

TwisTorr 74 FS PCB 24 Vdc Controller

**Model
X3510-64050**

**Manuale di istruzioni
Bedienungshandbuch
Notice de mode d'emploi
Manual de instrucciones
Manual de instruções
Bedrijfshandleiding
Instruktionsbog
Bruksanvisning**

**Instruksjon manual
Ohjekäsikirja
Felhasználói kézikönyv
Podrecznik instrukcji
Návod k použití
Návod na obsluhu
Priročnik za navodila
User Manual**

**87-901-056-01
07/2018**



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2018

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Manual Part Number

Publication Number: 87-901-056-01

Edition

Edition 07/2018

Printed in ITALY

Agilent Technologies Italia S.p.A.

Vacuum Products Division

Via F.lli Varian, 54

10040 Leini (TO)

ITALY

Warranty

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or

contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Trademarks

Windows and MS Windows are U.S. registered trademarks of Microsoft Corporation.

Safety Notices

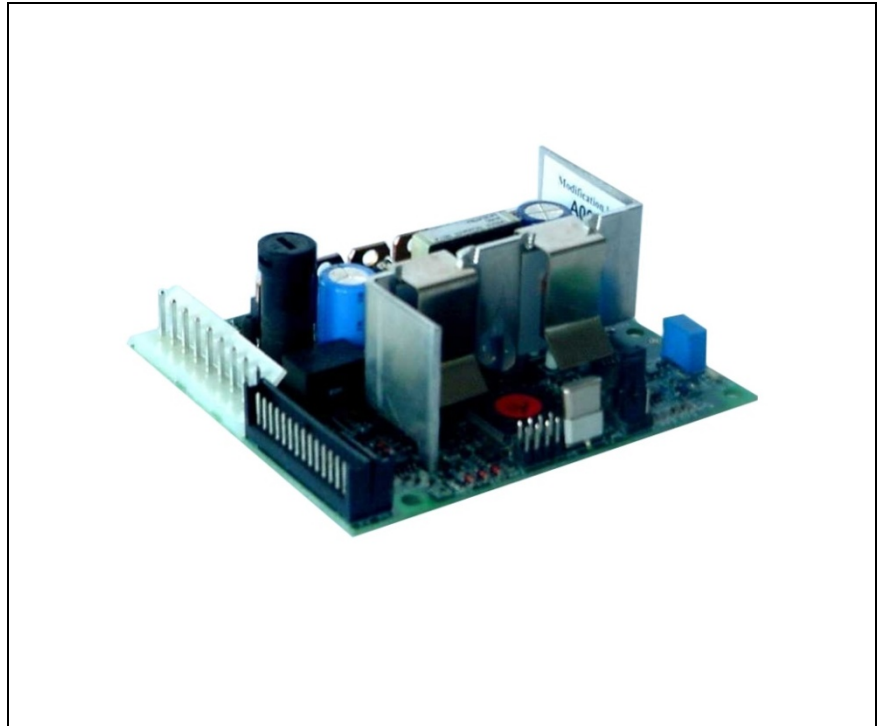
CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

TwisTorr 74 FS PCB 24 Vdc Controller



Contents

1	Istruzioni per l'uso	13
	Informazioni Generali	14
	Immagazzinamento	15
	Preparazione per l'installazione	16
	Installazione	18
	Uso	19
	Procedure di uso	19
	Manutenzione	20
	Smaltimento	21
2	Gebrauchsanleitung	23
	Allgemeines	24
	Lagerung	25
	Vor der Installation	26
	Installation	28
	Gebrauch	29
	Bedienung	29
	Wartung	30
	Entsorgung	31

Contents

3	Mode d'emploi	33
	Indications Générales	34
	Emmagasinage	35
	Préparation pour l'installation	36
	Installation	38
	Utilisation	39
	Procédures d'utilisation	39
	Entretien	40
	Mis au rebut	41
4	Manual de instrucciones	43
	Informaciones de carácter general	44
	Almacenamiento	45
	Preparación para la instalación	46
	Instalación	48
	Uso	49
	Procedimientos de uso	49
	Mantenimiento	50
	Eliminación	51
5	Manual de Instruções	53
	Informações gerais	54
	Armazenagem	55
	Preparação para a instalação	56

Instalação	58
Utilização	59
Procedimentos de uso	59
Manutenção	60
Eliminação	61

6 Bedrijfshandleiding 63

Algemene informatie	64
Opslag	65
Vorbereiding voor installatie	66
Installatie	68
Gebruik	69
Gebruiksprocedures	69
Onderhoud	70
Afvalverwerking	71

7 Istruktionsbog 73

Generel Information	74
Opbevaring	75
Forberedelser før installation	76
Installation	78
Anvendelse	79
Instruktion	79
Vedligeholdelse	80

Contents

	Bortskaffelse	81
8	Bruksanvisning	83
	Allmän information	84
	Förvaring	85
	Förberedelser för Installation	86
	Installation	88
	Användning	89
	Instruktioner för bruk	89
	Underhåll	90
	Bortskaffning	91
9	Instruksjon Manual	93
	Generell informasjon	94
	Lagring	95
	Forberede installasjonen	96
	Installasjon	98
	Bruk	99
	Instruksjoner for bruk	99
	Vedlikehold	100
	Eliminering	101
10	Ohjekäsikirja	103
	Yleisiä tietoja	104
	Varastointi	105

Valmistelut asennusta varten	106
Asennus	108
Käyttö	109
Käyttötoimenpiteet	109
Huolto	110
Hävittäminen	111
11 Felhasználói Kézikönyv	113
Általános információk	114
Tárolás	115
A telepítésre való előkészítés	116
Telepítés	118
Használat	119
Használati eljárások	119
Karbantartás	120
Megsemmisítés	121
12 Podrecznik Instrukcji	123
Informacje ogólne	124
Magazynowanie	125
Przygotowanie do instalacji	126
Instalacja	128
Użytkowanie	129
Procedury użytkowania	129

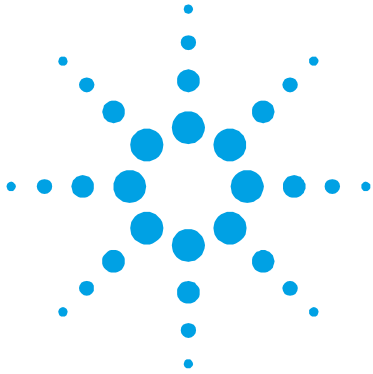
Contents

Konserwacja	130
Przetworstwo odpadów	131
13 Návod k Použití	133
Všeobecné informace	134
Uskladnění	135
Příprava k instalaci	136
Instalace	138
Použití	139
Používané procedury	139
Údržba	140
Likvidace	141
14 Návod na Obsluhu	143
Všeobecné informácie	144
Uskladňovanie	145
Príprava pre inštaláciu	146
Inštalácia	148
Použitie	149
Postup pri použití	149
Údržba	150
Likvidácia	151
15 Priročník za Navodila	153
Splošna navodila	154

Skladiščenje	155
Priprava za montažo	156
Montaža	158
Uporaba	159
Postopki za uporabo	159
Vzdrževanje	160
Odlaganje odpadkov	161
16 Instructions for Use	163
General Information	164
Storage	165
Preparation prior to Installation	166
Installation	168
Use	169
Operating Procedures	169
Maintenance	170
Disposal	171
17 Technical Information	173
General Description	174
TwisTorr 74 FS PCB 24 Vdc Controller Description	174
Controller Specifications	175
Controller Outline	178
Installation	179

Contents

Data Exchange Protocols	184
Example	194
Operation	199
Maintenance	201
Accessories and Spare Parts	202



1

Istruzioni per l'uso

Informazioni Generali	14
Immagazzinamento	15
Preparazione per l'installazione	16
Installazione	18
Uso	19
Procedure di uso	19
Accensione del Controller	19
Avvio della Pompa	19
Arresto della Pompa	19
Manutenzione	20
Smaltimento	21

Traduzione delle istruzioni originali



Informazioni Generali

Questa apparecchiatura è destinata ad uso professionale. L'utilizzatore deve leggere attentamente il presente manuale di istruzioni ed ogni altra informazione addizionale fornita dalla Agilent prima dell'utilizzo dell'apparecchiatura. La Agilent si ritiene sollevata da eventuali responsabilità dovute all'inosservanza totale o parziale delle istruzioni, ad uso improprio da parte di personale non addestrato, ad interventi non autorizzati o ad uso contrario alle normative nazionali specifiche. Il controller TwisTorr 74 FS PCB è un convertitore di frequenza, controllato da un microprocessore, realizzato con componenti a stato solido e con capacità di autodiagnostica e autoprotezione. Esso incorpora tutta la circuiteria necessaria per il funzionamento automatico delle pompe della serie TwisTorr 74 FS.

Il controller pilota le pompe della serie TwisTorr 74 FS (con un processo suddiviso in dieci passi) durante la fase di avvio controllando la tensione e la corrente in rapporto alla velocità raggiunta dalla pompa. Tramite connettori ausiliari sono disponibili i comandi per l'avvio e l'arresto della pompa da remoto, i segnali che indicano lo stato operativo della pompa, segnali di bloccaggio (per interruttori a pressione, interruttori di controllo del flusso dell'acqua, ecc.). Il controller deve essere alimentato con una tensione continua di 24 Vdc.

Nei paragrafi seguenti sono riportate tutte le informazioni necessarie a garantire la sicurezza dell'operatore durante l'utilizzo dell'apparecchiatura. Informazioni dettagliate sono fornite nell'appendice "Technical Information".

Questo manuale utilizza le seguenti convenzioni:

ATTENZIONE!

I messaggi di attenzione sono visualizzati prima di procedure che, se non osservate, potrebbero causare danni all'apparecchiatura.

AVVERTENZA!



I messaggi di avvertenza attirano l'attenzione dell'operatore su una procedura o una pratica specifica che, se non eseguita in modo corretto, potrebbe provocare gravi lesioni personali.

NOTA

Le note contengono informazioni importanti estrapolate dal testo.

Immagazzinamento

Durante il trasporto e l'immagazzinamento del controller devono essere soddisfatte le seguenti condizioni ambientali:

- temperatura: da -20 °C a +70 °C
- umidità relativa 0 – 95 % (non condensante)

Preparazione per l'installazione

Il controller viene fornito in un imballo protettivo speciale; se si presentano segni di danni, che potrebbero essersi verificati durante il trasporto, contattare l'ufficio vendite locale.

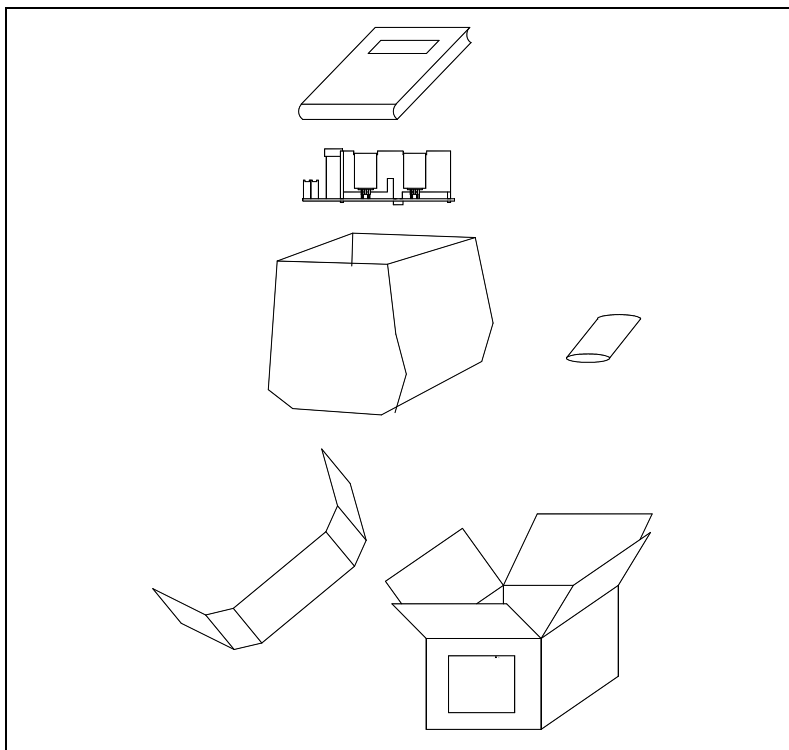


Figura 1 Imballo del controller

Durante l'operazione di disimballaggio, prestare particolare attenzione a non lasciar cadere il controller e a non sottoporlo ad urti. Non disperdere l'imballo nell'ambiente. Il materiale è completamente riciclabile e risponde alla direttiva CEE 85/399 per la tutela dell'ambiente. Prima di installare il controller occorre preparare alcuni cavi di interconnessione, quali ad esempio i cavi per le interconnessioni con l'alimentazione, quelli per le interconnessioni logiche e quello per il collegamento con la pompa. Nelle figure seguenti sono schematizzati i collegamenti da effettuare tra il controller e la pompa e tra il controller e l'alimentatore. Per ulteriori dettagli sulle suddette interconnessioni e le caratteristiche tecniche del trasformatore (disponibile come opzione) vedere l'appendice "Technical Information".

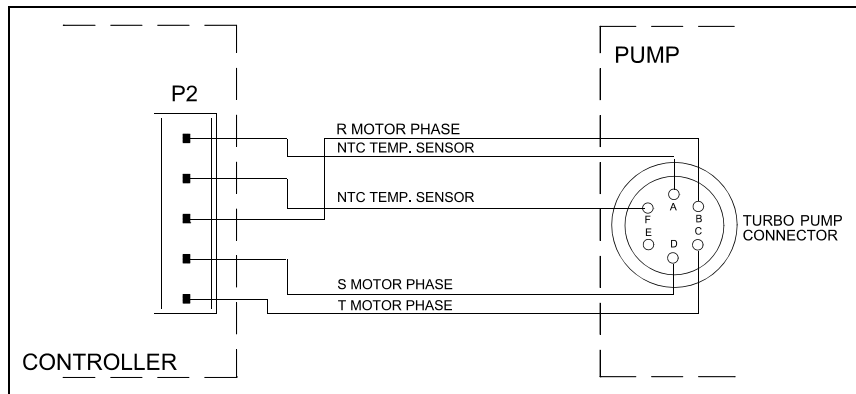


Figura 2 Cavo interconnessione tra Controller e pompa

Installazione

AVVERTENZA!



All'interno del controller si sviluppano tensioni che possono recare gravi danni. Prima di eseguire qualsiasi operazione di installazione o manutenzione del controller scollegarlo dalla alimentazione. All'interno del controller si sviluppano elevate temperature che possono recare gravi danni. Proteggere adeguatamente il controller nella sua installazione definitiva da contatti accidentali.

NOTA

Il controller installato nel sistema definitivo deve essere posizionato in modo tale che l'aria di raffreddamento possa circolare liberamente intorno all'apparato. Non installare e/o utilizzare il controller in ambienti esposti ad agenti atmosferici (pioggia, gelo, neve), polveri, gas aggressivi, in ambienti esplosivi o con elevato rischio di incendio.

Durante il funzionamento è necessario che siano rispettate le seguenti condizioni ambientali:

- temperatura: da 0 °C a +40 °C
- umidità relativa: 0 – 95 % (non condensante).

Uso

In questo paragrafo sono riportate le principali procedure operative. Per ulteriori dettagli e per procedure che coinvolgono collegamenti o particolari opzionali, fare riferimento al paragrafo "USE" dell'appendice "Technical Information". Prima di usare il controller effettuare tutti i collegamenti elettrici e pneumatici e fare riferimento al manuale della pompa collegata.

AVVERTENZA!

Fare riferimento al manuale relativo alla pompa in questione per utilizzarla in modo sicuro prima dell'avviamento.



Procedure di uso

Accensione del Controller

Per accendere il controller è sufficiente fornire la tensione di alimentazione al controller stesso.

Avvio della Pompa

Per avviare la pompa lasciare aperto il pin 12 del connettore J4, in modo tale che quando viene fornita la tensione di 24 Vdc al controller si ha l'avviamento della pompa.

Arresto della Pompa

Per arrestare la pompa occorre cortocircuitare i pin 12 e 15 del connettore J4.

Manutenzione

I controller della serie TwisTorr 74 FS non richiedono alcuna manutenzione. Qualsiasi intervento deve essere eseguito da personale autorizzato.

In caso di guasto è possibile usufruire del servizio di riparazione Agilent o del "Agilent advance exchange service", che permette di ottenere un controller ricondizionato in sostituzione di quello guasto.

AVVERTENZA!



Prima di effettuare qualsiasi intervento sul controller scollegare il cavo di alimentazione.

Qualora un controller dovesse essere rottamato, procedere alla sua eliminazione nel rispetto delle normative nazionali specifiche.

Smaltimento

Significato del logo "WEEE" presente sulle etichette. Il simbolo qui sotto riportato è applicato in ottemperanza alla direttiva CE denominata "WEEE". Questo simbolo (**valido solo per i paesi della Comunità Europea**) indica che il prodotto sul quale è applicato, NON deve essere smaltito insieme ai comuni rifiuti domestici o industriali, ma deve essere avviato ad un sistema di raccolta differenziata. Si invita pertanto l'utente finale a contattare il fornitore del dispositivo, sia esso la casa madre o un rivenditore, per avviare il processo di raccolta e smaltimento, dopo opportuna verifica dei termini e condizioni contrattuali di vendita.



Per maggiori informazioni riferirsi a:

<http://www.agilent.com/environment/product/index.shtml>

1 Istruzioni per l'uso

Smaltimento



2 Gebrauchsanleitung

Allgemeines	24
Lagerung	25
Vor der Installation	26
Installation	28
Gebrauch	29
Bedienung	29
Einschalten des Controllers	29
Pumpenstart	29
Pumpenstopp	29
Wartung	30
Entsorgung	31

Übersetzung der Originalanleitungen

Allgemeines

Dieser Apparat ist für Fachbetriebe bestimmt. Vor Gebrauch sollte der Benutzer dieses Handbuch sowie alle weiteren mitgelieferten Zusatzdokumentationen genau lesen. Bei Nichtbeachtung - auch teilweise - der enthaltenen Hinweise, unsachgemäßem Gebrauch durch ungeschultes Personal, nicht autorisierten Eingriffen und Missachtung der einheimischen, hier zur Geltung kommenden Bestimmungen übernimmt die Firma Agilent keinerlei Haftung. Der Controller der Serie TwisTorr 74 FS PCB ist ein mikroprozessorgesteuerter Frequenzwandler. Er ist aus Festkörperbauteilen gefertigt und verfügt über ein Selbstdiagnose- und ein Selbstschutzsystem. Er enthält alle für den automatischen Betrieb der Pumpenserie TwisTorr 74 FS erforderlichen Schaltungen. Der Controller steuert die Pumpen der Serie TwisTorr 74 FS (durch einen 10-Schritte-Prozess) in der Startphase, indem er die Spannung und die Stromstärke im Verhältnis zur Pumpengeschwindigkeit kontrolliert. Mittels Hilfsverbinder sind die Fernsteuerungen für Pumpenstart und -stopp, die Signale für die Betriebsanzeige der Pumpe und Sperrsignale (für Druckschalter, Wasserstrom-Kontrollschalter usw.) verfügbar. Der Controller soll mit einer Gleichstromspannung von 24 V versorgt werden. Für die Funktionskontrolle und die Visualisierung des Controllerzustandes ist ein entsprechendes Gerät (Hand Held Terminal) erhältlich; es besteht aus einer Steuertastatur und einem Display für die Visualisierung.

In den folgenden Abschnitten sind alle erforderlichen Informationen für die Sicherheit des Bedieners bei der Anwendung des Geräts aufgeführt. Detaillierte technische Informationen sind im Anhang "Technical Information" enthalten.

In dieser Gebrauchsanleitung werden Sicherheitshinweise folgendermaßen hervorgehoben:

VORSICHT! Die Vorsichtshinweise werden vor Vorgängen angegeben, die bei Nichtbeachtung Schäden an der Anlage verursachen könnten.

WARNUNG! Die Warnhinweise lenken die Aufmerksamkeit des Bedieners auf eine bestimmte Prozedur oder Praktik, die bei unkorrekter Ausführung schwere Verletzungen hervorrufen können.



HINWEIS Die Hinweise enthalten wichtige Informationen, die aus dem Text hervorgehoben werden.

Lagerung

Beim Transport und bei der Lagerung des Controllers müssen folgende klimatische Verhältnisse eingehalten werden:

- Temperatur: von -20 °C bis +70 °C
- Relative Luftfeuchtigkeit: 0 – 95 % (nicht kondensierend)

Vor der Installation

Der Controller wird mit einer speziellen Schutzverpackung geliefert. Eventuelle Transportschäden müssen der zuständigen örtlichen Verkaufsstelle gemeldet werden.

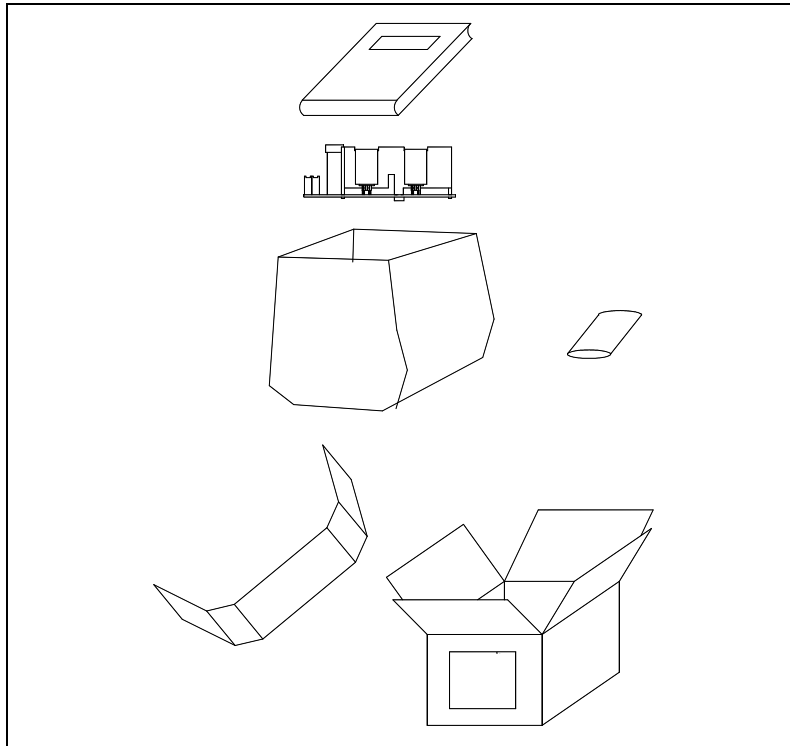


Abbildung 1 Verpackung des Controllers

Der Controller sollte vorsichtig ausgepackt werden, damit er nicht fällt oder Stößen ausgesetzt wird. Das Verpackungsmaterial muss korrekt entsorgt werden. Es ist vollständig recyclebar und entspricht der EG-Richtlinie 85/399 für Umweltschutz. Vor der Installation des Controllers müssen einige Verbindungskabel vorbereitet werden; wie zum Beispiel die Kabel für die Verbindung mit der Versorgung, die Kabel für die logischen Verbindungen und das Kabel für den Anschluss mit der Pumpe. Die nachstehenden Abbildungen schematisieren die Verbindungen, die zwischen dem Controller und der Pumpe sowie zwischen dem Controller und dem Netzgerät auszuführen sind. Weitere Einzelheiten zu den oben genannten Verbindungen und zu den Merkmalen des Versorgungstransformators (erhältlich als Optional), sind im Anhang "Technical Information" aufgeführt.

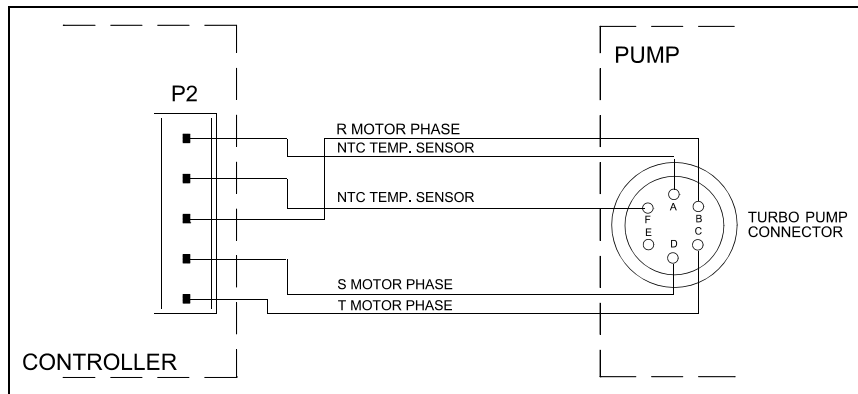


Abbildung 2 Verbindungskabel zwischen Controller und Pumpe

Installation

WARNUNG!



Im Innern des Controllers bilden sich Spannungen, die zu schweren Schäden führen können. Vor jedem Installations- oder Wartungseingriff den Controller von der Stromversorgung trennen. Im Innern des Controllers bilden sich erhöhte Temperaturen, die schwere Schäden verursachen können. Schützen Sie den Controller bei seiner definitiven Installation vor zufälligen Berührungen.

HINWEIS

Der im definitiven System installierte Controller muss so positioniert sein, dass die Kühlluft frei um das Gerät zirkulieren kann. Der Controller darf nicht in Umgebungen installiert u/o benutzt werden, die Witterungseinflüssen (Regen, Frost, Schnee), Staub und aggressiven Gasen ausgesetzt sind und in denen Explosions- und erhöhte Brandgefahr besteht.

Beim Betrieb müssen folgende Umgebungsbedingungen eingehalten werden:

- Temperatur: von 0 °C bis +40 °C;
- Relative Luftfeuchtigkeit: 0 – 95 % (nicht kondensierend).

Gebrauch

In diesem Kapitel sind die wichtigsten Betriebsvorgänge aufgeführt. Für weitere Hinweise bezüglich Anschlüsse und Montage des bestellbaren Zubehörs siehe Kapitel "USE" im Anhang zu "Technical Information". Vor Benutzung des Controllers sämtliche elektrischen und pneumatischen Anschlüsse ausführen und die Betriebsanleitung der angeschlossenen Pumpe durchlesen.

WARNUNG!



Bevor Sie die fragliche Pumpe in Betrieb setzen, beachten Sie zu Ihrer Sicherheit bitte die Hinweise im entsprechenden Handbuch.

Bedienung

Einschalten des Controllers

Zum Einschalten des Controllers genügt es, das Netzkabel an die Steckdose anzuschließen.

Pumpenstart

Zum Starten der Pumpe ist PIN 12 des Verbinders J4 offen zu lassen, so dass bei Zuschaltung der 24V-Gleichspannung am Controller der Pumpenstart erfolgt.

Pumpenstopp

Zum Stoppen der Pumpe sind die PIN 12 und 15 des Verbinders J4 kurzzuschließen.

Wartung

Die Controller der Serie TwisTorr 74 FS sind wartungsfrei. Eventuell erforderliche Eingriffe müssen von dazu befugtem Fachpersonal ausgeführt werden.

Bei einem Defekt kann der Agilent Reparaturdienst bzw. der "Agilent advanced exchange service" in Anspruch genommen werden, der für die Erneuerung des defekten Controllers sorgt.

WARNUNG!

Vor jedem Eingriff am Controller muss der Netzstecker gezogen werden.



Eine eventuelle Verschrottung hat unter Einhaltung der einschlägigen landesüblichen Vorschriften zu erfolgen.

Entsorgung

Bedeutung des "WEEE" Logos auf den Etiketten. Das folgende Symbol ist in Übereinstimmung mit der EU-Richtlinie WEEE (Waste Electrical and Electronic Equipment) angebracht. Dieses Symbol (**nur in den EU-Ländern gültig**) zeigt an, dass das betreffende Produkt nicht zusammen mit Haushaltsmüll entsorgt werden darf sondern einem speziellen Sammelsystem zugeführt werden muss. Der Endabnehmer sollte daher den Lieferanten des Geräts - d.h. die Muttergesellschaft oder den Wiederverkäufer - kontaktieren, um den Entsorgungsprozess zu starten, nachdem er die Verkaufsbedingungen geprüft hat.



