

TV 141 Navigator Vent Valve

**Model
969-9834**

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

**87-900-911-01 (F)
04/2011**



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: 87-900-911-01 (F)

Edition

Edition 04/2011

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.

TV 141 Navigator Vent Valve



Contents

1 Procedura per l'installazione 7

Generalità 8

Caratteristiche tecniche 9

Installazione 11

Smaltimento 13

2 Anleitung zur Installation 15

Allgemeines 16

Technische Daten 17

Installation 19

Entsorgung 21

3 Procédure pour l'installation 23

Généralités 24

Caractéristiques techniques 25

Installation 27

4 Installation procedure 31

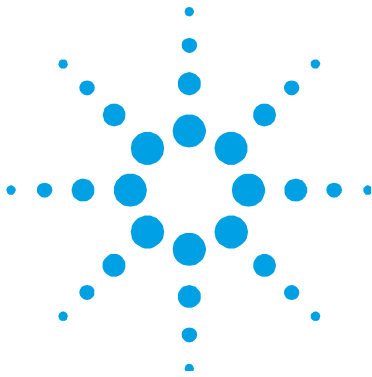
Overview 32

Technical Characteristics 33

Installation 35

Disposal 37

Contents



1 Procedura per l'installazione

Generalità	8
Caratteristiche tecniche	9
Unità di controllo	9
Cavi di connessione	9
Vent valve	10
Installazione	11
Smaltimento	13

Traduzione delle istruzioni originali



Generalità

Il kit “TV 141 Navigator Vent Valve” comprende un'unità di controllo ed una valvola, che realizzano un sistema completo per la ventilazione automatica della pompa nella fase di spegnimento o nel caso si verifichi una caduta di tensione. La valvola in condizioni di riposo (senza alimentazione) è normalmente aperta. L'attivazione avviene in modo elettromagnetico, mentre il fissaggio (Viton-sealed) viene realizzato tramite un raccordo di tipo M8 con relativo O-ring sul foro per alto vuoto della pompa. L'aria di ingresso nella valvola, viene filtrata tramite un opportuno filtro presente sull'ingresso aria della valvola stessa.

L'unità di controllo viene alimentata dal governo Turbo-V che non è predisposto per il montaggio su rack.

L'unità di controllo viene attivata con un ritardo prefissato di circa 0,8 secondi per evitare ventilazioni inopportune durante una caduta di tensione temporanea e per permettere la chiusura delle valvole di sistema prima della ventilazione.

Caratteristiche tecniche

Unità di controllo

Tensione di ingresso	24 Vdc \pm 10%
- potenza (max)	2,5 W
Tensione di uscita (max)	24 Vcc \pm 10%
- potenza (max)	1,2 W
Ritardo	circa 0,8 secondi
Temperatura operativa	da 0 a 40 °C
Temperatura di immagazzinamento	da -20 a 50 °C

Cavi di connessione

Ingresso	120 mm (4,72 inch)
Uscita dalla valvola	200 mm (7,87 inch)

1 Procedura per l'installazione

Caratteristiche tecniche

Vent valve

Stato valvola	Normalmente aperta (chiusa quando viene alimentata)
Connessione di vuoto	Raccordo M8
Filtro ingresso aria	Bronzo sinterizzato
Dimensione foro	1,2 mm (0,05")
Gamma di pressioni	da 10^{-6} mbar a 1 bar (da 10^{-7} Torr a 760 Torr)
Adattatore ingresso aria	\varnothing 6,35 mm (¼ inch)
Leak rate	$\leq 1 \times 10^{-7}$ mbar l/s
Vita	1 milione di cicli
Tensione di ingresso	24 Vdc \pm 10%
Potenza	2,5 W
Temperatura di bakeout	60 °C
Peso	140 g (0,3 lbs)

La figura seguente riporta le dimensioni di ingombro della valvola.

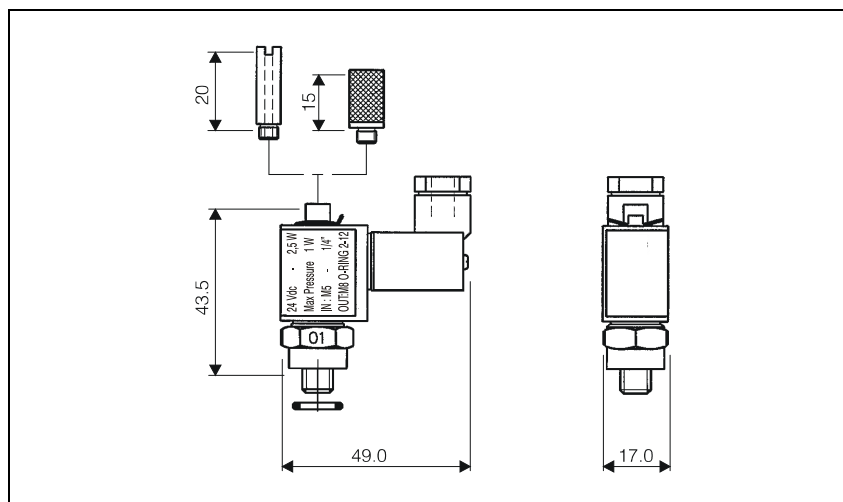


Figura 1 Dimensioni in mm

Installazione

In figura sono riportati i vari componenti presenti nel Kit TV 141 Navigator Vent Valve. Tali componenti sono forniti disassemblati; sarà quindi cura del cliente provvedere all'assemblaggio del Kit.

