

2 l/s Ion Pump

Models

**9190520, 9190521, 9190522,
9190523, 9190524**

Magnet 9190038

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

87-900-091-01 (F)

06/2019



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2019

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: 87-900-091-01 (F)

Edition

Edition 06/2019

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

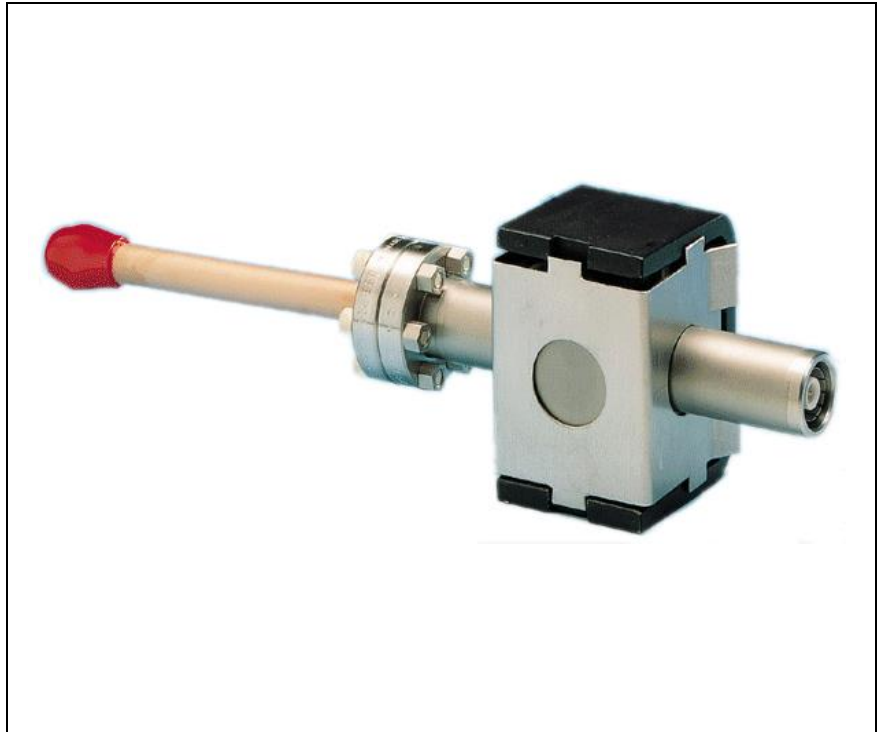
CAUTION

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.

2 l/s Ion Pump



Contents

1 Istruzioni per l'uso 7

Informazioni generali 8

Preparazione per l'installazione 9

Installazione 10

Uso 12

Manutenzione 14

Smaltimento 15

2 Gebrauchsanleitung 17

Allgemeine Hinweise 18

Vor der Installation 19

Installation 20

Gebrauch 22

Wartung 24

Entsorgung 25

3 Mode d'emploi 27

Indications Générales 28

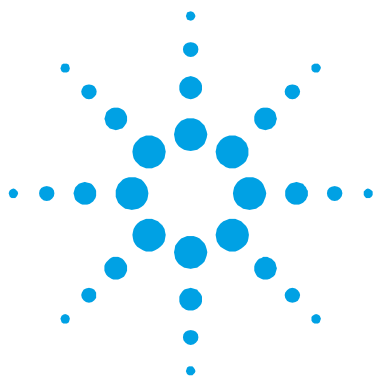
Préparation de l'installation 29

Installation 30

Utilisation 32

Contents

Maintenance	34
Mise au rebut	35
4 Instructions for Use	37
General Information	38
Preparation for Installation	39
Installation	40
Use	42
Maintenance	44
Disposal	45
5 Technical Information	47
Description of the Pump	48
Operating Notes	50
Technical Specification	51
Outline Drawing	53
Installation	56
Pump Installation	57
Control Unit Connection	58
Bakeout	59
Pump Troubleshooting	61
2 l/s Pump Replacement Parts and Accessories	62



1

Istruzioni per l'uso

Informazioni generali 8

Preparazione per l'installazione 9

Installazione 10

Uso 12

Procedure di uso 13

Manutenzione 14

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.