Weitere Informationen finden Sie unter:

<http://www.agilent.com/environment/product/index.shtml>

2 **Gebrauchsanleitung**

Entsorgung



3

Mode d'emploi

Indications Générales	34
Emmagasinage	35
Préparation pour l'installation	36
Installation	38
Utilisation	39
Procédures d'utilisation	39
Allumage du contrôleur	39
Mise en marche de la Pompe	39
Arrêt de la Pompe	39
Entretien	40
Mis au rebut	41

Traduction de la mode d'emploi originale



Indications Générales

Cet appareillage a été conçu en vue d'une utilisation professionnelle. Il est conseillé à l'utilisateur de lire attentivement cette notice d'instructions ainsi que toute autre indication supplémentaire fournie par Agilent, avant l'utilisation de l'appareil. Agilent décline par conséquent toute responsabilité en cas d'inobservation totale ou partielle des instructions données, d'utilisation incorrecte de la part d'un personnel non formé, d'opérations non autorisées ou d'un emploi contraire aux réglementations nationales spécifiques. Le contrôleur TwisTorr 74 FS PCB est un convertisseur de fréquence, contrôlé par un microprocesseur, réalisé avec des éléments à l'état solide et ayant des capacités d'autodiagnostic et d'autoprotection. Il incorpore l'ensemble de circuits nécessaire au fonctionnement automatique des pompes de la série TwisTorr 74 FS. Le contrôleur pilote les pompes de la série TwisTorr 74 FS (selon un processus subdivisé en dix pas) pendant la phase de mise en marche, en contrôlant la tension et le courant par rapport à la vitesse atteinte par la pompe. Des connecteurs auxiliaires permettent de disposer des commandes de mise en marche et d'arrêt de la pompe à distance, des signaux indiquant l'état opérationnel de la pompe, des signaux de blocage (pour interrupteurs à pression, interrupteurs de contrôle du flux de l'eau etc.) Le contrôleur doit être alimenté avec une tension continue de 24 Vdc.

Les paragraphes suivants donnent toutes les indications nécessaires à garantir la sécurité de l'opérateur pendant l'utilisation de l'appareillage.

Des renseignements plus détaillés se trouvent dans "Technical Information".

Cette notice utilise les signes conventionnels suivants:

ATTENTION!

Les messages d'attention apparaissent avant certaines procédures qui, si elles ne sont pas observées, pourraient endommager sérieusement l'appareillage.

AVERTISSEMENT!



Les messages d'avertissement attirent l'attention de l'opérateur sur une procédure ou une manoeuvre spéciale qui, si elle n'est pas effectuée correctement, risque de provoquer de graves lésions.

NOTE

Les notes contiennent des renseignements importants, isolés du texte.

Emmagasinage

Pendant le transport et l'emmagasinage des contrôleurs, il faudra veiller à respecter les conditions environnementales suivantes:

- température: de -20 °C à + 70 °C
- humidité relative: 0 – 95 % (non condensante).

Préparation pour l'installation

Le contrôleur est fourni dans un emballage de protection spécial; si l'on constate des marques de dommages pouvant s'être produits pendant le transport, contacter aussitôt le bureau de vente local.

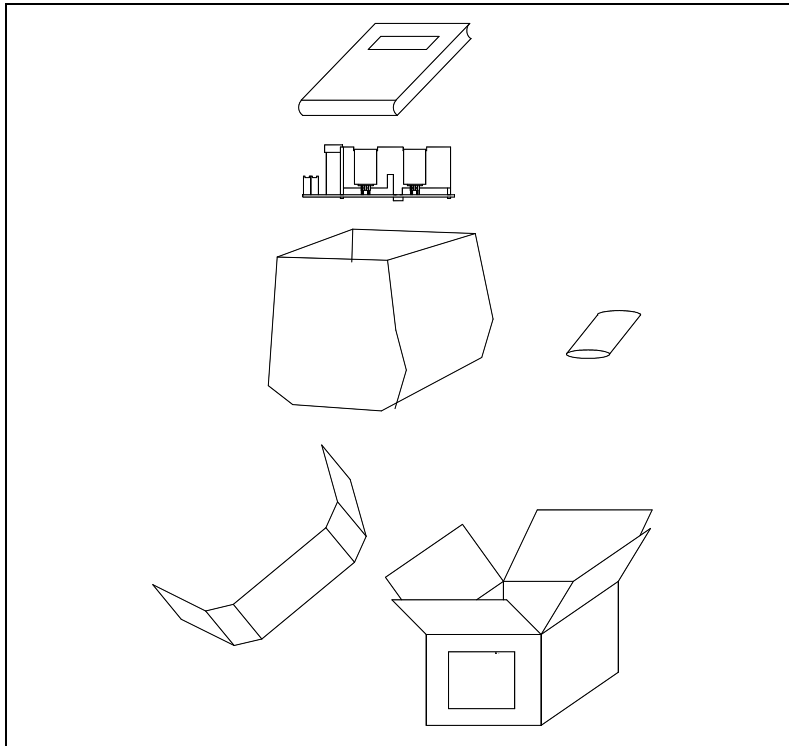


Figure 1 Emballage du contrôleur

Pendant l'opération d'ouverture de l'emballage, veiller tout particulièrement à ne pas laisser tomber le contrôleur et à ne lui faire subir aucun choc. Ne pas jeter l'emballage dans la nature. Le matériel est entièrement recyclable et il est conforme aux directives CEE 85/399 en matière de protection de l'environnement. Avant d'installer le contrôleur, il est nécessaire de préparer quelques câbles d'interconnexion, par exemple les câbles pour les interconnexions avec l'alimentation, ceux pour les interconnexions logiques et celui pour le branchement avec la pompe. Les schémas qui suivent indiquent les connexions à effectuer entre le contrôleur et la pompe et entre le contrôleur et l'alimentateur. Pour tous autres détails concernant les interconnexions susdites et les caractéristiques du transformateur d'alimentation (disponible en option), se reporter à l'appendice "Technical Information".

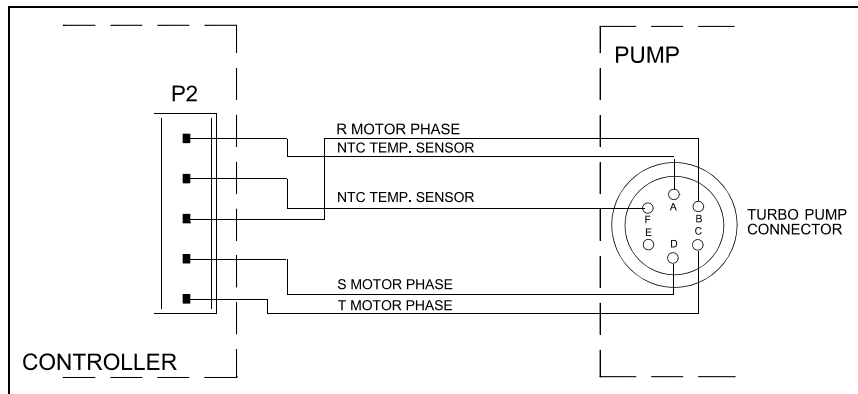


Figure 2 Câble d'interconnexion entre le contrôleur et la pompe

Installation

AVERTISSEMENT!



A l'intérieur du contrôleur se développent des tensions qui peuvent causer de graves dommages. Avant d'effectuer toute opération d'installation ou d'entretien du contrôleur, le débrancher de la prise d'alimentation. A l'intérieur du contrôleur se développent des températures élevées qui peuvent causer de graves dommages. Protéger de façon appropriée le contrôleur, dans son installation définitive, contre tous contacts accidentels.

NOTE

Le contrôleur installé dans le système définitif doit être positionné de façon à ce que l'air de refroidissement puisse circuler librement autour de l'appareil. Ne pas installer et/ou utiliser le contrôleur dans des milieux exposés à des agents atmosphériques (pluie, gel, neige), à des poussières, à des gaz de combat ainsi que dans des milieux explosifs ou à risque élevé d'incendie.

Pendant le fonctionnement, il est nécessaire de respecter les conditions environnementales suivantes:

- température: de 0 °C à + 40 °C
- humidité relative: de 0 % à 95 % (non condensante).

Utilisation

Ce paragraphe indique les principales procédures opérationnelles. Pour tous autres détails et pour les procédures concernant des connexions ou des éléments en option, se reporter au paragraphe "USE" de l'appendice "Technical Information". Avant d'utiliser le contrôleur, effectuer toutes les connexions électriques et pneumatiques et se référer à la notice de la pompe connectée.

AVERTISSEMENT!



Veillez-vous référer au manuel de la pompe spécifique pour l'utiliser en toute sécurité avant le démarrage.

Procédures d'utilisation

Allumage du contrôleur

Pour allumer le contrôleur, il suffit d'introduire le câble d'alimentation dans la prise du réseau.

Mise en marche de la Pompe

Pour mettre la pompe en marche il faut laisser ouvert le contact sur le pin 12 du connecteur J4 de façon que lorsque le contrôleur reçoit la tension de 24 Vdc, la pompe se met en marche .

Arrêt de la Pompe

Pour arrêter la pompe, il est nécessaire de court-circuiter les pins 12 et 15 du connecteur J4.

Entretien

Les contrôleurs de la série TwisTorr 74 FS n'exigent aucun entretien. Toute opération doit être effectuée par un personnel agréé. En cas de panne, il est possible de s'adresser au Service de réparation Agilent ou bien au "Agilent advance exchange service" qui permet d'obtenir un contrôleur régénéré à la place du contrôleur détraqué.

AVERTISSEMENT!



Avant d'effectuer toute opération sur le contrôleur, débrancher le câble d'alimentation.

Si un contrôleur doit être mis au rebut, procéder à son élimination dans le respect des normes nationales en vigueur.

Mis au rebut

Signification du logo "WEEE" figurant sur les étiquettes. Le symbole ci-dessous est appliqué conformément à la directive CE nommée "WEEE". Ce symbole (**uniquement valide pour les pays de la Communauté européenne**) indique que le produit sur lequel il est appliqué NE doit PAS être mis au rebut avec les ordures ménagères ou les déchets industriels ordinaires, mais passer par un système de collecte sélective. Après avoir vérifié les termes et conditions du contrat de vente, l'utilisateur final est donc prié de contacter le fournisseur du dispositif, maison mère ou revendeur, pour mettre en œuvre le processus de collecte et mise au rebut.

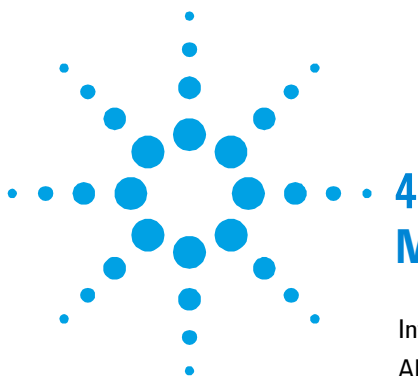


Pour en savoir plus, consulter :

<http://www.agilent.com/environment/product/index.shtml>

3 Mode d'emploi

Mis au rebut



4 Manual de instrucciones

Informaciones de carácter general	45
Almacenamiento	46
Preparación para la instalación	48
Instalación	49
Uso	49
Procedimientos de uso	49
Encendido del controlador	49
Puesta en marcha de la Bomba	49
Parada de la Bomba	49
Mantenimiento	50
Eliminación	51

Traducción de las instrucciones originales

Informaciones de carácter general

Este equipo se ha concebido para un uso profesional. El usuario deberá leer atentamente el presente manual de instrucciones y cualquier otra información suplementaria facilitada por Agilent antes de utilizar el equipo. Agilent se considera libre de cualquier responsabilidad debida al incumplimiento total o parcial de las instrucciones, al uso poco apropiado por parte de personal sin formación, a las operaciones no autorizadas o al uso que no cumpla con las normas nacionales específicas. El controlador TwisTorr 74 FS PCB es un convertidor de frecuencia, controlado por un microprocesador, realizado con componentes en estado sólido y con capacidad de autodiagnos y autoprotección. Éste incorpora todos los circuitos necesarios para el funcionamiento automático de las bombas de la serie TwisTorr 74 FS.

El controlador pilotea las bombas de la serie TwisTorr 74 FS (con un proceso dividido en diez pasos) durante la fase de puesta en marcha, controlando la tensión y la corriente en relación a la velocidad alcanzada por la bomba. Mediante conectores auxiliares están disponibles los mandos para la puesta en marcha y la parada de la bomba desde remoto, las señales que indican el estado operativo de la bomba y las señales de bloqueo (para interruptores a presión, interruptores de control del flujo del agua, etc.). El controlador debe ser alimentado con una tensión continua de 24 Vdc.

En los apartados siguientes se facilita toda la información necesaria para garantizar la seguridad del operador durante el uso del equipo. Una información más detallada se facilita en el Suplemento "Technical Information".

Este manual utiliza los siguientes símbolos convencionales:

¡ATENCIÓN!

Los mensajes de atención se visualizan antes de procedimientos que, al no respetarse, podrían provocar daños al equipo.

¡ADVERTENCIA!



Los mensajes de advertencia atraen la atención del operador sobre un procedimiento o una operación específica que, al no realizarse correctamente, podría provocar graves lesiones personales.

NOTA

Las notas contienen informaciones importantes extrapoladas del texto.

Almacenamiento

Durante el transporte y el almacenamiento del controlador se deberá cumplir con las condiciones ambientales siguientes:

- temperatura: de -20 °C a +70 °C
- humedad relativa: 0 – 95 % (no condensadora)

Preparación para la instalación

El controlador se suministra en un embalaje de protección especial; si se observan señales de daños, que podrían haberse producido durante el transporte, ponerse en contacto con la oficina de venta más cercana.

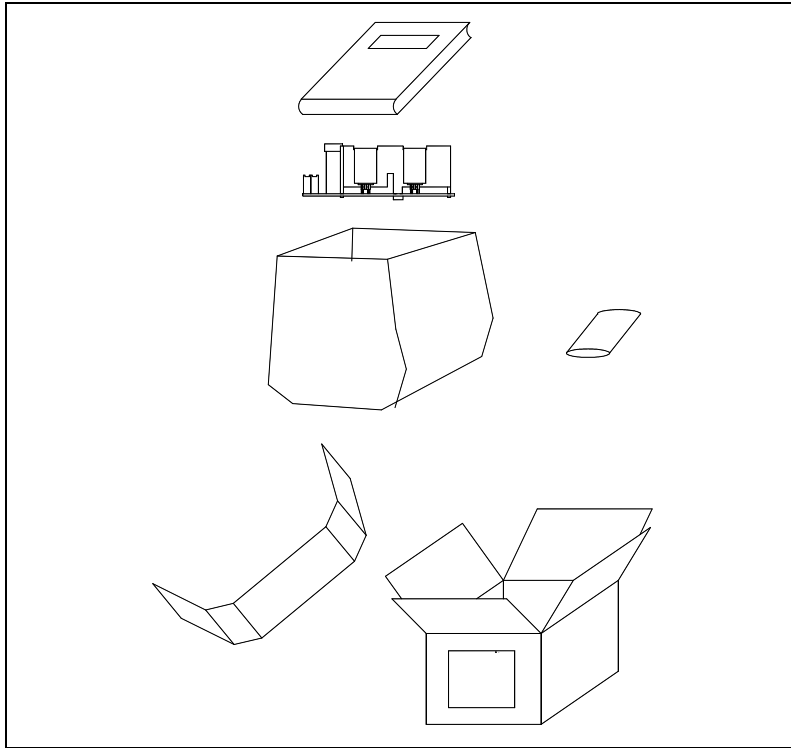


Figura 1 Embalaje del controlador

Durante la operación de desembalaje, prestar una atención especial a no dejar caer el controlador y evitarle golpes. No dispersar el embalaje en el medio ambiente. El material es totalmente reciclable y cumple con la directiva CEE 85/399 para la preservación del medio ambiente. Antes de instalar el controlador habrá que preparar algunos cables de interconexión, como por ejemplo los cables para las interconexiones con la alimentación, los de las interconexiones lógicas y el de acoplamiento con la bomba. En las figuras siguientes están esquematizados los acoplamientos que hay que realizar entre el controlador y la bomba y entre el controlador y el alimentador. Para más detalles sobre las interconexiones indicadas anteriormente y sobre las características técnicas del transformador (disponible como opción) véase el suplemento "Technical Information".

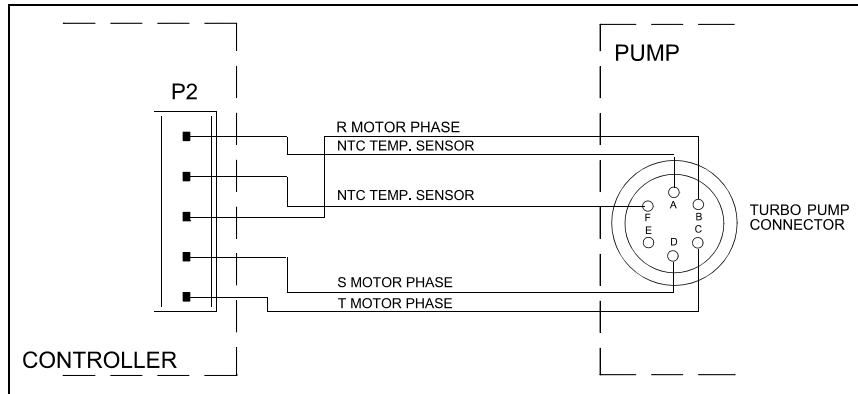


Figura 2 Cable de interconexión entre controlador y bomba

Instalación

¡ADVERTENCIA!



Dentro del controlador se desarrollan tensiones que pueden causar graves daños. Antes de efectuar cualquier operación de instalación o mantenimiento del controlador, desconectarlo de la alimentación. Dentro del controlador se desarrollan altas temperaturas que pueden causar graves daños. Proteger el controlador adecuadamente en su instalación definitiva contra contactos accidentales.

NOTA

El controlador instalado en el sistema definitivo ha de colocarse de manera que el aire de refrigeración pueda circular libremente alrededor del aparato. No instalar ni utilizar el controlador en ambientes expuestos a agentes atmosféricos (lluvia, hielo y nieve), polvos, gases agresivos, en ambientes explosivos o con alto riesgo de incendio.

Durante el funcionamiento es necesario que se respeten las siguientes condiciones ambientales:

- temperatura: de 0 °C a + 40 °C
- humedad relativa: 0 – 95 % (no condensadora).

Uso

En este apartado se indican los procedimientos operativos principales. Para más detalles y para procedimientos que impliquen conexiones o elementos opcionales, les remitimos al apartado “USE” del anexo “Technical Information”. Antes de usar el controlador efectuar todas las conexiones eléctricas y neumáticas y consultar el manual de la bomba conectada.

¡ADVERTENCIA!



Antes de arrancar la bomba consulte el correspondiente manual de instrucciones de la misma para operar de forma segura.

Procedimientos de uso

Encendido del controlador

Para encender el controlador es suficiente aprovisionar el controlador mismo con tensión de alimentación.

Puesta en marcha de la Bomba

Para poner en marcha la bomba dejar abierto el pin 12 del conector J4, de manera que al conectarse al controlador la tensión de 24 Vdc se verifique el arranque de la bomba.

Parada de la Bomba

Para detener la bomba es necesario cortocircuitar los pins 12 y 15 del conector J4.

Mantenimiento

Los controladores de la serie TwisTorr 74 FS no necesitan ningún mantenimiento. Toda operación ha de ser efectuada por personal autorizado. En caso de avería es posible utilizar el servicio de reparación Agilent o el “Agilent advance exchange service”, que permite obtener un controlador regenerado en vez del averiado.

¡ADVERTENCIA!



Antes de efectuar cualquier operación en el controlador se debe desenchufar el cable de alimentación.

En caso de que un controlador se tenga que desguazar, efectuar su eliminación respetando las normas nacionales específicas.

Eliminación

Significado del logotipo "WEEE" presente en las etiquetas. El símbolo que se indica a continuación, es aplicado en observancia de la directiva CE denominada "WEEE". Este símbolo (**válido sólo para los países miembros de la Comunidad Europea**) indica que el producto sobre el cual ha sido aplicado, NO debe ser eliminado junto con los residuos comunes sean éstos domésticos o industriales, y que, por el contrario, deberá ser sometido a un procedimiento de recogida diferenciada. Por lo tanto, se invita al usuario final, a ponerse en contacto con el proveedor del dispositivo, tanto si éste es la casa fabricante o un distribuidor, para poder proveer a la recogida y eliminación del producto, después de haber efectuado una verificación de los términos y condiciones contractuales de venta.



Para obtener más información, consulte:

<http://www.agilent.com/environment/product/index.shtml>

4 Manual de instrucciones

Eliminación



5

Manual de Instruções

Informações gerais	54
Armazenagem	55
Preparação para a instalação	56
Instalação	58
Utilização	59
Procedimentos de uso	59
Acendimento do Controller	59
Activação da bomba	59
Paragem da bomba	59
Manutenção	60
Eliminação	61

Tradução das instruções originais



Informações gerais

Esta aparelhagem destinase ao uso profissional. O utilizador deve ler atentamente o presente manual de instruções e todas as informações adicionais fornecidas pela Agilent antes de utilizar a aparelhagem. A Agilent não se responsabiliza pela inobservância total ou parcial das instruções, pelo uso indevido por parte de pessoas não treinadas, por operações não autorizadas ou pelo uso contrário às normas nacionais específicas. O controller TwisTorr 74 FS PCB é um conversor de frequência, controlado por um microprocessador, realizado com componentes em estado sólido e com capacidade de autodiagnóstico e autoprotecção. Incorpora todos os circuitos necessários para o funcionamento automático das bombas da série TwisTorr 74 FS. O controller comanda as bombas da série TwisTorr 74 FS (com um processo subdividido em dez passos) durante a fase de activação, controlando a tensão e a corrente em relação à velocidade atingida pela bomba. Através de conectores auxiliares, estão disponíveis os comandos para a activação e a paragem da bomba por controlo remoto, os sinais que indicam o estado operativo da bomba, os sinais de bloqueio (para interruptores de pressão, interruptores de controlo do fluxo de água, etc.). O controller deve ser alimentado por tensão contínua de 24 Vdc.

Nos parágrafos seguintes estão descritas todas as informações necessárias para garantir a segurança do operador durante o uso da aparelhagem. Informações pormenorizadas são fornecidas no apêndice "Technical Information".

Este manual utiliza as seguintes convenções:

CUIDADO!

As mensagens de cuidado são visualizadas antes de procedimentos que, se não observados, podem causar danos à aparelhagem.

ATENÇÃO!



As mensagens de atenção chamam a atenção do operador para um procedimento ou uma prática específica que, se não efectuada correctamente, pode provocar graves lesões pessoais.

NOTA

As notas contêm informações importantes destacadas do texto.

Armazenagem

Durante o transporte e a armazenagem do controller, devem ser satisfeitas as seguintes condições ambientais:

- temperatura: de -20 °C a + 70 °C
- humidade relativa: 0 – 95 % (não condensadora)

Preparação para a instalação

O controller é fornecido numa embalagem protectora especial; se se apresentarem sinais de danos, que poderiam verificarse durante o transporte, entrar em contacto com o escritório de vendas local.

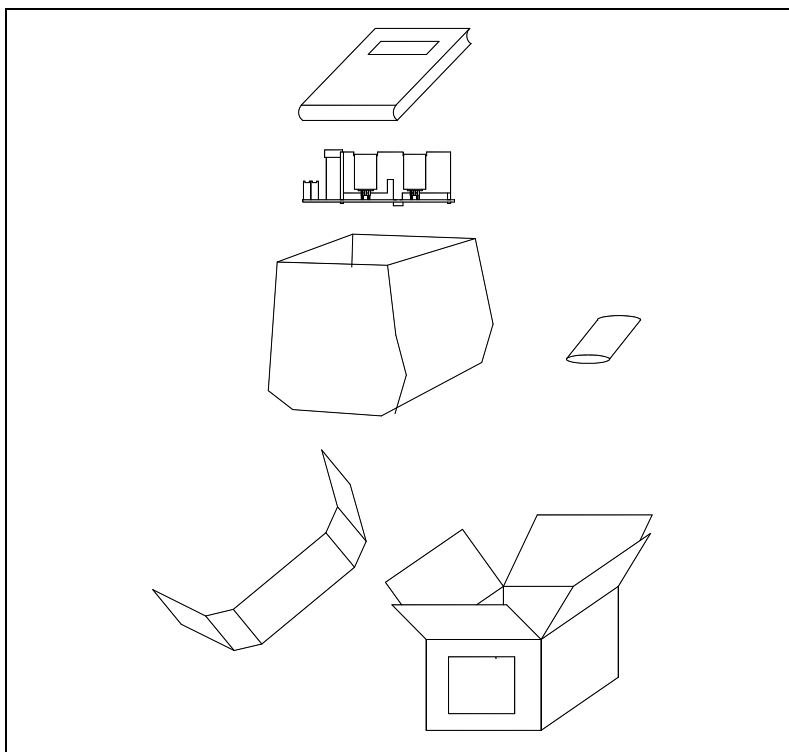


Figura 1 Embalagem do controller

Durante a remoção da embalagem, tomar muito cuidado para não deixar cair o controller e para não submetê-lo a choques. Não depositar a embalagem no meio ambiente. O material é completamente reciclável e responde à norma CEE 85/399 para a protecção do meio ambiente. Antes de instalar o controller, é necessário preparar alguns cabos de interconexão, como por exemplo os cabos para as interconexões com a alimentação, para as interconexões lógicas e para a ligação com a bomba. Nas figuras a seguir estão esquematizadas as ligações a efectuar entre o controller e a bomba, e entre o controller e o alimentador. Para maiores detalhes sobre as interconexões citadas acima ver o apêndice "Technical Information".

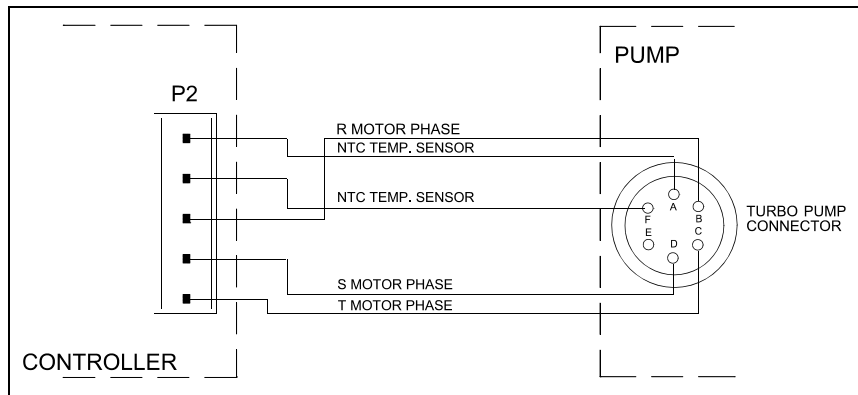


Figura 2 Cabo de Interconexão entre o Controller e a Bomba

Instalação

ATENÇÃO!



No interior do controller desenvolvem-se altas tensões que podem provocar graves danos ou a morte. Antes de efectuar qualquer operação de instalação ou manutenção do controller, desligar a tomada de alimentação. No interior do controller desenvolvem-se temperaturas elevadas que podem provocar graves danos. Proteger adequadamente o controller contra contactos acidentais durante a sua instalação definitiva.

NOTA

O controller instalado no sistema definitivo deve ser posicionado de modo que o ar de refrigeração possa circular livremente ao redor da aparelhagem. Não instalar e/ou utilizar o controller em ambientes expostos a agentes atmosféricos (chuva, gelo, neve), poeiras, gases agressivos ou em ambientes explosivos ou com elevado risco de incêndio.