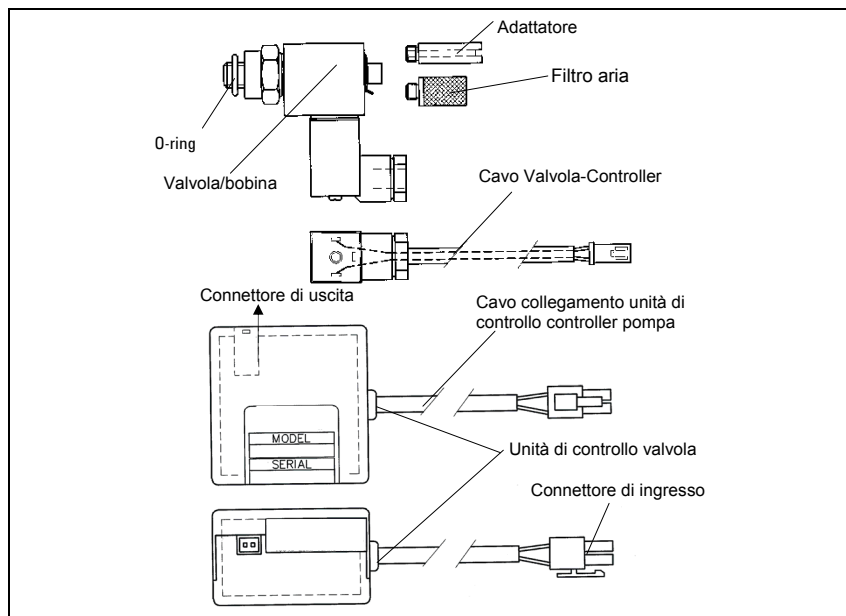


Figura 2 Componenti del Kit

ATTENZIONE! Durante la fase di assemblaggio del kit, fare attenzione a non svitare la ghiera ed il dado di fissaggio della bobina interna alla valvola.

1 Procedura per l'installazione

Installazione

Completato l'assemblaggio, procedere con l'installazione

Rimuovere il tappo a vite presente sulla pompa, quindi fissare la valvola alla pompa utilizzando una chiave esagonale da 16 mm, avendo cura di serrare il dado con una coppia pari a 2,5 Nm.

La figura seguente riporta un esempio d'installazione.

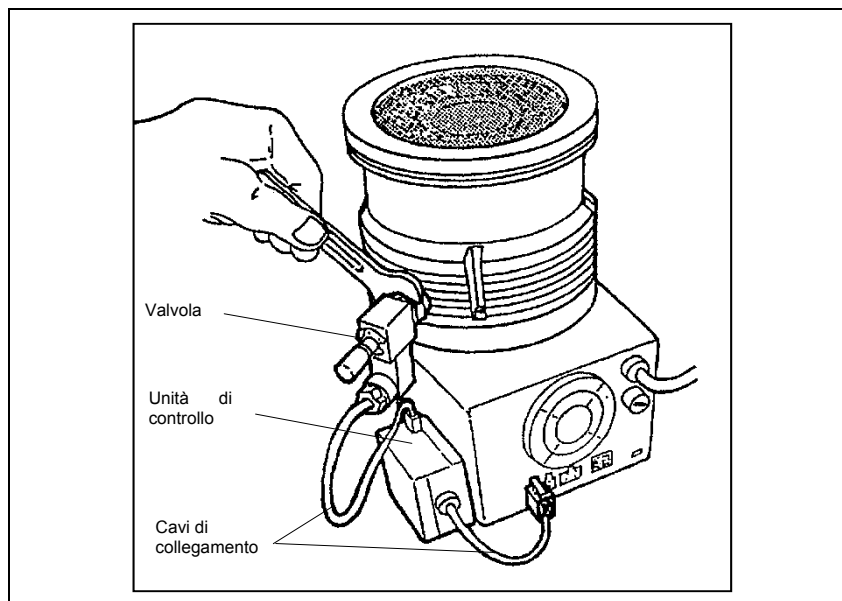


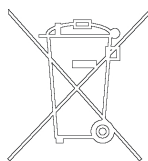
Figura 3 Installazione

ATTENZIONE! Per evitare danneggiamenti della pompa, non serrare il dado con una coppia maggiore a 2,5 Nm.

