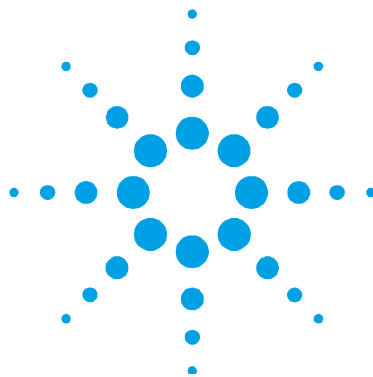




Cryopanel

Model
919-0180

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

87-900-078-01 (E)
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: 87-900-078-01 (E)

Edition

Edition 05/2011

Printed in ITALY

Agilent Technologies Italia S.p.A.

Vacuum Products Division

Via F.Ili 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

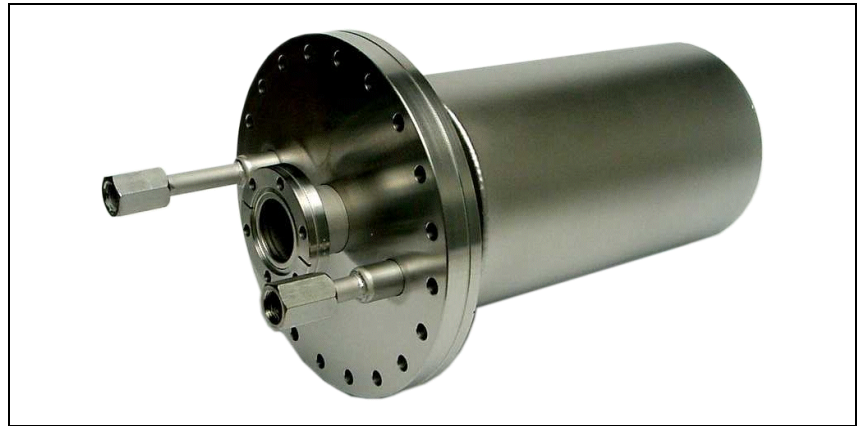
CAUTION

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

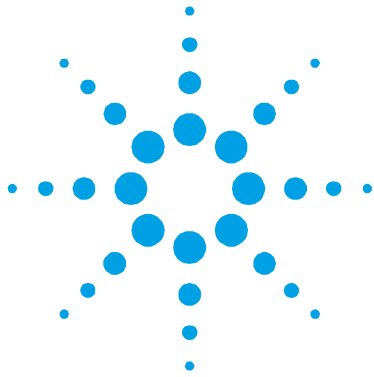
A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Cryopanel



Contents

1	Procedura per l'installazione	7
	Informazioni generali	8
	Generalità sul TSP Cryopanel Assy	9
	Smaltimento	15
2	Anleitung zur Installation	17
	Allgemeine Hinweise	18
	Allgemeine Angaben zum TSP Cryopanel	19
	Entsorgung	25
3	Procédure pour l'installation	27
	Indications Generales	28
	Generalites sur le TSP Cryopanel Assy	29
	Mise au rebut	35
4	Installation procedure	37
	General Information	38
	Overview on the TSP Cryopanel Assy	39
	Disposal	45



1

Procedura per l'installazione

Informazioni generali	8
Generalità sul TSP Cryopanel Assy	9
Specifiche Tecniche	10
Installazione del TSP Cryopanel	11
Installazione della Cartuccia TSP	12
Smaltimento	15

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.

Questo manuale utilizza le seguenti convenzioni:

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.

ATTENZIONE!

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

NOTA

Le note contengono informazioni importanti estrapolate dal testo.

Generalità sul TSP Cryopanel Assy

Il TSP Cryopanel è stato progettato per essere utilizzato in accoppiamento con la TSP (Titanium Sublimation Pump), tramite montaggio su una flangia 8" OD ConFlat®. Per il raffreddamento può essere utilizzata acqua o azoto liquido. Inoltre il sistema può funzionare in modalità UHV senza raffreddamento. Il Cryopanel può essere installato su pompe ioniche doubleended o sideported.

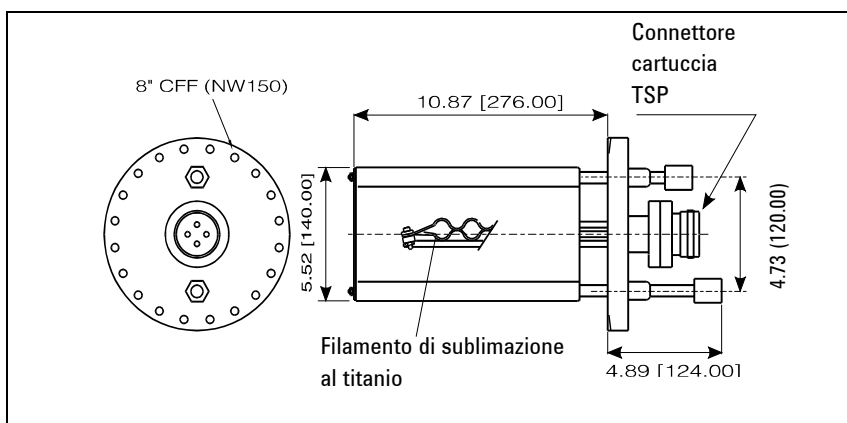


Figura 1 Ingombri TSP Cryopanel in mm [pollici]

Specifiche Tecniche

Tab. 1

	N ₂	H ₂	H ₂ O
Velocità di pompaggio a 20 °C con TSP e raffreddamento ad acqua (l/s)	516	1205	578
Velocità di pompaggio a -195 °C con TSP e raffreddamento ad LN ₂ (l/s)	555	1760	696*
Superficie di pompaggio interna (cm ²)	826		
Flangia principale	8" CF (NW 150)		
Vol. del serbatoio (litri)	1.8		
Raccordi di raffreddamento	3/8" Gas		
Flangia sorgente al titanio	2¾ CF		

* Tenendo presente che H₂O è un fluido condensabile, l'effetto del pompaggio sulla superficie di uscita del cryopanel è compreso. La velocità di pompaggio è 12.000 l/sec.