Durante o funcionamento é necessário que sejam respeitadas as seguintes condições ambientais:

- temperatura: de 0 °C a + 40 °C
- humidade relativa: 0 – 95 % (não condensadora)

Utilização

Neste parágrafo são descritos os principais procedimentos operativos. Para maiores detalhes e para procedimentos que envolvam ligações ou peças opcionais, consultar o parágrafo "USE" do apêndice "Technical Information". Antes de usar o controller, efectuar todas as ligações eléctricas e pneumáticas e consultar o manual da bomba ligada.

ATENÇÃO!

Consulte o respectivo manual da bomba para operá-la de forma segura antes de começar.



Procedimentos de uso

Acendimento do Controller

Para ligar o controller é suficiente fornecer a tensão de alimentação ao controller.

Activação da bomba

Para activar a bomba é necessário deixar o pin 12 do conector J4 aberto. Deste modo , quando o controller for alimentado com uma tensão de 24 Vdc, a bomba activa-se.

Paragem da bomba

Para parar a bomba é necessário pôr os pins 12 e 15 do conector J4 em curto-circuito.

Manutenção

O controller da série TwisTorr 74 FS não requer qualquer manutenção. Todas as operações devem ser efectuadas por pessoal autorizado. Em caso de defeito é possível utilizar o serviço de reparação Agilent ou o "Agilent advance exchange service", que permite obter um controller regenerado que substitua o controller com defeito.

ATENÇÃO!



Antes de efectuar qualquer operação no controller, desligar o cabo de alimentação.

Caso um controller deva ser destruído, proceder à sua eliminação respeitando as normas nacionais específicas.

Eliminação

Significado do logótipo "WEEE" presente nos rótulos. O símbolo abaixo indicado é aplicado de acordo com a directiva CE denominada "WEEE". Este **símbolo (válido apenas para os países da Comunidade Europeia)** indica que o produto no qual está aplicado NÃO deve ser eliminado juntamente com os resíduos domésticos ou industriais comuns, mas deve ser dirigido a um sistema de recolha diferenciada. Portanto, convidamos o utilizador final a contactar o fornecedor do dispositivo, seja este o fabricante ou um revendedor, para encaminhar o processo de recolha e eliminação, após a oportuna verificação dos termos e condições do contrato de venda.

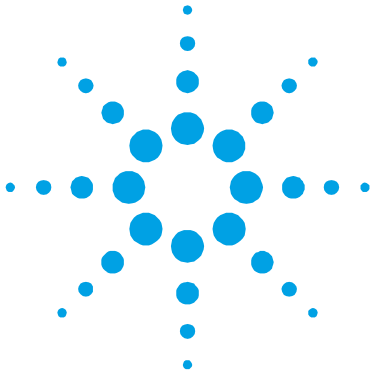


Para mais informações consulte:

<http://www.agilent.com/environment/product/index.shtml>

5 Manual de Instruções

Eliminação



6

Bedrijfshandleiding

Algemene informatie	64
Opslag	65
Vorbereiding voor installatie	66
Installatie	68
Gebruik	69
Gebruiksprocedures	69
Inschakelen van de controller	69
Starten van de pomp	69
Stoppen van de pomp	69
Onderhoud	70
Afvalverwerking	71

Vertaling van de originele instructies



Algemene informatie

Deze apparatuur is bestemd voor beroepsmatig gebruik. De gebruiker wordt verzocht aandachtig deze handleiding en alle overige door Agilent verstrekte informatie door te lezen alvorens het apparaat in gebruik te nemen. Agilent acht zich niet aansprakelijk voor de gevolgen van het niet of gedeeltelijk in acht nemen van de aanwijzingen, onoordeelkundig gebruik door niet hiervoor opgeleid personeel, reparaties waarvoor geen toestemming is verkregen of gebruik in strijd met de specifieke nationale wetgeving. De controller TwisTorr 74 FS PCB is een frequentieomzetter die gestuurd wordt door een microprocessor, is gemaakt van halfgeleiderelementen en is in staat om zelfdiagnose en zelfbescherming uit te voeren. De controller is van circuits voorzien die noodzakelijk zijn voor de automatische werking van de pompen van de serie TwisTorr 74 FS. De controller stuurt de pompen van de serie TwisTorr 74 FS (met een proces bestaande uit tien stappen) tijdens de startfase, en controleert hierbij de spanning en de stroom in verhouding tot de door de pomp bereikte snelheid. Via hulpconnectors zijn de sturingen voor het op afstand starten en stoppen van de pomp beschikbaar, de signalen die de bedrijfstoestand van de pomp aangeven, blokkeersignalen (voor drukschakelaars, regelschakelaars van de waterstroom, enz.). De controller moet met een gelijkspanning van 24 Vdc worden gevoed.

In de volgende paragrafen is alle informatie vermeld om de veiligheid van de operator tijdens het gebruik van de apparatuur te verzekeren. Gedetailleerde informatie is te vinden in de bijlage "Technical information".

Deze handleiding hanteert de volgende symbolen:

VOORZICHTIG!

Bij dit symbool staat tekst met procedures die, indien niet opgevolgd, schade aan apparatuur kunnen veroorzaken.

WAARSCHUWING!



Bij dit symbool staat tekst die de aandacht van de operator vestigt op een speciale procedure of methode die, indien niet correct uitgevoerd, ernstig lichamelijk letsel kan veroorzaken.

OPMERKING

De opmerkingen bevatten belangrijke informatie die uit de tekst is gelicht.

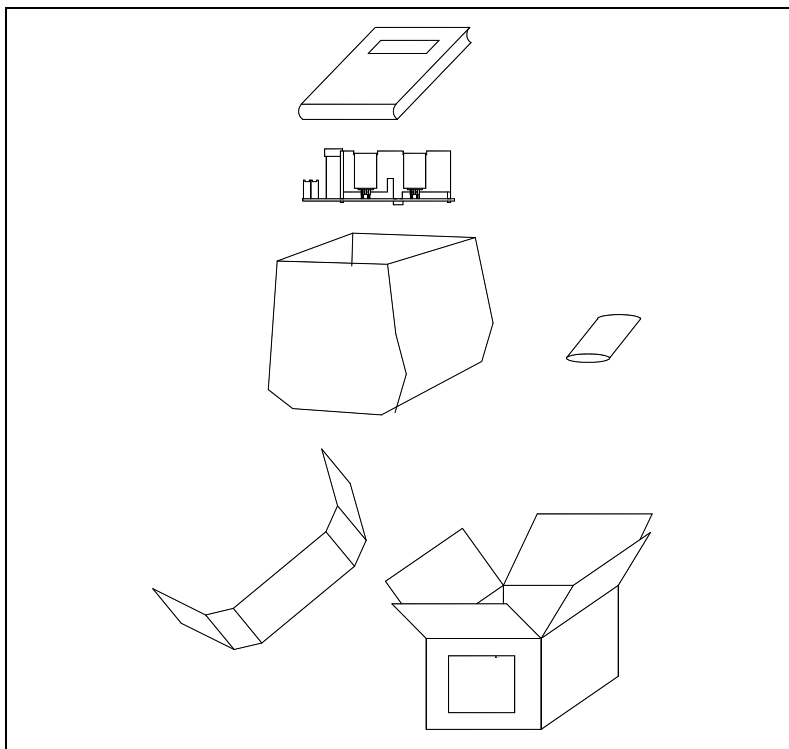
Opslag

Tijdens het transport en de opslag van de controllers moeten de volgende omgevingscondities aanwezig zijn:

- temperatuur: van -20 °C tot +70 °C
- relatieve vochtigheid: 0 – 95 % (niet condenserend)

Vorbereitung voor installatie

De controller wordt in een speciale beschermende verpakking geleverd; als er schade wordt geconstateerd die tijdens het transport veroorzaakt zou kunnen zijn, meteen contact opnemen met het plaatselijke verkoopkantoor.

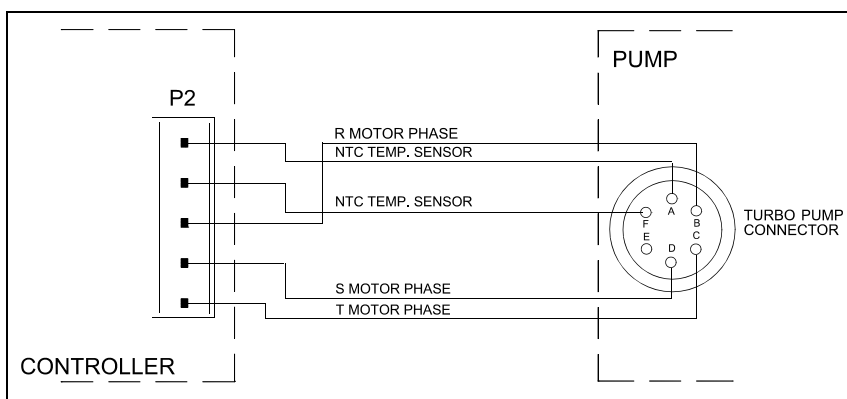


Figuur 1 Verpakking van de controller

Zorg er bij het uitpakken voor dat de controller niet kan vallen of stoten te verduren krijgt. Laat de verpakking niet ergens buiten achter. Het verpakkingsmateriaal is volledig recyclebaar en voldoet aan de EEG milieuriichtlijn 85/399.

Alvorens de controller te installeren, moeten enkele verbindingskabels worden voorbereid, zoals bijvoorbeeld kabels voor verbinding met de voeding, kabels voor de logicaverbindingen en de kabel voor aansluiting op de pomp. Op de volgende afbeeldingen zijn de verbindingen tussen controller en pomp en tussen controller en voedingsinrichting schematisch weergegeven.

Voor meer informatie over deze verbindingen en over de kenmerken van de voedingstransformator (leverbaar als extra) wordt verwezen naar de bijlage "Technical Information".



Figuur 2 Verbindingskabel tussen controller en pomp

Installatie

WAARSCHUWING!



In de controller ontwikkelen zich hoge spanningen die ernstige schade kunnen veroorzaken. Alvorens installatie- of onderhoudswerkzaamheden uit te voeren, de controller van de voeding afkoppelen. In de controller ontwikkelen zich hoge temperaturen die zware schade kunnen veroorzaken. Bij de definitieve installatie van de controller, deze op passende wijze tegen eventueel contact beschermen.

OPMERKING

De controller die definitief in het systeem geïnstalleerd wordt moet zodanig geplaatst worden dat de koellucht vrij rondom het apparaat kan circuleren. De controller mag niet geïnstalleerd en/of gebruikt worden in ruimten die blootgesteld zijn aan de weersomstandigheden (regen, vorst, sneeuw), stof, agressieve gassen, of in ruimten met explosiegevaar of zeer hoog brandgevaar.

Tijdens de werking moeten de volgende omgevings- condities aanwezig zijn:

- temperatuur: van 0 °C tot +40 °C
- relatieve vochtigheid: 0 – 95 % (niet condenserend).

Gebruik

In deze paragraaf worden de voornaamste bedieningswijzen uitgelegd. Voor meer informatie of procedures die aansluitingen of speciale opties betreffen wordt verwezen naar de paragraaf "USE" van de bijlage "Technical Information". Breng, alvorens de controller in gebruik te nemen, alle elektrische en pneumatische aansluitingen tot stand en raadpleeg hiervoor de handleiding van de aan te sluiten pomp.

WAARSCHUWING!



Raadpleeg de betreffende handleiding van de pomp, voor een veilige bediening, voorafgaand aan het starten.

Gebruiksprocedures

Inschakelen van de controller

Om de controller in te schakelen, de voedingskabel in de netcontactdoos inbrengen.

Starten van de pomp

Voor het starten van de pomp moet pin 12 van connector J4 open gelaten worden, zodat de pomp start wanneer 24 Vdc spanning aan de controller wordt geleverd.

Stoppen van de pomp

Voor het stoppen van de pomp moeten de pins 12 en 15 van connector J4 kortgesloten worden.

Onderhoud

De controllers van de serie TwisTorr 74 FS zijn onderhoudsvrij. Eventuele werkzaamheden moeten door bevoegd personeel worden uitgevoerd. In geval van storing is het mogelijk om de reparatiedienst van Agilent of de "Agilent advanced exchange service" in te schakelen: zo krijgt men een ruilcontroller ter vervanging van de defecte controller.

WAARSCHUWING!



Alvorens werkzaamheden aan de controller uit te voeren, de voedingskabel afkoppelen.

Mocht de controller gesloopt worden, ga dan overeenkomstig de specifieke nationale wetgeving te werk.

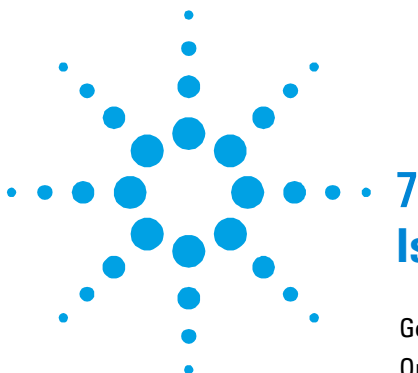
Afvalverwerking

Betekenis van het logo "WEEE" op de etiketten. Het onderstaande symbool wordt aangebracht in overeenstemming met de EG-richtlijn "WEEE". Dit symbool (**alleen geldig voor de landen van de Europese Gemeenschap**) geeft aan dat het product waarop het is aangebracht, NIET mag worden afgevoerd samen met normaal huisvuil of industrieel afval, maar gescheiden moet worden ingezameld. De eindgebruiker wordt dus verzocht contact op te nemen met de leverancier van het apparaat, zij het de fabrikant of een wederverkoper, om het proces van gescheiden inzameling en verwerking in gang te zetten, na de van toepassing zijnde termen en voorwaarden van het verkoopcontract te hebben gecontroleerd.



Voor meer informatie wordt verwezen naar:
<http://www.agilent.com/environment/product/index.shtml>

6 **Bedrijfshandleiding**
Afvalverwerking



7 Instruktionsbog

Generel Information	74
Opbevaring	75
Forberedelser før installation	76
Installation	78
Anvendelse	79
Instruktion	79
Start af styreenheden	79
Start af pumpen	79
Stop af pumpen	79
Vedligeholdelse	80
Bortskaffelse	81

Oversættelse af originalinstruktionerne



Generel Information

Dette materiel er beregnet til professionel anvendelse. Brugeren bør læse denne brugsanvisning og anden yderligere information fra Agilent, før udstyret anvendes. Agilent tager ikke ansvar for skader helt eller delvis som følge af tilsidesættelse af disse instruktioner, fejlagtig brug af personer uden tilstrækkelig kendskab, ukorrekt anvendelse af udstyret eller håndtering, der strider imod gældende lokale regler. Styreenheden i TwisTorr 74 FS PCB-serien er en mikroprocessorstyret frekvens-omformer, der består af komponenter med fast tilstand. Styreenheden er udstyret med selvdiagnose og selvbeskyttelsesfunktioner. Styreenheden omfatter alle midler, der kræves for automatisk drift af pumperne i TwisTorr 74 FS serien. Styreenheden kontrollerer pumperne i TwisTorr 74 FS serien (med en ti-trins-proces) i forbindelse med start. Spænding og strøm reguleres i forhold til pumpens opnåede hastighed. Hjælpekontakter forsyner kontrol til fjernstart og -stop af pumpen, signaler om pumpens tilstand, blokerings signaler (til tryk- og vandføringsafbrydere, osv.). Styreenheden skal forsynes med jævnspænding på 24 Vdc.

De følgende afsnit indeholder al information der behøves, for at garantere operatørens sikkerhed under anvendelsen. Detaljeret information findes i bilaget "Technical Information".

I brugsanvisningen anvendes følgende standard-rubrikker:

FORSIGTIG!

Denne advarselsmeddelelse vises før procedurer, der skal følges nøje for ikke at risikere maskinskader.

ADVARSEL!



Advarselsmeddelelserne informerer operatøren om, at en speciel procedure eller en vis type arbejde skal udføres præcist efter anvisningerne. I modsat fald er der risiko for svære personskader.

BEMÆRK

Dette gør opmærksom på vigtig information i teksten.

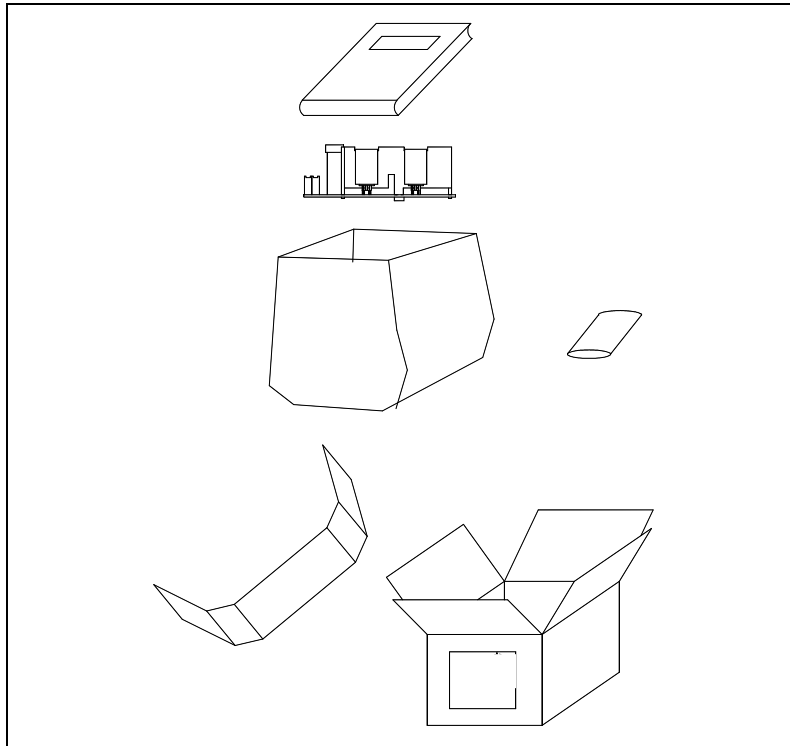
Opbevaring

Følgende krav til omgivelserforholdene gælder ved transport og opbevaring af styreenheden:

- temperatur: fra -20 °C til +70 °C
- relativ luftfugtighed: 0 – 95 % (ikke kondenserende)

Forberedelser før installation

Styreenheden leveres i en speciel beskyttende emballage. Kontakt den lokale forhandler, hvis emballagen viser tegn på skader, der kan være opstået under transporten.



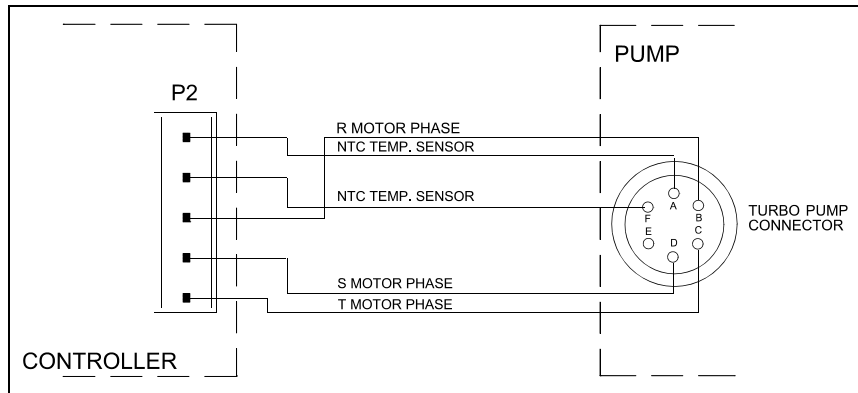
Figur 1 Styreenhedens emballage

Sørg for at styreenheden ikke tabes eller udsættes for stød ved udpakningen.

Smid ikke emballagen ud. Materialet kan genbruges 100 % og opfylder EU-direktiv 85/399 om miljøbeskyttelse.

Inden styreenheden installeres, skal man klargøre de nødvendige forbindelseskabler, d v s forbindelseskabler til strømforsyning og pumpe samt kabler til logiske signaler. Følgende billeder viser de nødvendige forbindelser mellem styreenhed og pumpe samt mellem styreenhed og strømkilde.

For yderligere information om forbindelser og strømkilde (ekstra tilbehør) henvises til bilag "Technical Information".



Figur 2 Forbindelseskabel mellem styreenheden og pumpe

Installation

ADVARSEL!



Spændinger frembragt i styreenheden kan nå høje værdier og forårsage stor skade. Frakobl altid strømkablet, inden der udføres installations- eller vedligeholdelsesarbejde på styreenheden. Temperaturen frembragt i styreenheden kan nå høje værdier og forårsage stor skade. I forbindelse med permanent installation skal styreenheden og transformatoren på egnet måde beskyttes mod utilsigtet tilslutning.

BEMÆRK

Styreenheden skal anbringes på en sådan måde, at luft kan cirkulere frit omkring apparatet. Installér og anvend ikke styreenheden i miljøer, der udsættes for påvirkninger fra atmosfæren (regn, sne, is), damp, aggressive gasser, og ligeledes ikke i eksplosivt eller brandfarligt miljø.

Følgende krav til omgivelsesforholdene gælder ved drift:

- temperatur: fra 0 °C til +40 °C
- relativ luftfugtighed: 0 – 95 % (ikke kondenserende)

Anvendelse

Dette afsnit beskriver de vigtigste driftsprocedurer. For en detaljeret beskrivelse samt procedurer, der involverer tilslutninger eller tilbehør, henvises til afsnittet "USE" i bilag "Technical Information". Inden styreenheden anvendes, bør samtlige elektriske og pneumatiske tilslutninger udføres. Læs brugsanvisningen før pumpen tilsluttes.

ADVARSEL!

For at operere den sikkert, henvises der til den relevantepumpemanual før start.

Instruktion

Start af styreenheden

Styreenheden startes ved at sætte strømkablet i vægudtaget.

Start af pumpen

Pumpen startes ved at lade stift 12 på konnektoren J4 være åben, således at pumpen startes, når kontrolenheden forsynes med spænding på 24 Vdc.

Stop af pumpen

Pumpen stopper ved at kortslutte stift 12 og stift 15 på konnektoren J4.

Vedligeholdelse

Styreenhederne i TwisTorr 74 FS-serien behøver ikke nogen vedligeholdelse. Ethvert indgreb på pumpen skal foretages af autoriseret personale. Hvis pumpen går i stykker, kan man benytte sig af Agilents reparations-service eller Agilent udvekslingsservice, hvor man kan få en repareret pumpe i bytte for den, der er gået i stykker.

ADVARSEL!

Inden der foretages noget som helst indgreb på styreenheden, skal strømmen først afbrydes.



Skrotning af pumpen skal foregå i overensstemmelse med det pågældende lands særlige love.

Bortskaffelse

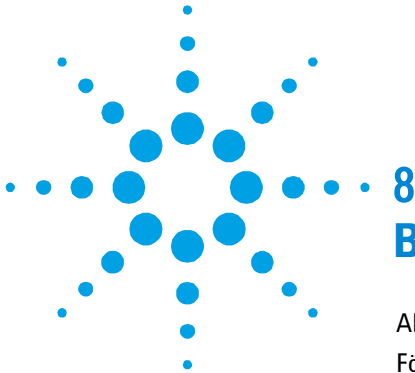
Betydningen af "WEEE" logoet på mærkaterne. Nedenstående symbol anvendes i overensstemmelse med det såkaldte EU-direktiv "WEEE". Symbolet (**kun gældende for EU-landene**) viser, at produktet, som det sidder på IKKE må bortskaffes sammen med affald fra private husholdninger eller industriel affald men skal indleveres på en godkendt affaldsstation. Vi opfordrer derfor slutbrugeren til at kontakte leverandøren af anordningen, enten fabrikken eller en forhandler, for igangsættelse af afhentnings- og bortskaffelsesprocessen efter nøje at have kontrolleret betingelserne i salgskontrakten.



For yderligere oplysninger henvises til:

<http://www.agilent.com/environment/product/index.shtml>

7 **Istruktionsbog**
Bortskaffelse



8

Bruksanvisning

Allmän information	84
Förvaring	85
Förberedelser för Installation	86
Installation	88
Användning	89
Instruktioner för bruk	89
Start av styrenheten	89
Start av pumpen	89
Stopp av pumpen	89
Underhåll	90
Bortskaffning	91

Översättning av originalinstruktionerna



Allmän information

Utrustningen är avsedd för yrkesmässig användning. Användaren bör läsa denna bruksanvisning, samt övrig dokumentation från Agilent före användning av utrustningen. Agilent tar inget ansvar för skador som helt eller delvis orsakats av åsidosättande av instruktionerna, olämplig användning av person utan tillräcklig kunskap, obehörigt bruk av utrustningen eller hantering som strider mot gällande lokala föreskrifter.

Styrenheten i TwisTorr 74 FS PCB-serien är en mikro-processorstyrd frekvensomvandlare som består av komponenter med fast tillstånd. Styrenheten är försedda med självdiagnos- och självskyddsfunktion. Styrenheten omfattar alla kretsar som behövs för automatisk drift av pumparna i TwisTorr 74 FS serien.

Styrenheten kontrollerar pumparna i TwisTorr 74 FS-serien (med en tiostegs-process) i samband med start. Spänning och ström regleras i förhållande till pumpens uppnådda hastighet. Hjälpkontakter erbjuder kontroller för fjärrstart och fjärrstopp av pumpen, signaler för pumpens tillstånd, blockeringssignaler (för tryckvakter, kontrollbrytare för vattenflöde osv). Styrenheten ska förses med likspänning på 24 Vdc.

De följande avsnitten innehåller all information som behövs för att garantera operatörens säkerhet under driften. Detaljerade uppgifter finns i bilagan "Technical information".

I bruksanvisningen används följande standard-rubriker:

OBSERVER!

Detta varningsmeddelande visas framför procedurer som måste följas exakt för att undvika skador på maskinen.

WARNING!



Varningsmeddelandena informerar operatören om att en speciell procedur eller en viss typ av arbete måste utföras exakt enligt anvisningarna. I annat fall finns risk för svåra personskador.

OBSERVERA

Detta visar på viktig information i texten.

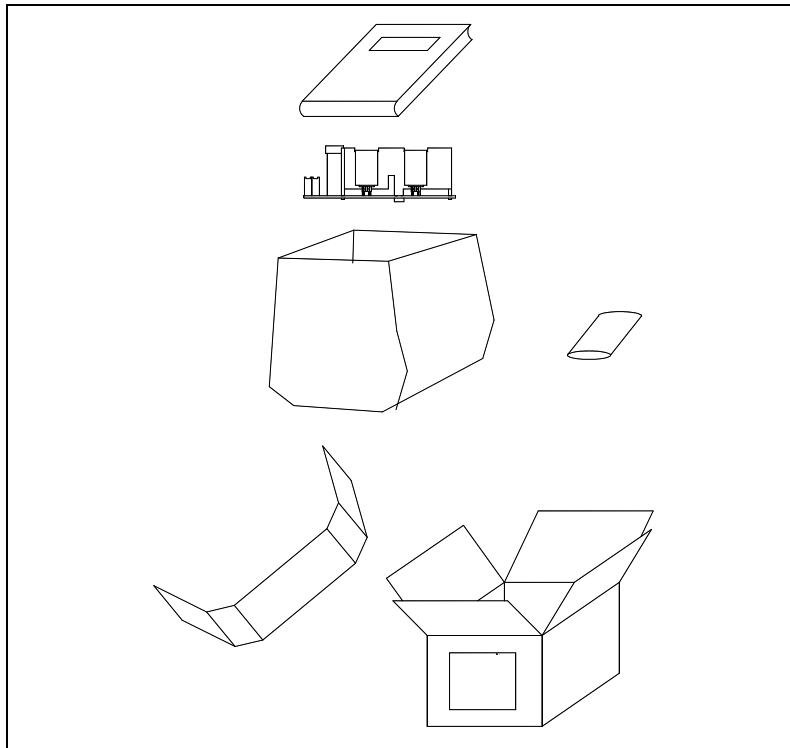
Förvaring

Följande krav på omgivningsförhållanden gäller vid transport och förvaring av styrenheten:

- temperatur: från -20 °C till +70 °C
- relativ luftfuktighet: 0 – 95 % (utan kondens)

Förberedelser för Installation

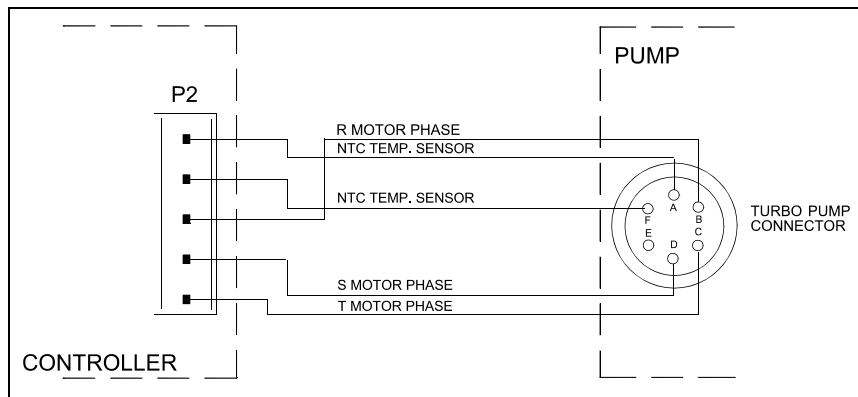
Styrenheten levereras i ett särskilt skyddande emballage. Kontakta det lokala försäljningskontoret om emballaget visar tecken på skador som kan ha uppstått under transporten. Se till att styrenheten inte tappas eller utsätts för stötar vid uppackningen.



