

Ion Pumps Controller

24 V

**Model
929-0198**

**Manuale di Istruzioni
Bedienungshandbuch
Notice de Mode D'Emploi
User Manual**

**87-900-127-01 (C)
05/2011**



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2011

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: 87-900-127-01 (C)

Edition

Edition 05/2011

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.

Ion Pumps Controller 24 V

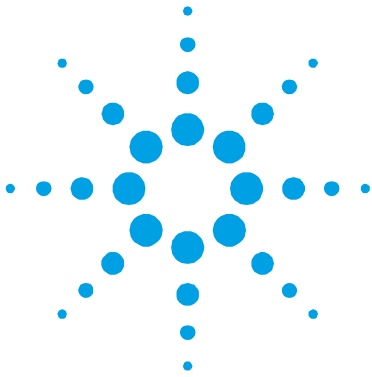


Contents

1	Procedura per l'installazione	7
	Descrizione del Controller Ion Pump 24 V	8
	Informazioni Generali	8
	Installazione	9
	Smaltimento	12
2	Anleitung zur Installation	13
	24 Volt Iongetterpumpen-Kontroller	14
	Allgemeines	14
	Installation	15
	Entsorgung	18
3	Procédure pour l'installation	19
	Description du controleur ion pump 24 V	20
	Informations Generales	20
	Installation	21
	Mise au rebut	24
4	Installation procedure	25
	Ion Pumps Controller 24V	26
	General Information	26
	Installation	27

Contents

	Disposal	30
5	Technical Information	31
	Controller Description	32
	Orderable Parts	36



1 Procedura per l'installazione

Descrizione del Controller Ion Pump 24 V8

Informazioni Generali 8

Installazione 9

Smaltimento 12

Traduzione delle istruzioni originali



Descrizione del Controller Ion Pump 24 V

AVVERTENZA!



L'alta tensione sviluppata nel controller può causare gravi danni fisici o il decesso. Prima di mettere in funzione l'unità, scollegare il cavo di alimentazione.

Per il collegamento del controller, utilizzare esclusivamente l'apposito cavo fornito.

ATTENZIONE!

Non installare o utilizzare il controller in ambienti esposti ad agenti atmosferici (pioggia, neve, ghiaccio), polvere, gas aggressivi, oppure in ambienti sottoposti al rischio di esplosione o incendio.

NOTA

Quando il controller viene acceso è udibile un rumore acuto di intensità e di frequenza variabile. Questo rumore è del tutto normale e non è indice di malfunzionamento.

Informazioni Generali

Questo controller è stato progettato per essere utilizzato da professionisti. Prima di utilizzare il controller, l'operatore deve leggere il presente Manuale d'Istruzioni e qualsiasi altro documento informativo fornito dalla Agilent. La Agilent non sarà tenuta responsabile di qualsiasi evento che possa verificarsi a causa del mancato rispetto, anche parziale, di dette istruzioni, nonché a causa dell'uso scorretto del controller da parte di personale non formato o non autorizzato o di qualsiasi operazione non conforme agli standard nazionali specifici.

Installazione

Durante il funzionamento, devono essere rispettate le seguenti condizioni ambientali:

- temperatura: da 0 °C a +45 °C
- umidità relativa: 0 – 95 % (in assenza di condensa)

Per collegare il controller alla pompa, utilizzare gli appositi cavi forniti. Per maggiori informazioni circa i collegamenti sopraccitati o di altro genere, fare riferimento all'appendice "Technical Information".

Attivazione/disattivazione HV mediante il pannello anteriore:

Agire sull'interruttore presente sul pannello anterior

Tab. 1

Tensione di input	
24 Vcc ± 10 %	OK
< 19 Vcc per più di 1 minuto	HV ON e Avvertimento batteria con accensione simultanea a luce lampeggiante di entrambi i LED con 1 lampeggio ogni 2 sec
< 15 Vdc per più di 2 minuti	HV OFF e assenza di lampeggiamento del LED

1 Procedura per l'installazione

Installazione

LED

Tab. 2 :

Led Pos: 1 lampeggio ogni 2 sec	HV POS ON: funzionamento normale
Led Neg: 1 lampeggio ogni 2 sec	HV NEG ON: funzionamento normale
Led Pos: 2 lampeggi ogni 2 sec	HV POS Corto Circuito: tensione di output < 100 V e corrente in uscita > 4 mA per più di 1 minuto
Led Neg: 2 lampeggi ogni 2 sec	HV NEG Corto Circuito: tensione di output < 100 V e corrente in uscita > 4 mA per più di 1 minuto
Led Pos: 3 lampeggi ogni 2 sec	HV POS fuori servizio: il valore della tensione di output è superiore del 120 % o inferiore dell'80 % rispetto al valore target, per più di 1 minuto
Led Neg: 3 lampeggi ogni 2 sec	HV NEG fuori servizio: il valore della tensione di output è superiore del 120 % o inferiore dell'80 % rispetto al valore target, per più di 1 minuto
Led Pos: 4 lampeggi ogni 2 sec	Cavo di Interlock non connesso
Led Neg: 4 lampeggi ogni 2 sec	Cavo di Interlock non connesso
1 lampeggio di entrambi i LED ogni 2 sec	La tensione di alimentazione è inferiore a 19 Vcc per più di 1 minuto.
Assenza di lampeggiamento di entrambi i LED	HV OFF: la tensione di alimentazione è inferiore a 15 Vcc per più di 2 minuti