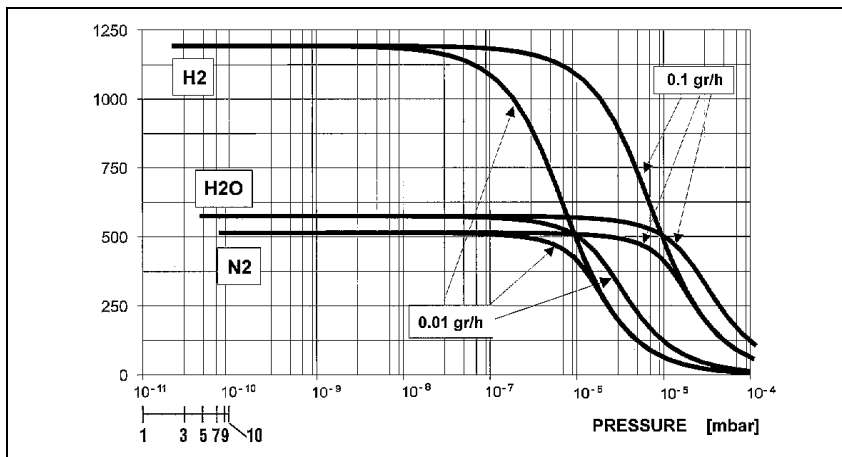


Figura 2 Velocità di pompaggio a 20 °C con TSP e raffreddamento ad acqua

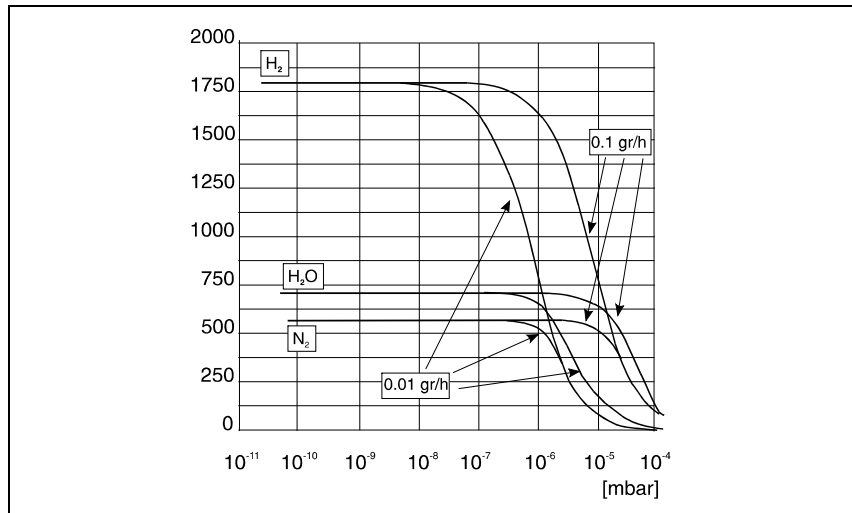


Figura 3 Velocità di pompaggio a -195 °C con TSP e raffreddamento ad azoto liquido

Installazione del TSP Cryopanel

AVVERTENZA!



Se per il raffreddamento viene utilizzato LN₂, prestare molta attenzione nel maneggiare il TSP Cryopanel durante le fasi d'installazione, in quanto il contatto del LN₂ con la pelle provoca gravi ustioni.

Per eseguire una corretta installazione del TSP Cryopanel, attenersi fedelmente alla procedura riportata di seguito.

- 1 Pulire accuratamente sia la flangia da 8" ConFlat® del Cryopanel, che la relativa flangia presente sulla pompa.
- 2 Posizionare i dadi ed avvitare ciascuno con una coppia da 6 – 11 Nm (4.5 - 8 ft.-lbs). Dopo aver avvitato un dado, avvitare sempre quello dalla parte opposta rispetto al centro della flangia.

1 Procedura per l'installazione

Generalità sul TSP Cryopanel Assy

- 3 Ripetere questa sequenza di avvitamento per altri due cicli.
- 4 Continuare ad avvitare i dadi fino a che i lati delle flange si accostano e fino a rilevare un incremento di coppia.
- 5 Fissare i condotti del fluido di raffreddamento ai relativi attacchi presenti sul TSP Cryopanel, avendo cura di stringere in modo adeguato i dadi di bloccaggio.

Installazione della Cartuccia TSP

Normalmente la cartuccia TSP viene fornita con una flangia OD 2 $\frac{3}{4}$ " ConFlat® con tre filamenti al titanio-molibdeno; ciascuno bakeable (riscaldabile) a 400 °C. La sublimazione massima si ottiene con l'erogazione di 300 W di potenza.

Per il fissaggio della cartuccia sul Cryopanel, procedere come descritto nel paragrafo precedente, facendo attenzione a fissare la flangia in modo tale da non creare interferenze tra il connettore del cavo proveniente dall'unità di controllo ed i condotti di raffreddamento.

NOTA

Durante la fase d'installazione, occorre rispettare il verso di circolazione del fluido di raffreddamento, come riportato negli esempi seguenti.

ATTENZIONE!

Prestare attenzione in fase di rimozione ai liquidi di raffreddamento dal Cryopanel prima del baking e tenere aperto il collegamento del tubo di raffreddamento durante il baking.

Le figure riportano alcuni esempi di posizionamento del TSP Cryopanel.

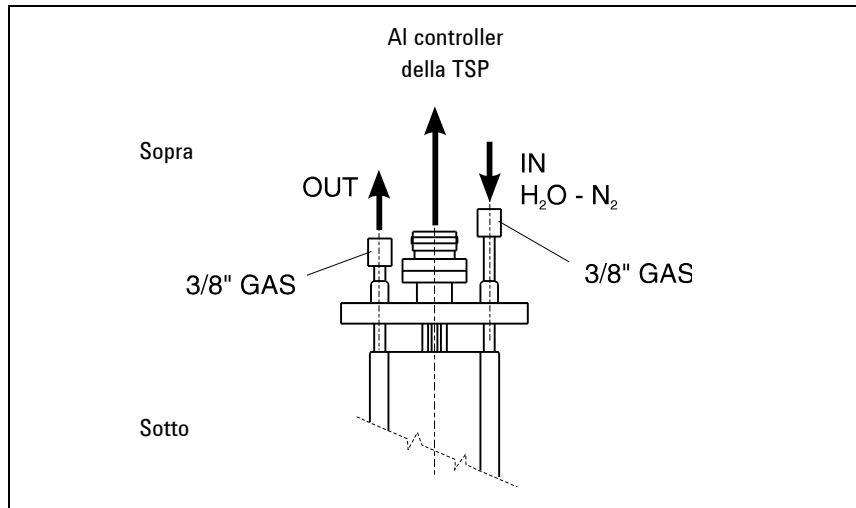


Figura 4 Esempio di posizionamento verticale

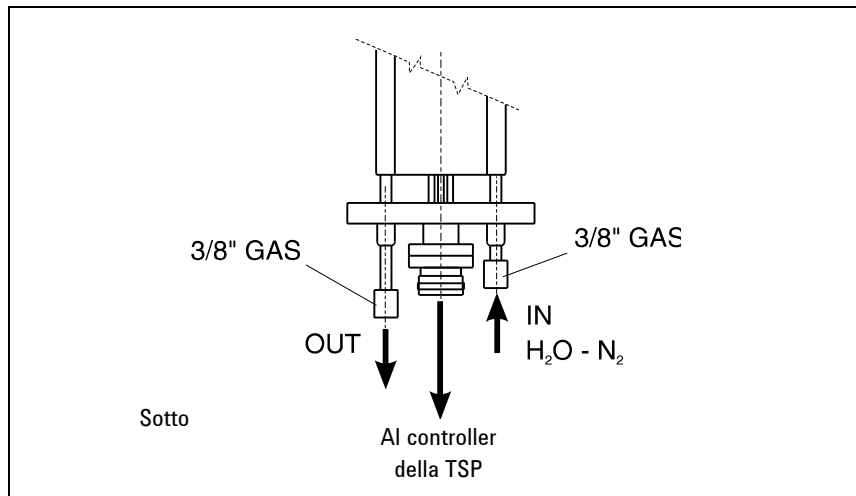


Figura 5 Esempio di posizionamento verticale

1 Procedura per l'installazione

Generalità sul TSP Cryopanel Assy

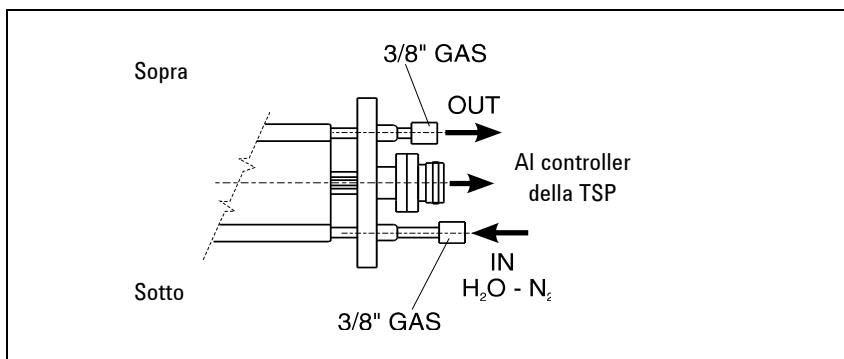


Figura 6 Esempio di posizionamento orizzontale

- 1 Collegare il cavo proveniente dal Titanium Sublimation Controller al relativo connettore presente al centro del TSP Cryopanel.