Le pompe della serie 2l/s sono pompe ioniche utilizzate comunemente per applicazioni di ultra alto vuoto, grazie alla loro pulizia, capacità di pompare qualsiasi tipo di gas, e del loro funzionamento senza vibrazioni e necessità di manutenzione.

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

Questo manuale utilizza le seguenti convenzioni:

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.

Preparazione per l'installazione

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

Durante l'operazione di disimballaggio, prestare particolare attenzione a non lasciar cadere la pompa e a non sottoporla ad urti o vibrazioni.

Non disperdere l'imballo nell'ambiente. Il materiale è completamente riciclabile e risponde alla direttiva CEE 85/399 per la tutela dell'ambiente.

ATTENZIONE!

Onde evitare problemi di degassamento, non toccare con le mani nude i componenti destinati ad essere esposti al vuoto. Utilizzare sempre i guanti o altra protezione adeguata.

NOTA

La pompa non può essere danneggiata rimanendo semplicemente esposta all'atmosfera. Si consiglia comunque di mantenerla chiusa fino al momento dell'installazione sul sistema onde evitare eventuale inquinamento da polvere.

1 Istruzioni per l'uso Installazione

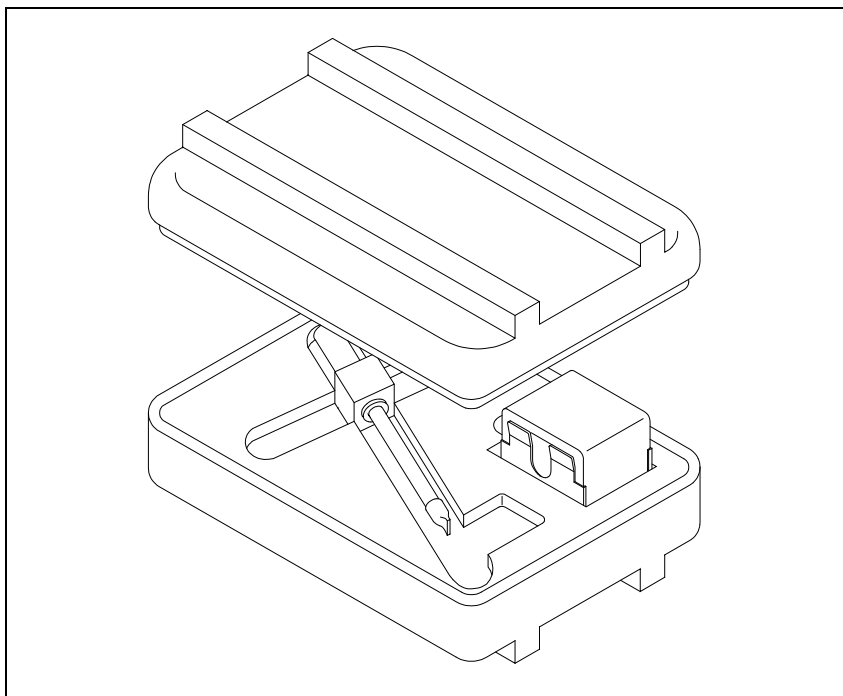


Figura 1

Installazione

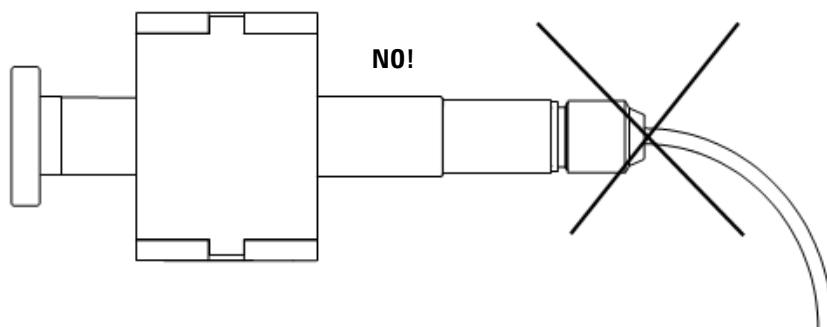
Non installare e/o utilizzare la pompa in ambienti esposti ad agenti atmosferici (pioggia, gelo, neve), polveri, gas aggressivi, in ambienti esplosivi o con elevato rischio di incendio. Durante il funzionamento, per ottenere le specifiche tecniche dichiarate, la temperatura ambiente deve essere compresa tra 0 °C e +85 °C.

ATTENZIONE! La pompa deve essere tenuta sigillata finchè non è pronta per essere collegata al sistema.

AVVERTENZA!



Per evitare lesioni alla persona, non collegare l'alta tensione alla pompa prima che sia installata nel sistema e che tutte le flange di ingresso siano adeguatamente collegate o chiuse.



ATTENZIONE!

Il rispetto delle normative di sicurezza nell'uso delle pompe è garantito solo con l'uso delle unità di controllo e dei cavi Agilent.