Dopo aver completato l'installazione meccanica, collegare il cavo di collegamento dalla valvola all'unità di controllo, ed il cavo di collegamento dall'unità di controllo dell'elettrovalvola al controller della pompa. Per fissare la scatola del controller, utilizzare il velcro in dotazione.

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.



1 **Procedura per l'installazione**
Smaltimento



2

Anleitung zur Installation

Allgemeines	16
Technische Daten	17
Steuereinheit	17
Anschlußkabel	17
Vent valve	18
Installation	19
Entsorgung	21

Übersetzung der Originalanleitungen



Allgemeines

Die Baugruppe "TV 141 Navigator Vent Valve" umfaßt eine Steuereinheit und ein Ventil, die ein komplettes System zur automatischen Belüftung der Pumpe beim Ausschalten oder bei Spannungsabfall darstellen. In Ruhestellung (ohne Versorgung) ist dieses Ventil normalerweise geöffnet. Es wird elektromagnetisch betätigt, während die Befestigung (Viton-sealed) durch einen Anschluß Typ M8 mit entsprechendem O-Ring an der Bohrung für Hochvakuum der Pumpe erfolgt. Die Eintrittsluft in das Ventil wird durch einen entsprechenden Filter am Lufteinlauf des Ventils gefiltert.

Die Steuereinheit wird durch eine Turbo-V-Steuerung versorgt, die nicht zur Montage auf dem Baugruppenträger vorgesehen ist.

Die Steuereinheit wird mit einer auf ca. 0,8 Sekunden voreingestellten Verzögerung eingeschaltet, um eine unerwünschte Belüftung bei zeitweisem Spannungsabfall zu verhindern und um das Schließen der Ventile des Systems vor der Belüftung zu ermöglichen.

Technische Daten

Steuereinheit

Eingangsspannung	24 Vdc \pm 10%
- Leistung (max)	2,5 W
Ausgangsspannung (max)	24 Vcc \pm 10%
- Leistung (max)	1,2 W
Verzögerung	ca. 0,8 Sekunden
Betriebstemperatur	von 0 bis 40 °C
Lagerungstemperatur	von -20 bis 50 °C

Anschlußkabel

Eingang	120 mm (4,72 inch)
Ausgang vom Ventil	200 mm (7,87 inch)

Vent valve

Ventilzustand	Normalerweise offen (bei Versorgung geschlossen)
Vakuumbuleitung	M8-Anschluß
Lufteingangsfiter	Sinterbronze
Bohrungsdurchmesser	1,2 mm (0,05")
Druckwerte	von 10^{-6} mbar bis 1 bar (von 10^{-7} Torr bis 760 Torr)
Adapter Lufttritt	\varnothing 6,35 mm (1/4 inch)
Leak rate	$\leq 1 \times 10^{-7}$ mbar l/s
Lebensdauer	1 Million Zyklen
Eingangsspannung	24 Vdc \pm 10%
Leistung	2,5 W
Bakeout-Temperatur	60 °C
Gewicht	140 g (0,3 lbs)

Aufnachstehender Abbildung sind die Abmessungen des Ventils angegeben.

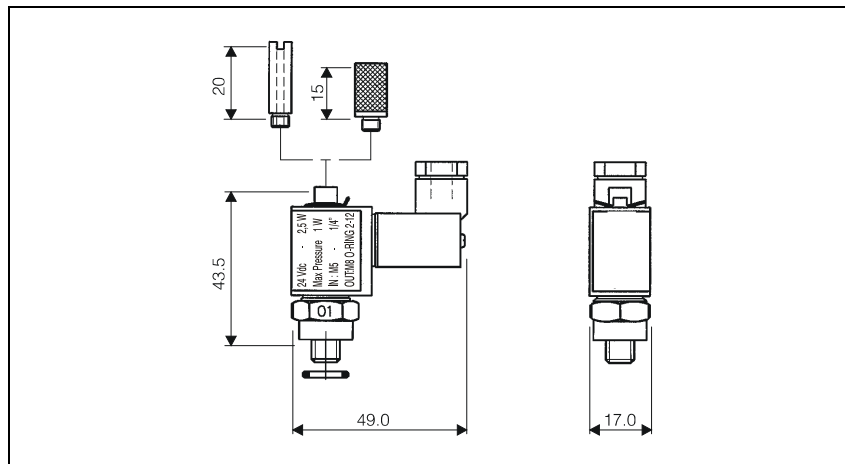


Abbildung 1 Abmessungen in mm

Installation

In der Abbildung sind die verschiedenen Bauteile der Baugruppe TV 141 Navigator Vent Valve dargestellt. Diese Bauteile werden lose geliefert und müssen daher vom Kunden zusammengesetzt werden.