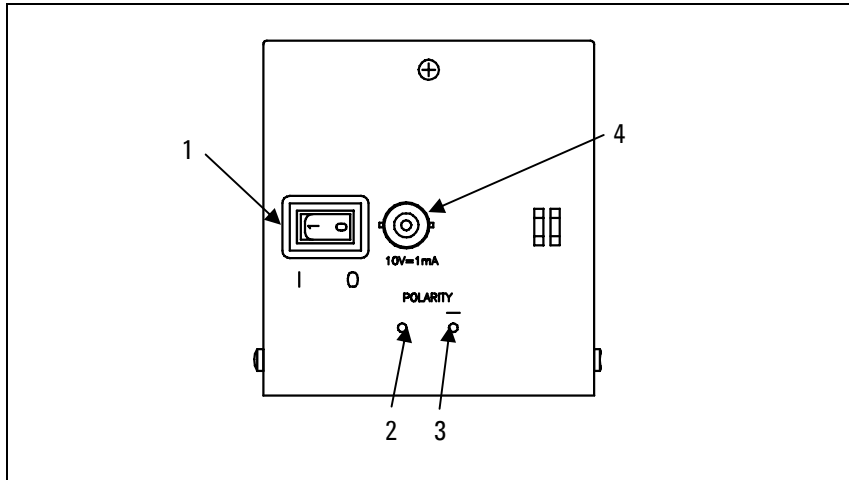


Figura 1 Pannello anteriore

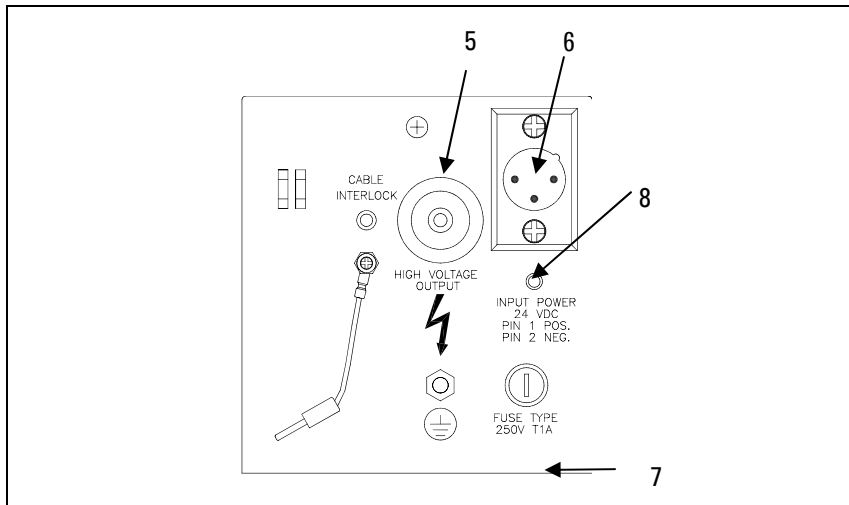


Figura 2 Pannello posteriore

1 Procedura per l'installazione

Smaltimento

Tab. 3

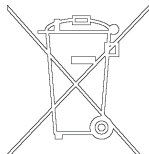
RIF.	SEGNALE	DESCRIZIONE
1	ON/OFF	Accende l'unità e l'Alta Tensione
2	Positive LED	1 lampeggio ogni 2 secondi quando l'unità e HV pos. sono ON
3	Negative LED	1 lampeggio ogni 2 secondi quando l'unità e HV neg. sono ON
4	Recorder Output Connector	Da 0 a 10 V lineari 1 μ A = 10 mV e 1 mA = 10 V
5	HV Connector	Connettore Alta Tensione
6	DC INPUT	24 V \pm 10 %
7	FUSE	Supporto fusibile T1A SLOW

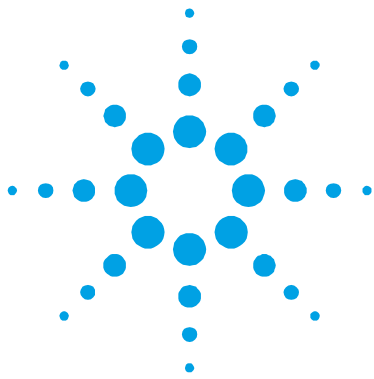
Smaltimento

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.





2

Anleitung zur Installation

24 Volt Ionengetterpumpen-Kontroller	14
Allgemeines	14
Installation	15
Entsorgung	18

Übersetzung der Originalanleitungen



24 Volt Ionengetterpumpen-Kontroller

WARNUNG!



Die im Kontroller erzeugte Hochspannung kann ernsthafte Verletzungen oder Tod verursachen. Vor Wartungsarbeiten am Gerät unbedingt das Stromkabel ziehen.

Um den Kontroller in Betrieb zu nehmen, das mitgelieferte Stromkabel verwenden.

VORSICHT!

Den Kontroller nicht in Umgebungen installieren oder benutzen, die atmosphärischen Bedingungen (Regen, Schnee, Eis, etc.), Staub, aggressiven Gasen, Explosionsgefahr oder solchen mit hoher Feuergefahr ausgesetzt sind.

HINWEIS

Die Geräusche in verschiedener Lautstärke und Frequenz, die bei angeschaltetem Gerät hörbar sind, sind völlig normal und bedeuten keine Fehlfunktion.

Allgemeines

Der Kontroller ist für Fachbetriebe bestimmt. Vor Gebrauch sollte der Benutzer dieses Handbuch sowie alle weiteren von der Agilent mitgelieferten Zusatzdokumente sorgfältig lesen. Bei - auch nur teilweiser - Nichtbeachtung der enthaltenen Hinweise, unsachgemäßem Gebrauch durch ungeschultes Personal, nicht autorisierten Eingriffen und Missachtung der einheimischen, hier zur Geltung kommenden Bestimmungen übernimmt die Agilent keinerlei Haftung.

Installation

Beim Gebrauch müssen die folgenden Umgebungsbedingungen eingehalten werden:

- Temperatur: 0 °C bis + 45 °C
- rel. Luftfeuchtigkeit: 0 - 95 % (nicht kondensierend)

Zum Anschluss des Kontrollers an die Pumpe die mitgelieferten Kabel benutzen. Für Einzelheiten zu den erwähnten und anderen Anschlüsse den Anhang "Technical Information" zu Rate ziehen.

Hochspannung An/Aus auf dem vorderen Bedienfeld:

Benutzen Sie den Schalter auf dem vorderen Bedienfeld

Tab. 1

Eingangsspannung	
24 V Gleichspannung ± 10 %	OK
< 19 V Gleichspannung für mehr als eine Minute	Hochspannung liegt an und eine Batterie-Warnung mit beiden LEDs wird angezeigt (blinken 1 mal alle 2 Sekunden)
< 15 V Gleichspannung für mehr als 2 Minuten	Hochspannung ausgeschaltet und keine blinkenden LEDs

2 Anleitung zur Installation

Installation

LED-Anzeigen:

Tab. 2

Positiv LED: Blink 1x alle 2 Sekunden	POSITIVE HOCHSPANNUNG AN: Normalbetrieb
Negativ LED: Blink 1x alle 2 Sekunden	NEGATIVE HOCHSPANNUNG AN: Normalbetrieb
Positiv LED: Blink 2x alle 2 Sekunden	Kurzschluß positive Hochspannung: Ausgangsspannung <100 V und Ausgangsstrom > 4mA für länger als 1 Minute
Negativ LED: Blink 2x alle 2 Sekunden	Kurzschluß negative Hochspannung: Ausgangsspannung <100 V und Ausgangsstrom > 4mA für länger als 1 Minute
Positiv LED: Blink 3x alle 2 Sekunden	POS.HV- Fehler: Der Wert der Ausgangsspannung ist länger als 1 Minute größer als 120 % oder kleiner als 80 % des Zielwertes
Negativ LED: Blink 3x alle 2 Sekunden	NEG.HV- Fehler: Der Wert der Ausgangsspannung ist länger als 1 Minute größer als 120 % oder kleiner als 80 % des Zielwertes
Positiv LED: Blink 4x alle 2 Sekunden	Interlock-Kabel nicht angeschlossen
Positiv LED: Blink 4x alle 2 Sekunden	Interlock-Kabel nicht angeschlossen
Beide LED's blinken 1x alle 2 Sekunden	Die Versorgungsspannung liegt für länger als 1 Minute unter 19 Volt Gleichspannung
Keine der beiden LED's blinkt	Hochspannung AUS: Die Versorgungsspannung liegt für länger als 2 Minuten unter 15 V Gleichspannung