ATTENZIONE!

Il passante per la connessione del cavo di alta tensione è previsto per l'uso con il cavo Agilent P/N 9290706.

Solo in questo caso è disponibile la funzione "H.V. Cable Safety Interlock" che fa sì che l'alta tensione venga spenta quando il relativo cavo viene scollegato dalla pompa.

1 Istruzioni per l'uso

Uso

La pompa può essere installata in qualsiasi posizione. Per convenienza normalmente viene montata in posizione verticale con la flangia di ingresso in alto, o in posizione orizzontale.

Le pompe possono anche essere mantenute sospese in ogni posizione tramite la loro flangia di ingresso.

Per informazioni dettagliate sull'installazione della pompa, vedere l'appendice "Technical Information".

Uso

Tutte le istruzioni per il corretto funzionamento della pompa sono contenute nel manuale dell'unità di controllo.

Leggere attentamente tale manuale prima dell'utilizzo.

Si raccomanda di portare la pompa ad una pressione di 1×10^{-4} mbar in modo da ottenere un avvio più rapido. A questo scopo è meglio non utilizzare una pompa meccanica sigillata con olio, ma se deve essere utilizzata una tale pompa, si raccomanda di inserire una trappola lungo la linea di vuoto per ridurre la pressione causata dai vapori di acqua e di olio provenienti dalla pompa meccanica. Cercare di ridurre al minimo il tempo in cui la pompa meccanica è aperta verso il sistema e verso la pompa ionica, poiché i suoi vapori si diffondono nel sistema a pressioni inferiori a 1×10^{-1} mbar causando una contaminazione dello stesso. Nei sistemi in cui l'olio deve essere completamente assente, è meglio utilizzare pompe Turbo per la suddetta operazione.

Depositi igroscopici e l'assorbimento dell'idrogeno all'interno del composto di titanio possono provocare l'allungamento del tempo di avvio con l'invecchiamento della pompa. Durante l'esposizione all'aria, il deposito del composto di titanio assorbe vapore acqueo; in conseguenza di ciò, al successivo avvio, il riscaldamento della pompa provoca il rilascio del vapore e di parte dell'idrogeno pompato precedentemente, con l'effetto dell'allungamento del tempo di avvio.