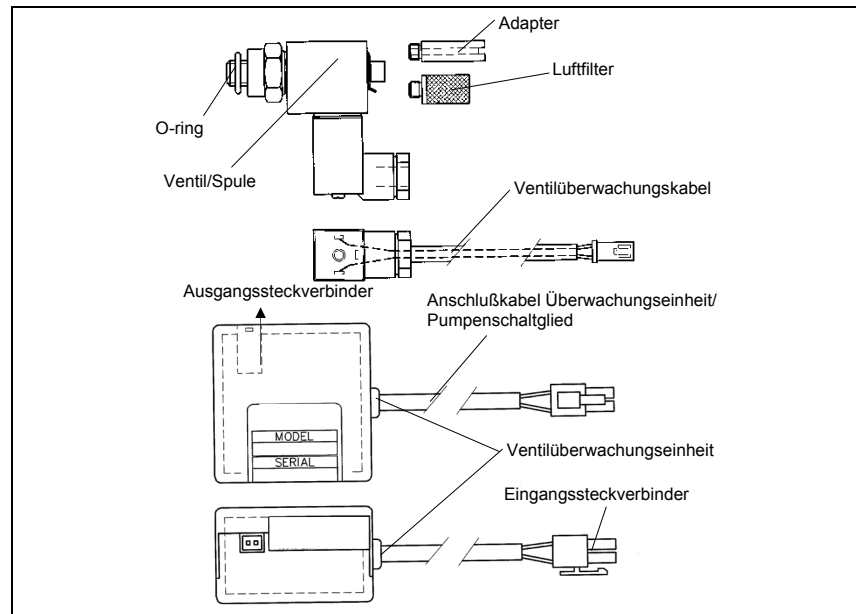


Abbildung 2 Bauteile der Baugruppe

VORSICHT!

Beim Zusammenbau der Baugruppe dürfen der Gewinding und die Befestigungsmutter der Spule im Ventilinneren nicht losgeschraubt werden.

2 Anleitung zur Installation

Installation

Nach dem Zusammenbau, Installation vornehmen.
Verschlußschraube an der Pumpe entfernen, danach Ventil unter Verwendung eines Imbusschlüssels zu 16 mm an der Pumpe befestigen, indem darauf zu achten ist, daß die Mutter mit einem Anziehmoment von 2,5 Nm festgezogen wird.

Nachfolgende Abbildung stellt ein Installationsbeispiel dar.

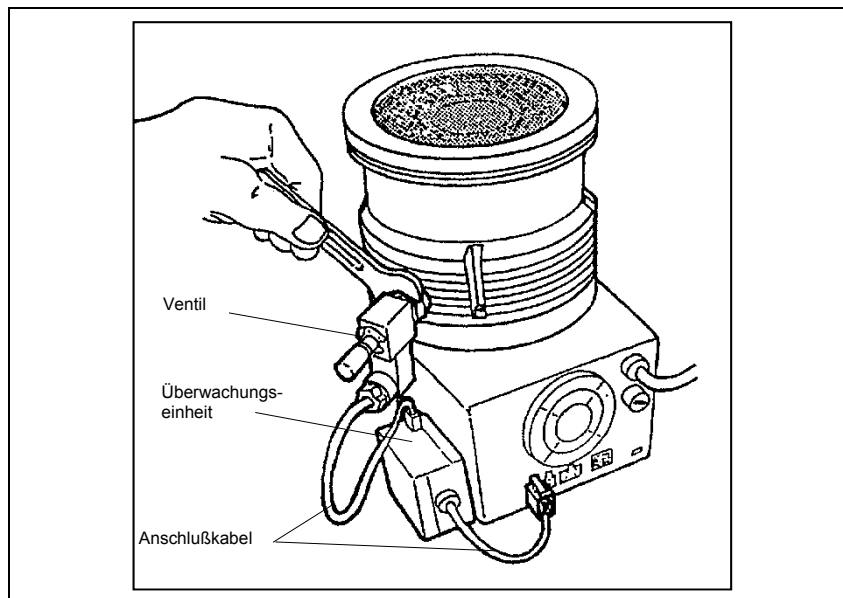


Abbildung 3 Installation

VORSICHT!

Um Beschädigungen der Pumpe zu vermeiden, die Mutter nicht mit einem Anziehmoment höher als 2,5 Nm festziehen.

Nach Abschluß der mechanischen Installation, Verbindungskabel vom Ventil zur Überwachungseinheit, sowie Verbindungskabel von der Überwachungseinheit des Elektroventils zum Schaltglied der Pumpe anschließen. Zum Befestigen des Schaltgehäuses, beigefügtes Klettband verwenden.

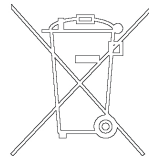
Entsorgung

Bedeutung des "WEEE" Logos auf den Etiketten.