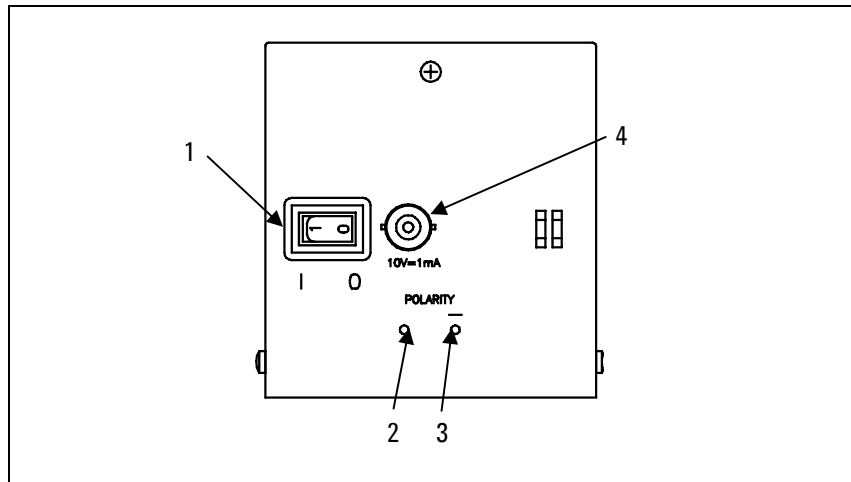


Abbildung 1 Vorderseite

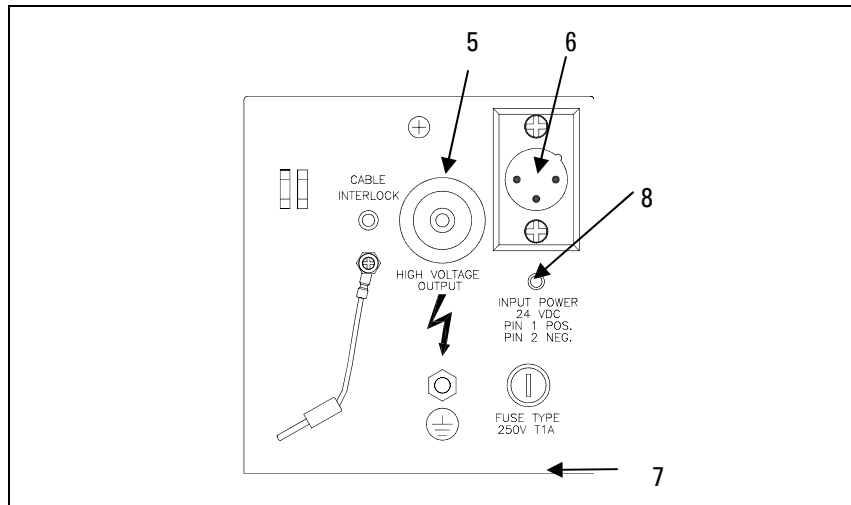


Abbildung 2 Rückseite

2 Anleitung zur Installation

Entsorgung

Tab. 3

REF.	SIGNAL	BESCHREIBUNG
1	EIN- / AUS-Schalter	Gerät und Hochspannung werden gleichzeitig ein bzw. ausgeschaltet
2	Positiv-LED	Blinkt 1x alle 2 Sekunden, wenn das Gerät und die positive Hochspannung eingeschaltet sind
3	Negativ-LED	Blinkt 1x alle 2 Sekunden, wenn das Gerät und die negative Hochspannung eingeschaltet sind
4	Recorder Ausgangs-Verbindung	0 bis 10 V linear $1\mu\text{A} = 10\text{mV}$ und $1\text{mA} = 10\text{V}$
5	Hochspannungsausgang	Zum Anschluß der Ionengetterpumpe mittels eines geeigneten Hochspannungskabels
6	GLEICHSPANNUNG-EINGANG	$24\text{V} \pm 10\%$
7	FUSE(Sicherung)	Sicherungshalter für T1A TRÄGE

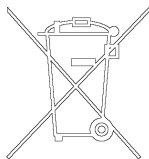
Entsorgung

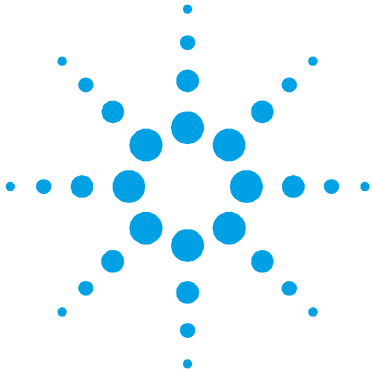
Bedeutung des "WEEE" Logos auf den Etiketten.

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.





3

Procédure pour l'installation

Description du controleur ion pump 24 V 20

Informations Generales 20

Installation 21

Mise au rebut 24

Traduction de la mode d'emploi originale



Description du contrôleur ion pump 24 V

AVERTISSEMENT!



La Haute Tension développée dans le contrôleur peut provoquer de graves lésions physiques voire la mort. Avant de mettre en fonction l'unité, débrancher le câble de l'alimentation.

Pour le branchement du contrôleur, utiliser exclusivement le câble spécial en fourniture.

ATTENTION!

Ne pas installer ou utiliser le contrôleur dans des lieux exposés aux agents atmosphériques (pluie, neige, gelée), poussières, gaz agressifs ou bien dans des lieux à risque d'explosion ou d'incendie.

NOTE

Lorsque le contrôleur est allumé, on entend un son aigu avec une intensité et une fréquence variable. Ce bruit est tout à fait normal et n'est pas signe de dysfonctionnement.

Informations Generales

Ce contrôleur a été projeté pour être utilisé par des professionnels. Avant de l'utiliser, l'opérateur doit lire le Manuel d'Instructions et tout autre document fourni par Agilent. Agilent décline toute responsabilité pour les incidents ayant lieu à cause du non respect, même partiel, de ces instructions ainsi que de l'utilisation incorrecte du contrôleur de la part du personnel non formé ou non autorisé ou de toute autre opération non conforme aux standards nationaux spécifiques.