Procedure di uso

Osservare la seguente procedura per l'uso della pompa:

1. Tramite una pompa di pre-vuoto portare il sistema ad una pressione di 1×10^{-4} mbar od inferiore.
2. Collegare l'unità di controllo ad una apposita fonte di alimentazione ed accenderla.
3. Lasciare la valvola di pre-vuoto aperta durante l'avvio della pompa ionica finchè non si raggiunge una adeguata pressione di avvio. Se la tensione della pompa ionica scende dopo la chiusura della valvola, riaprirla per un pre-pompaggio aggiuntivo. Appena la pressione diminuisce, la tensione cresce nuovamente e la valvola di pre-vuoto deve essere chiusa.
4. Quando si porta la pompa alla pressione atmosferica, usare azoto secco in modo da evitare l'assorbimento di vapore acqueo da parte delle pareti della pompa.

AVVERTENZA!



Quando la pompa viene utilizzata per il pompaggio di gas tossici, infiammabili o radioattivi, seguire le appropriate procedure tipiche di ciascun gas. Non usare la pompa in presenza di gas esplosivi.

ATTENZIONE!

Non avvicinare dispositivi elettronici alla pompa. Il campo magnetico attorno ad essa può provocare dei malfunzionamenti dei dispositivi stessi.

Manutenzione

Le pompe della serie 2 l/s non richiedono alcuna manutenzione.
Qualsiasi intervento deve essere eseguito da personale autorizzato.

AVVERTENZA!



Prima di effettuare qualsiasi intervento sulla pompa scollegarla dall'alta tensione.

Qualora una pompa dovesse essere rottamata, procedere alla sua eliminazione nel rispetto delle normative nazionali specifiche.

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.

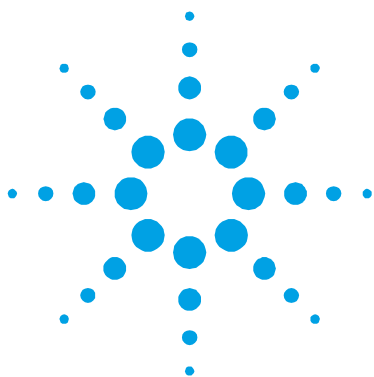
Si invita pertanto l'utente finale a contattare il fornitore del dispositivo, sia esso la casa madre o un rivenditore, per avviare il processo di raccolta e smaltimento, dopo opportuna verifica dei termini e condizioni contrattuali di vendita.



Per maggiori informazioni riferirsi a:

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

1 Istruzioni per l'uso Smaltimento



2

Gebrauchsanleitung

Allgemeine Hinweise	18
Vor der Installation	19
Installation	20
Gebrauch	22
Bedienungsschritte	23
Wartung	24
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.

Die Pumpen der Serie 2 l/s sind Ionenpumpen, die aufgrund ihrer Reinheit, ihrer Fähigkeit, alle Arten von Gas zu pumpen, und ihres vibrations- und wartungsfreien Betriebes, allgemein für Ultrahochvakuumanwendungen zum Einsatz kommen.

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

In dieser Gebrauchsanleitung werden Sicherheitshinweise folgendermaßen hervorgehoben:

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.

Vor der Installation

Die Pumpe wird in einer speziellen Schutzverpackung geliefert. Eventuelle Transportschäden sind der zuständigen örtlichen Verkaufsstelle zu melden.

Beim Auspacken vorsichtig vorgehen, damit die Pumpe nicht fällt oder Stößen oder Vibrationen ausgesetzt wird.

Das Verpackungsmaterial ist ordnungsgemäß zu entsorgen. Es ist vollständig recyclebar und entspricht der EG-Richtlinie 85/399 für den Umweltschutz.

VORSICHT!

Um Entgasungsprobleme zu vermeiden, dürfen die Komponenten, die mit dem Vakuum in Kontakt kommen, nicht mit bloßen Händen berührt werden. Stets Handschuhe oder einen anderen geeigneten Schutz tragen.

HINWEIS

Die Pumpe kann, wenn sie einfach der Atmosphäre ausgesetzt ist, nicht beschädigt werden. Sie sollte jedoch bis zur Installation an der Anlage geschlossen bleiben, um Verunreinigungen durch Staub zu vermeiden.