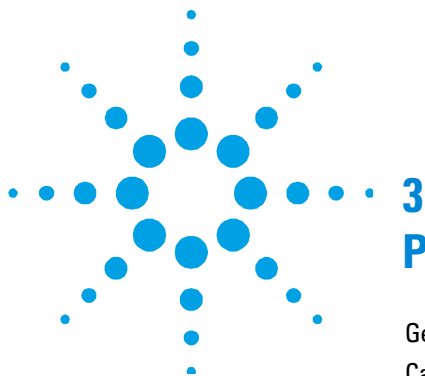
Das folgende Symbol ist in Übereinstimmung mit der EURichtlinie WEEE (Waste Electrical and Electronic Equipment) angebracht.

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.



2 **Anleitung zur Installation** **Entsorgung**



3 Procédure pour l'installation

Généralités	24
Caractéristiques techniques	25
Unité de commande	25
Câbles de connexion	25
Vent valve	26
Installation	27
Mise au Rebut	29

Traduction de la mode d'emploi originale



Généralités

Le kit "TV 141 Navigator Vent Valve" est composé d'une unité de commande et d'une valve qui forment un dispositif complet pour la ventilation automatique de la pompe pendant la phase d'extinction ou en cas de chute de tension. La valve en état de repos (non alimentée) est normalement ouverte. L'activation est réalisée d'une manière électromagnétique, alors que la fixation (Viton-sealed) est réalisée au moyen d'un adaptateur de type M8 doté d'un joint torique situé sur l'orifice pour vide poussé de la pompe. L'air en entrée de la valve est filtré au moyen d'un filtre monté sur l'entrée d'air de la valve même.

Le dispositif de commande Turbo-V, qui n'est pas conçu pour le montage sur support, alimente l'unité de commande du système.

L'unité de commande est activée avec un retard prédéfini de 0,8 secondes environ, pour éviter toute ventilation inopportune en cas de chute de tension temporaire et pour permettre la fermeture des valves de système avant la ventilation.

Caractéristiques techniques

Unité de commande

Tension en entrée	24 Vdc \pm 10%
- puissance max.	2,5 W
Tension de sortie max.	24 Vcc \pm 10%
- puissance max.	1,2 W
Retard	0,8 secondes environ
Température de service	0 à 40 °C
Température de stockage	-20 à 50 °C

Câbles de connexion

Entrée	120 mm (4,72 inch)
Sortie de la valve	200 mm (7,87 inch)

3 Procédure pour l'installation

Caractéristiques techniques

Vent valve

Etat de la valve	Normalement ouverte (fermée lorsqu'elle est alimentée)
Connexion de vide	Adaptateur M8
Filtre entrée d'air	Bronze fritté
Dimension orifice	1,2 mm (0,05")
Plage de pression	10 ⁻⁶ mbar à 1 bar (10 ⁻⁷ Torr à 760 Torr)
Adaptateur entrée d'air	Ø 6,35 mm (¼ inch)
Leak rate	≤ 1x10 ⁻⁷ mbar l/s
Durée de vie	1 million de cycles
Tension d'entrée	24 Vdc ± 10%
Puissance	2,5 W
Température de "bakeout"	60 °C
Poids	140 g (0,3 lbs)

Les dimensions d'encombrement de la valve sont indiquées sur la figure ci-après.

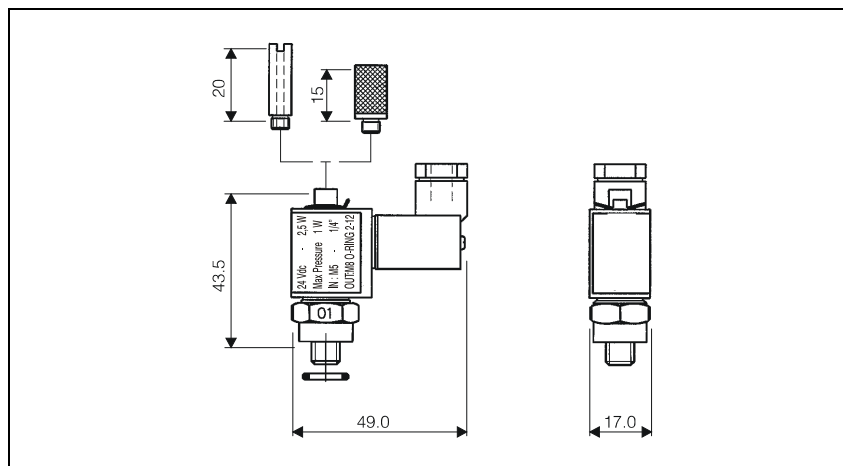


Figure 1 Dimensions en mm

Installation

Les éléments composant le kit TV 141 Navigator Vent Valve sont illustrés sur la figure ci-après. Ces éléments sont livrés désassemblés; par conséquent le client devra effectuer l'assemblage du kit.