Installation

Durant le fonctionnement, les conditions environnementales suivantes doivent être suivies:

- température: de 0 °C à +45 °C
- humidité relative: 0 – 95 %
(en absence de vapeur d'eau condensée)

Pour brancher le contrôleur à la pompe, utiliser les câbles appropriés en fourniture. Pour de plus amples renseignements concernant les branchements indiqués ci-dessus, se référer au supplément "Technical Information".

Activation/désactivation HV au moyen du panneau antérieur:

Agir sur l'interrupteur présent sur le panneau antérieur

Tab. 1

Tension de d'entrée	
24 Vcc ± 10 %	OK
< 19 Vcc pendant plus d' 1 minute	HT ON et Avertissement batterie avec allumage simultané avec lumière clignotante des deux LED (voyants) avec 1 clignotement toutes les 2 secondes
< 15 Vdc pendant plus de 2 minutes	HT OFF et absence de clignotement du LED

3 Procédure pour l'installation

Installation

LED:

Tab. 2

Led Pos: 1 clignotement toutes les 2 secondes	HT POS ON: fonctionnement normal
Led Nég: 1 clignotement toutes les 2 secondes	HT NEG ON: fonctionnement normal
Led Pos: 2 clignotements toutes les 2 secondes	HT POS Court-Circuit: tension de sortie < 100 V et courant en sortie > 4 mA pendant plus d'1 minute
Led Nég: 2 clignotements toutes les 2 secondes	HT NEG Court-Circuit: tension de sortie < 100 V et courant en sortie > 4 mA pendant plus d'1 minute
Led Pos: 3 lampeggi ogni 2 sec	HT POS hors service: la valeur de la tension de sortie est supérieure de 120 % ou inférieure de 80 % par rapport à la valeur ciblée pendant plus d' 1 minute
Led Neg: 3 clignotements toutes les 2 secondes	HT NEG hors service: la valeur de la tension de sortie est supérieure de 120 % ou inférieure de 80 % par rapport à la valeur ciblée pendant plus d' 1 minute
Led Pos: 4 3 clignotements toutes les 2 sec	Câble d'interlock non connecté
Led Neg: 4 3 clignotements toutes les 2 sec	Câble d'interlock non connecté
1 clignotement des deux LED toutes les 2 secondes	La tension d'alimentation est inférieure à 19 Vcc pendant plus de 1 minute.
Absence de clignotement des deux LED	HT OFF: la tension d'alimentation est inférieure à 15 Vcc pendant plus de 2 minutes

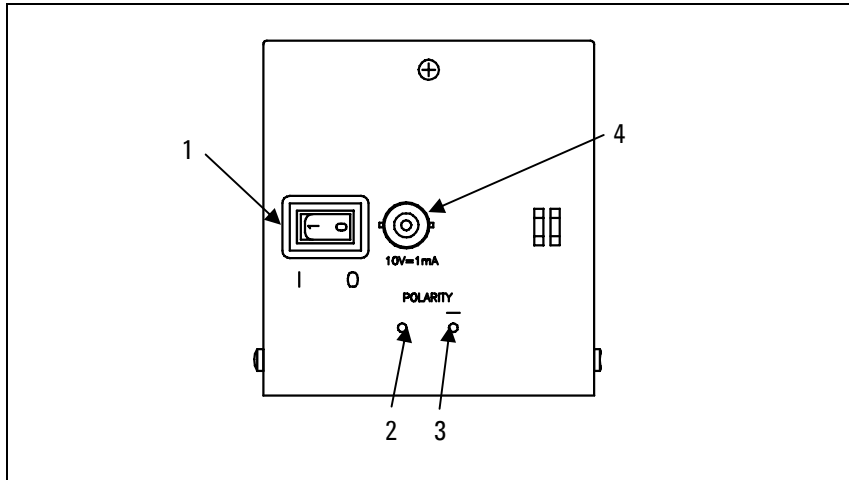


Figure 1 Panneau antérieur

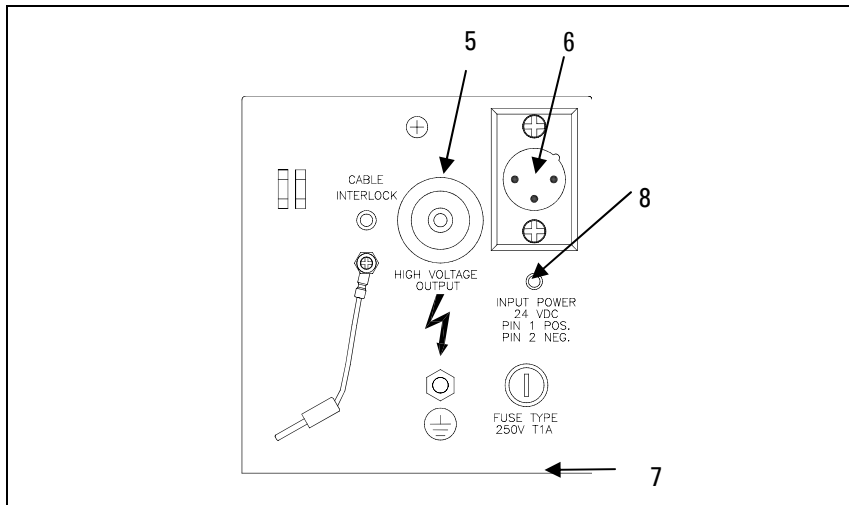


Figure 2 Panneau postérieur

3 Procédure pour l'installation

Mise au rebut

Tab. 3

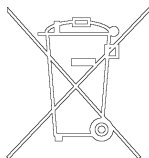
RIF.	SIGNAL	DESCRIPTION
1	ON/OFF	Il allume l'unité et l'HT
2	LED positif	1 clignotement toutes les 2 secondes lorsque l'unité et l'HT post. sont sur ON
3	LED négatif	1 clignotement toutes les 2 secondes lorsque l'unité et l'HT nég. sont sur ON
4	Recorder Output Connector	De 0 à 10 V linéaires $1\mu\text{A} = 10\text{ mV}$ et $1\text{mA} = 10\text{ V}$
5	Connecteur HV	Connecteur HT
6	DC INPUT	$24\text{ V} \pm 10\%$
7	FUSE (FUSIBLE)	SUPPORT FUSIBLE T1A SLOW

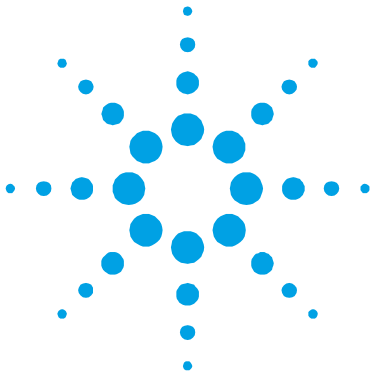
Mise au rebut

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.