AVVERTENZA!



L'elevata corrente che transita nel cavo di collegamento all'unità di controllo potrebbe causare gravi infortuni o pericolo di morte. Prima di collegare il cavo al connettore della cartuccia TSP, oppure prima di scollegarlo, assicurarsi di aver spento l'unità di controllo.

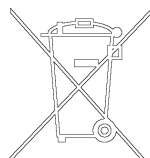
Smaltimento

Significato del logo "WEEE" presente sulle etichette.

Il simbolo qui sotto riportato applicato in ottemperanza alla direttiva CE denominata "WEEE".

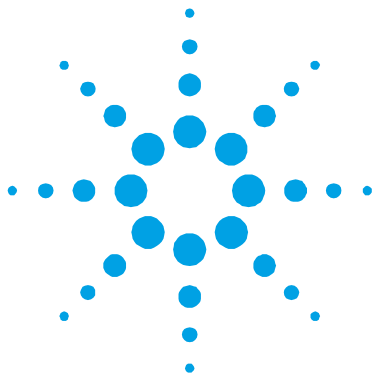
Questo simbolo (**valido solo per i paesi della Comunità Europea**) indica che il prodotto sul quale è applicato, NON deve essere smaltito insieme ai comuni rifiuti domestici o industriali, ma deve essere avviato ad un sistema di raccolta differenziata.

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

Allgemeine Hinweise	18
Allgemeine Angaben zum TSP Cryopanel	19
Technische Daten	20
Installation des TSP Cryopanel	21
Installation des TSP-Einsatzes	22
Entsorgung	25

Übersetzung der Originalanleitungen



Allgemeine Hinweise

Dieses Gerät ist für den professionellen Gebrauch bestimmt. Vor dem Gebrauch soll der Benutzer dieses Handbuch sowie alle weiteren von Agilent mitgelieferten Zusatzdokumentationen genau lesen. Bei vollständiger bzw. teilweiser Nichtbeachtung der enthaltenen Hinweise, unsachgemäßem Gebrauch durch ungeschultes Personal, nicht autorisierten Eingriffen und Mißachtung der nationalen Bestimmungen übernimmt Firma Agilent keinerlei Haftung.

In dieser Gebrauchsanleitung werden Sicherheitshinweise folgendermaßen hervorgehoben:

WARNUNG!



Die Warnhinweise lenken die Aufmerksamkeit des Bedieners auf eine spezielle Prozedur oder Praktik, die bei unkorrekter Ausführung schwere Verletzungen hervorrufen könnte.

VORSICHT!

Die Vorsichtshinweise vor bestimmten Prozeduren machen den Bediener darauf aufmerksam, daß bei Nichteinhaltung Schäden am Gerät entstehen können.

HINWEIS

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

Allgemeine Angaben zum TSP Cryopanel

Der TSP Cryopanel wurde entwickelt, um auf einen Flansch 8" OD ConFlat® montiert und in Verbindung mit TSP (Titanium Sublimation Pump) verwendet zu werden. Das Cryopanel ist für den Betrieb im UHV vorgesehen und kann sowohl mit Wasser oder flüssigen Stickstoff gekühlt, als auch ungekühlt eingesetzt werden. Das Cryopanel kann an Ionengetterpumpen mit zusätzlichen Flansch am Boden (doubleended) oder an der Seite (side port) installiert werden.

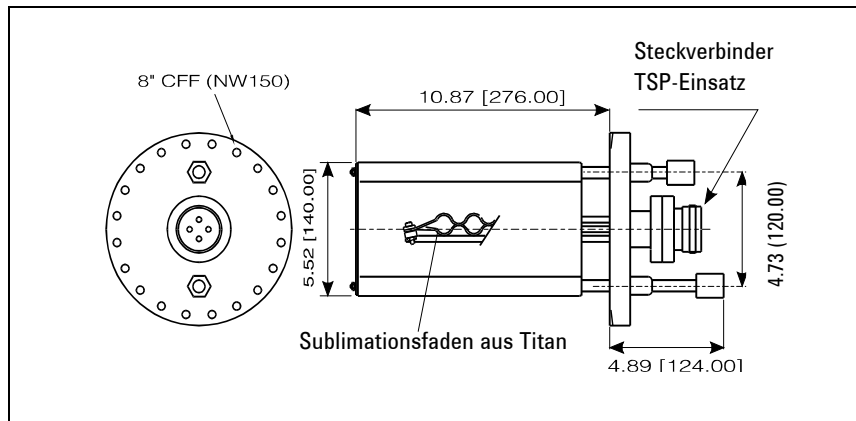


Abbildung 1 Abmessungen TSP Cryopanel in mm [Zoll]

Technische Daten

Tab. 1

	N ₂	H ₂	H ₂ O
Pumpgeschwindigkeit bei 20 °C mit TSP und Wasserkühlung (l/s)	516	1205	578
Pumpgeschwindigkeit bei -195 °C mit TSP und Kühlung mit LN ₂ (l/s)	555	1760	696*
Interne Pumpoberfläche (cm ²)	826		
Hauptflansch	8" CF (NW 150)		
Behälterinhalt (Liter)	1.8		
Kühlanschlüsse	3/8" Gas		
Flansch der Titanquelle	2 3/4 CF		

* Da H₂O eine kondensierbare Flüssigkeit ist, berücksichtigt dieser Wert die Pumpwirkung auf die Ausgangsfläche des Cryopanel. Die Pumpgeschwindigkeit beträgt 12.000 l/sec.

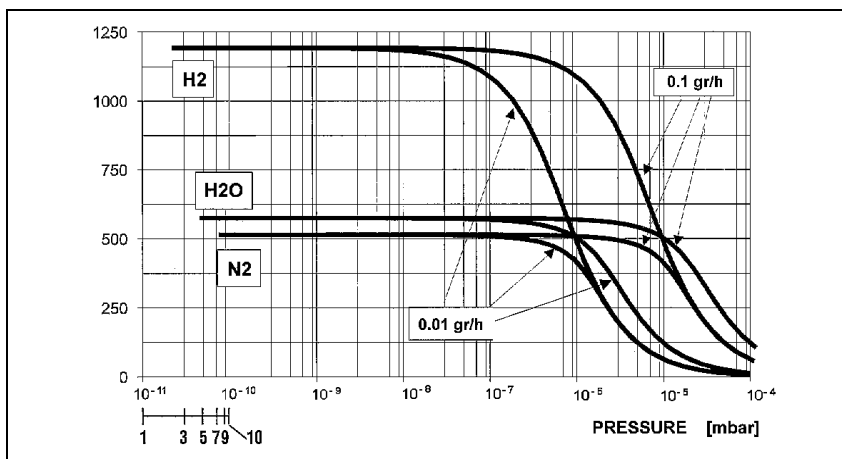


Abbildung 2 Pumpgeschwindigkeit bei 20 °C mit TSP und Wasserkühlung

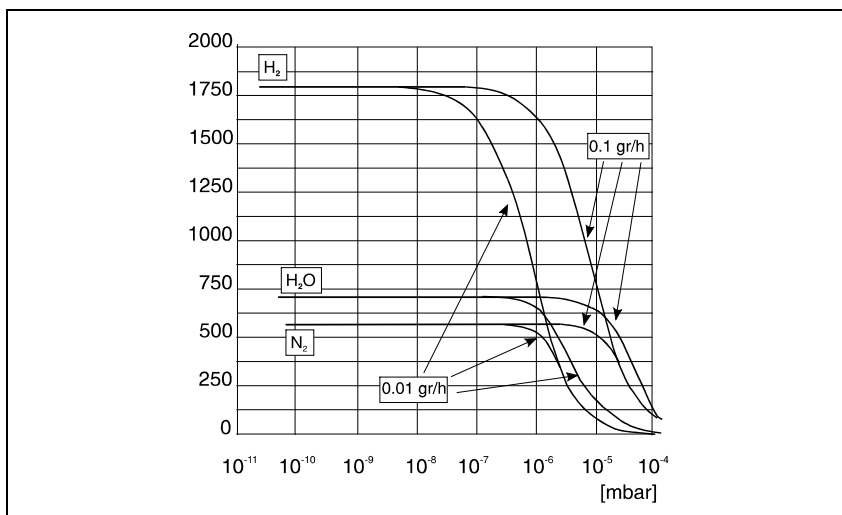


Abbildung 3 Pumpgeschwindigkeit bei -195 °C mit TSP und Kühlung mit flüssigem Stickstoff

Installation des TSP Cryopanel

WARNUNG!