Figur 1 Styrenhetensförpackning

Kasta inte packmaterialet i soporna. Materialet är återvinningsbart till 100 % och uppfyller EU-direktiv 85/399 om miljöskydd. Innan styrenheten installeras ska du förbereda de nödvändiga anslutningskablarna, det vill säga anslutningskablar till matarspänning och pump samt kablar för logiska signaler. Följande bilder visar de nödvändiga anslutningarna mellan styrenhet och pump samt mellan styrenhet och strömkälla.

För ytterligare information om anslutningar och strömkälla (tillval) hänvisas till bilaga "Technical Information".



Figur 2 Anslutningskabel mellan styrenhet och pump

Installation

WARNING!



Spänningen inuti styrenheten kan nå höga värden och förorsaka allvarliga skador. Frånkoppla alltid strömkabeln innan något installations- eller underhålls-moment utförs på styrenheten. Temperaturen inuti styrenheten kan nå höga värden och förorsaka allvarliga skador. I samband med permanent installation ska styrenhet och transformator lämpligen skyddas mot oavsiktlig kontakt.

OBSERVERA

Styrenheten ska installeras permanent på ett sådant ställe, att kylluften kan cirkulera fritt kring apparaten. Installera och använd inte styrenheten i miljöer som utsätts för påverkan från atmosfären (regn, snö, is), damm, aggressiva gaser, och inte heller i explosiv eller brandfarlig miljö.

Följande krav på omgivningsförhållanden gäller vid drift:

- temperatur: från 0 °C till +40 °C
- relativ luftfuktighet: 0 – 95 % (utan kondens)

Användning

Detta avsnitt beskriver de viktigaste driftmomenten. För en detaljerad beskrivning samt beträffande moment som involverar anslutningar eller tillbehör hänvisas till avsnittet "USE" i bilaga "Technical Information". Innan styrenheten används bör samtliga elektriska och pneumatiska anslutningar utföras. Läs bruksanvisningen för den anslutna pumpen.

WARNING!



Referera till den relevanta pumpmanualen innan start för att använda den på ett säkert sätt.

Instruktioner för bruk

Start av styrenheten

Styrenheten startas enkelt genom att strömkabeln sätts i vägguttaget.

Start av pumpen

Pumpen startas genom att stift 12 öppnas på kontakt J4 så att pumpen startar när spänningen 24 Vdc når styrenheten.

Stopp av pumpen

Pumpen stoppas genom att stift 12 och 15 kortsluts på kontakt J4.

Underhåll

Styrenheterna i TwisTorr 74 FS-serien är underhållsfria. Allt servicearbete måste utföras av auktoriserad personal.

Om styrenheten havererar, kontakta Agilent reparation- sverkstad eller Agilent utbytesservice, som kan ersätta styrenheten med en renoverad styrenhet.

WARNING!



Innan något arbete utförs på styrenheten måste dess strömförsörjning brytas.

Skrotning av pumpen skall ske enligt gällande lagstiftning.

Bortskaffning

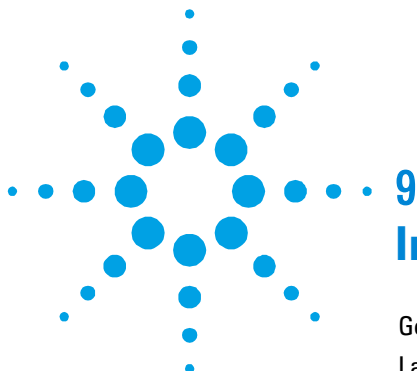
Betydelse av logotypen "WEEE" på etiketterna. Symbolen som visas nedan har tillämpats i enlighet med CD-direktivet som har betecknats som "WEEE". Den här symbolen (**gäller endast i de länder som tillhör den Europeiska Unionen**) indikerar att produkten på vilken symbolen har applicerats INTE får skaffas bort tillsammans med vanliga hushålls- eller industriavfall, men att däremot ett differentierat uppsamlingsystem måste upprättas. Vi rekommenderar därför att slutanvändaren tar kontakt med leverantören av anordningen, oberoende om det handlar om moderföretaget eller återförsäljaren, för att kunna starta uppsamlings- och bortskaffningsprocessen, detta efter lämplig kontroll av kontraktsevenliga tidsgränser och försäljningsvillkor.



För mer information, se:

<http://www.agilent.com/environment/product/index.shtml>

8 Bruksanvisning
Bortskaffning



9

Instruksjon Manual

Generell informasjon	94
Lagring	95
Forberede installasjonen	96
Installasjon	98
Bruk	99
Instruksjoner for bruk	99
Starte styreenheten	99
Starte pumpen	99
Stoppe pumpen	99
Vedlikehold	100
Eliminering	101

Oversetting av den opprinnelige samsvarserklæringen



Generell informasjon

Dette utstyret er beregnet til bruk av profesjonelle brukere. Brukeren bør lese denne brukerveiledningen og all annen informasjon fra Agilent før utstyret tas i bruk.

Agilent kan ikke holdes ansvarlig for hendelser som skjer på grunn av manglende oppfølging, selv delvis, av disse instruksjonene, feilaktig bruk av utrenet personell, ikke godkjente endringer av utstyret eller handlinger som på noen måte er i strid med nasjonale bestemmelser.

Styreenheten i TwisTorr 74 FS PCB serien er en mikroprosessorstyrt frekvensomvender, som består av komponenter med fast tilstand. Styreenheten har funksjoner for selvdiagnose og selvbeskyttelse.

Styreenheten omfatter alle kretser som er nødvendige for automatisk drift av pumpene i TwisTorr 74 FS serien.

Styreenheten kontrollerer pumpene i TwisTorr 74 FS serien (med en titrinnspesess) ved oppstart. Spenning og strømstyrke justeres i forhold til pumpens oppnådde hastighet. Hjelpekontakter gir muligheter for fjernstyrt start og stopp av pumpen, signaler for pumpens tilstand, blokkeringssignaler (for trykksensorer, kontrollbrytere for vannstrøm osv). Styreenheten skal forsynes med jevnspenning på 24 VDC.

De følgende avsnittene inneholder all informasjon som er nødvendig for å sikre brukeren når utstyret er i bruk. For mer detaljert bruk vises det til tillegget "Technical Information".

Denne veiledningen bruker følgende standard- protokoll:

FORSIKTIG!

Denne advarselen vises foran fremgangsmåter som, dersom de ikke følges, kan føre til at utstyret skades.

ADVARSEL!



Disse meldingene skal tiltrekke seg brukerens oppmerksomhet til en spesiell fremgangsmåte eller praksis som, hvis den ikke følges, kan medføre alvorlige skader.

MERK

Merknadene inneholder viktig informasjon som er hentet fra teksten.

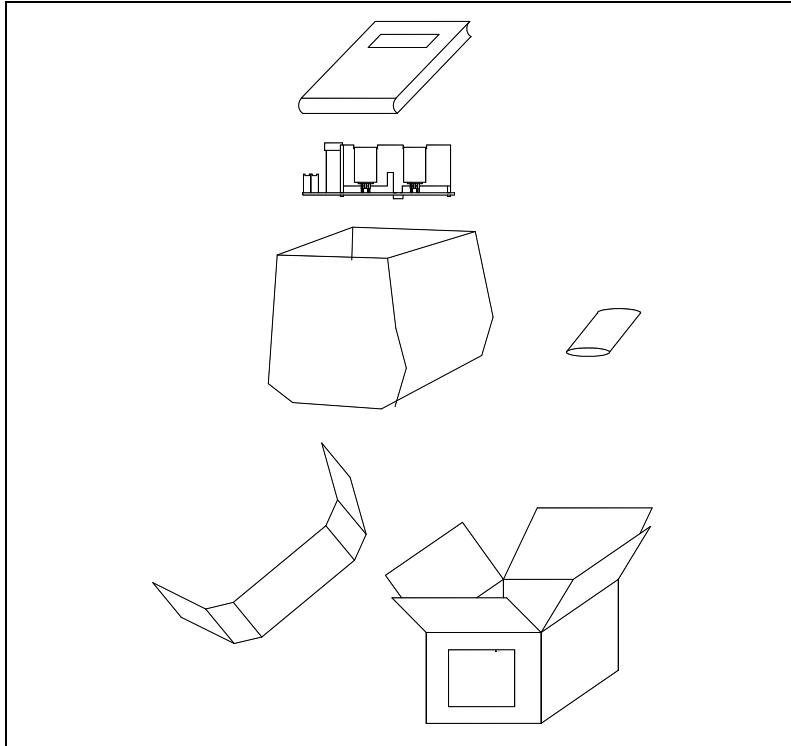
Lagring

Når styreenhetene transporteres eller lagres, må følgende forhold være oppfylt:

- temperatur: fra 20 °C til +70 °C
- relativ fuktighet: 0 – 95 % (uten kondens)

Forberede installasjonen

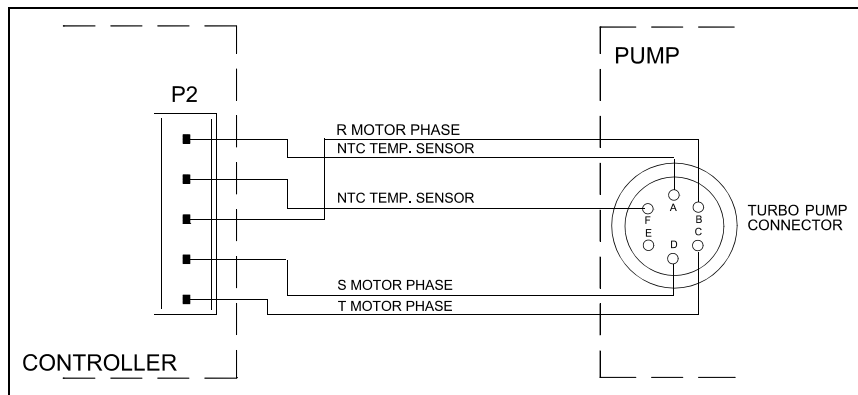
Styreenheten leveres i en spesiell beskyttelsesemballasje. Viser denne tegn på skader som kan ha oppstått under transporten, må du ta kontakt med det lokale salgskontoret.



Figur 1 Styreenhetens emballasje

Når styreenheten pakkes ut, må du passe på at den ikke slippes ned eller utsettes for noen form for støt. Emballasjen må ikke kastes på en ulovlig måte. Alle materialer er 100 % resirkulerbare og er i samsvar med EU-direktiv 85/399 om miljøbeskyttelse.

Før styreenheten installeres skal du klargjøre de nødvendige forbindelseskablene, det vil si forbindelseskabler til matespenning og pumpe samt kabler for logiske signaler. Følgende bilder viser de nødvendige forbindelsene mellom styreenhet og pumpe samt mellom styreenhet og strømkilde. For ytterligere informasjon om forbindelser og strømkilde (ekstrautstyr) vises det til vedlegget "Technical Information".



Figur 2 Forbindelseskabel mellom styreenheten og pumpen

Installasjon

ADVARSEL!



Spenningen inne i styreenheten kan nå høye verdier og kan føre til alvorlige skader. Kople alltid strømkabelen fra strømmettet før alle installasjons- eller vedlikeholds-arbeider som utføres på styreenheten. Temperaturen inne i styreenheten kan nå høye verdier og kan føre til alvorlige skader. I forbindelse med permanent installasjon skal styreenhet og transformator beskyttes mot utilsiktet kontakt på en passende måte.

MERK

Styreenheten skal installeres på slik måte, at kjøleluften kan sirkulere fritt rundt apparatet. Ikke installer eller bruk styreenheten i miljøer som utsettes for regn, snø eller is, støv, aggressive gasser, eksplosjonsfarlige miljøer eller i miljøer med stor brannfare.

Under bruk må følgende forhold respekteres:

- temperatur: fra 0 °C til + 40 °C
- relativ fuktighet: 0 – 95 % (uten kondens)

Bruk

Dette avsnittet beskriver de viktigste driftsmomentene. For en detaljert beskrivelse samt moment som omfatter tilkoplinger eller ekstrautstyr vises det til avsnittet "Use" i vedlegget "Technical Information". Før styreenheten tas i bruk bør samtlige elektriske og pneumatiske tilkoplinger gjøres. Les brukerveiledningen for pumpen som er tilkopleet.

ADVARSEL!

Se i den relevante pumpehåndboken for sikker bruk før du starter.



Instruksjoner for bruk

Starte styreenheten

Styreenheten startes ved å sette strømkabelen i veggkontakten.

Starte pumpen

Pumpen startes ved å la stift 12 på konnektoren J4 være åpen, slik at pumpen starter når styreenheten får spenningen på 24 Vdc.

Stoppe pumpen

Pumpen stopper når stiftene 12 og 14 på konnektoren J4 kortsluttes.

Vedlikehold

TwisTorr 74 FS seriens styreenheter er vedlikeholdsfrie. Alt arbeid på styreenheten må kun utføres av autorisert personell. Dersom styreenheten stanser, må du ta kontakt med Agilents reparasjonsservice eller med Agilents avanserte bytteservice, som kan tilby overhalte styreenheter til erstatning for den ødelagte styreenheten.

ADVARSEL!

Før noe arbeid utføres på styreenheten, må den frakoples strømmettet.



Dersom en styreenhet skal kasseres, må dette skje i henhold til nasjonale bestemmelser.

Eliminering

Betydelsen av symbolet på logo "WEEE" på etikettene. Symbolet nedenunder som finnes, er anvendt i henhold til EC-direktiv kalt "WEEE". Dette symbolet (**som bare gjelder for land i Det europeiske fellesselskap**), viser at produktet som det sitter på, IKKE må behandles som vanlig hus-industriavfall, men må legges i kildesortering. Derfor oppfordrer man den sluttelige brukeren av anordningen å henvende seg til leverandøren av anordningen, som kan være et firma eller en forhandler, som sørger for oppsamling og eliminering etter å ha kontrollert avtal og betingelser i kjøpekontrakten.



For mer informasjon se:

<http://www.agilent.com/environment/product/index.shtml>

9 Instruksjon Manual

Eliminering



10 Ohjekäsikirja

Yleisiä tietoja	104
Varastointi	105
Valmistelut asennusta varten	106
Asennus	108
Käyttö	109
Käyttötoimenpiteet	109
Valvojan päälle pano	109
Pumpun käynnistys	109
Pumpun pysäyttäminen	109
Huolto	110
Hävittäminen	111

Alkuperäisen ohjeiden käännös



Yleisiä tietoja

Tämä laite on tarkoitettu ammattimaiseen käyttöön. Ennen laitteen käyttöönottoa käyttäjän tulee lukea huolellisesti mukana seuraava käyttöohje sekä kaikki muut Agilentin toimittamat lisätiedot. Agilent ei vastaa seurauksista, jotka johtuvat laitteen käyttöohjeiden täydellisestä tai osittaisesta laiminlyömisestä, ammattitaidottomien henkilöiden suorittamasta laitteen virheellisestä käytöstä, valtuuttamattomista toimenpiteistä tai maakohtaisten säädösten ja normien vastaisesta käytöstä. Sarjan TwisTorr 74 FS PCB valvoja on mikroprosessorin valvoma kiinteistä materiaaleista tehtyjä taajuudenmuuntimia, jotka kykenevät itsemäärittelyyn ja itsesuojaukseen. Ne yhdistävät kaikki sähköpiirit, jotka ovat välttämättömiä TwisTorr 74 FS-sarjan pumpun automaattiselle toiminnalle. Valvoja ajaa TwisTorr 74 FS-sarjan pumppuja (kymmenportaisessa järjestelmässä) käynnistysvaiheessa valvoen jännitettä ja sähkövirtaa suhteessa pumpunsaavuttamaan nopeuteen.

Apuliittimiä käyttäen on mahdollista käyttää kauko-ohjattua pumpun käynnistystä ja pysähdystä, signaaleja, jotka ilmaisevat pumpun toimintatilan, pysäytyssignaaleja (painokatkaisimille, veden virtauksen säätelykatkaisijalle jne.). Valvojan syöttöjännitteen tulee olla 24 Vdc.

Seuraavilla sivuilla on luettavissa tarpeelliset tiedot laitteen käyttäjän turvallisuuden takaamiseksi laitteen käytön aikana. Yksityiskohtaiset tiedot löytyvät liitteestä "Technical Information".

Tämä käsikirja käyttää seuraavia merkintöjä:

HUOMIO!

Huomio-merkinnät varoittavat toiminnoista, joiden laiminlyönti voi johtaa laitteen vaurioitumiseen.

VAROITUS!



Vaara-merkinnät saavat käyttäjän huomion kiinnittymään erityisiin toimintotapoihin, joiden seuraamatta jättäminen voi aiheuttaa vakavia henkilövaurioita.

HUOM

Huomiot sisältävät tärkeätä tekstistä otettua tietoa.

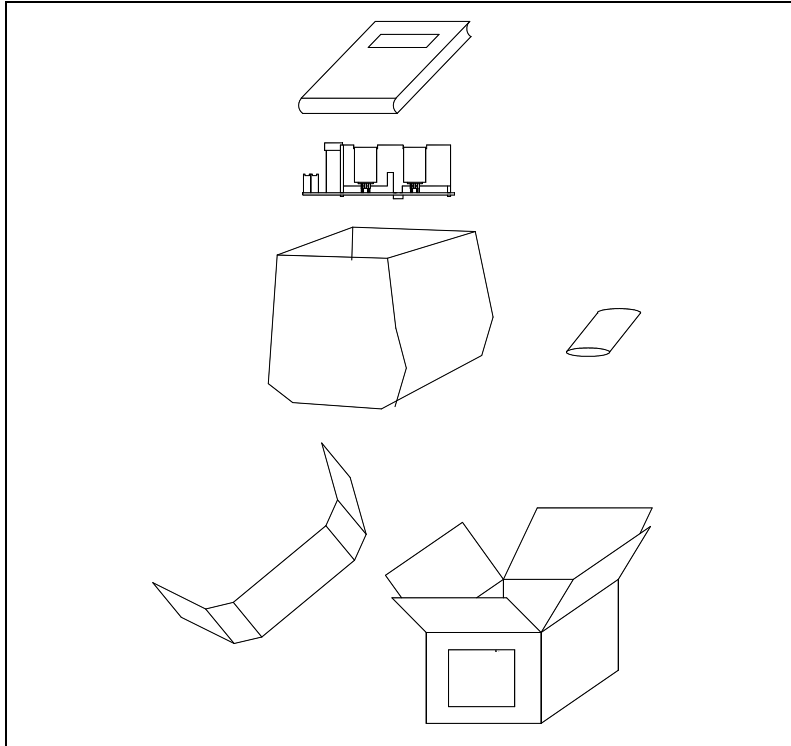
Varastointi

Valvojan kuljetuksen ja varastoinnin aikana tulevat seuraavat ympäristövaatimukset olla täytettyinä:

- lämpötila: -20 °C ja +70 °C asteen välillä
- suhteellinen kosteus: 0 – 95 % (ilman lauhdetta)

Valmistelut asennusta varten

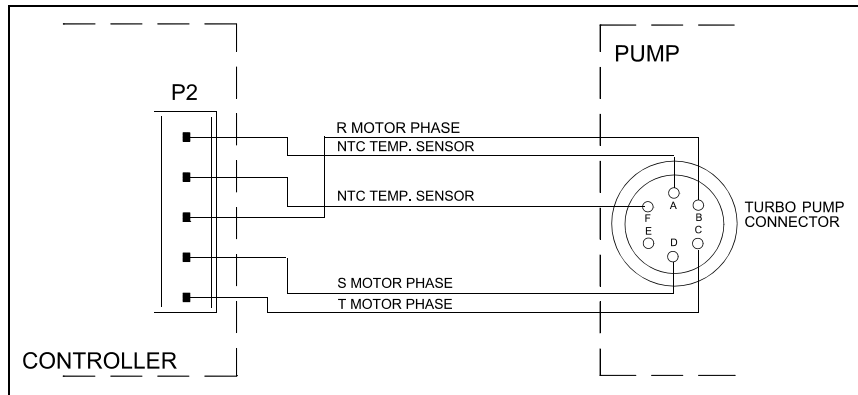
Valvoja toimitetaan erityisessä suojaavassa pakkauksessa. Mikäli havaitsette mahdollisesti kuljetuksen aikana sattuneita vaurioita, ottakaa yhteys paikalliseen myyntitoimistoon.



Kuva 1 Valvojan pakkaus

Pakkauksen purkamisen yhteydessä huolehtikaa, että valvoja ei pääse putoamaan ja välttäkää sen joutumista iskujen kohteeksi. Älkää jättäkö pakkausta ympäristöön. Materiaali voidaan kokonaisuudessaan kierrättää ja se vastaa EY:n 85/399 direktiiviä ympäristön suojelusta.

Ennen kuin asennatte valvojan, pitää valmistaa liitoskaapeleita, esimerkiksi sähkövirtakaapeleita, loogisten signaleiden kaapeleita ja pumpun liitoskaapeli. Seuraavissa piirroksissa näkyvät valvojan ja pumpun välillä ja valvojan ja sähkön syötön välillä suoritettavat liitännät. Liitännän ja (lisä-) muuntajan lähimmät tiedot löytyvät liitteestä "Technical Information".



Kuva 2 Valvojan ja pumpun välinen liitoskaapeli

Asennus

VAROITUS!



Valvojan sisällä syntyy korkeajännitettä, joka voi aiheuttaa vakavia vammoja. Ennen minkätahansa valvojan huolto- tai asennustoimenpiteen suorittamista, irroittakaa valvoja sähköverkosta. Valvojan sisällä syntyy ylikuumentumista, joka voi aiheuttaa vakavia vammoja. Suojatkaa valvoja ja muuntaja tahattomalta kosketukselta, kun ne ovat jo asennetut.

HUOM

Lopulliseen järjestelmään asennettu valvoja on asetettava siten, että jäähdytysilma pystyy kiertämään vapaasti laitteen sisällä. Älä asenna ja/tai käytä valvojaa ilmaston tekijöille (sade, jää, lumi), pölylle, syövyttävillä kaasuilla, räjähdyksille tai tulelle alttiissa paikoissa.

Toiminnan aikana tulee noudattaa seuraavia ympäristönoloja koskevia sääntöjä:

- lämpötila: 0 °C ja +40 °C välillä
- suhteellinen kosteus: 0 – 95 % välillä (ilman lauhdetta)

Käyttö

Tähän kappaleeseen on kirjattu tärkeimmät käyttötoimenpiteet. Tarkempia lisätietoja sekä kytkentöjä, että valinnaisia lisälaitteita koskevien toimenpiteiden suorittamista käsittäviä tietoja löydätte kappaleesta "Käyttö", joka on "Technical Information" -kappaleen liitteenä. Ennen valvojan käyttöä suorittaakaa kaikki sähkökytkennät seuraten kytkettävän pumpun käyttöohjeita.

VAROITUS!



Tutustu pumpun käyttöohjeeseen ennen aloittamista pumpun turvallisesti käyttämiseksi.

Käyttötoimenpiteet

Valvojan päälle pano

Valvoja käynnistyy asettamalla virtakaapeli pistorasiaan.

Pumpun käynnistys

Jättäkää pumpun käynnistämiseksi liittimen J4 pistike 12 auki siten, että pumppu käynnistyy valvojan 24 Vdc jännitteen päälle panon yhteydessä.

Pumpun pysäyttäminen

Pumppu pysähtyy, kun liittimen J4 pistikkeet 12 ja 15 asetetaan oikosulkuun.

Huolto

TwisTorr 74 FS sarjan valvojat eivät kaipaakaan minkäänlaista huoltoa. Mahdolliset valvojan tehtävät toimenpiteet tulee jättää aina valtuutetun henkilön tehtäviksi. Toimintahäiriön sattuessa on mahdollista käyttää Agilentin korjauspalvelua tai "Agilent advance exchange service" -palvelua, jolloin on mahdollista vaihtaa rikkiöntunut valvoja ladattuun valvojaan.

VAROITUS!



Ennen minkätähansa valvojan tehtävän toimenpiteen suorittamista irroittakaa sähkökaapeli pistorasiasta.

Mikäli valvoja täytyy romuttaa, toimikaa sen hävittämisessä kansallisten säädösten ja normien määrävällä tavalla.

Hävittäminen

Pakkausmerkinnöissä olevan WEEE-logon merkitys. Alla näkyvä merkki on lisätty pakkaukseen EY:n ns. WEEE-direktiivin mukaisesti. Merkki (**koskee ainoastaan Euroopan Unionin jäsenmaita**) tarkoittaa, että tuotetta EI saa hävittää tavallisen kotitalous- tai teollisuusjätteen mukana, vaan se on toimitettava erilliseen keräyspisteeseen. Loppukäyttäjää kehoitetaan sen vuoksi ottamaan keräys- ja hävittämisprosessia varten yhteyttä laitteen toimittajaan, olipa se sitten laitteen valmistaja tai jälleenmyyjä, tarkastettuaan ensin kaupan sopimusehdot.

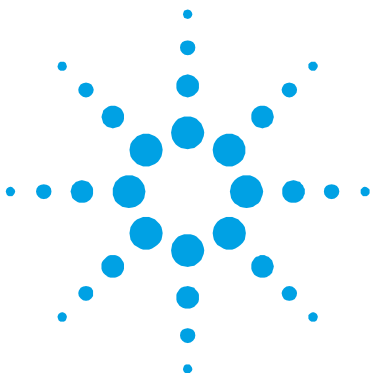


Katso lisätietoja kohdasta:

<http://www.agilent.com/environment/product/index.shtml>

10 Ohjekäsikirja

Hävittäminen



11 Felhasználói Kézikönyv

Általános információk	114
Tárolás	115
A telepítésre való előkészítés	116
Telepítés	118
Használat	119
Használati eljárások	119
A vezérlő bekapcsolása	119
A szivattyú beindítása	119
A szivattyú leállítása	119
Karbantartás	120
Megsemmisítés	121

Az eredeti utasítás fordítása



Általános információk

Ez a berendezés hivatásos használatra tervezett. A felhasználónak, a berendezés használatát megelőzően, gondosan el kell olvasnia a felhasználói kézikönyvet és minden egyéb, a Agilent által szolgáltatott információt. A Agilent elhárít magától minden esetleges felelősséget, mely a felhasználói utasítások teljes vagy részleges figyelmen kívül hagyásához, a nem kiképzett személyzet által történő nem megfelelő használathoz, meghatalmazás nélküli beavatkozásokhoz, vagy a specifikus nemzeti irányelvekkel ellentétes használathoz kötődik.