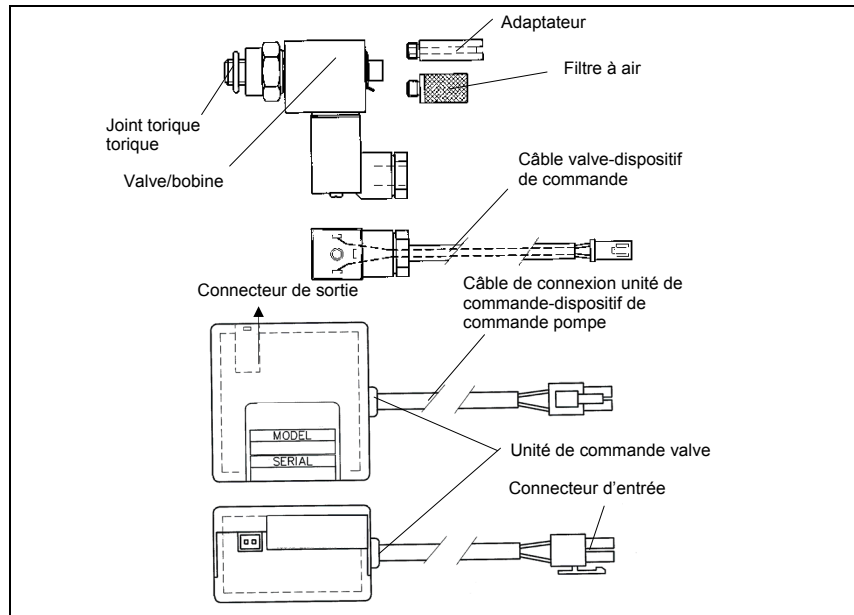


Figure 2 Eléments du kit

ATTENTION!

En cours d'assemblage du kit faire attention à ne pas dévisser la douille et l'écrou de fixation de la bobine interne de la valve.

3 Procédure pour l'installation

Installation

Après avoir terminé l'assemblage, installer le kit.

Déposer le bouchon à vis situé sur la pompe, puis fixer la valve à la pompe au moyen d'une clé à six pans de 16 mm et serrer l'écrou au couple de 2,5 Nm.

La figure ci-après illustre un exemple d'installation.

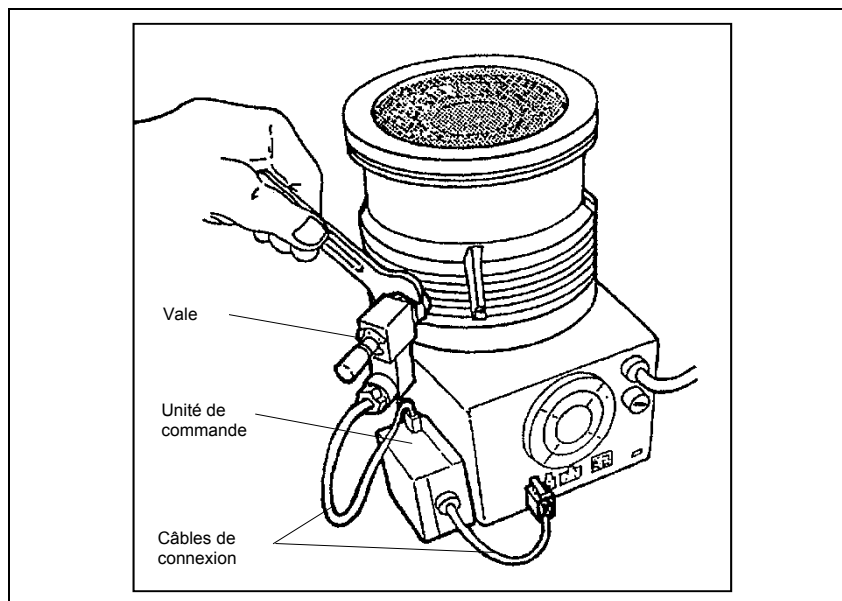


Figure 3 Installation

ATTENTION!

Pour éviter tout endommagement de la pompe ne pas serrer l'écrou à un couple supérieur à 2,5 Nm.

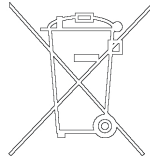
Après avoir terminé l'installation mécanique, brancher le câble de connexion de la valve sur l'unité de commande et le câble de connexion de l'unité de commande de l'électrovanne sur le dispositif de commande de la pompe. Pour fixer le boîtier du dispositif de commande utiliser le velcro livré avec le kit.

Mise au Rebut

Signification du logo "WEEE" figurant sur les étiquettes.