Wird zur Kühlung LN₂ verwendet, ist bei der Installation des TSP Cryopanel besondere Vorsicht geboten, da LN₂ schwere Verbrennungen verursacht, wenn es mit der Haut in Berührung kommt.

Zur richtigen Installation des TSP Cryopanel sind die nachstehende Anleitungen genau zu befolgen:

- 1 Den 8" ConFlat® Flansch des Cryopanel und den pumpseitigen Flansch sorgfältig reinigen.

2 Anleitung zur Installation

Allgemeine Angaben zum TSP Cryopanel

- 2 Die Muttern einsetzen und mit einem Drehmoment von jeweils 6 – 11 Nm (4.5 - 8 ft.-lbs) anschrauben. Sobald eine Mutter eingeschraubt ist, muß danach immer die der Flanschmitte gegenüberliegende Mutter eingeschraubt werden.
- 3 Diesen Vorgang zweimal wiederholen.
- 4 Die Muttern soweit anziehen, bis sich die Flansche berühren und eine Erhöhung des Drehmoments festzustellen ist.
- 5 Die Leitungen der Kühlflüssigkeit mit den Anschlüssen am TSP Cryopanel verbinden und die Befestigungsmuttern entsprechend anziehen.

Installation des TSP-Einsatzes

Normalerweise wird der TSP-Einsatz mit einem Flansch OD 2¾" ConFlat® mit drei Titan-Molybdän-Fäden geliefert, die bis 400 °C aufheizbar (bakeable) sind. Die größte Sublimation erfolgt bei 300 W Leistung.

Die Befestigung des Einsatzes am Cryopanel ist entsprechend der im vorherigen Abschnitt enthaltenen Beschreibung vorzunehmen. Dabei ist da-rauf zu achten, dass der Flansch so angebracht wird, dass sich der Stecker des von der Steuereinheit kommenden Kabels und die Kühlleitungen sich nicht gegenseitig behindern.

HINWEIS

Bei der Installation muß die Strömungsrichtung der Kühlflüssigkeit beachtet werden, wie in den folgenden Beispielen dargestellt.

VORSICHT!

Es ist darauf zu achten, daß evtl. vorhandene Kühlflüssigkeiten vor dem Ausheizen aus dem Cryopanel entfernt werden und Kühlleitungen während des Ausheizvorgangs nicht verschlossen sind.

Die folgenden Abbildungen enthalten einige Beispiele für die Anordnung des TSP Cryopanel.

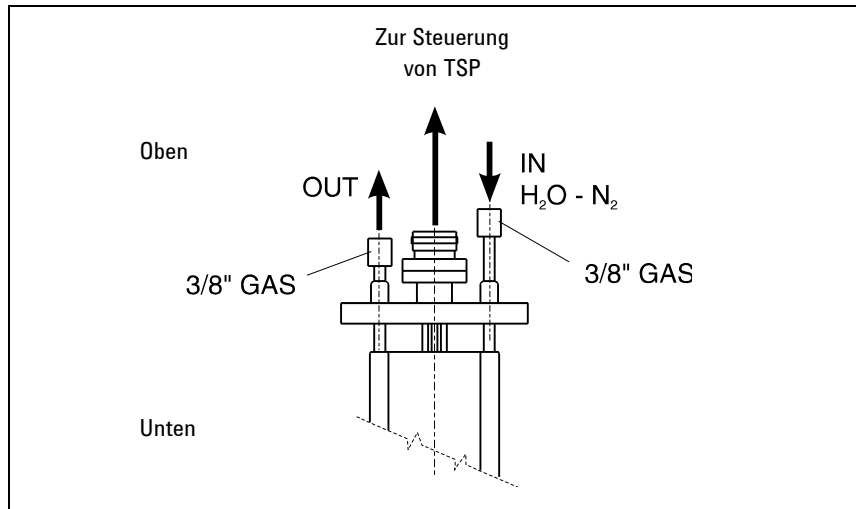


Abbildung 4 Beispiel für senkrechte Anordnung

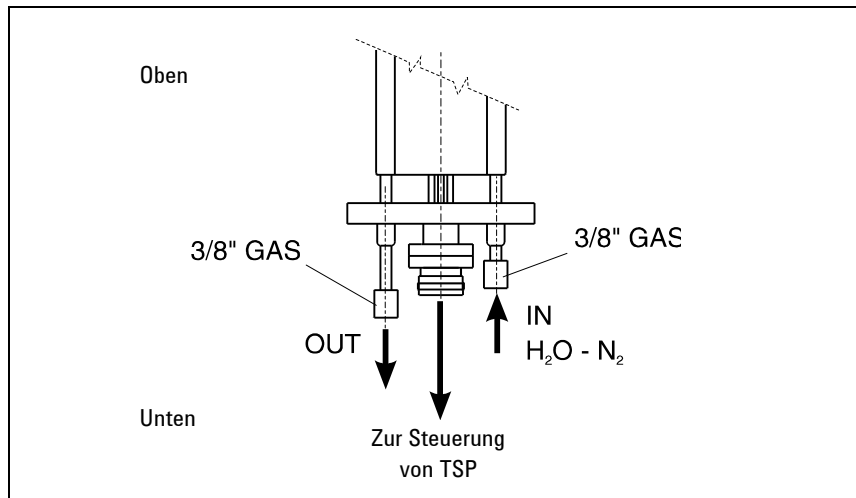


Abbildung 5 Beispiel für senkrechte Anordnung

2 Anleitung zur Installation

Allgemeine Angaben zum TSP Cryopanel

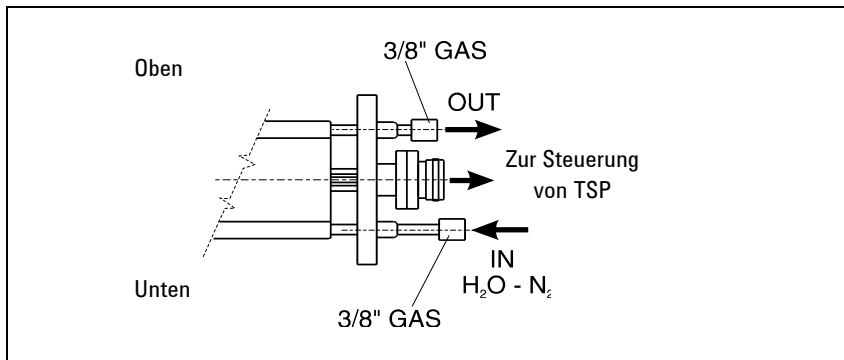


Abbildung 6 Beispiel für waagrechte Anordnung

- 1 Das vom Titanium Sublimation Controller kommende Kabel an den entsprechenden Steckverbinder auf dem TSP Cryopanel anschließen.