2 Gebrauchsanleitung Installation

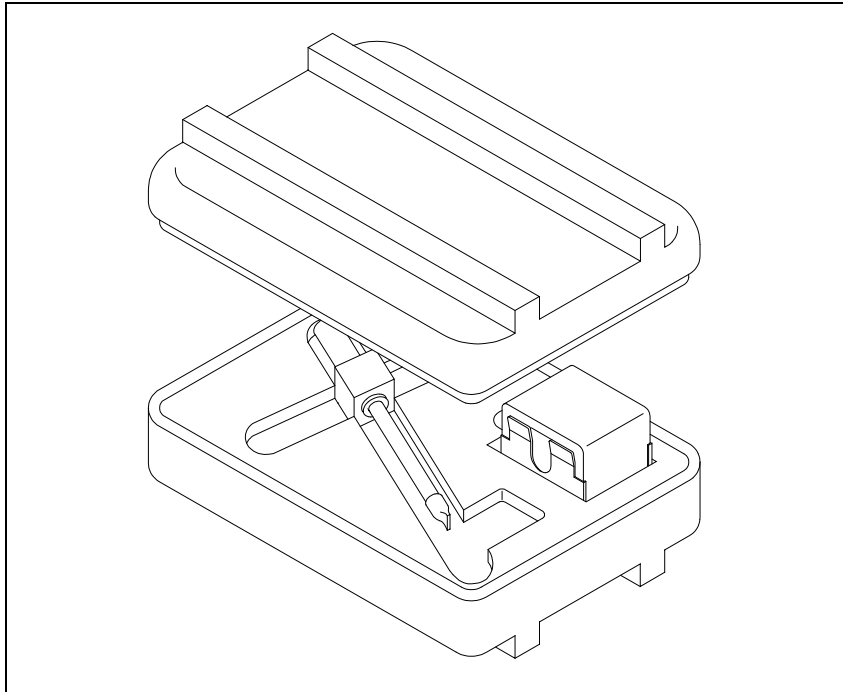


Abbildung 1

Installation

Die Pumpe darf nicht in Umgebungen installiert und/oder benutzt werden, die ungeschützt vor Witterungsbedingungen (Regen, Frost, Schnee), Staub und aggressiven Gasen sind, und in denen Explosions- und erhöhte Brandgefahr besteht.

Während des Betriebes soll die Umgebungstemperatur zwischen 0 °C und +85 °C betragen, um die angegebenen technischen Merkmale zu gewährleisten.

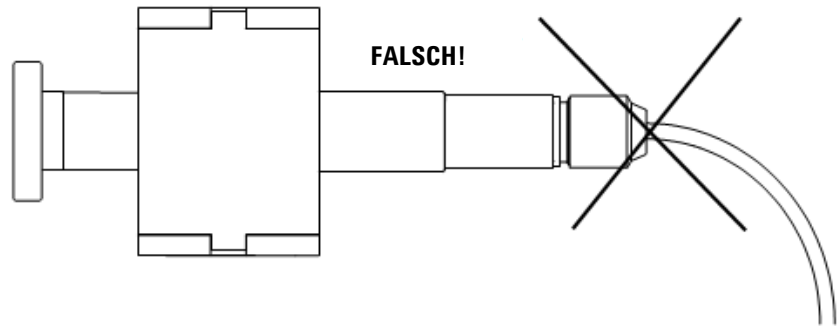
VORSICHT!

Die Pumpe soll versiegelt bis sie für den Anschluß an das System bereit ist.

WARNUNG!



Um Personenschäden zu vermeiden, darf die Hochspannungsleitung der Pumpe erst angeschlossen werden, wenn die Pumpe im System installiert ist und alle Eintrittsflansche entsprechend angeschlossen oder geschlossen sind.



VORSICHT!

Die Einhaltung der Sicherheitsvorschriften im Gebrauch der Pumpen wird nur durch den Einsatz der Kontroll- und Verkabelungssysteme Agilent garantiert.