A TwisTorr 74 FS PCB sorozatú vezérlők mikroprocesszoros vezérlésű, szilárdtest-elemekkel realizált frekvencia-átalakítók, öndiagnosztikai és saját védelmi képességekkel. A TwisTorr 74 FS sorozatú szivattyúk automatikus működtetéséhez szükséges valamennyi áramkört. A vezérlők TwisTorr 74 FS sorozatú szivattyúkat irányítanak (egy tíz lépésből álló folyamat révén), az indítási fázisban a feszültséget és az áramot a szivattyú által elért sebesség függvényében szabályozva. Segédcsatlakozók révén rendelkezésre állnak a szivattyú indítási és leállítási távparancsai, a szivattyú működési állapotát mutató jelek, az elő-vákuum szivattyú indításához és leállításához szükséges parancsok, valamint a reteszelő jelek (nyomáskapcsolók, vízáram-érzékelő kapcsolók stb. számára). A vezérlőt 24 Vdc egyenfeszültséggel kell táplálni.

Az elkövetkező szakaszok tartalmazzák mindazokat az ismereteket, melyek a kezelő biztonságát hivatottak garantálni a berendezés használata során. Részletes információk találhatóak a "Technical Information" című mellékletben.

E kézikönyv az alábbi megegyezésszerű jelöléseket alkalmazza:

FIGYELEM!

A figyelmeztető üzenetek azon eljárások előtt kerülnek megjelenítésre, melyek figyelmen kívül hagyása a berendezés károsodásához vezethet.

VESZÉLY!



A veszélyt jelző üzenetek olyan eljárásra vagy specifikus tevékenységre hívják fel a kezelő figyelmét, melyek, ha nem megfelelő módon kerülnek végrehajtásra, súlyos személyi sérüléseket okozhatnak.

MEGJEGYZÉS

A megjegyzések fontos, a szövegből kivonatolt információkat tartalmaznak.

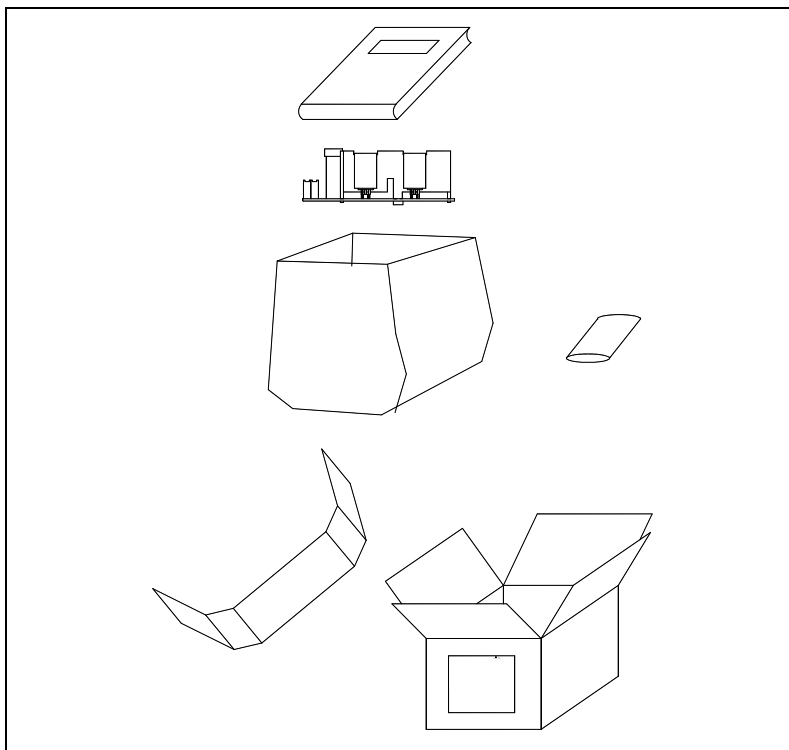
Tárolás

A vezérlők szállítása és tárolása során az alábbi környezeti feltételeket kell biztosítani:

- hőmérséklet: -20 °C és +70 °C között
- relatív nedvességtartalom: 0 – 95 % (nem lecsapódó).

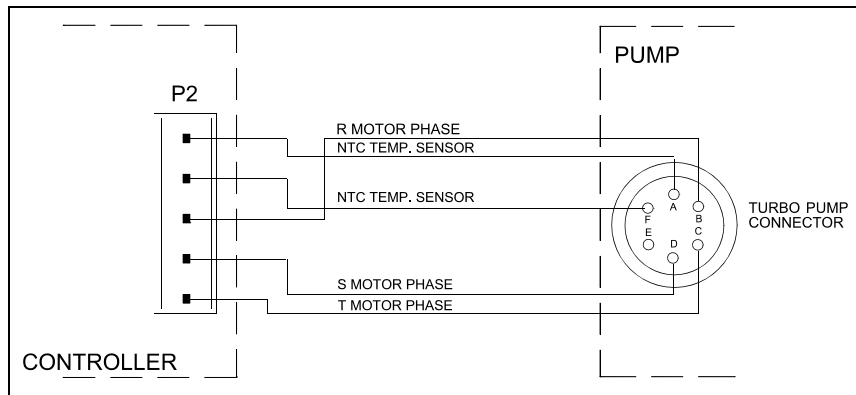
A telepítésre való előkészítés

A vezérlő speciális védőcsomagolásban kerül leszállításra; ha ez károsodás jeleit mutatja, melyek a szállítás során keletkezettek, kapcsolatba kell lépni a helyi forgalmazó irodával.



Ábra 1 A vezérlő csomagolása

A kicsomagolás során különleges figyelmet kell fordítani arra, hogy a vezérlő le ne essék, és ne szenvedjen ütődéseket. A csomagolóanyagot nem szabad szétszórni a környezetben. Az anyag teljes mértékben újrafelhasználható, s megfelel a környezetvédelemmel kapcsolatos EGK 85/399-es irányelvnek. A vezérlő telepítését megelőzően, elő kell készíteni néhány összekötőkábelt; mint például a táplálással való összekötéshez, a logikai összeköttetésekhez és a szivattyúval való összekötéshez. Az alábbi ábrák a vezérlő és a szivattyú, valamint a vezérlő és a táphálózat között létesítendő összeköttetéseket vázolja fel. A fent említett összeköttetésekkel kapcsolatos további részletek és a transzformátor (mely opcióként áll rendelkezésre) műszaki jellemzői a "Technical Information" című mellékletben található.



Ábra 2 Vezérlő és szivattyú közötti összekötő kábel

Telepítés

VESZÉLY!



A vezérlő belsejében nagyfeszültségek alakulnak ki, melyek súlyos károkat vagy halált okozhatnak. A vezérlő bármely telepítési vagy karbantartási műveletének megkezdése előtt ki kell húzni a hálózati csatlakozót. A vezérlő belsejében magas hőmérséklet képződhet, mely súlyos károkat okozhat. A vezérlőt, végleges telepítéskor, megfelelően védeni kell, a véletlenszerű érintések elkerülése érdekében.

MEGJEGYZÉS

A végleges rendszerben telepített vezérlő elhelyezésének biztosítania kell a hűtőlevegő szabad áramlását a berendezés körül. A vezérlőt nem szabad légköri hatásoknak (eső, fagy, hó) kitett, poros, maró hatású gázt tartalmazó, robbanás- vagy fokozottan tűzveszélyes környezetben telepíteni és/vagy használni.

Üzem közben az alábbi környezeti feltételeket kell biztosítani:

- hőmérséklet: 0 °C és +40 °C között
- relatív nedvességtartalom: 0 – 95 % (nem lecsapódó).

Használat

Ebben a szakaszban a fontosabb működési eljárások kerülnek ismertetésre. További részletek és az összekötéseket vagy az opcionális elemeket érintő eljárások találhatóak a “Technical Information” melléklet “USE” című szakaszában. A vezérlő használata előtt valamennyi elektromos és pneumatikus összeköttetést létre kell hozni, és figyelembe kell venni a csatlakoztatott szivattyú kézikönyvét.

VESZÉLY!



Kérjük, beindítás előtt gondosan olvassa el a szóban forgó szivattyú kézikönyvét, a biztonságos használat érdekében.

Használati eljárások

A vezérlő bekapcsolása

A vezérlő bekapcsolásához elegendő számára a tápfeszültség biztosítása.

A szivattyú beindítása

A szivattyú beindításához a J4 csatlakozó 12-es pin-jét nyitva kell hagyni; ily módon, amikor a 24 Vdc feszültség a vezérlőre rákapcsolásra kerül, a szivattyú beindul.

A szivattyú leállítása

A szivattyú leállítása a J4 csatlakozó 12-es és 15-ös pin-jének rövidre zárásával történhet.

Karbantartás

A TwisTorr 74 FS sorozatú vezérlők semmiféle karbantartást sem igényelnek. Valamennyi beavatkozást meghatalmazott személyzet hajthat csak végre. Meghibásodás esetén igénybe vehető a Agilent javítási szolgáltatása, vagy a "Agilent advanced exchange service", mely a meghibásodott helyett, egy javított vezérlőt biztosít.

VESZÉLY!



A vezérlőn való bármely beavatkozás végrehajtása előtt ki kell húzni a hálózati csatlakozót.

Amikor egy vezérlőt le kell selejtezni, megsemmisítésekor a tárgykörbe tartozó nemzeti normáknak megfelelően kell eljárni.

Megsemmisítés

A címkén jelenlévő "WEEE" logo jelentése.

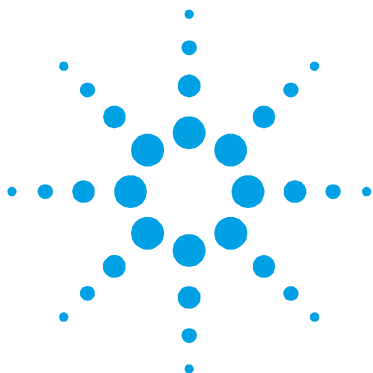
Az alább látható szimbólum az EK "WEEE" elnevezésű irányelvvel összhangban kerül alkalmazásra. Ez a szimbólum **(mely csak az Európai Közösség országaiban érvényes)**, azt jelzi, hogy a termék, melyen megtalálható, NEM kerülhet közöségi háztartási vagy ipari hulladékkal együtt megsemmisítésre, hanem azt egy szelektív hulladékgyűjtő rendszerbe kell továbbítani. A végfelhasználónak, a gyűjtési és megsemmisítési eljárás beindítása céljából, az adásvételi szerződés határidőinek és feltételeinek alapos áttanulmányozása után, fel kell vennie a kapcsolatot a berendezés szállítójával, legyen ez a gyártó vagy egy viszonteladó.



További információkért lásd:

<http://www.agilent.com/environment/product/index.shtml>

11 Felhasználói Kézikönyv
Megsemmisítés



12 Podrecznik Instrukcji

Informacje ogólne	124
Magazynowanie	125
Przygotowanie do instalacji	126
Instalacja	128
Użytkowanie	129
Procedury użytkowania	129
Zaswiecenie kontrolera	129
Uruchomienie pompy	129
Zatrzymanie pompy	129
Konserwacja	130
Przetworstwo odpadów	131

Tłumaczenie instrukcji oryginalnej



Informacje ogólne

Urządzenie to jest przeznaczone do użytku zawodowego. Przed przystąpieniem do korzystania użytkownik powinien wnikliwie przesledzić zarówno ten podrecznik zawierający instrukcje jak również każda inna dostarczona przez Agilent informację. Agilent uchyła się od jakiegokolwiek odpowiedzialności w przypadku całkowitego lub częściowego nie przestrzegania swoich instrukcji, przeznaczenia urządzenia niezgodnie z jego przeznaczeniem i jego obsługi przez personel uprzednio nie przeszkolony, naprawy urządzenia bez odpowiedniej autoryzacji lub przeznaczenia niezgodnego z obowiązującymi, specyficznymi normami krajowymi. Kontroler z serii TwisTorr 74 FS PCB jest przetwornikiem częstotliwości kontrolowanym przez mikroprocesor zrealizowanym z komponentów w stanie stałym i posiadającym zdolności samodiagnostyczne i samoochronne.

Kontroler ten pilotuje pompe z serii TwisTorr 74 FS (przy pomocy procesu podzielonego na dziesięć etapów) Podczas fazy rozruchu kontroluje napięcie i prąd elektryczny w stosunku do szybkości pompy. Do jego zadania należy połączenie całego obwodu niezbędne do prawidłowego automatycznego funkcjonowania pompy z serii TwisTorr 74 FS. Poprzez dodatkowy przetwornik są dostępne sterowniki sterujące uruchomieniem i zatrzymywaniem pompy na odległość, sygnały wskazujące stan operacyjny pompy, i sterowniki sterujące uruchomieniem i zatrzymywaniem pompy wstępnej próżni, sygnały blokujące (dla przerywaczy ciśnieniowych, przerywaczy kontrolujących przepływ wody itp). Kontroler powinien być zasilany napięciem stałym 24 Vdc.

W następujących paragrafach zawarte są wszystkie niezbędne informacje dotyczące gwarancji bezpieczeństwa personelu obsługującego urządzenie.

Bardziej szczegółowe informacje są zawarte w dodatku do "Technical Information".

Podrecznik ten opiera sie na nastepujacych konwencjach:

UWAGA!

Sygnaly wzrokowe ukazujace sie przed procedura jezeli zostana zlekcewazone moga spowodowac uszkodzenia urzadzenia.

ZAGROZENIE!



Sygnaly ostrzegajace o zagrozeniu powinny zwrocic uwage operatora na odpowiednia specyficzna procedure ktora jezeli nie jest przestrzegana w sposob prawidlowy moze przyczynic sie do powaznych obrazen osoby obslugujacej.

PRZYPIS

Przypisy zawieraja najwazniejsze informacje wyciagniete z tekstu.

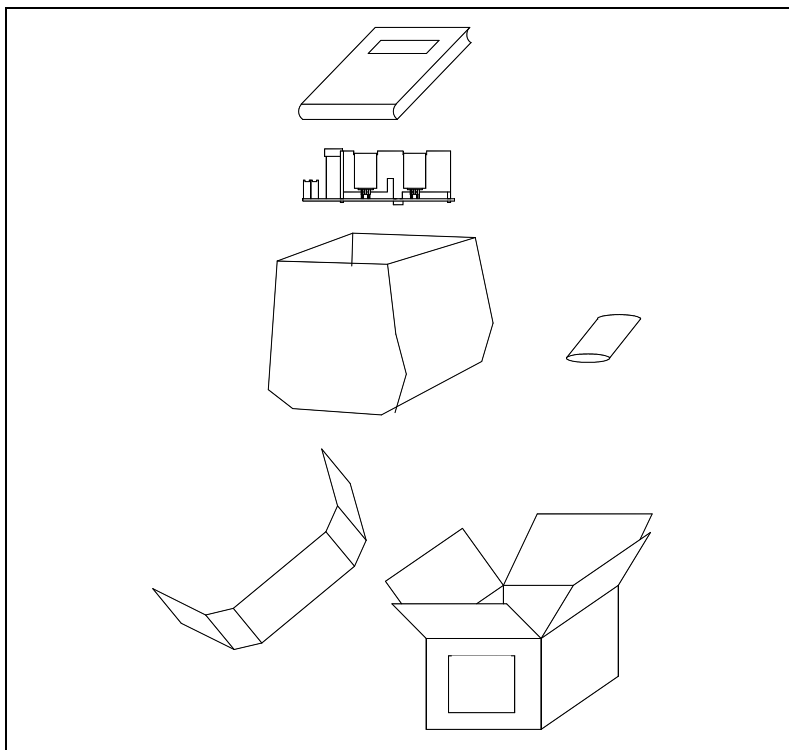
Magazynowanie

Kontrolery podczas transportu i przechowywania w magazynie powinny miec nastepujace warunki otoczenia:

- temperatura: od -20 °C do +70 °C
- wilgotnosc wzgledna: od 0 – 95 % (nie kondensujaca).

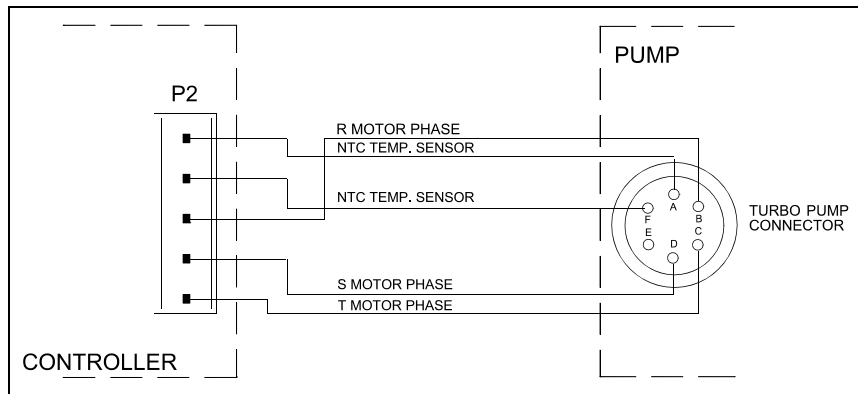
Przygotowanie do instalacji

Kontroler zostaje dostarczony w specjalnym, ochronnym opakowaniu; jeżeli opakowanie prezentuje jakiegokolwiek uszkodzenia które mogłyby powstać podczas transportu wówczas należy niezwłocznie powiadomić najbliższy, lokalny punkt sprzedaży.



Postać 1 Opakowanie kontrolera

Podczas operacji rozpakowywania należy zachować specjalną ostrożność, aby nie spowodować upadku kontrolera lub nie narazić na uderzenia. Nie wyrzucać opakowania, ponieważ jest ono całkowicie recyklowalne i odpowiada normom CEE 855/399 dotyczącym ochrony środowiska. Przed zainstalowaniem kontrolera należy przygotować niektóre kable służące do połączenia kilku elementów, takie jak na przykład kabel zasilający każdy z elementów połączonych z sobą, kable służące do wspólnych połączeń logicznych i te do połączenia z pompą. Schematy zamieszczone poniżej przedstawiają połączenia, które należy wykonać pomiędzy kontrolerem i pompą oraz pomiędzy kontrolerem i zasilaniem. Dla uzyskania bardziej szczegółowych wyjaśnień na temat tych połączeń, jak i danych technicznych transformatora (do dyspozycji jako opcja) należy konsultować dodatek do "Technical Information".



Postać 2 Kabel łączący kontroler z pompą

Instalacja

ZAGROZENIE!



Wewnątrz kontrolera powstają wysokie napięcia, które mogą spowodować poważne obrażenia lub śmiertelne wypadki. Przed wykonaniem jakiegokolwiek operacji instalacji lub konserwacji należy odłączyć kontroler od źródła zasilania. Wewnątrz kontrolera powstają również bardzo wysokie temperatury, które mogą spowodować poważne obrażenia. Podczas definitywnej instalacji kontrolera chronić go odpowiednio od przypadkowych kontaktów.

PRZYPIS

Kontroler zainstalowany w systemie definitywnie powinien być tak usytuowany, aby wokół niego mogło swobodnie krążyć powietrze chłodzące. Nie należy instalować i/lub użytkować kontroler w miejscach narażonych na działanie warunków atmosferycznych (deszcz, mroz, śnieg), pyłów, gazów agresywnych, w środowiskach wybuchowych lub o dużym zagrożeniu pożarowym.

Podczas pracy kontrolera należy zapewnić następujące warunki otoczenia:

- temperatura: od 0 °C do +40 °C
- wilgotność względna: 0 – 95 % (nie kondensująca).

Uzytkowanie

Paragraf ten zawiera podstawowe procedury operatywne. Dla uzyskania dodatkowych bardziej szczegolowych informacji lub dostepu do procedur na temat polaczen lub akcesoriow skonsultowac paragraf "USE" znajdujacy sie w dodatku do "Technical Information". Przed przystapieniem do uzytkowania kontrolera wykonac wszystkie polaczenia elektryczne i pneumatyczne podlaczenie pompy konsultowac z odnosa instrukcja.

ZAGROZENIE!



Aby obslugiwac pompe w bezpieczny sposob, nalezy zapoznac sie z odpowiednia instrukcja obslugi.

Procedury uzytkowania

Zaswiecenie kontrolera

Aby zaswiecic kontroler wystarczy dostarczyc do niego napiecie zasilania.

Uruchomienie pompy

Zeby uruchomic pompe nalezy zostawic otwarty pin 12 lacznika J4 w ten sposob dostarczone kontrolerowi napiecie 24 Vdc pozwoli na uruchomienie pompy.

Zatrzymanie pompy

Do zatrzymania pompy nalezy zewrzec pin 12 i 15 lacznika J4.

Konservacja

Kontroler z serii TwisTorr 74 FS nie wymaga żadnej konserwacji. Jakakolwiek interwencja musi być przeprowadzona wyłącznie przez osoby autoryzowane.

.W przypadku uszkodzenia możliwe jest korzystanie z serwisów naprawczych Agilent lub “Agilent advance exchange service”, w którym można otrzymać kontroler odnowiony zastępując go tym uszkodzonym.

ZAGROZENIE!



Przed wykonaniem jakiegokolwiek interwencji na kontrolerze należy odłączyć kabel zasilający.

W przypadku decyzji złomowania należy przeprowadzić operacje respektując krajowe, specyficzne normy.

Przetworstwo odpadów

Objasnienie znajdującego się na etykiecie znaczenia logo “WEEE”.

Użyty poniżej symbol jest zgodny z wymogiem zarządzenia “WEEE” Unii Europejskiej. Symbol ten (**prawomocny tylko w krajach Unii Europejskiej**) oznacza, że wyrob który nim został oznaczony NIE może być przetworzony jako odpad razem z innymi domowymi lub przemysłowymi natomiast musi być składowany w miejscu przeznaczonym dla odpadów do przerobki zroznicowanej .

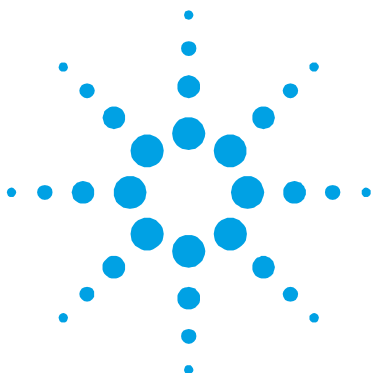
Dlatego też poleca się użytkownikowi, po uprzedniej weryfikacji terminu i warunków zawartych w kontrakcie sprzedaży, nawiązać kontakt z dostawcą lub sprzedawcą urządzenia w celu uruchomienia procesu zbiórki i przerobu.



Aby uzyskać więcej informacji, patrz:

<http://www.agilent.com/environment/product/index.shtml>

12 Podrecznik Instrukcji
Przetworstwo odpadów



13 Návod k Použití

Všeobecné informace	134
Uskladnění	135
Příprava k instalaci	136
Instalace	138
Použití	139
Používané procedury	139
Zapalování Kontrolérů	139
Spuštění čerpadla	139
Zastavení čerpadla	139
Údržba	140
Likvidace	141

Překlad originálního návodu



Všeobecné informace

Tento přístroj je určen pro profesionální potřebu. Uživatel si musí před použitím tuto příručku velmi pozorně přečíst, ale i každou další aktualizaci, kterou podnik bude předávat. V důsledku nedostatečného i částečně nesprávného použití, nebo v případě chyb nevyškoleného personálu či neoprávněných zásahů a neodpovídajících specifickým státním normám, Agilent nepřijímá a nenese žádnou zodpovědnost.

Kontrolér TwisTorr 74 FS PCB je frekvenčním konvertorem, je kontrolovaný mikroprocesorem a skládá se z pevných komponentů s autodiagnostickou a samoochrannou schopností.

Obsahuje veškeré obvody nutné pro automatický provoz čerpadel série TwisTorr 74 FS. Kontrolér řídí čerpadla série TwisTorr 74 FS (s procesem rozděleným na deset kroků), během spouštěcí fáze kontrolují napětí a proud ve vztahu k rychlosti, kterou čerpadlo dosáhne. Dálkové příkazy ke spuštění a zastavení čerpadla jsou vykonávány prostřednictvím výpomocného konektoru, signály, které poukazují na operativní stav čerpadla, příkazy ke spuštění a zastavení pre-prázdného čerpadla, blokovací signály (pro tlakové vypínače, kontrolní vypínače vodního toku, atd.). Kontrolér musí být napájený stejnosměrným 24 Vdc napětím.

V následujících odstavcích jsou uvedeny veškeré nutné informace, týkající se operátora a bezpečnostních podmínek během provozu přístroje. Detailní informace se nacházejí v dodatku "Technical Information".

Tato příručka používá následující běžná označení:

POZOR!

Tyto zprávy bývají vizualizovány před jakoukoliv procedurou, je třeba se jimi řídit, aby nedošlo k poškození přístroje.

NEBEZPEČÍ!



Odkazy, které upozorňují na nebezpečí a donutí operátora, aby se řídil dle procedury nebo specifické praktiky, která nebude-li řádně vykonána, může ohrozit vážným způsobem jeho zdraví.

POZNÁMKA

Poznámky, vyjmuté z textu obsahují velmi důležité informace.

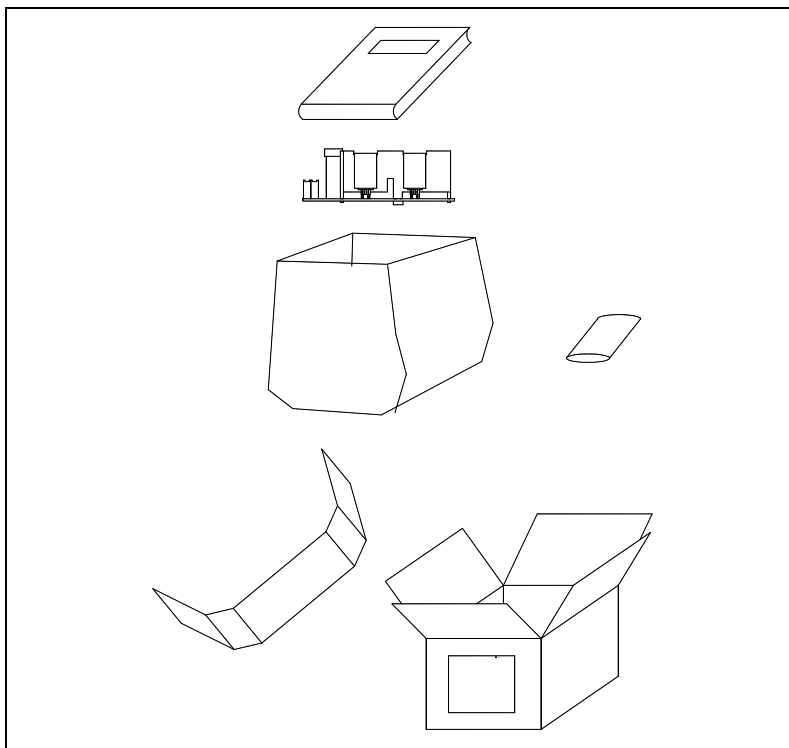
Uskladnění

Během dopravy a skladování kontrolérů, se musí dbát na následující podmínky prostředí:

- teplota: od -20 °C do +70 °C
- relativní vlhkost: 0 – 95 % (nesmí kondenzovat).

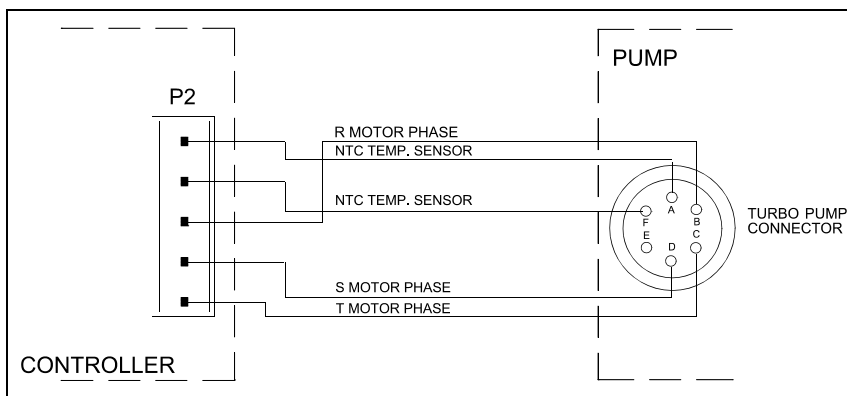
Příprava k instalaci

Kontrolér je dodáván v speciálním ochranném balení; objeví-li se následkem dopravy nějaké poškození, je třeba se okamžitě obrátit na místní prodejní servis.