WARNUNG!



Durch den starken Strom, der durch das Anschlusskabel der Steuereinheit fließt, besteht große Unfall- bzw. Todesgefahr. Bevor das Kabel an den Steckverbinder des TSP-Einsatzes angeschlossen wird und bevor es von diesem abgezogen wird, ist unbedingt sicherzustellen, dass die Steuereinheit ausgeschaltet ist.

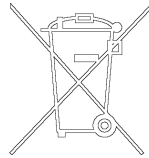
Entsorgung

Bedeutung des "WEEE" Logos auf den Etiketten.

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

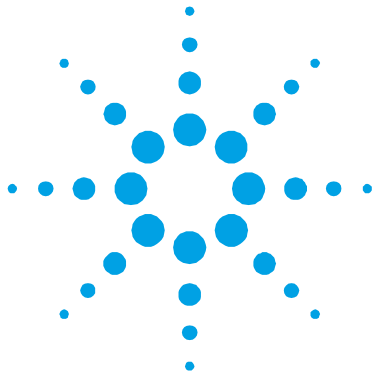
Dieses Symbol (**nur in den EU-Ländern gültig**) zeigt an, dass das betreffende Produkt nicht zusammen mit Haushaltsmüll entsorgt werden darf sondern einem speziellen Sammelsystem zugeführt werden muss.

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.



2 Anleitung zur Installation

Entsorgung



3

Procédure pour l'installation

Indications Generales 28

Generalites sur le TSP Cryopanel Assy 29

Caractéristiques Techniques 30

Installation du TSP Cryopanel 31

Installation de la Cartouche TSP 32

Mise au rebut 35

Traduction de la mode d'emploi originale



Indications Generales

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 d'utiliser l'appareil. Agilent décline toute responsabilité en cas de non respect total ou partiel des instructions fournies, d'opérations non autorisées, d'utilisation impropre par du personnel non formé ou contraires aux réglementations nationales spécifiques.

Cette notice utilise les signes conventionnels suivants:

AVERTISSEMENT!



Les messages d'avertissement attirent l'attention de l'opérateur sur une procédure ou une manœuvre spéciale dont la mauvaise exécution risque de provoquer de graves lésions.

ATTENTION!

Les messages d'attention apparaissent avant certaines procédures dont le non-respect peut endommager sérieusement l'appareillage.

NOTE

Les notes contiennent des renseignements importants, extrapolés du texte.

Generalites sur le TSP Cryopanel Assy

Le TSP Cryopanel a été conçu pour être utilisé conjointement avec la TSP (Titanium Sublimation Pump) par un montage sur une bride 8" OD Con-Flat®. On peut utiliser de l'eau ou de l'azote liquide pour le refroidissement. En outre, le système peut fonctionner en modalité UHV sans refroidissement. Le Cryopanel peut être installé sur des pompes ioniques doubleended ou sideported.

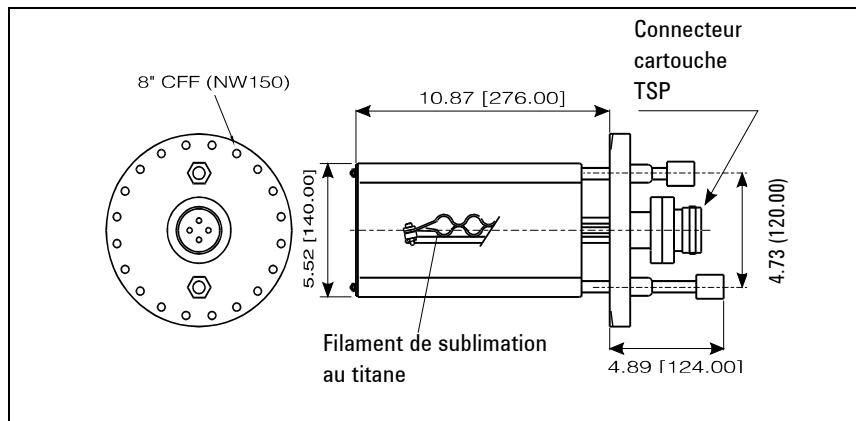


Figure 1 Encombrements TSP Cryopanel en mm [pouces]

Caractéristiques Techniques

Tab. 1

	N ₂	H ₂	H ₂ O
Vitesse de pompage à 20 °C avec TSP et refroidissement à l'eau (l/s)	516	1205	578
Vitesse de pompage à -195 °C avec TSP et refroidissement au LN ₂ (l/s)	555	1760	696*
Surface de pompage interne (cm ²)	826		
Bride principale	8" CF (NW 150)		
Vol. du réservoir (litres)	1.8		
Raccords de refroidissement	3/8" Gas		
Bride source au titane	2 3/4 CF		

* Compte tenu du fait que l'H₂O est un fluide condensable, l'effet du pompage sur la surface de sortie du cryopanel est compris. La vitesse de pompage est de 12.000 l/sec.

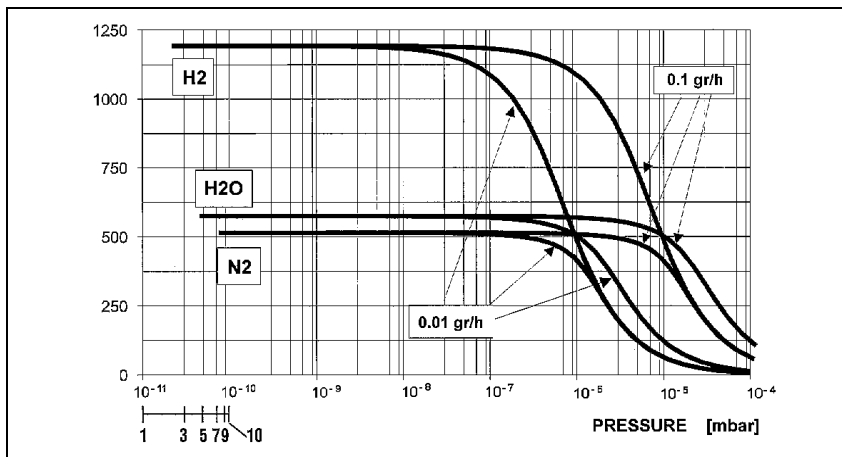


Figure 2 Vitesse de pompage à 20 °C avec TSP et refroidissement à l'eau

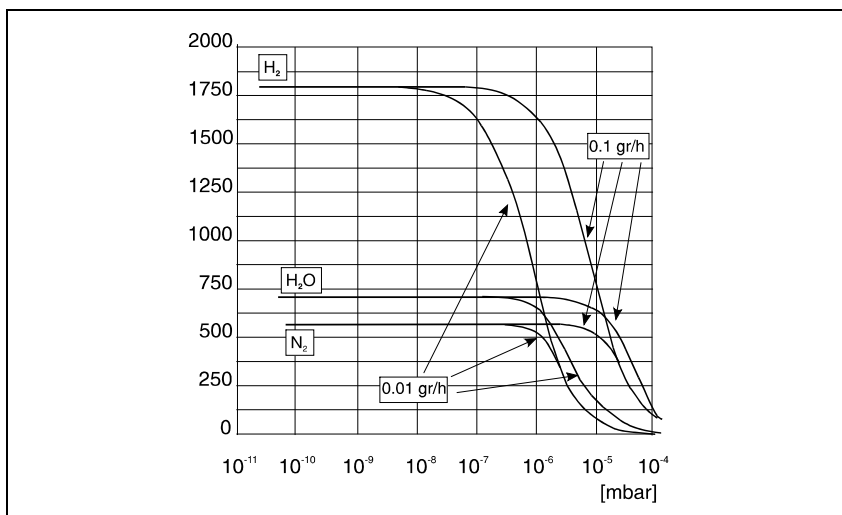


Figure 3 Vitesse de pompage à -195 °C avec TSP et refroidissement à l'azote liquide

Installation du TSP Cryopanel

AVERTISSEMENT!