4 Installation Procedure

Ion Pumps Controller 24V	26
General Information	26
Installation	27
Disposal	30

Original Instructions



Ion Pumps Controller 24V

WARNING!

High voltage developed in the controller can cause severe injury or death. Before servicing the unit, disconnect the power cable.



To connect the controller use the specified cable supplied.

CAUTION!

Do not install or use the controller in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

NOTE

The acute noise of variable intensity and frequency heard when the controller is switched on is absolutely normal and does not indicate malfunctioning.

General Information

This controller is destined for use by professionals. The user should read this instruction manual and any other additional information supplied by Agilent before operating the controller. Agilent will not be held responsible for any events occurring due to non-compliance, even partial, with these instructions, improper use by untrained persons, non-authorized interference with the controller or any action contrary to that provided for by specific national standards.

Installation

During operation, the following environmental conditions must be respected:

- temperature: from 0 °C to +45 °C
- relative humidity: 0 – 95 % (non-condensing)

To connect the controller to the pump use the specific cables supplied. See the appendix “Technical Information” for detailed information about the above mentioned and the other connections.

HV On/Off by front panel:

Move the switch on Front Panel

Tab. 1

Input Voltage	
24 Vdc ± 10 %	OK
< 19 Vdc for more of 1 minute	HV On and battery Warning with both LED Blink simultaneously 1 blink every 2 sec
< 15 Vdc for more of 2 minutes	HV OFF and no blink of LED

4 Installation Procedure

Installation

Led Indicator:

Tab. 2

Pos Led: 1 Blink every 2 sec	HV POS ON: Normal Operation
Neg Led: 1 Blink every 2 sec	HV NEG ON: Normal Operation
Pos Led: 2 Blink every 2 sec	HV POS Short Circuit: Output Voltage < 100 V and Output Current > 4 mA for more of 1 minute
Neg Led: 2 Blink every 2 sec	HV NEG Short Circuit: Output Voltage < 100 V and Output Current > 4 mA for more of 1 minute
Pos Led: 3 Blink every 2 sec	HV POS Out of Order: The value of Output Voltage is major of 120 % or minor of 80 % of the Target Value, for more of 1 minute
Neg Led: 3 Blink every 2 sec	HV NEG Out of Order: The value of Output Voltage is major of 120 % or minor of 80 % of the Target Value, for more of 1 minute
Led Pos: 4 Blink every 2 sec	Interlock cable not connected
Led Neg: 4 Blink every 2 sec	Interlock cable not connected
1 Blink of both LED every 2 sec	The supply voltage is under 19 Vdc for more of 1 minute.
No Blink of both LED	HV OFF: The supply voltage is under 15 Vdc for more of 2 minutes

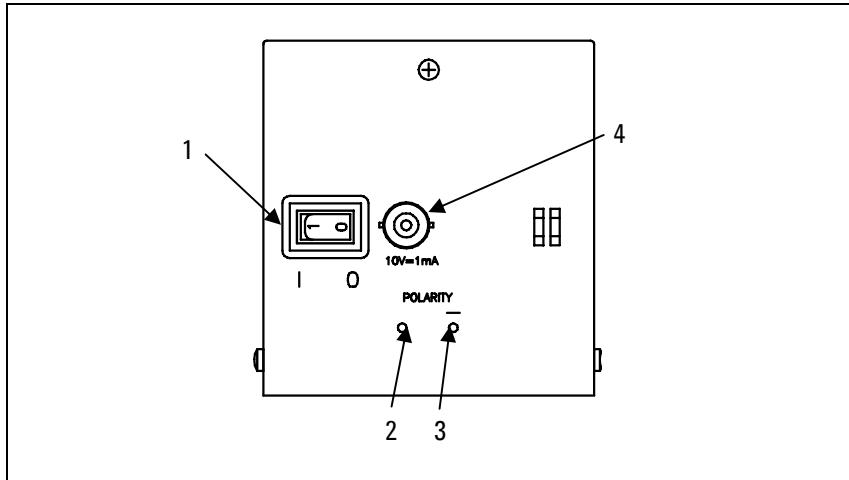


Figure 1 Front Panel

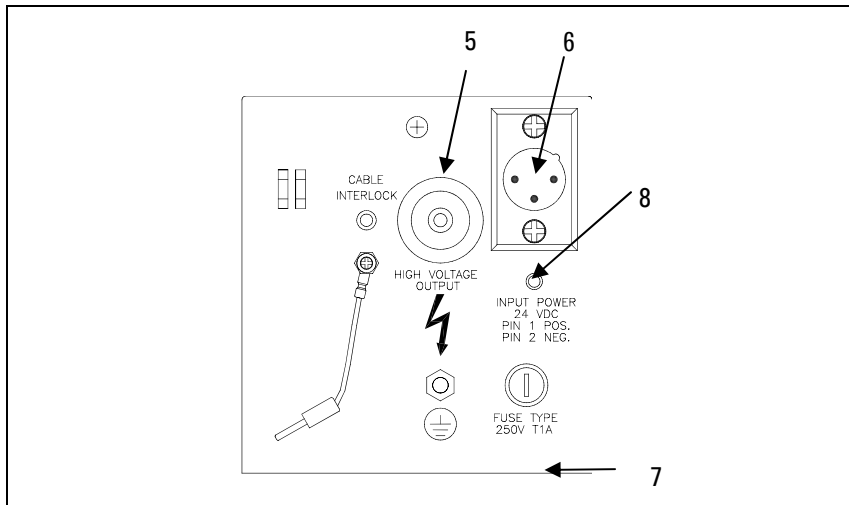


Figure 2 Rear Panel

4 Installation Procedure

Disposal

Tab. 3

REF.	NAME	DESCRIPTION
1	ON/OFF	It switches On the Unit and the High Voltage
2	Positive LED	1 blink every 2 seconds when the unit and Positive HV are ON
3	Negative LED	1 blink every 2 seconds when the unit and negative HV are ON
4	Recorder Output Connector	0 to 10 V linear 1 μ A = 10 mV and 1 mA = 10 V
5	HV Connector	HV Connector
6	DC INPUT	24 V \pm 10 %
7	FUSE	Fuse holder T1A SLOW

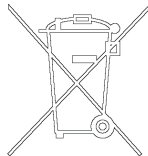
Disposal

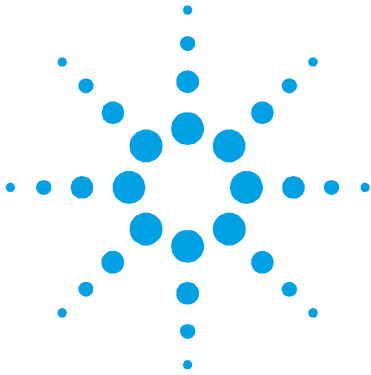
Meaning of the "WEEE" logo found in labels

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.