Postava 1 Obal kontroléru

Během rozbalování se obzvláště musí dávat pozor, aby kontrolér nespadol nebo aby neutrpěl nárazy. Balicí materiál se nesmí odkládat v přírodě. Tento materiál se může kompletně recyklovat, neboť se shoduje se směrnicemi EU, CEE 85/399, které se týkají životního prostředí. Před instalací kontroléru je třeba připravit kabely na vzájemná propojení, jako například vzájemné propojení s napájením, vzájemná logická propojení a kabely na čerpadlové spojení. Následující obrázky zobrazují potřebné propojení mezi kontrolérem a čerpadlem a mezi kontrolérem a napáječem. Pro další detaily výše uvedených vzájemných spojení a technických charakteristik transformátoru (k dispozici jako opce) viz dodatek "Technical Information".



Postava 2 Propojovací kabel mezi Kontrolérem a čerpadlem

Instalace

NEBEZPEČÍ!



Uvnitř kontroléru dochází k vysokým napětím, která jsou velmi nebezpečná i smrtící. Před provedením jakékoliv instalační či údržbové operace s kontrolérem, je třeba ho odpojit od elektrického příkonu. Uvnitř kontroléru vznikají velmi vysoké teploty, které mohou být velice škodlivé. Kontrolér je třeba adekvátně ochránit proti nahodilým kontaktům.

POZNÁMKA

Kontrolér nainstalovaný v definitivním systému musí být umístěn tak, aby kolem přístroje mohl volně proudit ochlazovací vzduch. Kontroléry se nesmí instalovat v místech vystavených atmosferickým vlivům (déšť, mráz, sníh), prachu, agresivním plynům, na výbušných místech nebo na místech kde existuje velké riziko požárů.

Během provozu se musí dbát na následující podmínky prostředí:

- teplota: od 0 °C do +40 °C
- relativní vlhkost: 0 – 95 % (nesmí kondenzovat).

Použití

V tomto odstavci jsou uvedeny hlavní operativní postupy. Pro další detaily i pro procedury týkající se spojení nebo dalších opčních detailů, je třeba se obracet na odstavec "USE" v příloze "Technical Information". Před použitím kontroléru se musí vykonat veškerá elektrická a pneumatická spojení, při čemž je třeba se obracet na příslušnou příručku napojeného čerpadla.

NEBEZPEČÍ!



Před spuštěním se pro bezpečný provoz obraťte na příslušný manuál k pumpě.

Používané procedury

Zapalování Kontrolérů

Zapálení kontrolérů se děje tak, že se přivede elektrické napětí do kontroléru.

Spuštění čerpadla

Pro spuštění čerpadla je třeba nechat otevřený pin 12 konektoru J4 tak, aby se, až kontrolér dostane napětí 24 Vdc, čerpadlo spustilo.

Zastavení čerpadla

Na zastavení čerpadla je třeba uskutečnit zkrat pinu 12 a 15 konektoru J4.

Údržba

Kontrolér série TwisTorr 74 FS nevyžaduje žádnou údržbu. Jakýkoliv zásah musí být vykonáván jen kompetentním personálem. V případě poruchy se může využít opravářského servisu Agilent nebo "Agilent advanced exchange service", kde poškozený kontrolér bude vyměněn za nový.

NEBEZPEČÍ!



Nejdříve, před jakýmkoliv zásahem na kontroléru, se musí odpojit elektrický kabel.

Budeli zapotřebí kontrolér likvidovat, musí se postupovat dle platných specifických státních norem.

Likvidace

Význam loga "WEEE" **nacházejícího se na štítku**. Níže uvedený symbol odpovídá směrnici CE pojmenovaným "WEEE". Tento symbol (**platný jen pro státy Evropské Unie**) určuje, že výrobek, který je takto označen **NESMÍ** být likvidován společně s ostatními domácími nebo průmyslovými odpady, ale je nutno ho předat do patřičných provozních sběrů, kde musí být tříděn a likvidován odděleně od městského odpadu. Uživateli, který má výrobek likvidovat doporučujeme, aby se nakontaktoval přímo s výrobcem nebo prodejcem, který se po patřičném prověření termínů a smluvních podmínek postará o kompletní likvidaci uvedeného výrobku.

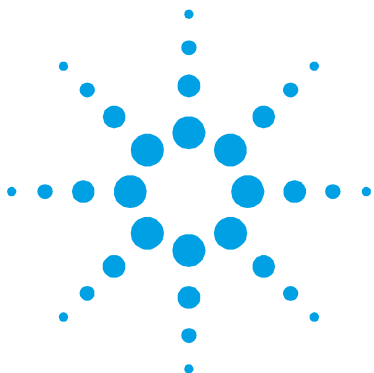


Více informací naleznete na:

<http://www.agilent.com/environment/product/index.shtml>

13 **Návod k Použití**

Likvidace



14 Návod na Obsluhu

Všeobecné informácie	144
Uskladňovanie	145
Príprava pre inštaláciu	146
Inštalácia	148
Použitie	149
Postup pri použití	149
Zapnutie Controller	149
Spustenie čerpadla	149
Zastavenie čerpadla	149
Údržba	150
Likvidácia	151

Preklad originálneho návodu



Všeobecné informácie

Tento prístroj slúži k profesionálnemu použitiu. Užívateľ si pred použitím tohto prístroja musí pozorne prečítať tento návod k použitiu a jeho ďalšie doplnky, dodané podnikom Agilent. Podnik Agilent nepreberá na seba žiadnu zodpovednosť za úplné, alebo čiastočné nedodržiavanie návodu, za nesprávne používanie prístroja nekompetentnými osobami, za nepovolené zásahy a za používanie, ktoré nie je v zhode s príslušnými národnými normami.

“Controllers” série TwisTorr 74 FS PCB je menič kmitočtu, kontrolovaný mikroprocesorom, vyrobený prvkami pevnej fázy a so schopnosťou samodiagnostiky a samokontroly.

Obsahuje všetky obvody, potrebné pre automatickú prevádzku čerpadiel série TwisTorr 74 FS. “Controllers” riadi čerpadlá série TwisTorr 74 FS (s procesom, rozdeleným do desať rozpätí) počas štartovacej fázy, kontrolujú napätie a prúd vo vzťahu s rýchlosťou, dosiahnutou čerpadlom. Pomocná spojka umožní diaľkové povely pre štart a vypnutie čerpadla, signály, ktoré ukazujú operatívny stav čerpadla, povely pre štart a vypnutie pred-prázdňového čerpadla blokacie signály (pre tlakové tlačidlá, tlačidlá pre kontrolu toku vody, atď.). Controller musí byť napájaný rovnakosmerným 24 Vdc napätím.

Ďalej sú popísané všetky potrebné informácie, ktorých úlohou je zabezpečiť bezpečnosť pracovníka počas používania prístroja.

Podrobné informácie sú uvedené v dodatku "Technical Information".

Tento návod používa nasledujúce signály:

POZOR!

Signály “pozor” sú pred tými postupmi, ktoré, v prípade ich nedodržania, môžu poškodiť prístroj.

NEBEZPEČÍ!



Signál nebezpečie upozorňuje pracovníka na určitý pracovný postup, ktorý, keď nie je prevedený správne, môže zapríčiniť vážne osobné poranenia.

POZNÁMKA

Poznámky, vybrané z textu obsahujú veľmi dôležité informácie.

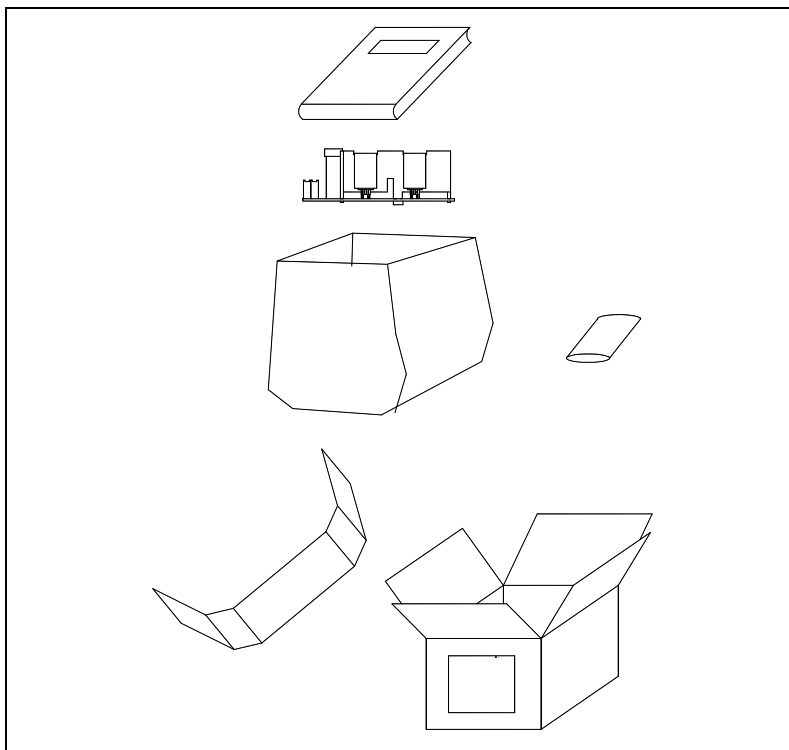
Uskladňovanie

Počas prepravy a uskladňovania “Controllers” musia byť dodržané tieto podmienky prostredia:

- teplota: od -20 °C do +70 °C
- relatívna vlhkosť: 0 – 95 % (bez kondenzátu).

Príprava pre inštaláciu

“Controller” je dodaný v špeciálnom ochrannom obale; keby ste objavili škodu, spôsobenú prepravou, skontaktujte sa s kanceláriou miestneho predaja.

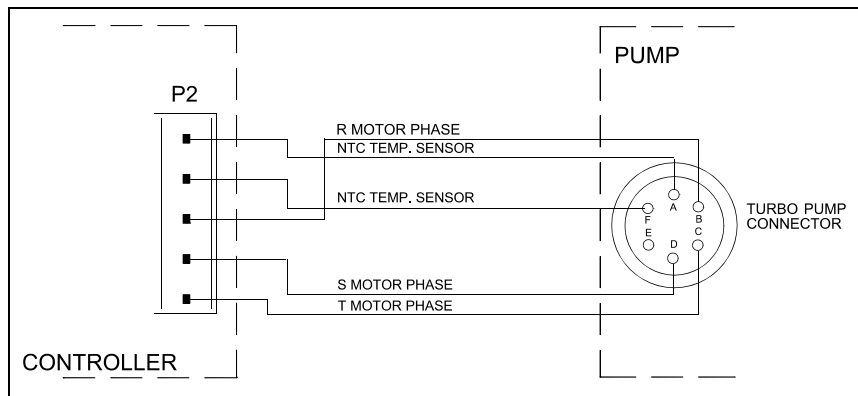


Postava 1 Obal controller

Pri rozbal'ovaní dávajte pozor, aby vám "Controller" nespadol, alebo niekde nenarazil. Obal nezhadzujte. Je vyrobený z recyklovateľného materiálu a zodpovedá smernici CEE 85/399 pre ochranu životného prostredia.

Pred inštaláciou controller je treba pripraviť káble na vzájomné prepojenie, ako napríklad vzájomné prepojenie s napájaním, vzájomné logické prepojenia a káble na čerpadlové spojenie. Nasledujúce obrázky zobrazujú potrebné prepojenia medzi controllerom a čerpadlom a medzi controllerom a napájaním.

Pre ďalšie detaily, týkajúce sa vzájomných spojení a technických charakteristík transformátora (k dispozícii ako opcia) viď dodatok "Technical Information".



Postava 2 Kábel prepojenia medzi Controllerom a čerpadlom

Inštalácia

NEBEZPEČÍ!



Vo vnútri controller dochádza k vysokému napätiu, ktoré môže spôsobiť veľké škody alebo smrť. Pred prevedením akékoľvek inštalácie alebo údržby controller, je treba tento odpojiť od elektrického príkonu. Vo vnútri controller vznikajú vysoké teploty, ktoré môžu byť veľmi škodlivé. Controller je treba adekvátne ochrániť voči náhodným kontaktom.

POZNÁMKA

Controller nainštalovaný v definitívnom systéme musí byť umiestnený tak, aby okolo prístroja mohol voľne prúdiť chladný vzduch. Controllers sa nesmú inštalovať v miestach, vystavených atmosferickým vplyvom (dážď, mráz, sneh), prachu, agresívnym plynom, na výbušných miestach, alebo na miestach, kde existuje veľké riziko požiarov.

Počas chodu je dôležité, aby boli rešpektované tieto podmienky prostredia:

- teplota: od 0 °C do +40 °C
- relatívna vlhkosť: 0 – 95 % (nekondenzovaná).

Použitie

V tomto odstavci sú uvedené hlavné operatívne postupy. Pre ostatné detaily a pre postupy, ktoré sa vzťahujú na napojenia, alebo na zvolené príslušenstvo, postupujte podľa časti "USE" v dodatku "Technical Information". Všetky elektrické a pneumatické napojenia treba previesť pre použitím controller. Pritom treba konzultovať príslušnú príručku napojeného čerpadla.

NEBEZPEČÍ!



Pred začiatkom manipulácie si z dôvodu zaistenia bezpečného použitia dôkladne preštudujte priložený manuál.

Postup pri použití

Zapnutie Controller

Na zapnutie controller stačí priviesť elektrické napätie do controller.

Spustenie čerpadla

Pre spustenie čerpadla nechajte otvorený pin 12 spojky J4 tak, že keď controller dostane napätie 24 Vdc, čerpadlo sa spustí.

Zastavenie čerpadla

Na zastavenie čerpadla treba skratovať pin 12 a 15 spojky J4.

Údržba

“Controllers” série TwisTorr 74 FS nevyžadujú žiadnu údržbu. Akýkoľvek zásah musí previesť poverený personál. V prípade vady je možné využiť opravnú službu Agilent alebo “Agilent advanced exchange service”, ktorá vymení vadný controller.

NEBEZPEČÍ!



Pred akýmkoľvek zásahom na “controller” vytiahnite šnúru z elektrickej siete.

V prípade, že “controller” musí ísť do zberu, riadte sa platnými príslušnými národnými normami.

Likvidácia

Význam loga "WEEE" nachádzajúceho sa na štítkoch. Aplikovanie doluo značeného symbolu dodržiava smernicu EÚ s názvom "WEEE". Tento symbol (**platný iba pre štáty Európskej Únie**) znamená, že výrobok s týmto štítkom NEMIE byť odstránený spolu s bežným domácim alebo priemyselným odpadom, ale sa musí odstrániť ako delený odpad. Vyzývame preto konečného užívateľa, aby sa skontaktoval s dodávateľom prístroja, či už je to výrobca alebo predajca za účelom jeho likvidácie podľa zmluvných podmienok predaja.

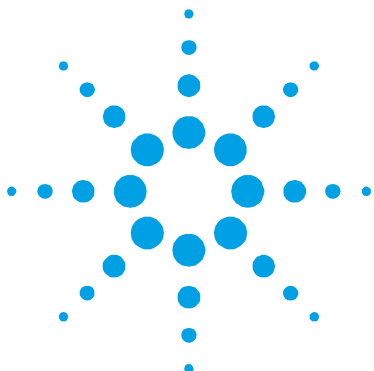


Viac informácií nájdete na:

<http://www.agilent.com/environment/product/index.shtml>

14 **Návod na Obsluhu**

Likvidácia



15 Priročnik za Navodila

Splošna navodila	154
Skladiščenje	155
Priprava za montažo	156
Montaža	158
Uporaba	159
Postopki za uporabo	159
Vkllopitev Controllera	159
Zagon Črpalke	159
Zaustavitev Črpalke	159
Vzdrževanje	160
Odlaganje odpadkov	161

Prevod navodil v izvirniku



Splošna navodila

Naprava je namenjena samo za strokovno rabo. Uporabnik mora temeljito prebrati priročnik in pri tem upoštevati vsako dodatno informacijo, ki jo nudi Agilent. Agilent ne odgovarja v slučaju celotnega ali delnega nespoštovanja navodil, pri nepravilni uporabi s strani nestrokovnega osebja, v slučaju nedovoljenih posegov ali pri zanemarjenju specifičnih nacionalnih meril.

Controller serije TwisTorr 74 FS PCB je frekvenčni pretvornik, na katerega nadzor ima mikrokrmilnik, sestavljen iz solidnih delov za katere so značilne lastnosti avtodiagnoze in avtozaščite. Controller vsebuje celotno električno vezje, ki je potrebno za avtomatično delovanje črpalke serije TwisTorr 74 FS. Controller krmi črpalke serije TwisTorr 74 FS (s 10-faznim postopkom) med pogonom in nadzoruje napetost ter električni tok v sorazmerju s hitrostjo, ki jo doseže črpalka.

Preko pomočnih konektorjev so na razpolago komande za pogon in zaustavitev črpalke (ki se regulira s daljinskim upravljanjem), komande ki pokažejo operativno stanje črpalke, komande za zagon in zaustavitev črpalke predhodnega črpanja, komande za blokiranje (za stikala na pritisk, nadzorna stikala vodnega toka, itn.). Controlleru dovodite trajno napetost od 24 Vdc. V naslednjih odstavkih so navedene vse informacije potrebne za zagotavljanje sigurnosti uporabnika med rokovanjem s to napravo. Podrobnejše informacije dobite v priponki "Technical Information".

V tem priročniku so varnostne informacije razvrščene v dva razreda:

POZOR!

Znak 'pozor' se pojavi pred postopke, ki če jih uporabnik zanemari, lahko napravo poškoduje.

SVARILO!



Znak svarila opozarja uporabnika, da mora pri določenem postopku ali pri posebnem delovanju paziti. Kršitev svarila lahko privede do lažjih ali hudih telesnih poškodb.

OPOMBA

V opombah so vsebovane vse najbolj pomembne informacije iz besedila.

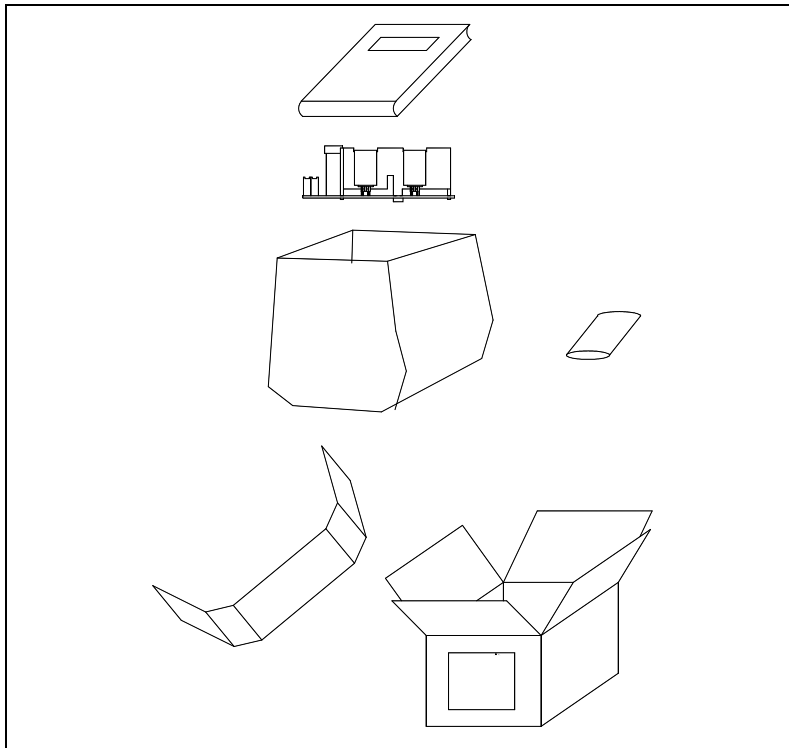
Skladiščenje

Med prevozom in skladiščenjem controllera morate zagotoviti naslednje pogoje okolice:

- temperatura: od -20 °C do +70 °C
- relativna vlaga 0 – 95 % (ne-kondenzna)

Priprava za montažo

Controller dobite v posebni zaščitni embalaži; če so na embalaži znaki poškodbe, ki naj bi nastale med transportom obrnite se krejavnemu prodajalnemu biroju.

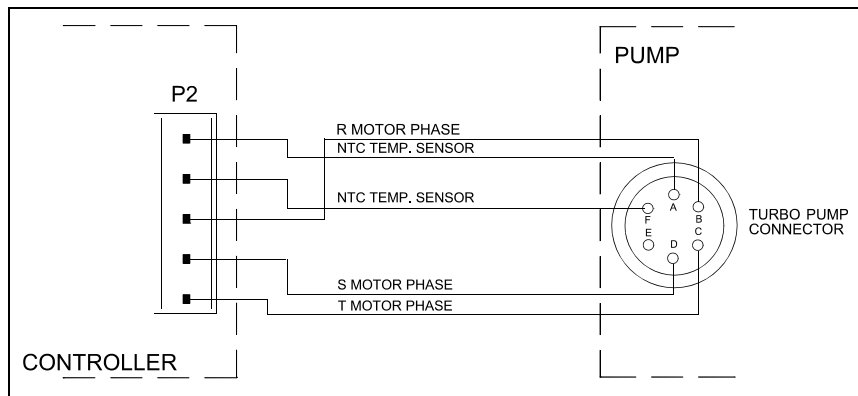


Slika 1 Embalaža controllera

Pri odstranitvi embalaže morate še posebno paziti, da controller ne pade na tla ali da ne podlegne udarcem. Ne zavržite embalaže v okolico. Material lahko v celoti recikirate in ustreza direktivi CEE 85/399 za zaščito okolja. Preden bi vgradili controller, morate pripraviti nekaj interkonektorskih kablov, kot na primer interkonektorske kable za napoj, interkonektorske logične kable in kabel za priključitev črpalke.

Naslednje slike prikazujejo priključke in vezave, ki jih morate izpeljati med controllerjem in črpalko ter med controllerjem in napajalnikom.

Za podrobnejše informacije o raznih prikljopitvah in o tehničnih značilnosti transformatorja (mogoč kot opcija) si oglejte priponko "Technical Information".



Slika 2 Interkonektorski kabel med Controllerjem in črpalko

Montaža

SVARILO!



Znotraj controllera se razvija visoka napetost, ki lahko resno poškoduje ali celo povzroči smrt. Preden bi nadaljevali s katerokoli dejavnostjo montaže ali vzdrževanja, izklopite controller iz vtičnice. V notranjosti controllera se razvijajo visoke temperature, ki lahko povzročijo hude telesne poškodbe.

OPOMBA

Ko controller vgradite v dokončni sistem, mora imeti taki položaj, da lahko zrak za hlajenje prosto kroži skozi napravo. Controller ne smete vgraditi in/ali uporabiti v okolih, ki so ogrožene s klimatskimi dejavniki (dež, led, sneg), ki so izpostavljene prahu, agresivnim plinom ali v okolih, kjer obstaja nevarnost vnetja ali požara.

Med delovanjem controllera morate zagotoviti naslednje pogoje okolice:

- temperatura: od 0 °C do +40 °C
- relativna vlaga: 0 – 95 % (ne-kondenzna).

Uporaba

V tem odsatvku so navedeni glavni operativni postopki. Za podrobnejše informacije in postopke, ki se nanašajo na priključitve ali posamezne opcije ogledajte si paragraf "USE" v dodatku "Technical Information".

Predno vkločite controller, morate izvesti vse električne in pnevmatske priključke; pri tem se nanajšate na priročnik priložene črpalke.

SVARILO!



Pred zagonom si pogledajte ustrezen priročnik za varno upravljanje in delovanje črpalke.

Postopki za uporabo

Vklopitev Controllera

Controllera vklopite tako, da controlleru priključite napajalno napetost.

Zagon Črpalke

Za zagon črpalke mora biti pin 12 konektorja J4 odprt, tako da ko Controller dobi napetost od 24 Vdc lahko začne zagon črpalke.

Zaustavitev Črpalke

Za zaustavitev črpalke morata biti pin 12 in 15 konektorja J4 v kratkem stiku.

Vzdrževanje

Controllerji serije TwisTorr 74 FS ne potrebujejo vzdrževanja. Samo pooblaščen osebje lahko rokuje s to napravo. V slučaju okvare se lahko obrnite servisu Agilent ali "Agilent advance exchange service", ki Vam stavlja na razpolago popravljeni controller.

SVARILO!

Pred katerikoli posegom, izključite napajalni kabel iz controllera.



Controller zavrzite kot industrijski odpadek v skladu z državnimi predpisi.

Odlaganje odpadkov

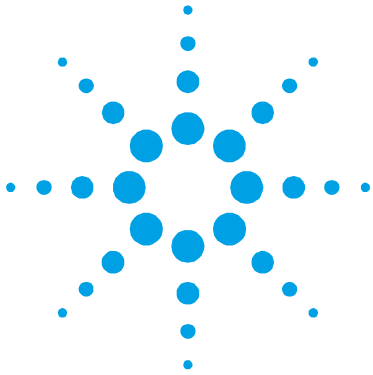
Pomen znamke "WEEE" na etiketah. Spodaj navedeni simbol je v skladu z direktivo ES znano pod imenom "WEEE". Ta simbol (**ki velja samo v državah Evropske Skupnosti**) pomeni, da izdelek NE SMETE ODSTRANITI skupaj z ostalimi komunalnimi ali pa industrijskimi odpadki, temveč morate poskrbeti za njihovo primerno ločevanje. Zato pozivamo uporabnike, da se ali pri prodajalnem centru ali pa pri prodajalcu seznanijo o postopku ločevanja in odstranitve odpadkov, šele nato, ko se je seznanil s pogoji in z merili kupopordajne pogodbe.



Za dodatne informacije se obrnite na:

<http://www.agilent.com/environment/product/index.shtml>

15 Priročnik za Navodila
Odlaganje odpadkov



16 Instructions for Use

General Information	164
Storage	165
Preparation prior to Installation	166
Installation	168
Use	169
Operating Procedures	169
Powering on the Controller	169
Starting the Pump	169
Stopping the Pump	169
Maintenance	170
Disposal	171

Original Instructions



General Information

This equipment is intended for professional use. Prior to using the equipment, the operator must carefully read this instruction manual and any other additional information provided by Agilent. Agilent declines all responsibility which may become as the result of the total or partial misinterpretation of the information provided in the documentation, by the improper use of the equipment by untrained personnel or by the use of the equipment by unauthorized personnel or not in compliance with specific national norms.

The TwisTorr 74 FS PCB controller is a microprocessor-controlled frequency converter made up of solid state components and capable of autodiagnostic and auto-protection functionality. It incorporates the circuitry needed for the automatic operation of TwisTorr 74 FS series pumps.

The TwisTorr 74 FS series pumps are driven by the controller (in a 10-step procedure) during the start-up phase that also checks the voltage and current according to the speed reached by the pump. Auxiliary connectors render the commands needed to start and stop the pump from remote, the signals that indicate the operating state of the pump and the locking commands (for pressure switches, water flow control switches, etc.) available. The controller must be provided with 24 Vdc.

In the next sections you will find all the information needed to guarantee operator safety during the use of the equipment. Detailed information is provided in the Appendix entitled "Technical Information".

The following conventions are adopted in this manual:

CAUTION!

Caution messages are provided before a procedure that could damage the equipment is not executed in the correct way.

WARNING!



Warning messages call the operator's attention to a specific procedure or operation that could cause serious personal injury if not executed in the correct way.

NOTE

Notes provide important information extracted from the text.

Storage

The controller must be transported or stored within the following environmental conditions:

- Temperature: -20 °C to +70 °C
- Relative humidity 0 to 95 % (non-condensing)

Preparation prior to Installation

The controller comes in a special protective packaging; if you notice any sign of damage that could have been caused during transportation, notify our local sales office immediately.