Si l'on utilise du LN₂ pour le refroidissement, faire très attention en manipulant le TSP Cryopanel pendant les phases d'installation, car le contact du LN₂ avec la peau provoque de graves brûlures.

Pour installer correctement le TSP Cryopanel, suivre fidèlement la procédure indiquée cidessous:

- 1 Nettoyer soigneusement la bride de 8" ConFlat® du Cryopanel ainsi que la bride correspondante sur la pompe.
- 2 Positionner les écrous et les visser chacun à un couple de 6 – 11 Nm (4.5 - 8 ft.-lbs). Après avoir vissé un écrou, toujours visser celui à l'opposé par rapport au centre de la bride.

3 Procédure pour l'installation

Generalites sur le TSP Cryopanel Assy

- 3 Répéter cette séquence de vissage pendant encore deux cycles.
- 4 Continuer à visser les écrous jusqu'à ce que les côtés des brides entrent en contact et jusqu'à avoir une augmentation du couple.
- 5 Fixer les tubulures du fluide de refroidissement aux fixations correspondantes sur le TSP Cryopanel, en prenant soin de serrer les écrous de blocage de manière appropriée.

Installation de la Cartouche TSP

Normalement, la cartouche TSP est fournie avec une bride OD 2 $\frac{3}{4}$ " ConFlat® avec trois filaments au titane - molybdène; chacun étant bakable (chauffable) à 400 °C. La sublimation maximum est obtenue en fournissant une puissance de 300 W.

Pour fixer la cartouche sur le Cryopanel, procéder comme le décrit le paragraphe précédent, en prenant garde à fixer la bride de manière à ne pas créer d'interférences entre le connecteur du câble provenant du système de commande et les tubulures de refroidissement.

NOTE

Pendant la phase d'installation, il faut respecter le sens de circulation du fluide de refroidissement, ainsi que l'illustrent les exemples suivants.

ATTENTION!

Veuillez retirer les liquides de refroidissement du Cryopanel avant de cuire au four et laissez le raccord du tuyau de refroidissement ouvert pendant la cuisson.

Les figures illustrent quelques exemples de positionnement du TSP Cryopanel.

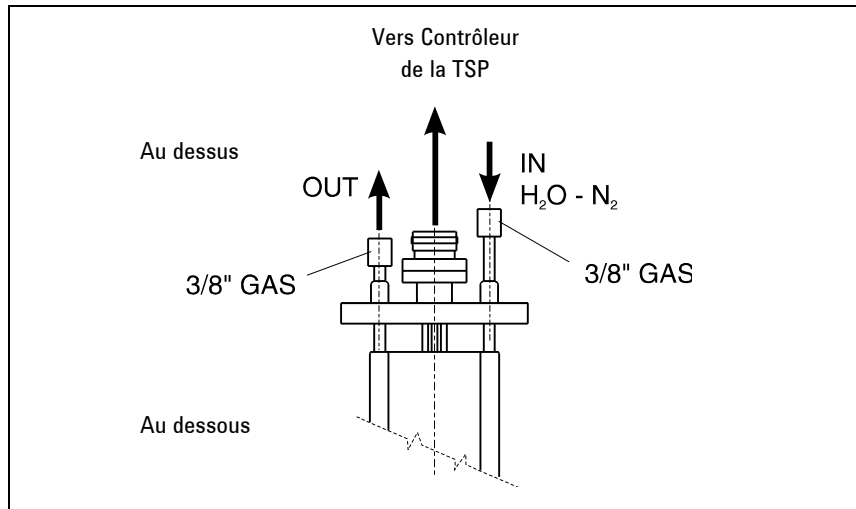


Figure 4 Exemple de positionnement vertical

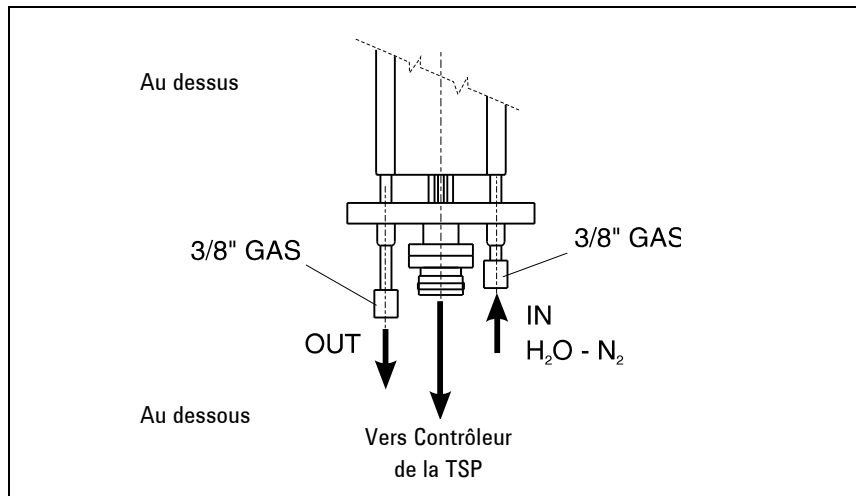


Figure 5 Exemple de positionnement vertical

3 Procédure pour l'installation

Generalites sur le TSP Cryopanel Assy

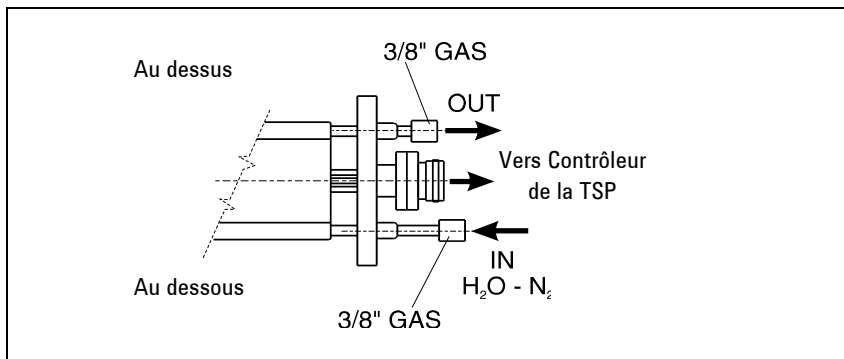


Figure 6 Exemple de positionnement horizontal

- 1 Brancher le câble provenant du Titanium Sublimation Controller au connecteur correspondant sur le centre du TSP Cryopanel.

AVERTISSEMENT!



Le courant élevé qui passe dans le câble de connexion au système de commande pourrait provoquer de graves accidents et mettre les personnes en danger de mort. Avant de brancher les câbles au connecteur de la cartouche TSP, et de même avant de le débrancher, vérifier que l'on a éteint le système de commande.

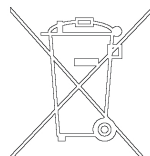
Mise au rebut

Signification du logo "WEEE" imprimé sur les étiquettes.

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

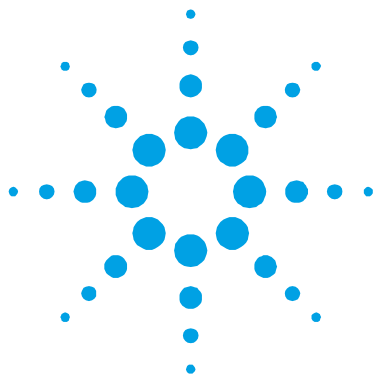
Ce symbole (**uniquement valide pour les pays de la Communauté européenne**) indique que le produit sur lequel il est appliqué NE doit PAS être mis au rebut avec les ordures ménagères ou les déchets industriels ordinaires, mais passer par un système de collecte sélective.