5 Technical Information

Controller Description	32
Controller Specifications	33
Output Polarity Selection	35
Polarity Change Procedure	35
Orderable Parts	36

Original Instructions



Controller Description

Agilent's controller is an ion pumps controller.

It can drive 1 Ion Pump. The controller is designed to give the HV to the Ion Pumps when it is connected to 24 Vdc (Max. Output Voltage = ± 3000 Vdc).

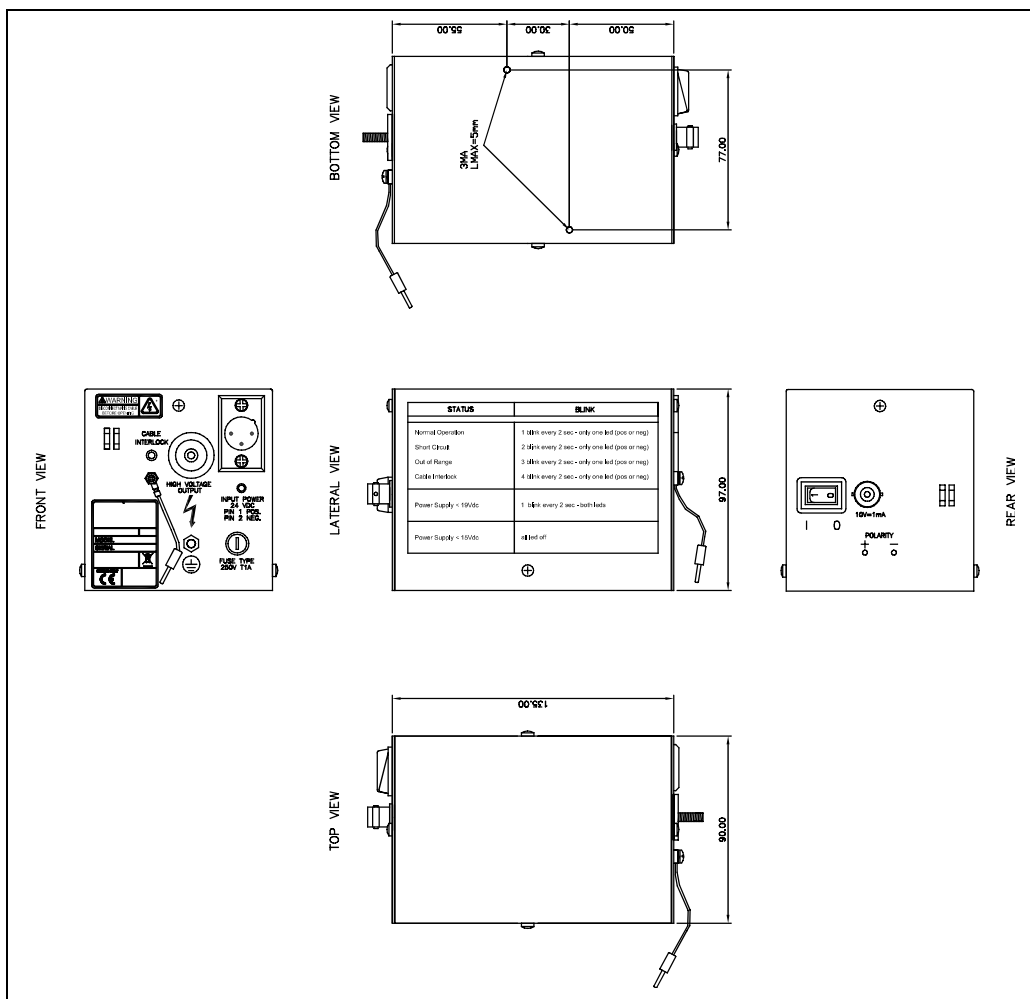


Figure 4 Controller Outline

Controller Specifications

Tab. 4

Input:	
Voltage	24 Vdc +/- 10 %
Max Power	7 W
Operating conditions: Max Altitude:	
	3000 m
HV-OUTPUT:	
Voltage	Default Value = $\pm 3000\text{ V} \pm 5\%$ (mean Value)
Ripple VPP MAX @ 3Kv	300 V
Current max	4 mA $\pm 10\%$
Maximum Output Power	4 W
Current Recorder Output signal	0 to 10 V linear 1 $\mu\text{A} = 10\text{ mV}$ and 1 mA = 10 V
Operating Temperature	0 °C to 45 °C
Storage Temperature	-10 °C to 75 °C
Compliance to Norms	EN 61010 EN 61326-1
Line fuse	T1A slow
High Voltage Connection	1 High Voltage Fischer connectors