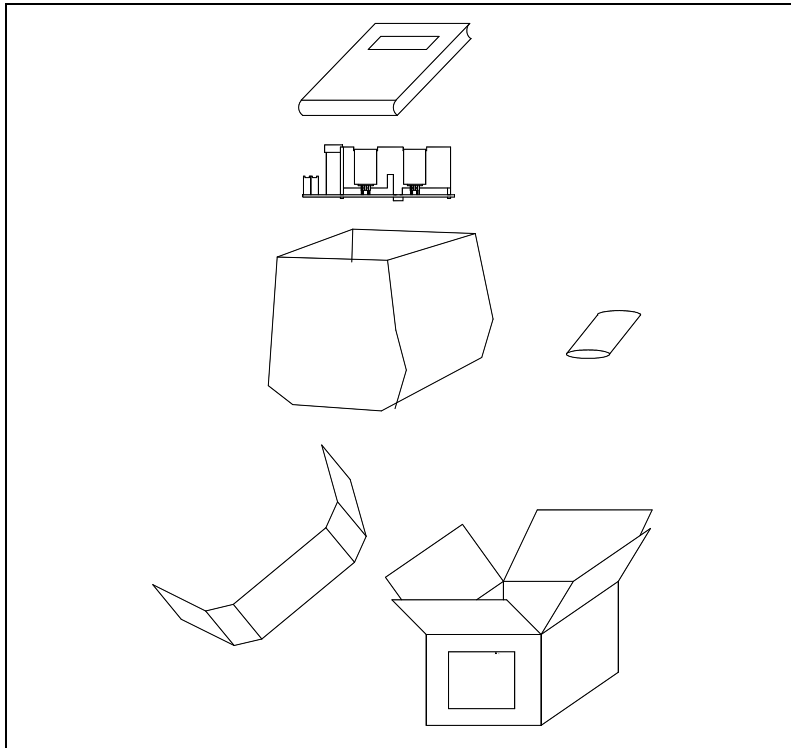


Figure 1 Controller Packaging

During the unpacking procedure, be careful to avoid dropping the controller or hitting it against anything. Do not discard the packaging material in the environment. The packaging material is completely recyclable and complies with the EEC environmental protection directive 85/399. Before installing the controller you need to prepare certain interconnection cables such as, for example, the power supply interconnection cables, the logic interconnection cables and the pump interconnection cable. The following figures provide the connection diagrams between the controller and pump and between the controller and power supply. For further information on these interconnections and for the technical characteristics of the transformer (available as option), refer to the Appendix entitled "Technical Information".

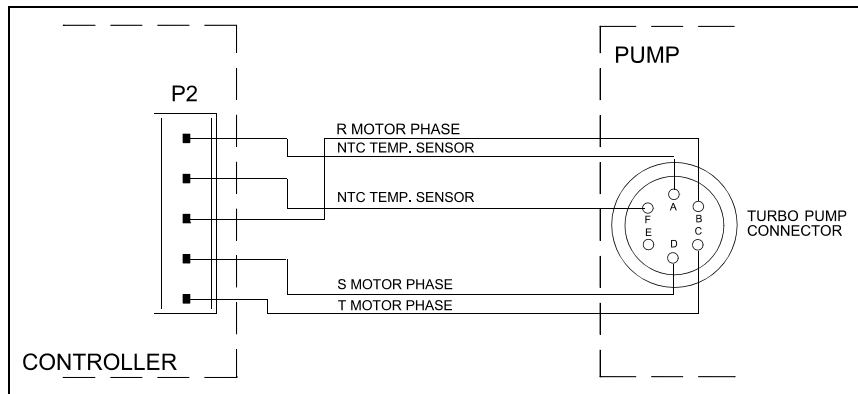


Figure 2 Controller – Pump Interconnection Cable

Installation

WARNING!



The controller develops internal voltages that could cause serious injury. Unplug the controller from the electrical wall outlet before beginning with any type of installation or maintenance procedure. In addition, the controller develops internal temperatures that could also cause serious injury. Adequately protect the controller in its final point of installation so as to avoid accidental contacts.

NOTE

When definitively installed in the system, the controller must be positioned so that the cooling air can circulate freely around the equipment. Do not install and/or use the controller in environments exposed to the atmosphere (rain, ice, and snow), to dust, to aggressive gases and in explosive areas where there is a high risk of fire.

The controller must be used within the following environmental conditions:

- Temperature: 0 °C to +40 °C
- Relative humidity: 0 to 95 % (non-condensing).

Use

This section provides all the major operating procedures. For further details and for the procedures that involve connections or specific options, refer to the section entitled "USE" in the Appendix "Technical Information".

Perform all electrical and pneumatic connections before using the controller, with reference to the manual of the connected pump.

WARNING!

Refer to the relevant pump manual to operate it in a safe way before starting.



Operating Procedures

Powering on the Controller

To power on the controller, simply provide electrical power to it.

Starting the Pump

To start the pump, leave pin 12 of connector J4 open so that the pump starts-up when 24 Vdc is provided to the controller.

Stopping the Pump

To stop the pump, short-circuit pins 12 and 15 of connector J4.

Maintenance

The TwisTorr 74 FS series controllers do not require any form of maintenance. Any intervention must be performed by authorized personnel.

In case of failure you can use the Agilent customer support service or the "Agilent advance exchange service", the latter consisting of replacing a faulty controller with a refurbished one.

WARNING!



Unplug the power cord from the electrical socket before performing any type of intervention on the controller.

If a controller needs to be scraped, proceed to do so in compliance with the specific national laws.

Disposal

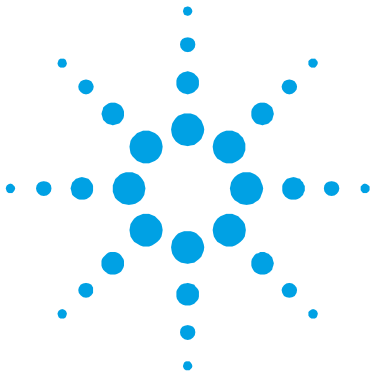
Meaning of the "WEEE" logo found in labels. The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive. This symbol (**valid only in countries of the European Community**) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system. The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale



For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

16 Instructions for Use
Disposal



17 Technical Information

General Description	174
TwisTorr 74 FS PCB 24 Vdc Controller Description	174
Controller Specifications	175
Controller Connector and Jumper	176
Jumper Configurations	177
Run-up Time Adjustment	177
Controller Outline	178
Installation	179
Data Exchange Protocols	184
RS-232 Protocol 1 Communication Description	184
Communication Protocol	185
Error Code Table	190
RS-232 Protocol 2 Communication Descriptions	190
Transmission Channel Characteristics	190
Message Structure	191
Example	194
Operation	199
Low Speed Activation/Deactivation	200
Maintenance	201
Accessories and Spare Parts	202



General Description

The TwisTorr 74 FS PCB controller is a microprocessor-controlled, solid-state, frequency converter with self-diagnosis and protection features.

TwisTorr 74 FS PCB 24 Vdc Controller Description

The controller is driven by a single chip microcomputer and is composed of a PCB that includes all the circuitry necessary for its operation. The microcomputer generates the variable output voltage according to the software and the gas load condition of the pump. Moreover, it manages signals from sensors, input/output connection information, and gives output for a fully automatic operation. The controller can be operated via remote signals through an RS-232 connection. The controller can be operated in local mode through suitable switches connected between the input pins of J4 connector and ground (see section J4 connector for a detailed description of the input signal present on J4 connector).

Controller Specifications

Tab. 1

Input:	
- Voltage	20.4 to 32.2 Vdc with 2 Vpp max ripple
- Power	128 W max.
Output:	
- Voltage	42 Vac nominal $\pm 10\%$, 3-phase
- Frequency	1167 Hz, $\pm 2\%$
- Power	70 W maximum
Fuse	T 4A
Operating temperature	0 °C to + 40 °C
Storage temperature	-20 °C to + 70 °C
Weight	0.5 Kg (1.1 lbs)

Controller Connector and Jumper

The following figure shows the controller and indicates its connectors and jumpers.

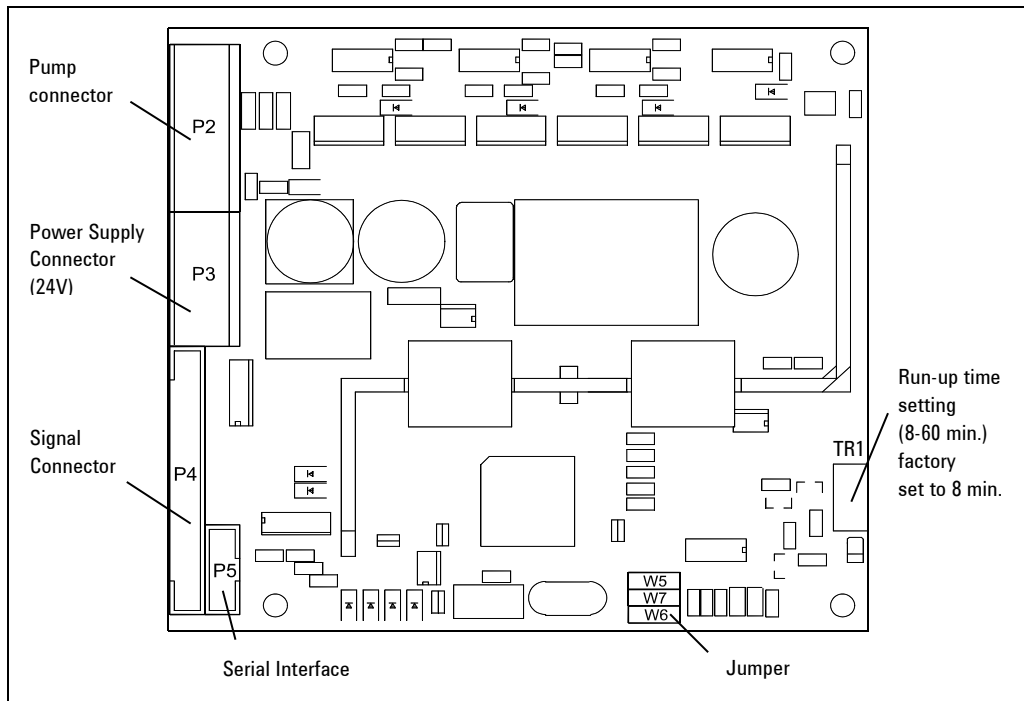


Figure 3 Connector and Jumper

Controller connector are named with prefix Px; in the following pages the connector counterpart are named as Jx to indicate the use of the appropriate connector housing and pins.

Jumper Configurations

The following figures show the configuration and meaning of the individual jumpers on the board.

Jumper W7 is used to select the data exchange rate, as shown in the figure.

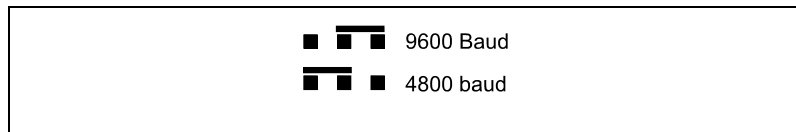


Figure 4 W7 Jumper

Jumper W6 is used to set the Soft Start, as shown in the figure.

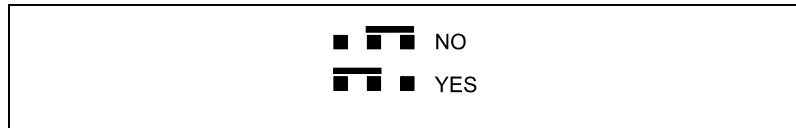


Figure 5 W6 Jumper

The W5 jumper is just for factory setting and must not be moved.

Run-up Time Adjustment

TR1 can be used to change the run-up time between 8 and 60 min. The controller is factory set to 8 min. If the gas load is large and the pump cannot reach the nominal speed in the set time, the pump is stopped and the controller gives Error 6 (Normal not reached at run-up time). In this case, the run-up time setting can be increased up to 60 min. by rotating TR1. The run-up time can be read through the serial communication at parameter 00700 (if protocol 1 is used) or at window 127 (if protocol 2 is used).

Controller Outline

The outline dimension for the controller is shown in the following figure.

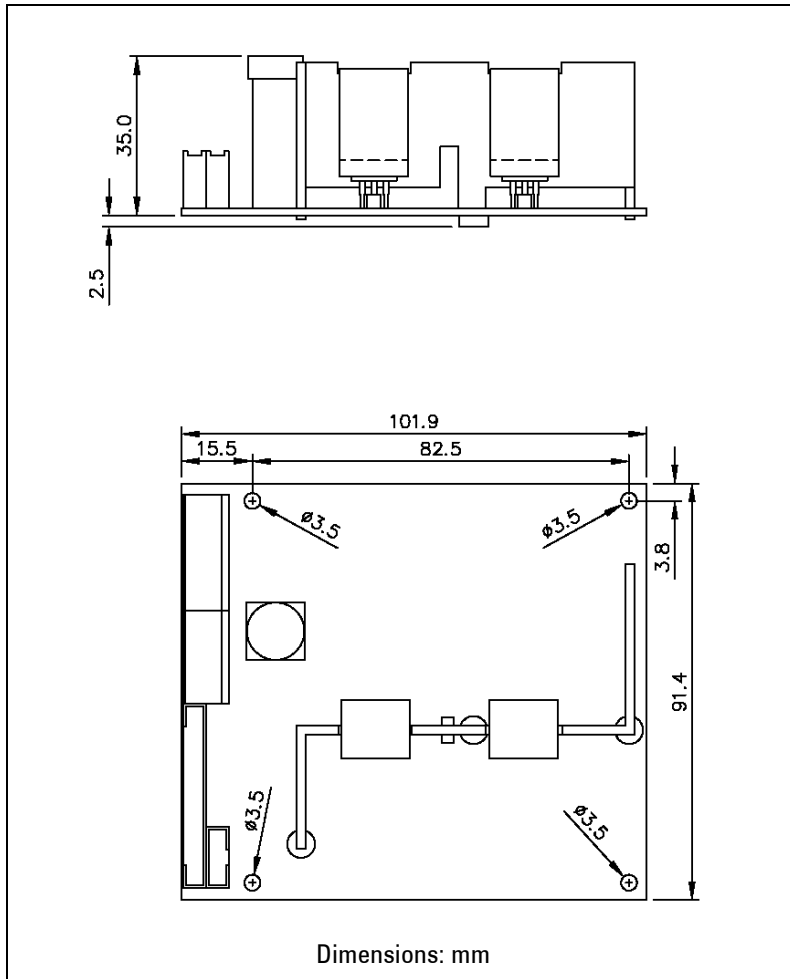


Figure 6 Controller Outline

Installation

Inspect the controller for any shipping damage. The controller is designed to be installed into a rack. All the connections are fitted to J2, J3, J4 and J5 connectors. Should the controller be connected to a host computer via the RS-232 interface, a suitable cable must be prepared. In the following paragraphs are detailed the input/output signals.

NOTE

The PCB installed into the customer system must be positioned so that cold air (forced or natural convection) can flow through the PCB components.

J3 Connector

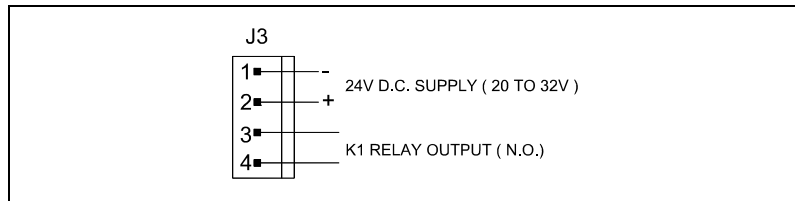


Figure 7 J3 – Power Connector

The signals of J3 connector are the following:

- **Pins 1/2** Power supply 24. Vdc (20 to 32 Vdc) (pin 2 positive).
- **Pins 4/3** K1 relay contacts. The relay switches when the output frequency overrides the set threshold (1117 Hz). The threshold overriding is indicated also on pin 6 of J4 connector.

J2 Connector

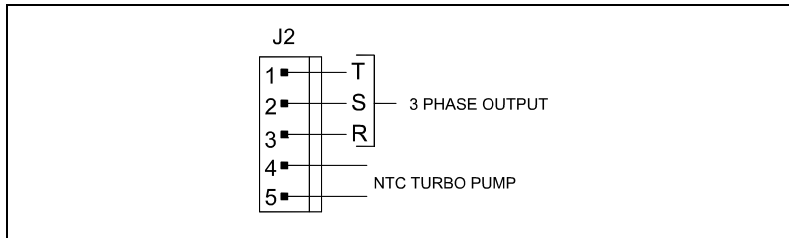


Figure 8 J2 – Connector to pump

The signals of J2 connector are the following:

- **Pin 1** 42 Vac 3-phase output to pump motor stator (phase T).
- **Pin 2** 42 Vac 3-phase output to pump motor stator (phase S).
- **Pin 3** 42 Vac 3-phase output to pump motor stator (phase R).
- **Pins 5/4** Pump temperature sensor.

J4 Connector

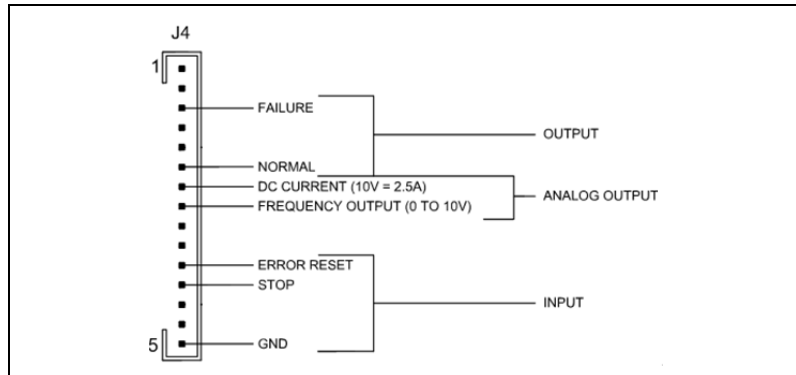


Figure 9 J4 – Signal Connector

The signals of J4 connector are the following:

- **Pin 3** 24V 60mA open collector Output failure. When a fault condition is detected, the pump is stopped and this output is ON.
- **Pin 6** 24V 60mA open collector Output When the output frequency is less than 1117 Hz, this output is ON.
- **Pin 7** Analog output. A voltage from 0 to 10 Vdc proportional to the output pump supplied current (from 0 to 2.5 A).
- **Pin 8** Analog output. A voltage from 0 to 10 Vdc proportional to the output pump motor driving frequency (from 0 to 1167 Hz).

17 Technical Information

Installation

- **Pin 11** Error reset command. Should an operation fault be detected (Pin 3 of J4 is ON), the pump is stopped. To reset the fault you must connect pin 11 to ground (pin 15 of J4) for at least one sec.
- **Pin 12** ON/OFF pump. To stop the pump you must connect pin 12 to ground (pin 15 of J4); to start the pump you must remove the contact between pin 12 and ground.
- **Pin 15** Ground

The following figure shows a typical contact logic input connection and the related simplified circuit of the controller.

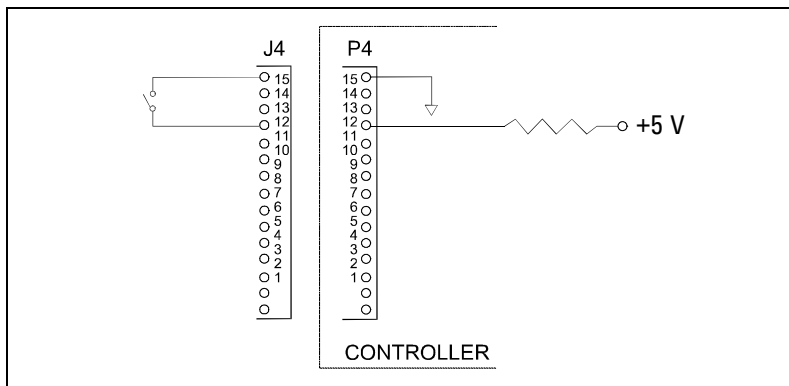


Figure 10 Typical Logic Input

The following figure shows a typical logic output connection, but any other device may be connected, and the related simplified circuit of the controller.

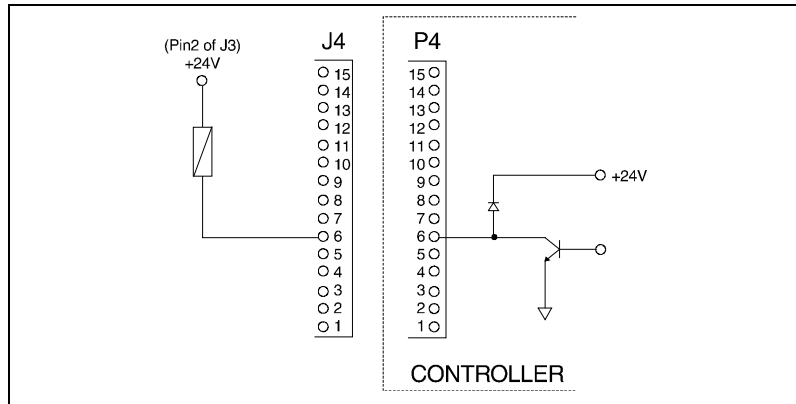


Figure 11 Typical Logic Output

J5 RS-232 Interface Connector

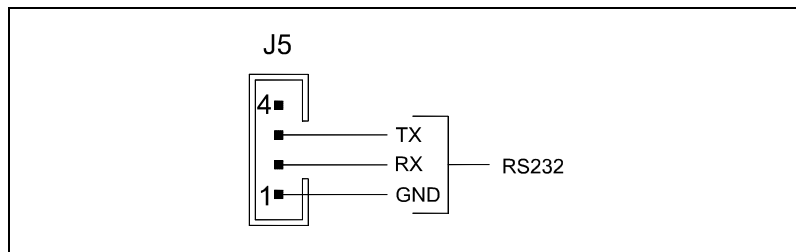


Figure 12 J2 – Interface Connector

The signal of J5, the RS-232 interface connector, are the following:

- **Pin 1** Ground.
- **Pin 2** RXD. Received data.
- **Pin 3** TXD. Transmitted data.
- **Pin 4** 24 Vdc Output (power supply to HHT, if necessary)

Data Exchange Protocols

This controller can manage two different serial communication protocols (1 or 2) on its RS-232 port. During the first communication frame after power on, the controller automatically recognizes the type of protocol used by the host to which it will interface.

NOTE

The identified protocol is kept active until the controller is powered off.

RS-232 Protocol 1 Communication Description

Transmission Channel Characteristics

Tab. 2

levels:	RS-232
baud rate:	9600/4800 programmable by W7 jumper on the board
character length:	8 bits
parity:	none
stop bit:	1 bit
protocol:	master (PC) / slave (converter)

Communication Protocol

The communication protocol is master-slave type, where:

- Host = Master
- Controller = slave

NOTE To write data through the serial interface, the controller must be set to "REMOTE OFF" operation, otherwise only read data command will be available.

The exchanged frame format is composed of ASCII values. The frame composition is:

slave address			Parameter number					data field length	data					check sum		CR			

where:

- **Slave Address:** made of three characters containing the controller address requested by the master. In RS-232 version the address is "001";
- **Parameter Number:** made of five characters, it points out to the controller about which parameter the master wants to operate, and about the kind of operation (read/write). It consists of:

1	2	3	4	5
0 = parameter reading 1 = parameter writing	always 0	000 to 299 = commands 300 to 699 = status 700 to 999 = parameter		

17 Technical Information

Data Exchange Protocols

- **Data Field Length:** made of 2 characters, it points out the next field length in characters. In this protocol the possible combinations are:

02 = parameter reading request

06 = parameter, writing
- **Data:** variable length, it could be 2 or 6 characters according to the previous field content, it can be:
"=?": in the case of parameter reading request
"000000" or "111111": in the case of YES/NO command or parameter writing
a six numeric character string 0 to 9
pointing out the value of a configuration parameter or an analog measure
a six character string pointing out an error
- **Checksum:** made of 3 characters, it is the sum of the string ASCII values up to the first character of the checksum with a 8 bit cutting-off converted to decimal number.
- **CR:** made of 1 character < 0xD>

All the fields are aligned to the right with the "0" character used as pad character.

When the controller receives a frame with the correct address value (001 in this case) it answers as detailed in the following:

- for a parameter request: a frame having the same format described before, but the first character of the "Parameter Number" field set to 1.

Example for a low speed status request:

Master to Controller:
0010000202= ?097< 0xD >

Controller to Master:
0011000206111111016< 0xD >
where " 111111" is for low speed status ON

- for a parameter writing: a frame including the parameter value that has been set out.

Example for a low speed status ON writing:

Master to Controller:
0011000206111111016< 0xD >

Controller to Master:
0011000206111111016< 0xD >

If the controller receives wrong frames it answers as following:

- in the case of frame with length greater than 40 characters:

			N	A	K	CR
Controller address						

- in case of frame with wrong checksum:

			N	A	K	CR
Controller address						

- in case of frame containing parameter number not present:

		1	0				0	6	N	O	-	D	E	F				
control. address		param. address													check-sum	CR		

- in case of writing values out of limits:

		1	0				0	6	-	R	A	N	G	E				
control. address		param. address													check-sum	CR		

17 Technical Information
Data Exchange Protocols

- in the case of an attempt of writing a logic parameter with a value not equal to "000000" or "111111", or when the parameter is a read only one:

			1	0				0	6	-	L	O	G	I	C				
control. address			param. address												check-sum	CR			

- in case of an excessive time between two characters (greater than 1 second):

			N	A	K	CR
Controller address						

Information available on the serial interface are listed in the following table where the column "TYPE" points out a logic variable with "D" (allowed values "000000" or "111111") and a numeric variable with "N".

Variables with the first character of "Parameter Number" equal to X can be either written or read. Variables with that character equal to "0" can be only read, equal to "1" can be only written.

Tab. 3

PARAMETER NUMBER	TYPE	DESCRIPTION
X0002	D	Low speed ON/OFF
X0003	D	Pump ON/OFF
X0008	D	Remote ON/OFF
10009	D	Error acknowledgment: error status clearing
00302	D	K1 status
00303	N	Error code: error code displaying (see the following table)

PARAMETER NUMBER	TYPE	DESCRIPTION
00306	D	Normal YES/NO
00307	D	Starting YES/NO
00309	N	Output frequency (Hz)
00310	N	Pump current * 100 (A)
00311	N	Temperature (°C)
00312	N	Software version The controller sends "QExxxx" where "xxxx" is the CRC16 of EPROM
00313	N	Power (W)
00314	N	Pump life (hours)
00315	N	Number of cycles
00316	N	Last cycle time (minutes)
00317	N	Software version. The controller sends "PAxxxx" where "xxxx" is the CRC16 of EEPROM
00700	N	Run up time (minutes)
X0701	N	Threshold (Hz)
X0702	D	Soft-start YES/NO
X0703	D	Reduced output power (YES) Full output power (NO)

Error Code Table

Tab. 4

CODE	DESCRIPTION
00E000	No error
00E001	Output current > 15 A
00E002	Pump not connected
00E003	Output current > 1.5 A for more than 15 sec.
00E004	Bearing temperature > 60 °C
00E005	Heat exchanger temperature > 60 °C
00E006	Normal not reached within run up time
00E009	Vp voltage 8V lower than nominal for more than 5 sec.
00E010	Soft start ramp not ended within the expected time

RS-232 Protocol 2 Communication Descriptions

Transmission Channel Characteristics

Tab. 5

levels:	RS-232
baud rate:	9600/4800 programmable by W7 jumper on the board
character length:	8 bits
parity:	none
stop bit:	1 bit
protocol:	master (PC) / slave (converter)

Message Structure

The communication protocol is a MASTER/SLAVE type where:

- Host = MASTER
- Controller = SLAVE

The communication is performed in the following way:

1. the host (MASTER) send a MESSAGE to the controller (SLAVE);
2. the controller answer with a MESSAGE to the host.