VORSICHT!

Der Durchgangsisolator für den Anschluß des Hochspannungskabels ist für den Gebrauch des Kabels Agilent P/N 9290706. Nur in diesem Fall steht die Funktion "H.V. Cable Safety Interlock" zur Verfügung, die für die Ausschaltung der Hochspannung sorgt, sobald das entsprechende Kabel von der Pumpe gelöst wird.

Die Pumpe kann in einer beliebigen Position installiert werden. Aus praktischen Gründen wird sie in der Regel senkrecht mit nach oben gerichtetem Eintrittsflansch oder waagerecht eingebaut.

Die Pumpen können auch in einer beliebigen Position an ihrem Eintrittsflansch hängend eingebaut werden.

Detaillierte Informationen zur Installation der Pumpe sind dem Anhang "Technical Information" zu entnehmen.

Gebrauch

Sämtliche Hinweise für den korrekten Betrieb der Pumpen sind im Handbuch der Steuereinheit enthalten.

Dieses Handbuch ist vor der Inbetriebnahme genau durchzulesen.

Es wird empfohlen, die Pumpe auf einen Druck von 1×10^{-4} mbar zu bringen, um einen rascheren Anlauf zu gewährleisten. Zu diesem Zweck empfiehlt es sich, keine ölversiegelte mechanische Pumpe zu verwenden. Falls hingegen eine derartige Pumpe benutzt werden muß, soll an der Vakuumleitung eine Falle eingesetzt werden, um den durch die Wasser- und Öldämpfe aus der mechanischen Pumpe erzeugten Druck zu verringern. Es ist zu versuchen, die Zeit auf ein Mindestmaß zu reduzieren, während der die mechanische Pumpe zum System und zur Ionenpumpe geöffnet ist, da sich ihre Dämpfe bei Drücken unter 1×10^{-1} mbar im System verbreiten und dessen Verunreinigung verursachen. Bei Systemen, die vollkommen ölfrei sein müssen, erweisen sich für den obengenannten Vorgang Turbopumpen als geeigneter.

Hygroskopische Ablagerungen und die Absorption von Wasserstoff in der Titanverbindung können die Anlaufzeit verlängern und eine kürzere Standzeit der Pumpe verursachen. Die Ablagerungen der Titanverbindungen absorbieren Wasserdampf, wenn sie der Luft ausgesetzt werden. Dadurch bewirkt beim anschließenden Anlaufvorgang die Aufheizung der Pumpe die Abgabe des Dampfes und eines Teils des zuvor gepumpten Wasserstoffs, so daß sich die Anlaufzeit verlängert.

Bedienungsschritte

Beachten Sie das folgende Verfahren für die Verwendung der Pumpe:

1. Mittels einer Vorvakuumpumpe das System auf einen Druck von 1×10^{-4} mbar oder einen niedrigeren Druck bringen.
2. Die Steuereinheit an eine entsprechende Versorgungsquelle anschließen und einschalten.
3. Während des Anlaufes der Ionenpumpe soll das Vorvakuumventil geöffnet bleiben, bis ein angemessener Anlaufdruck erreicht ist. Wenn die Spannung der Ionenpumpe nach Schließung des Ventils abfällt, ist das Ventil für eine zusätzliche Vorvakuumpumpe zu öffnen. Sobald der Druck sinkt, steigt die Spannung erneut an und ist das Vorvakuumventil zu schließen.
4. Wenn die Pumpe auf atmosphärischen Druck gebracht wird, sollte Trockenstickstoff verwendet werden, damit die Wände der Pumpen keinen Wasserdampf aufnehmen.

WARNUNG!



Wenn die Pumpe zur Förderung von giftigen, leicht entflammaren oder radioaktiven Gasen benutzt wird, sind die für das jeweilige Gas vorgeschriebenen Vorgänge zu befolgen. Die Pumpe nie bei Vorhandensein von explosivem Gas benutzen.