5 Technical Information

Controller Description

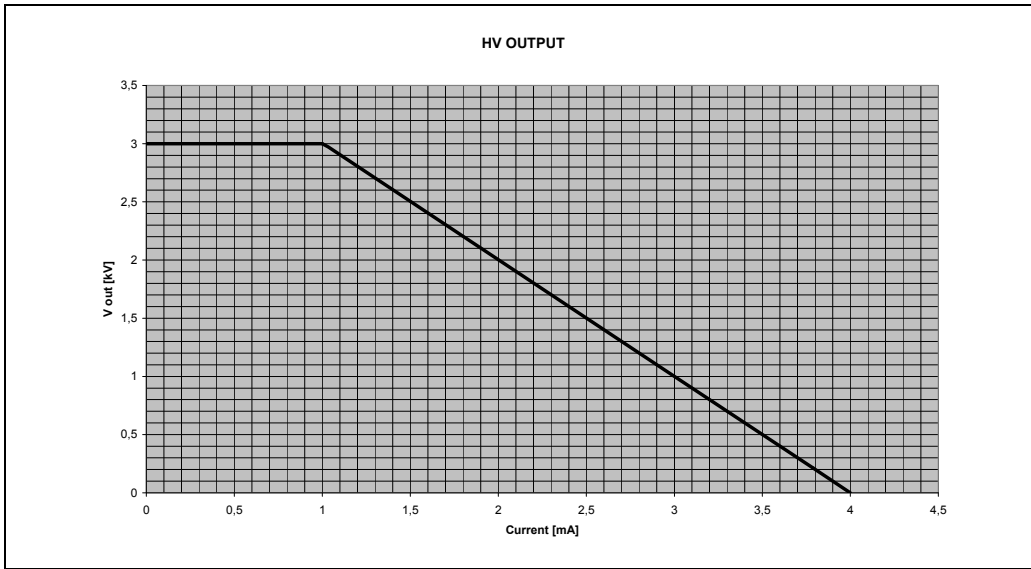


Figure 5

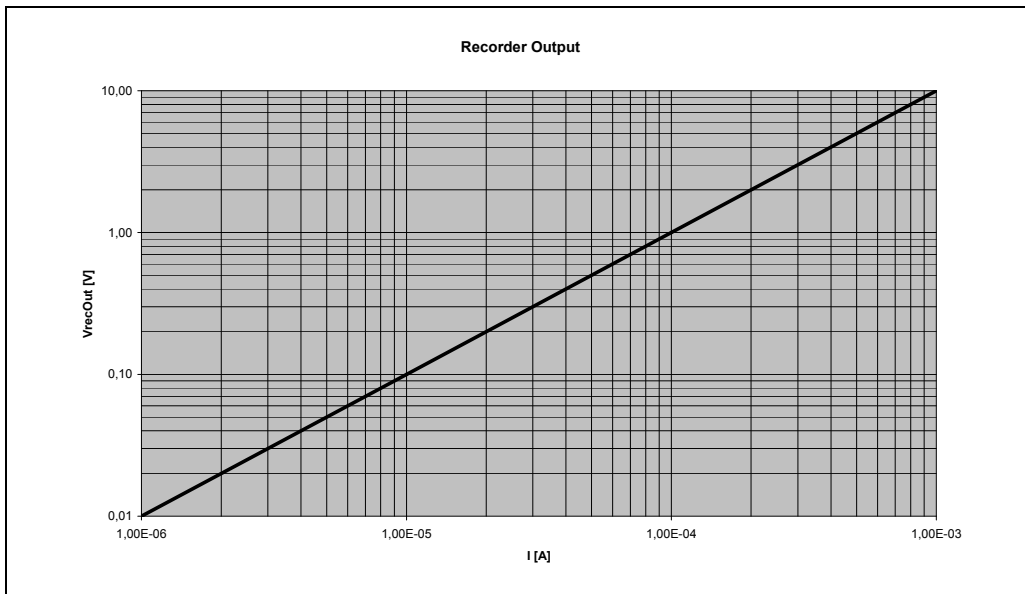


Figure 6

Tab. 5

Recorder Output Resolution:	1 μ A
Recorder Output Sensibility:	1 μ A
Recorder Output Rload >	100 k Ω
Recorder Output Ripple \max_{pp} <	10 mVpp

Output Polarity Selection

The output polarity is selected by the high voltage multiplier card.

Polarity Change Procedure

1. Switch off the power and unplug the power cord from the controller rear panel socket.
2. Undo the 4 screws and then remove the cover
3. Disconnect the H.V. white wires (1) from the multiplier board, remove the board and rotate it 180 ° horizontally; reinstall the board into the socket and reconnect the H.V. white wires.
4. Install the cover and screws.
5. Connect the power cord then switch on the unit and check the output polarity on the front panel indication.

5 Technical Information

Orderable Parts

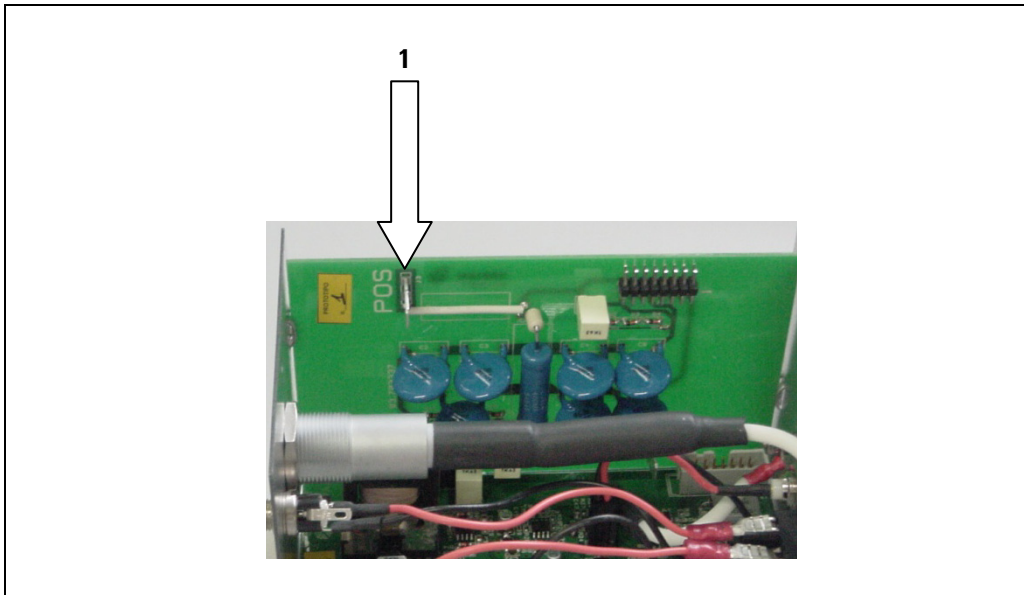


Figure 7

Orderable Parts

Tab. 6

DESCRIPTION	PART NUMBER
IPC-24 V	929-0198
HV Bakeable cable, radiation resistant, 4 m (13ft) long	929-0705
HV Bakeable cable, radiation resistant, 7 m (23ft) long	929-0707



Agilent Technologies

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,

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.

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 Leinì (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.):		
CORRECTIVE ACTION PLAN / ACTUATION (by AGILENT VPD)		DATE _____
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.

- 1) Complete the attached Request for Return form and send it to Agilent Technologies (see below), taking particular care to identify all products that have pumped or been exposed to any toxic or hazardous materials.
- 2) 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).
- 3) **Important steps for the shipment of returning product:**
 - Remove all accessories from the core product (e.g. inlet screens, vent valves).
 - Prior to shipment, 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.
 - Agilent Technologies is not responsible for returning customer provided packaging or containers.
 - **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.
- 4) Return only products for which the RA was issued.
- 5) **Product being returned under a RA must be received within 15 business days.**
- 6) **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.
- 7) Return shipments must comply with all applicable **Shipping Regulations** (IATA, DOT, 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
vpt-customer@agilent.com

NORTH AMERICA:
Fax: 1 781 860 9252
Toll Free: 800 882 7426, Option 3
vpl-ra@agilent.com

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



Please read important policy information on Page 3 that applies to all returns.