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.



3 Procédure pour l'installation

Mise au rebut



4 Installation procedure

General Information	38
Overview on the TSP Cryopanel Assy	39
Technical Specifications	40
TSP Cryopanel Installation	41
TSP Cartridge Installation	42
Disposal	45

Original Instructions



General Information

This equipment is destined for use by professionals. The user should read this instruction manual and any other additional information supplied by Agilent before operating the equipment. 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 equipment or any action contrary to that provided for by specific national standards.

This manual uses the following standard protocol:

WARNING!



The warning messages are for attracting the attention of the operator to a particular procedure or practice which, if not followed correctly, could lead to serious injury.

CAUTION!

The caution messages are displayed before procedures which, if not followed, could cause damage to the equipment.

NOTE

The notes contain important information taken from the text.

Overview on the TSP Cryopanel Assy

The TSP Cryopanel has been designed for use with the TSP (Titanium Sublimation Pump) and is mounted on an 8" OD ConFlat® flange. It can operate with water or liquid nitrogen cooling, or un-cooled if used at UHV. The Cryopanel can be mounted on double-ended or side ported ion pumps.

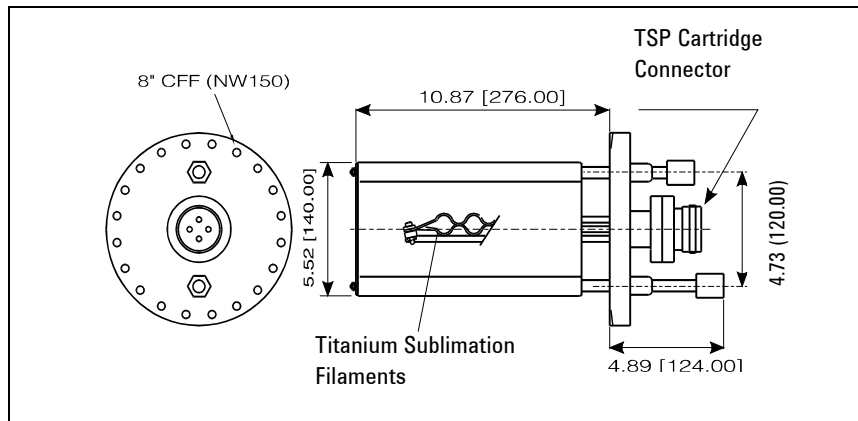


Figure 1 TSP Cryopanel Outline Drawing in mm [Inches]

Technical Specifications

Tab. 1

	N ₂	H ₂	H ₂ O
Pumping speed at 20 °C with TSP and water cooled (l/s)	516	1205	578
Pumping speed at -195 °C with TSP and LN ₂ cooled (l/s)	555	1760	696*
Inner pumping surface (cm ²)	826		
Main flange	8" CF (NW 150)		
Reservoir volume (liters)	1.8		
Cooling connection	3/8" Gas		
Titanium source flange	2¾ CF		

* It must be considered that H₂O is a condensable gas and therefore the pumping effect of the outlet surface of the Cryopanel should be included. This pumping speed is = 12.000 l/s.

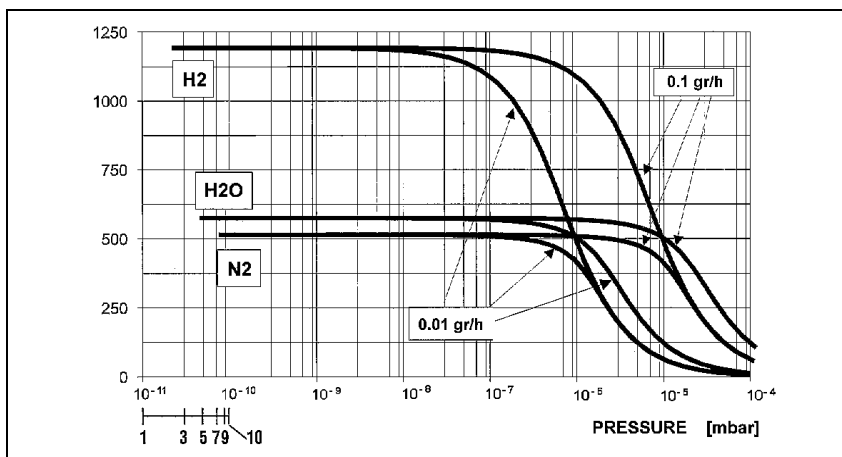


Figure 2 Pumping Speed at 20 °C with TSP and Water Cooling

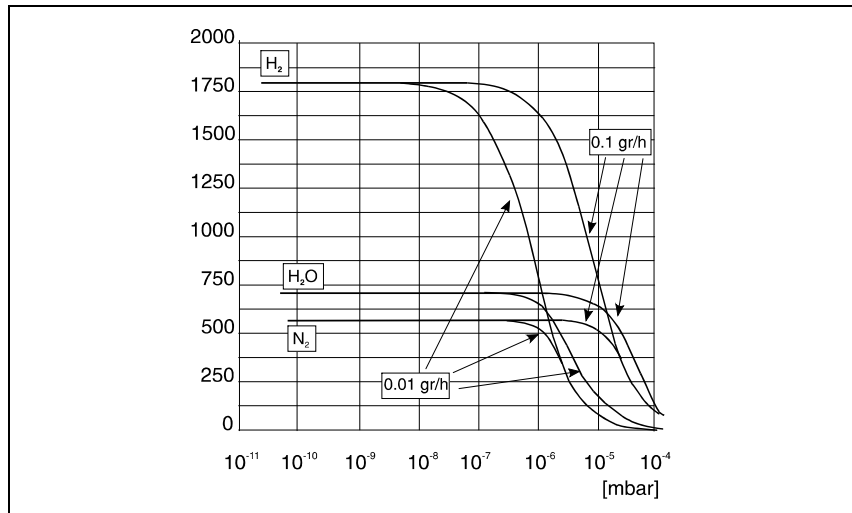


Figure 3 Pumping Speed at -195 °C with TSP and Liquid Nitrogen Cooling

TSP Cryopanel Installation

WARNING!



If LN₂ is used for cooling, be very careful when handling the TSP Cryopanel during installation. Contact with LN₂ causes severe burns to the skin.

To correctly install the TSP Cryopanel, carefully follow the procedure indicated below.

- 1 Adequately clean the Cryopanel's 8" ConFlat® flange and the flange present on the pump.
- 2 Attach the nuts and tighten each one turn to 6 – 11 Nm (4.5 - 8 ft.-lbs) of torque. After tightening a nut, always tighten the opposite nut with respect to the center of the flange.
- 3 Repeat this tightening sequence for another two cycles.

4 Installation procedure

Overview on the TSP Cryopanel Assy

- 4 Continue tightening the bolts until the flange faces meet and a pronounced increase in torque is felt.
- 5 Secure the cooling tubes for the liquid to the respective fittings on the TSP Cryopanel making sure to adequately tighten the connecting nuts.

TSP Cartridge Installation

The TSP cartridge is usually provided with an OD 2¾" ConFlat® flange with three titanium-molybdenum filaments, each bakeable at 400 °C. Maximum sublimation is obtained with a power of 300 W.

To secure the cartridge on the Cryopanel, proceed as explained in the previous section, making sure to position the flange in a way to avoid interference between the connector of the cable from the control unit and the cooling tubes.