VORSICHT!

Keine elektronischen Geräte in die Nähe der Pumpe bringen. Das darum befindliche Magnetfeld kann zu Funktionsstörungen der Geräte führen.

Wartung

Die Pumpen der Serie 2 l/s erfordern keine Wartung. Sämtliche Eingriffe dürfen nur von autorisiertem Personal vorgenommen werden.

WARNUNG!



Vor Eingriffen an der Pumpe ist diese von der Hochspannungsquelle zu trennen.

Bei eventueller Verschrottung einer Pumpe ist diese entsprechend der einschlägigen nationalen Vorschriften zu entsorgen.

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.

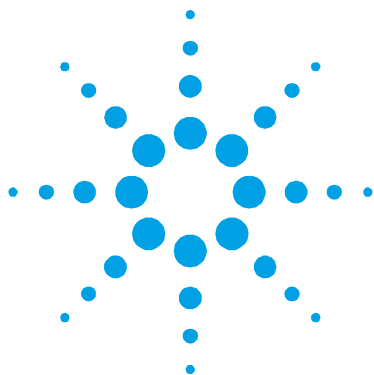
Der Endabnehmer sollte daher den Lieferanten des Geräts - d.h. die Muttergesellschaft oder den Wiederverkäufer - kontaktieren, um den Entsorgungsprozess zu starten, nachdem er die Verkaufsbedingungen geprüft hat.



Für weitere Informationen:

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

2 Gebrauchsanleitung Entsorgung



3

Mode d'emploi

Indications Générales 28

Préparation de l'installation 29

Installation 30

Utilisation 32

Procédure d'utilisation 33

Maintenance 34

Mise au rebut 35

Traduction de la mode d'emploi originale



Indications Générales

Cet appareillage a été conçu en vue d'une utilisation professionnelle. Il est conseillé à l'utilisateur de lire attentivement cette notice d'instructions ainsi que toute autre indication supplémentaire fournie par Agilent avant 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.

Grâce à leur propreté, à leur capacité de pomper tous les types de gaz, à leur fonctionnement sans vibrations et à l'absence d'entretien, les pompes de la série 2 l/s sont des pompes ioniques généralement utilisées pour des applications de vide ultra poussé.

Les paragraphes suivants fournissent toutes les indications nécessaires à garantir la sécurité de l'opérateur pendant l'utilisation de l'appareillage. Des renseignements plus détaillés se trouvent dans l'appendice "Technical Information".

Cette notice utilise les signes conventionnels suivants:

AVERTISSEMENT!



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

ATTENTION!

Les messages d'attention apparaissent avant certaines procédures dont le non respect peut provoquer des dommages irréversibles à l'appareillage.

NOTE

Les notes contiennent des renseignements importants, extrapolés du texte.

Préparation de l'installation

La pompe est fournie dans un emballage de protection spécial; si l'on constate des signes d'endommagement imputables au transport, contacter aussitôt le revendeur local.

Pendant l'opération de déballage, veiller tout particulièrement à ne pas laisser tomber la pompe et à ne lui faire subir aucun choc ni aucune vibration.

Ne pas abandonner l'emballage dans la nature. Le matériel est entièrement recyclable et il est conforme à la directive CEE 85/399 en matière de protection de l'environnement.

ATTENTION!

En vue d'éviter tout problème de dégazage, ne pas toucher à mains nues les éléments devant être exposés au vide. Mettre toujours des gants ou toute autre protection appropriée.

NOTE

La pompe ne peut être endommagée si elle reste simplement exposée à l'atmosphère. Il est quoi qu'il en soit conseillé de ne pas la retirer de son emballage avant le moment de l'installation, afin d'éviter toute pollution due à la poussière.