1) CUSTOMER INFORMATION

Form with fields for Company Name, Contact Name, Tel, Email, Fax, Customer Ship To, Customer Bill To, and VAT/USA/Canada tax information.

2) PRODUCT IDENTIFICATION

Table with 4 columns: Product Description, Agilent P/N, Agilent S/N, Original Purchasing Reference.

3) TYPE OF RETURN (Choose one from each row and supply Purchase Order if requesting a billable service)

- 3A. [] Non-Billable [] Billable -> New PO # (hard copy must be submitted with this form):
3B. [] Exchange [] Repair [] Upgrade [] Consignment/Demo [] Calibration [] Evaluation [] Return for Credit

4) HEALTH and SAFETY CERTIFICATION

Health and Safety Certification section containing warnings, equipment listing instructions, hazard checkboxes (Toxic, Corrosive, etc.), and signature fields.

5) FAILURE INFORMATION:

Form with fields for Failure Mode, Detailed Description of Malfunction, and Application (system and model).

Final agreement section: I understand and agree to the terms of Section 6, Page 3/3. Includes Print Name, Authorized Signature, and Date fields.



Vacuum Products Division
Request for Return Form
(Health and Safety Certification)

Please use these Failure Mode to describe the concern about the product on Page 2.

TURBO PUMPS and TURBO CONTROLLERS

Table with 3 columns: APPARENT DEFECT/MALFUNCTION, POSITION, and PARAMETERS. Includes sub-sections like OPERATING TIME.

ION PUMPS/CONTROLLERS

Table listing failure modes for Ion Pumps/Controllers such as Bad feedthrough, Vacuum leak, and Error code on display.

VALVES/COMPONENTS

Table listing failure modes for Valves/Components such as Main seal leak, Solenoid failure, and Damaged sealing area.

LEAK DETECTORS

Table listing failure modes for Leak Detectors such as Cannot calibrate, Vacuum system unstable, and Failed to start.

INSTRUMENTS

Table listing failure modes for Instruments such as Gauge tube not working, Communication failure, and Error code on display.

SCROLL AND ROTARY VANE PUMPS

Table listing failure modes for Scroll and Rotary Vane Pumps such as Pump doesn't start, Doesn't reach vacuum, and Pump seized.

DIFFUSION PUMPS

Table listing failure modes for Diffusion Pumps such as Heater failure, Doesn't reach vacuum, and Vacuum leak.

Section 6) ADDITIONAL TERMS

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.

- 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.
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.
A Special Cleaning fee will apply to all exposed products per Section 4 of this document.
If requesting a calibration service, units must be functionally capable of being calibrated.

Sales and Service Offices

United States

**Agilent Technologies
Vacuum Products Division**
121 Hartwell Avenue
Lexington, MA 02421 - USA
Tel.: +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426

Benelux

**Agilent Technologies Netherlands B.V.
Vacuum Products Division**
Herculesweg 8
4338 PL Middelburg
The Netherlands
Tel.: +31 118 671570
Fax: +31 118 671569
Toll-Free: 00 800 234 234 00

Canada

**Central coordination through: Agilent Technologies
Vacuum Products Division**
121 Hartwell Avenue
Lexington, MA 02421 - USA
Tel.: +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426

China

**Agilent Technologies (China) Co. Ltd
Vacuum Products Division**
No.3, Wang Jing Bei Lu,
Chao Yang District,
Beijing, 100102
China
Tel.: +86 (10) 6439 7718
Toll-Free: 800 820 6556

France

**Agilent Technologies France
Vacuum Products Division**
7 Avenue des Tropiques
Z.A. de Courtaboeuf - B.P. 12
91941 Les Ulis cedex - France
Tel.: +33 (0) 1 69 86 38 84
Fax: +33 (0) 1 69 86 29 88
Toll free: 00 800 234 234 00

Germany and Austria

**Agilent Technologies
Vacuum Products Division**
Alsfelder Strasse 6 Postfach 11
14 35
64289 Darmstadt – Germany
Tel.: +49 (0) 6151 703 353
Fax: +49 (0) 6151 703 302
Toll free: 00 800 234 234 00

India

**Agilent Technologies India Pvt. Ltd.
Vacuum Product Division**
G01. Prime corporate Park,
230/231, Sahar Road, Opp. Blue Dart Centre,
Andheri (East), Mumbai – 400 099.India
Tel: +91 22 30648287/8200
Fax: +91 22 30648250
Toll Free: 1800 113037

Italy

**Agilent Technologies Italia S.p.A.
Vacuum Products Division**
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

Japan

**Agilent Technologies Japan, Ltd.
Vacuum Products Division**
8th Floor Sumitomo Shibaura Building
4-16-36 Shibaura Minato-ku Tokyo 108-0023 - Japan
Tel.: +81 3 5232 1253
Fax: +81 3 5232 1710
Toll-Free: 0120 655 040

Korea

**Agilent Technologies Korea, Ltd.
Vacuum Products Division**
Shinsa 2nd Bldg. 2F 966-5 Daechi-dong
Kangnam-gu, Seoul
Korea 135-280
Tel.: +82 2 3452 2452
Fax: +82 2 3452 2451
Toll-Free: 080 222 2452

Mexico

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

Singapore

**Agilent Technologies Singapore Pte. Ltd,
Vacuum Products Division**
Agilent Technologies Building,
1 Yishun Avenue 7,
Singapore 768923
Tel : (65) 6215 8045
Fax : (65) 6754 0574

Southeast Asia

**Agilent Technologies Sales Sdn Bhd
Vacuum Products Division**
Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya,
Selangor, Malaysia
Tel : +603 7712 6106
Fax: +603 6733 8121

Taiwan

**Agilent Technologies Taiwan Limited
Vacuum Products Division (3F)**
20 Kao-Shuang Rd.,
Pin-Chen City, 324
Taoyuan Hsien , Taiwan, R.O.C.
Tel. +886 34959281
Toll Free: 0800 051 342

UK and Ireland

**Agilent Technologies UK, Ltd.
Vacuum Products Division**
6 Mead Road Oxford Industrial Park
Yarnton, Oxford OX5 1QU – UK
Tel.: +44 (0) 1865 291570
Fax: +44 (0) 1865 291571
Toll free: 00 800 234 234 00

Other Countries

**Agilent Technologies Italia S.p.A.
Vacuum Products Division**
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

Customer Support & Service

NORTH AMERICA:

Toll Free: 800 882 7426, Option 3
vpl-ra@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>

Worldwide Web Site, Catalog and Order

On-line:

www.agilent.com
Representative in most countries
12/10

© Agilent Technologies, Inc. 2011

Printed in ITALY

05/2011

Publication Number: 87-900-127-01 (C)



Agilent Technologies