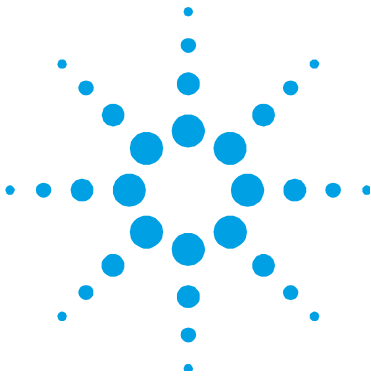
Le symbole ci-dessous est appliqué conformément à la directive CE nommée "WEEE".

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 oeuvre le processus de collecte et mise au rebut.



3 Procédure pour l'installation

Installation



4 Installation procedure

Overview	32
Technical Characteristics	33
Control Unit	33
Connection Cables	33
Vent Valve	34
Installation	35
Disposal	37

Original Instructions



4 Installation procedure

Overview

Overview

The “TV 141 Navigator Vent Valve” kit, consisting of a control unit and valve, is a complete unit for automatic pump venting when the pump is switched off or during a power failure. The valve is normally open during power off conditions, is electromechanically activated and secured (Viton-sealed) by means of an M8 thread with related O-ring on the high vacuum port. The air input into the valve is filtered by means of a specific filter located on the valve’s own air inlet.

The control unit is powered by the Turbo-V controller which is not suitable for rack mounting.

The control unit is activated with a preset delay of approximately 0.8 seconds to avoid undesired venting during a temporary power failure and to allow the closure of the system valves before venting.

Technical Characteristics

Control Unit

Input voltage	24 Vdc \pm 10%
- power (max)	2.5 W
Output voltage (max)	24 Vcc \pm 10%
- power (max)	1,2 W
Delay	Approx. 0.8 seconds
Operating temperature	0 to 40 °C
Storage temperature	-20 to 50 °C

Connection Cables

Input	120 mm (4.72 inches)
Output from valve	200 mm (7.87 inches)

Vent Valve

Valve status	Normally open (closed when power is applied)
Vacuum connection	M8 thread
Air intake filter	Sintered bronze
Hole dimension	1.2 mm (0.05 inch)
Pressure ranges	10 ⁻⁶ mbar to 1 bar (10 ⁻⁷ Torr to 760 Torr)
Air intake adapter	Ø 6.35 mm (¼ inch)
Leak rate	≤ 1x10 ⁻⁷ mbar l/s
Life span	1 million cycles
Input voltage	24 Vdc ± 10%
Power	2.5 W
Bakeout temperature	60 °C
Weight	140 g (0.3 lbs)

The following figure shows the valve's overall dimensions.

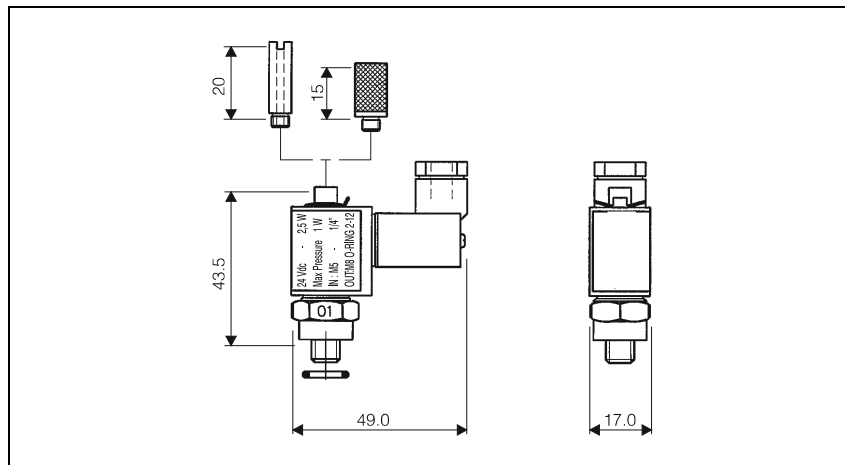


Figure 1 Dimensions in mm

Installation

The following figure shows the various components of the TV 141 Navigator Vent Valve kit. These components come disassembled and it is therefore up to the customer to assemble the kit.