The MESSAGE is a string with the following format:

<STX>+<ADDR>+<WIN>+<COM>+<DATA>+<ETX>+<CRC>

where:

NOTE

When a data is indicated between two quotes ('...') it means that the indicated data is the corresponding ASCII character.

-
- <STX> (Start of transmission) = one character containing 0x02
 - <ADDR> (Unit address) = one character containing 0x80 (for RS-232)
<ADDR> (Unit address) = one character containing 0x80 + device number (0 to 31) (for RS-485)
 - <WIN> (Window) = a string of 3 numeric characters indicating the window number (from '000' to '999'); for the meaning of each window see the relevant paragraph.
 - <COM> (Command) = one character containing 0x30 to read the window, 0x31 to write into the window.

17 Technical Information

Data Exchange Protocols

- For host messages this field is present only in case of writing containing logic, numeric or alphanumeric data types according to the following table.
- For controller messages this field is always present containing the required data (logic, numeric or alphanumeric) in case of reading or the relevant protocol answer in case of writing acknowledge or protocol errors.

Tab. 6

	Length	Characters Permitted
Logic (L)	1	'0'= off '1'= on
Numeric (N)	6	'-', '.', '0' ... '9' right justified with '0'
Alphanumeric (A)	10	from blank to ' _ ' (ASCII)

- <ETX> (End of transmission) = one character containing 0x03
- <CRC> = two characters containing XOR of all characters subsequent to <STX> and including the <ETX> terminator. The value is hexadecimal coded and indicated by two ASCII character.

Only the addressed SLAVE will respond.

NOTE

Using the RS-485 interface, the message structure remains identical to the one used for the RS-232 interface, the only difference being that the value assigned to the ADDRESS <ADDR>

The controller can answers with the following response types:

Tab. 7 Response type

Type	Length	Value	Description
Logic	1 byte	-	Logic value
Numeric	6 bytes	-	Numeric value
Alphanumeric	10 bytes	-	Alphanumeric value
ACK	1 byte	(0x6)	The command execution has been successfully completed
NACK	1 byte	(0x15)	The command execution has been failed
Unknown Window	1 byte	(0x32)	The specified window in the command is not a valid window
Data Type Error	1 byte	(0x33)	The data type specified in the command (Logic, Numeric or Alphanumeric) is not accorded with the specified Window
Out of Range	1 byte	(0x34)	The value expressed during a write command is out of the range value of the specified window
Win Disabled	1 byte	(0x35)	The specified window is Read Only or temporarily disabled (for example you can't write the Soft Start when the Pump is running)

17 Technical Information

Example

Example

Command: START

Source: PC

Destination: Inverter

02	80	30	30	30	31	31	03	42	33
STX	ADDR	WINDOW			WR	ON	ETX	CRC	

Source: Inverter

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

Command: STOP

Source: PC

Destination: Inverter

02	80	30	30	30	31	30	03	42	32
STX	ADDR	WINDOW			WR	OFF	ETX	CRC	

Source: Inverter

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

Command: SOFT-START (ON)

Source: PC

Destination: Inverter

02	80	31	30	30	31	31	03	42	32
STX	ADDR	WINDOW			WR	ON	ETX	CRC	

Source: Inverter

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

Command: SOFT-START (OFF)

Source: PC

Destination: Inverter

02	80	31	30	30	31	30	03	42	33
STX	ADDR	WINDOW			WR	OFF	ETX	CRC	

Source: Inverter

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

17 Technical Information

Example

Command: CURRENT

Source: PC

Destination: Inverter

02	80	32	30	30	30	03	38	31
STX	ADDR	WINDOW	RD	ETX	CRC			

Source: Inverter

Destination: PC

02	80	32	30	30	30	30	30	30	2E	30	30	03	39	44
STX	ADD	WINDOW	RD	000.00				ETX	CRC					

Command: FREQUENCY

Source: PC

Destination: Inverter

02	80	32	30	33	30	03	38	32
STX	ADDR	WINDOW	RD	ETX	CRC			

Source: Inverter

Destination: PC

02	80	32	30	33	30	30	30	30	30	34	32	03	38	34
STX	ADD	WINDOW	RD	000042				ETX	CRC					

Command: ERR-CODE

Source: PC

Destination: Inverter

02	80	32	30	36	30	03	38	37
STX	ADDR	WINDOW	RD	ETX	CRC			

Source: Inverter

Destination: PC

02	80	32	30	36	30	30	30	30	30	30	30	03	38	37
STX	ADD	WINDOW	RD	000000				ETX	CRC					

Serial Communication Windows

Tab. 8

WIN	TYPE	R	W	Description
000	L	X	X	START/STOP (1= START ; 0= STOP)
008	L	X	X	REMOTE/SERIAL Configuration (1= Remote ; 0= Serial)
100	L	X	X	SOFT START YES/NO (1= YES ; 0= NO) Default = 0
101	N	X	X	SET POINT TYPE (0 =Frequency; 1=Current; 2=Time) Default = 0
102	N	X	X	SET POINT VALUE [Hz, Ma, SEC] Default = 1000 Hz
103	N	X	X	SET POINT MASK TIME [s] Default = 0
104	L	X	X	OUTPUT STATE (0= OFF if Val > threshold; 1 = ON if Val > threshold) Default = 0
105	N	X	X	HYSTERESIS IN THRESHOLD % [0-100] Default = 2
107	L	X	X	ACTIVE STOP (0=NO; 1=YES) Default = 1
108	N	X	X	BAUD RATE (3-4) [4800-9600] Default = 4
109	L		X	PUMP LIFE RESET [Write "1" to Reset]
111	N	X	X	SET ANALOG OUTPUT TYPE [0=Power; 1=Current] Default = 0
120	N	X	X	SET ROTATIONAL FREQUENCY [Hz] $1100 \text{ Hz} \leq F_{\text{imp}} \leq F_{\text{MAX}}$
121	N	X	X	MAX SETTABLE ROTATIONAL FREQUENCY [Hz] $F < 1167 \text{ Hz}$
127	N	X		RUN UP TIME (8 – 60 minute)
130	N	X		RAMP CURRENT [mA]
200	N	X		CURRENT [mA]
201	N	X		VOLTAGE [V]
202	N	X		POWER [W]
203	N	X		DRIVING FREQUENCY [Hz]

17 Technical Information

Example

WIN	TYPE	R	W	Description
204	N	X		PUMP TEMPERATURE [°C]
205	N	X		STATUS [0=stop; 1=interlock; 2=ramp; 3=regulation; 4=brake; 5=normal; 6=failure]
206	N	X		ERROR CODE:
211	N	X		CONTROLLER SENSOR TEMPERATURE READING [208= 25 °C - 128= 60 °C]
216	N	X		AMBIENT SENSOR TEMPERATURE READING
300	N	X		CYCLE TIME [min]
301	N	X		CYCLE NUMBER
302	N	X		PUMP LIFE [h]
319	A	X		Controller Model
320	A	X		Base Pump Model Number (8 characters)
321	A	X		Modified Standard Model Number (4 characters)
323	A	X		Controller Serial Number (5 characters)
325	A	X		Electrical Modification Level (10 characters)
400	A	X		CRC PROGRAM LISTING [QE7xxxx]
401	A	X		CRC BOOTLOADER [BL1xxxx]
402	A	X		CRC PARAMETER LISTING [PA7xxxx]
404	A	X		CRC FILE PARAMETER STRUCTURE
406	A	X		PROGRAM LISTING CODE & REVISION
407	A	X		PARAMETER LISTING CODE & REVISION

WIN = Window

R = Read

W = Write

T = Type

L = Logical

N = Numeric

A = Alphanumeric

Operation

Make all vacuum manifold and electrical connections and refer to Turbo pump instruction manual prior to operating the Turbo controller.

The controller operates completely automatically after the remote start command is given.

Switching On/Off the Pump

To switch on the pump is necessary to remove the connection between pin 12 of J4 connector and pin 15 (ground) of the same connector. To switch off the pump is necessary to connect pin 12 of J4 connector to pin I5 (ground) of the same connector.

Low Speed Activation/Deactivation

To activate the Low Speed it is necessary to connect pin 13 of J4 connector to pin 15 (ground) of the same connector.

To deactivate the Low Speed status it is necessary to disconnect pin 13 of J4 connector from pin 15 (ground) of the same connector.

The low speed status is available on pin 5 of the same connector.

The low speed frequency is 1100 Hz.

Error reset

If the controller identifies an error, the pump is switched off. To start again the pump the error cause has been removed and the error status has to be reset connecting pin 11 of J4 connector to pin 15 (ground) of the same connector for at least 1 second.

Maintenance

Replacement controllers are available on an advance exchange basis through Agilent. If necessary, information is provided to aid the operator in determining malfunctions and corrective steps to be taken.

WARNING!



Voltages developed in the unit are dangerous and may be fatal. Service must be performed by authorized personnel only.

Error Messages

The controller will self-diagnose some error causes. The controller shows the error occurred by means of a diagnostic LED located on the board, and through the serial communication port.

The LED blinks in a coded mode:

it flashes a number of time equal to the error code (see the following table) and then stays off for a while; this sequence has been repeated until the error status has been reset.

LED Error Code Table

Tab. 9

LED BLINKING NUMBER	DESCRIPTION
0	No error
1	Output current > 15 A
2	Not connected pump
3	Output current > 1.5 A for 15 sec.
4	Bearing temperature > 60 °C
5	Heat exchanger temperature > 60 °C
6	Normal not reached at run up time
9	V _p voltage 8V lower than nominal for 5 sec.
10	Soft start ramp not ended within the expected time

Accessories and Spare Parts

Tab. 10

DESCRIPTION	PART NUMBER
Pump Cable 1 m long with connector	9699869



Vacuum Products Division

Dear Customer,

Thank you for purchasing an Agilent vacuum product. At Agilent Vacuum Products Division we make every effort to ensure that you will be satisfied with the product and/or service you have purchased.

As part of our Continuous Improvement effort, we ask that you report to us any problem you may have had with the purchase or operation of our products. On the back side you find a Corrective Action request form that you may fill out in the first part and return to us.

This form is intended to supplement normal lines of communications and to resolve problems that existing systems are not addressing in an adequate or timely manner.

Upon receipt of your Corrective Action Request we will determine the Root Cause of the problem and take the necessary actions to eliminate it. You will be contacted by one of our employees who will review the problem with you and update you, with the second part of the same form, on our actions.

Your business is very important to us. Please, take the time and let us know how we can improve.

Sincerely,

A handwritten signature in black ink, appearing to read "Giampaolo LEVI".

Giampaolo LEVI

*Vice President and General Manager
Agilent Vacuum Products Division*

Note: Fax or mail the Customer Request for Action (see backside page) to Agilent Vacuum Products Division (Torino) – Quality Assurance or to your nearest Agilent representative for onward transmission to the same address.

United States

Agilent Technologies
121 Hartwell Avenue
Lexington, MA 02421 - USA
Ph. +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426
vppl-customer@agilent.com

Netherlands

Agilent Technologies Netherlands B.V.
Customer Contact Center
Laan van Langerhuize 1, toren A-8
1186 DS Amstelveen
Tel. +31 020 547 2600
Fax +31 020 654 5748
customer@agilent.com

Belgium

Agilent Technologies Belgium S.A./N.V.
Customer Contact Center
Pegasus Park
De Kleetlaan 12A bus 12
B-1831 Diegem
Tel. +32 2 404 92 22
Fax +32 2 626 46 30
customer@agilent.com

Brazil

Agilent Technologies Brasil
Avenida Marcos Penteados de Ulhoa
Rodrigues, 939 - 6° andar
Castelo Branco Office Park
Torre Jacarandá - Tamboré
Barueri, Sao Paulo CEP: 06460-040
Toll free: 0800 728 1405

China

Agilent Technologies (China) Co. Ltd
No.3, Wang Jing Bei Lu, Chao Yang District
Beijing, 100102, China
Tel: +86 (0)10 64397888
Fax: +86 (0)10 64392765
Toll free: 800 820 6778 (ph); 400 820 6778 (cell.ph.)
vacuum.cnmarketing@agilent.com
vpk-customer@agilent.com

France

Agilent Technologies
Parc Technopolis - Z.A. de Courtaboeuf
3, avenue du Canada - CS 90263
91978 Les Ulis cedex, France
Tel: +33 (0) 1 64 53 61 15
Fax: +33 (0) 1 64 53 50 01
vpf.sales@agilent.com

Southeast Asia

Agilent Technologies Sales Sdn Bhd
Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya, Selangor, Malaysia
Ph. +603 7712 6181
Fax: +603 7727 1239
Toll free: 1 800 880 805
vps-customer@agilent.com

India (Sales)

Agilent Technologies India Pvt. Ltd.
Unit Nos 110- 116, & Part of 101 & 109
First Floor, Splendor Forum,
Plot No.-3 , District Centre, Jasola
New Delhi-110025
Ph. +91 11 4623 7100
Fax: +91 4623 7105
Toll Free: 1 800 180 1517

Italy

Agilent Technologies Italia S.p.A.
Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel: +39 011 9979 111
Fax: +39 011 9979 350
Toll free: 00 800 234 234 00
vppl-customer@agilent.com

Japan

Agilent Technologies Japan, Ltd.
9-1 Takakura-cho Hachioji-city,
Tokyo, Japan
Tel.: +81-3-5232-1253
Fax: +81-120-565-154
Toll-Free: +81-120-477-111
jp-vvt-sales.pdl-ext@agilent.com

Singapore

Agilent Technologies Singapore Pte. Ltd.
1 Yishun Avenue 7,
Singapore 768923
Tel : (65) 6215 8045
Fax : (65) 6754 0574
Toll free: 1 800 2762622
vps-customer@agilent.com

Korea

Agilent Technologies Korea, Ltd.
Ilshin Building 4F
Yongsan-gu Hannam-daero
Seoul Korea 04418
Tel: +82 (0)2 2194 9449
Fax: +82 (0)2 2194 9853
Toll free: 080 222 2452
vpk-customer@agilent.com

UK and Ireland

Agilent Technologies LDA UK Limited
Lakeside Cheadle Royal Business Park
Cheadle, Cheshire SK8 3GR,
United Kingdom
Ph. +44 01865291570
Fax +44 01865291571
Toll free: 00 800 234 234 00
Toll free fax: 00 800 345 345 00
vpt-customer@agilent.com

India (Service)

Agilent Technologies India Pvt. Ltd.
C-Block, RMZ Centennial Plot Number-
8A, 8B, 8C, 8D,
Doddanakundi Industrial Area, ITPL
Road,
Mahadevapura Post, Bangalore- 560048
Ph. +91 80 40614000
Fax: +91 80 40148991

Taiwan

Agilent Technologies Taiwan Limited
No. 20 Gao-shuang Road, Ping-zhen Dist
Tao-Yuan City
32450 Taiwan, R.O.C.
Tel: +886 3 4959004
Toll free: 0800 018 768
vpw-customer@agilent.com

Germany and Austria

**Agilent Technologies
Sales & Services GmbH & Co. KG**
Lyoner Str. 20
60 528 Frankfurt am Main
GERMANY
Tel: +49 69 6773 43 2230
Fax: +49 69 6773 43 2250

Mexico

Agilent Technologies
Concepcion Beistegui No 109
Col Del Valle
C.P. 03100 – Mexico, D.F.
Tel.: +52 5 523 9465
Fax: +52 5 523 9472

Other Countries

Agilent Technologies Italia S.p.A.
Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel.: +39 011 997 9111
Fax: +39 011 997 9350
Toll-Free: 00 800 234 234 00
vpt-customer@agilent.com

Customer Support & Service

NORTH AMERICA:
Toll Free: 800 882 7426
vppl-ra@agilent.com
Lexington-service@agilent.com

EUROPE:

Toll Free: 00 800 234 234 00
vpt-customer@agilent.com

PACIFIC RIM:

please visit our website for individual
office information
<http://www.agilent.com/chem/vacuum>

Worldwide Web Site, Catalog and Order

On-line:
www.agilent.com/chem/vacuum
Representatives in most countries

© Agilent Technologies, Inc. 2016

Printed in ITALY

07/2018

Publication Number: 87-901-056-01



Agilent Technologies

CUSTOMER REQUEST FOR CORRECTIVE / PREVENTIVE / IMPROVEMENT ACTION

TO: AGILENT VACUUM PRODUCTS DIVISION TORINO – QUALITY ASSURANCE FAX

N°: XXXX-011-9979350

ADDRESS: AGILENT TECHNOLOGIES ITALIA S.p.A. – Vacuum Products Division –
Via F.lli Varian, 54 – 10040 Leini (TO) – Italy

E-MAIL: vpd-qualityassurance_pdl-ext@agilent.com

NAME	COMPANY	FUNCTION
ADDRESS:		
TEL. N° :		FAX N° :
E-MAIL:		
PROBLEM / SUGGESTION :		
REFERENCE INFORMATION (model n°, serial n°, ordering information, time to failure after installation, etc.):		
		DATE
CORRECTIVE ACTION PLAN / ACTUATION (by AGILENT VPD)		LOG N°

XXX = Code for dialing Italy from your country (es. 01139 from USA; 00139 from Japan, etc.)

Vacuum Products Division Instructions for returning products

Dear Customer,

Please follow these instructions whenever one of our products needs to be returned.

Complete the attached **Request for Return form** and send it to Agilent Technologies (see below), taking particular care to include the completed **Health and Safety** declaration Section. No work can be started on your unit until we receive a completed copy of this form.

After evaluating the information, Agilent Technologies will provide you with a **Return Authorization (RA) number** via email or fax, as requested. Note: Depending on the type of return, a Purchase Order may be required at the time **the Request for Return is submitted**. We will quote any necessary services (evaluation, repair, special cleaning, eg).

Product preparation

- Remove all accessories from the core product (e.g. inlet screens, vent valves).
- Prior to shipment and if applicable for your product, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
- If ordering an Advance Exchange product, please use the packaging from the Advance Exchange to return the defective product.
- Seal the product in a plastic bag, and package product carefully to avoid damage in transit. You are responsible for loss or damage in transit.
- Include a copy of the Health and Safety Declaration in the shipping documentation on the outside of the shipping box of your returning product.
- Clearly label package with RA number. Using the shipping label provided will ensure the proper address and RA number are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will be returned.
- Return only products for which the RA was issued.

Shipping

- Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information. Customer is responsible for freight charges on returning product.
- Return shipments must comply with all applicable Shipping Regulations (IATA, DOT, ADR, etc.) and carrier requirements.

RETURN THE COMPLETED **REQUEST FOR RETURN** FORM TO YOUR NEAREST LOCATION:

EUROPE:

Fax: 00 39 011 9979 330
Fax Free: 00 800 345 345 00
Toll Free: 00 800 234 234 00

NORTH AMERICA:

Fax: 1 781 860 9252
Toll Free: 800 882 7426, Option 3

PACIFIC RIM:

Please visit our website for individual office information

vpt-customer@agilent.com

vpl-ra@agilent.com

<http://www.agilent.com>




TERMS AND CONDITIONS

Please read the terms and conditions below as they apply to all returns and are in addition to the Agilent Technologies Vacuum Product Division – Products and Services Terms of Sale.

- Unless otherwise pre-negotiated, customer is responsible for the freight charges for the returning product. Return shipments must comply with all applicable **Shipping Regulations** (IATA, DOT, etc.) and carrier requirements.
- Agilent Technologies is not responsible for returning customer provided packaging or containers.
- Customers receiving an Advance Exchange product agree to return the defective, rebuildable part to Agilent Technologies **within 15 business days**. Failure to do so, or returning a non-rebuildable part (crashed), will result in an invoice for the non-returned/non-rebuildable part.
- Returns for credit toward the purchase of new or refurbished Products are subject to prior Agilent approval and may incur a restocking fee. Please reference the original purchase order number.
- Units returned for evaluation will be evaluated, and a quote for repair will be issued. If you choose to have the unit repaired, the cost of the evaluation will be deducted from the final repair pricing. A Purchase Order for the final repair price should be issued within 3 weeks of quotation date. Units without a Purchase Order for repair will be returned to the customer, and the evaluation fee will be invoiced.
- Products returned that have not been drained from oil will be disposed.
- A Special Cleaning fee will apply to all exposed products.
- If requesting a calibration service, units must be functionally capable of being calibrated.

Customer information		
Company :		Contact Name:
Address:	Tel:	Fax:
		Email:

Equipment			
Product description	Agilent PartNo	Agilent Serial No	Original Purchasing Reference
Failure description		Type of process (for which the equipment was used)	

Type of return
<input type="checkbox"/> Non Billable <input type="checkbox"/> Billable  New PO # (hard copy must be submitted with this form): _____ <input type="checkbox"/> Exchange <input type="checkbox"/> Repair <input type="checkbox"/> Upgrade <input type="checkbox"/> Demo <input type="checkbox"/> Calibration <input type="checkbox"/> Evaluation <input type="checkbox"/> Return for Credit

Health and safety	Substances (please refer to MSDS forms)																																																																																				
The product has been exposed to the following substances: (by selecting 'YES' you MUST complete the table to the right)	* Agilent will not accept delivery of any product that is exposed to radioactive, biological, explosive substances or dioxins, PCB's without written evidence of decontamination.																																																																																				
	<table border="1"> <thead> <tr> <th>Trade name</th> <th>Chemical name</th> <th>Chemical Symbol</th> <th>CAS Number</th> </tr> </thead> <tbody> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> <tr><td> </td><td> </td><td> </td><td> </td></tr> </tbody> </table>	Trade name	Chemical name	Chemical Symbol	CAS Number																																																																																
Trade name	Chemical name	Chemical Symbol	CAS Number																																																																																		
Toxic	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Harmful	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Corrosive	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Reactive	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Flammable	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Explosive (*)	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Radioactive (*)	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Biological (*)	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Oxidizing	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Sensitizer	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				
Other dangerous substances	<input type="checkbox"/> YES <input type="checkbox"/> NO																																																																																				

Goods preparation
If you have replied YES to one of the above questions. Has the product been purged? <input type="checkbox"/> YES <input type="checkbox"/> NO
If yes, which cleaning agent/method:
Has the product been drained from oil? <input type="checkbox"/> YES <input type="checkbox"/> NOT APPLICABLE
I confirm to place this declaration on the outside of the shipping box. <input type="checkbox"/>

**I declare that the above information is true and complete to the best of my knowledge and belief.
I understand and agree to the terms and conditions on page 2 of this document.**

Name:	Authorized Signature:
Position:	
Date:	

NOTE: If a product is received at Agilent which is contaminated with a toxic or hazardous material that was not disclosed, **the customer will be held responsible** for all costs incurred to ensure the safe handling of the product, and is liable for any harm or injury to Agilent employees as well as to any third party occurring as a result of exposure to toxic or hazardous materials present in the product.

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies

121 Hartwell Avenue
Lexington, MA 02421 - USA
Ph. +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426
vpl-customer@agilent.com

Netherlands

Agilent Technologies Netherlands B.V.

Customer Contact Center
Laan van Langerhuize 1, toren A-8
1186 DS Amstelveen
Tel. +31 020 547 2600
Fax +31 020 654 5748
customer@agilent.com

Belgium

Agilent Technologies Belgium S.A./N.V.

Customer Contact Center
Pegasus Park
De Kleetlaan 12A bus 12
B-1831 Diegem
Tel. +32 2 404 92 22
Fax +32 2 626 46 30
customer@agilent.com

Brazil

Agilent Technologies Brasil

Avenida Marcos Penteado de Ulhoa
Rodrigues, 939 - 6° andar
Castelo Branco Office Park
Torre Jacarandá - Tamboré
Barueri, Sao Paulo CEP: 06460-040
Toll free: 0800 728 1405

China

Agilent Technologies (China) Co. Ltd

No.3, Wang Jing Bei Lu, Chao Yang District
Beijing, 100102, China
Tel: +86 (0)10 64397888
Fax: +86 (0)10 64392765
Toll free: 400 8206778 (mobile)
Toll free: 800 8206778 (landline)
vacuum.cnmarketing@agilent.com
vpc-customerservice@agilent.com

France

Agilent Technologies
Parc Technopolis - Z.A. de Courtaboeuf
3, avenue du Canada - CS 90263
91978 Les Ulis cedex, France
Tel: +33 (0) 1 64 53 61 15
Fax: +33 (0) 1 64 53 50 01
vpf.sales@agilent.com

Southeast Asia

Agilent Technologies Sales Sdn Bhd

Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya, Selangor, Malaysia
Ph. +603 7712 6181
Fax: +603 7727 1239
Toll free: 1 800 880 805
vps-customerservice@agilent.com

India (Sales)

Agilent Technologies India Pvt. Ltd.

Unit Nos 110- 116, & Part of 101 & 109
First Floor, Splendor Forum,
Plot No.-3, District Centre, Jasola
New Delhi-110025
Ph. +91 11 4623 7100
Fax: +91 4623 7105
Toll Free: 1 800 180 1517

Italy

Agilent Technologies Italia S.p.A.

Via F.Ili Varian, 54
10040 Leini, (Torino) - Italy
Tel: +39 011 9979 111
Fax: +39 011 9979 350
Toll free: 00 800 234 234 00
vpt-customerservice@agilent.com

Japan

Agilent Technologies Japan, Ltd.

9-1 Takakura-cho Hachioji-city,
Tokyo, Japan
Tel.: +81- 3-5232-1253
Fax: +81-120-565-154
Toll-Free: +81-120-477-111
jp-vvt-sales.pdl-ext@agilent.com

Singapore

Agilent Technologies Singapore Pte. Ltd.

1 Yishun Avenue 7,
Singapore 768923
Tel : (65) 6215 8045
Fax : (65) 6754 0574
Toll free: 1 800 2762622
vps-customerservice@agilent.com

Korea

Agilent Technologies Korea, Ltd.

Ilshin Building 4F
Yongsan-gu Hannam-daero
Seoul Korea 04418
Tel: +82 (0)2 2194 9449
Fax: +82 (0)2 2194 9853
Toll free: 080 222 2452
vpc-customerservice@agilent.com

UK and Ireland

Agilent Technologies LDA UK Limited

Lakeside Cheadle Royal Business Park
Cheadle, Cheshire SK8 3GR,
United Kingdom
Ph. +44 01865291570
Fax +44 01865291571
Toll free: 00 800 234 234 00
Toll free fax: 00 800 345 345 00
vpt-customer@agilent.com

© Agilent Technologies, Inc. 2018
07/2018

India (Service)

Agilent Technologies India Pvt. Ltd.
C-Block, RMZ Centennial Plot Number- 8A, 8B, 8C,
8D,
Doddanakundi Industrial Area, ITPL Road,
Mahadevapura Post, Bangalore- 560048
Ph. +91 80 40614000
Fax: +91 80 40148991

Taiwan

Agilent Technologies Taiwan Limited

No. 20 Gao-shuang Road, Ping-zhen Dist
Tao-Yuan City
32450 Taiwan, R.O.C.
Tel: +886 3 4959004
Toll free: 0800 018 768
vpw-customerservice@agilent.com

Germany and Austria

Agilent Technologies Sales & Services GmbH & Co. KG

Lyoner Str. 20
60 528 Frankfurt am Main
GERMANY
Tel: +49 69 6773 43 2230
Fax: +49 69 6773 43 2250

Mexico

Agilent Technologies

Concepcion Beistegui No 109
Col Del Valle
C.P. 03100 – Mexico, D.F.
Tel.: +52 5 523 9465
Fax: +52 5 523 9472

Other Countries

Agilent Technologies Italia S.p.A.

Via F.Ili Varian, 54
10040 Leini, (Torino) - Italy
Tel.: +39 011 997 9111
Fax: +39 011 997 9350
Toll-Free: 00 800 234 234 00
vpt-customer@agilent.com

Customer Support & Service

NORTH AMERICA:

Toll Free: 800 882 7426
vpl-ra@agilent.com
Lexington-service@agilent.com

EUROPE:

Toll Free: 00 800 234 234 00
vpt-customer@agilent.com

PACIFIC RIM:

please visit our website for individual office
information
<http://www.agilent.com/chem/vacuum>

Worldwide Web Site, Catalog and Order On-line:

www.agilent.com/chem/vacuum
Representatives in most countries