NOTE

During installation consider the direction in which the cooling liquid flows, as shown in the following examples.

CAUTION!

Take care to remove cooling liquids from the Cryopanel prior baking and keep the cooling tube connection open during baking.

The following figures show some examples for how to position the TSP Cryopanel.

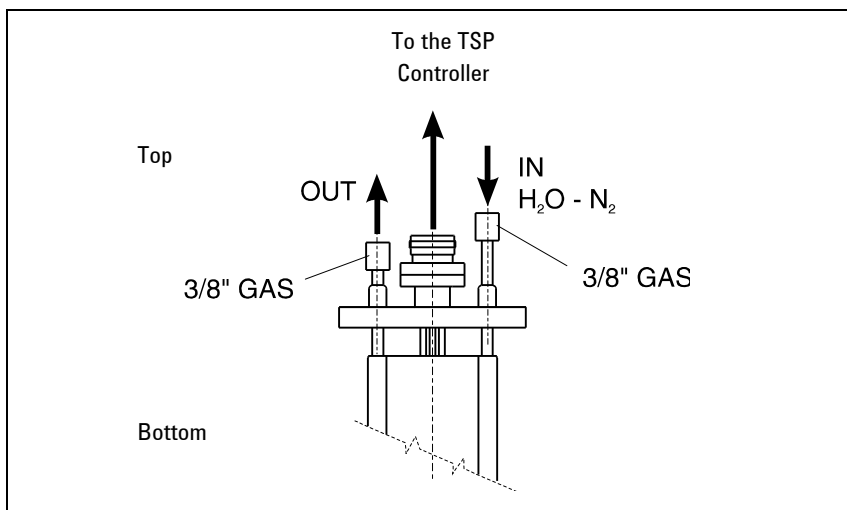


Figure 4 Example of a Vertical Positioning

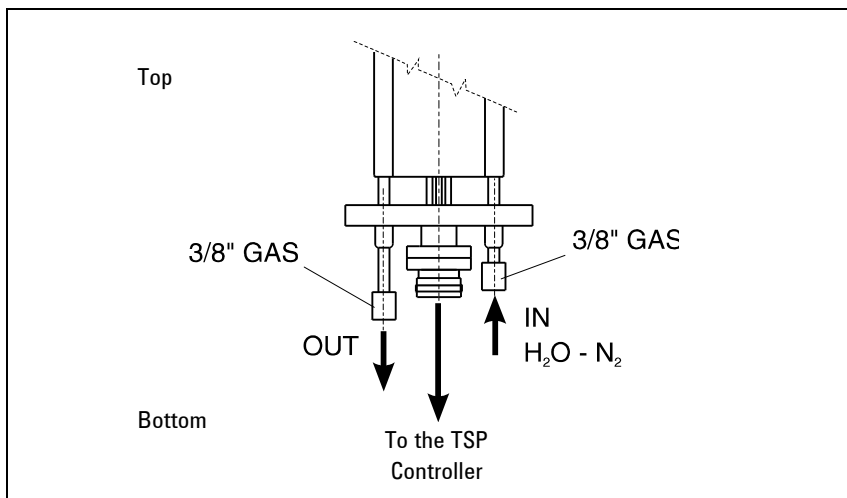


Figure 5 Example of a Vertical Positioning

4 Installation procedure

Overview on the TSP Cryopanel Assy

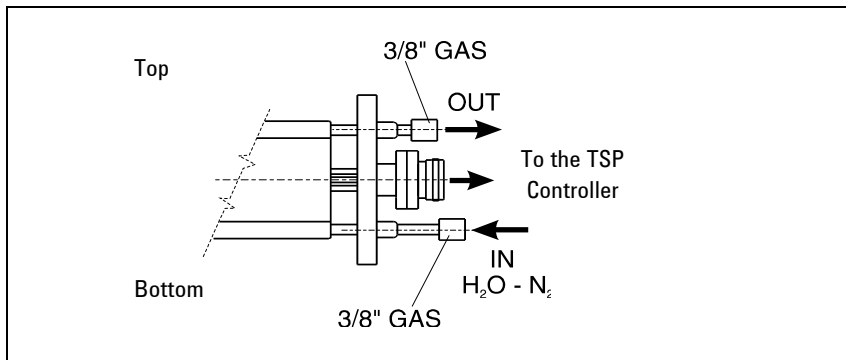


Figure 6 Example of a Horizontal Positioning

- 1 Attach the cable from the Titanium Sublimation Controller to the connector in the center of the TSP Cryopanel.

WARNING!



The high voltage present in the control unit connection cable could cause severe injury or death. Be sure to power off the control unit before attaching the cable to the TSP cartridge connector, or before disconnecting it.

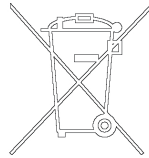
Disposal

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>



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

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

1) CUSTOMER INFORMATION

Company Name:		Contact Name:
Tel:	Email:	Fax:
Customer Ship To:		Customer Bill To:
Europe only: VAT reg. Number:		USA/Canada only: ☐ Taxable ☐ Non-taxable

2) PRODUCT IDENTIFICATION

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

AGILENT TECHNOLOGIES CANNOT ACCEPT ANY PRODUCTS CONTAMINATED WITH BIOLOGICAL OR EXPLOSIVE HAZARDS, RADIOACTIVE MATERIAL, OR MERCURY AT ITS FACILITY.

Call Agilent Technologies to discuss alternatives if this requirement presents a problem.

The equipment listed above (check one):

- ☐ **HAS NOT** pumped or been exposed to any toxic or hazardous materials. OR
- ☐ **HAS** pumped or been exposed to the following toxic or hazardous materials. If this box is checked, the following information must also be filled out. Check boxes for all materials to which product(s) pumped or was exposed:

☐ Toxic ☐ Corrosive ☐ Reactive ☐ Flammable ☐ Explosive ☐ Biological ☐ Radioactive

List all toxic/hazardous materials. Include product name, chemical name, and chemical symbol or formula:

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.

Print Name: _____ **Authorized Signature:** _____ **Date:** _____

5) FAILURE INFORMATION:

Failure Mode (REQUIRED FIELD. See next page for suggestions of failure terms):
Detailed Description of Malfunction: (Please provide the error message)
Application (system and model):

I understand and agree to the terms of Section 6, Page 3/3.

Print Name: _____ **Authorized Signature:** _____ **Date:** _____



**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

APPARENT DEFECT/MALFUNCTION	POSITION	PARAMETERS
- Does not start - Does not spin freely - Does not reach full speed - Mechanical Contact - Cooling defective	- Noise - Vibrations - Leak - Overtemperature - Clogging	- Vertical - Horizontal - Upside-down - Other:
		Power: Rotational Speed: Current: Inlet Pressure: Temp 1: Foreline Pressure: Temp 2: Purge flow: OPERATING TIME:

ION PUMPS/CONTROLLERS

- Bad feedthrough - Vacuum leak - Error code on display	- Poor vacuum - High voltage problem - Other
---	--

VALVES/COMPONENTS

- Main seal leak - Solenoid failure - Damaged sealing area	- Bellows leak - Damaged flange - Other
--	---

LEAK DETECTORS

- Cannot calibrate - Vacuum system unstable - Failed to start	- No zero/high background - Cannot reach test mode - Other
---	--

INSTRUMENTS

- Gauge tube not working - Communication failure - Error code on display	- Display problem - Degass not working - Other
--	--

SCROLL AND ROTARY VANE PUMPS

- Pump doesn't start - Doesn't reach vacuum - Pump seized	- Noisy pump (describe) - Over temperature - Other
---	--

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
05/2011
Publication Number: 87-900-078-01 (E)



Agilent Technologies