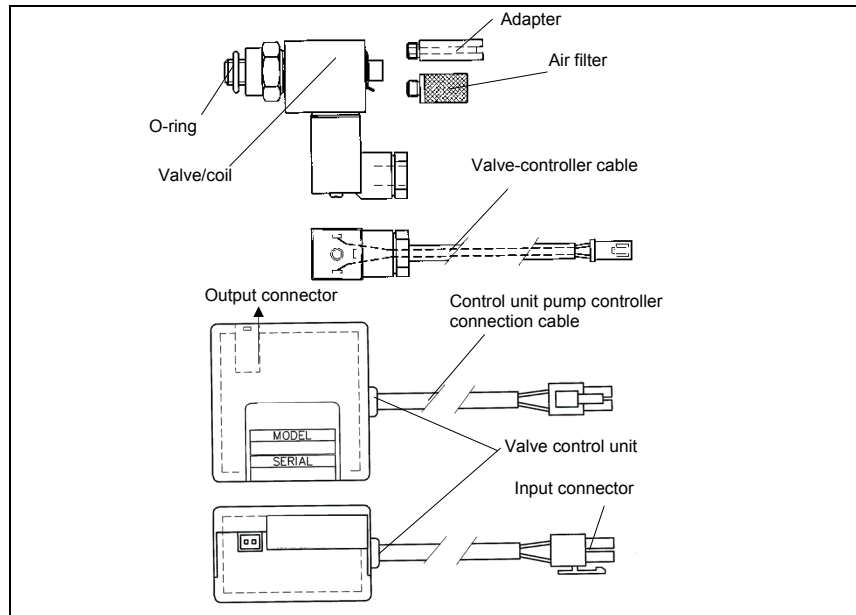


Figure 2 Kit Components

CAUTION!

When assembling the kit be careful not to unscrew the coil securing ring and nut inside the valve.

4 Installation procedure

Installation

Once the kit is assembled, install it on the pump.

Remove the screw cap from the pump, then secure the valve to the pump using a 16 mm wrench making sure to tighten the nut with a torque equivalent to 2.5 Nm.

The following figure shows an example of installation.

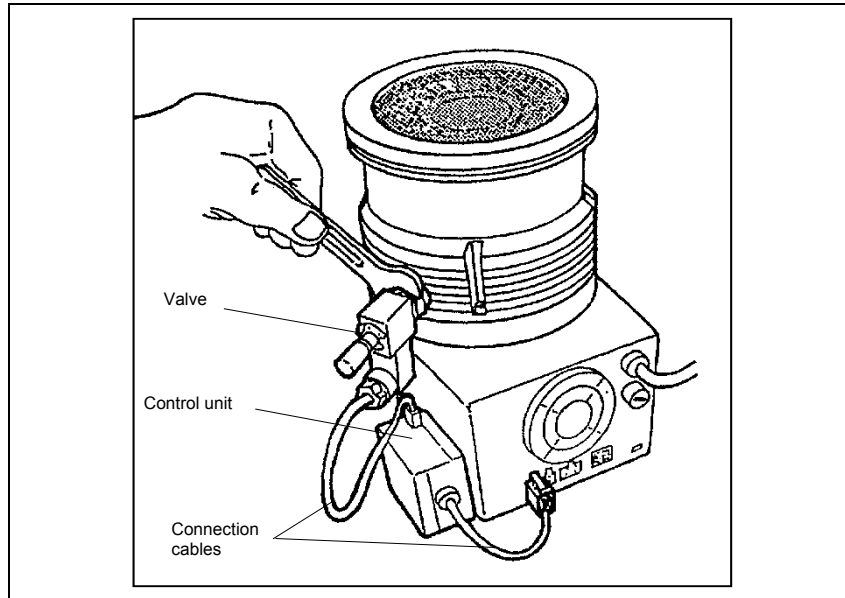


Figure 3 Installation

CAUTION!

To avoid damaging the pump, do not tighten the nut with a torque greater than 2.5 Nm.

Upon completion of the mechanical installation, attach the connection cable between the valve to the control unit, and the connection cable between the electrovalve control unit to the pump's controller. Use the velcrose provided to secure the controller box.

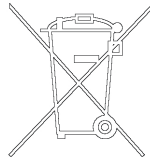
Disposal

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.



4 Installation procedure

Disposal



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.): _____ _____ _____ 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.

- 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. Lists various failure modes like 'Does not start', 'Noise', 'Vertical', 'Horizontal', etc.

ION PUMPS/CONTROLLERS

Table listing failure modes for Ion Pumps/Controllers: Bad feedthrough, Vacuum leak, Error code on display, Poor vacuum, High voltage problem, Other.

VALVES/COMPONENTS

Table listing failure modes for Valves/Components: Main seal leak, Solenoid failure, Damaged sealing area, Bellows leak, Damaged flange, Other.

LEAK DETECTORS

Table listing failure modes for Leak Detectors: Cannot calibrate, Vacuum system unstable, Failed to start, No zero/high background, Cannot reach test mode, Other.

INSTRUMENTS

Table listing failure modes for Instruments: Gauge tube not working, Communication failure, Error code on display, Display problem, Degas not working, Other.

SCROLL AND ROTARY VANE PUMPS

Table listing failure modes for Scroll and Rotary Vane Pumps: Pump doesn't start, Doesn't reach vacuum, Pump seized, Noisy pump (describe), Over temperature, Other.

DIFFUSION PUMPS

Table listing failure modes for Diffusion Pumps: Heater failure, Doesn't reach vacuum, Vacuum leak, Electrical problem, Cooling coil damage, Other.

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

04/2011

Publication Number: 87-900-911-01 (F)



Agilent Technologies