3 Mode d'emploi

Installation

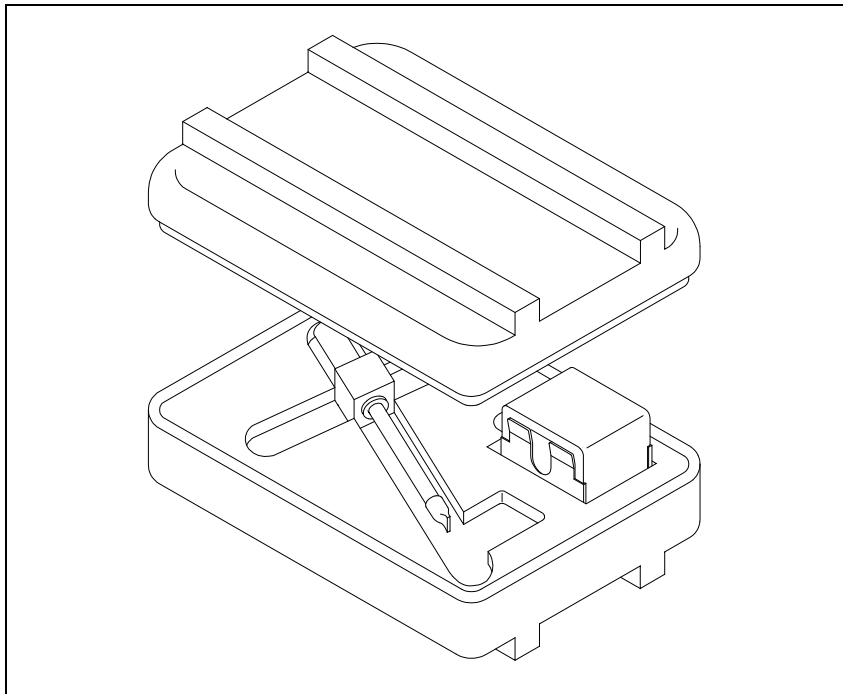


Figure 1

Installation

Ne pas installer et/ou utiliser la pompe dans des milieux exposés à des agents atmosphériques (pluie, gel, neige), à des poussières, à des gaz de combat ainsi que dans des milieux explosifs ou à fort risque d'incendie. Pendant le fonctionnement, pour respecter les spécifications techniques déclarées la température ambiante doit être comprise entre 0 °C et +85 °C;

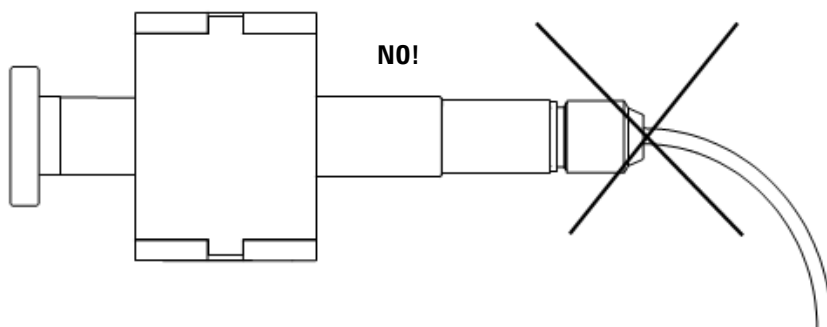
ATTENTION!

La pompe doit être conservée scellée jusqu'à ce qu'elle soit prête à être branchée au système.

AVERTISSEMENT!



Pour éviter toute lésion aux personnes, ne pas brancher la haute tension à la pompe avant que celle-ci soit installée dans le système et avant que toutes les brides d'entrées soient correctement assemblées ou fermées.



ATTENTION!

Lors de l'utilisation des pompes, le respect des normes de sécurité n'est garanti que par l'usage des unités de contrôle et des câbles Agilent.

ATTENTION!

Le passage pour la connexion du câble de haute tension est conçu pour l'utilisation du câble Agilent P/N 9290706. Dans ce cas uniquement la fonction "H.V. Cable safety Interlock" est opérationnelle. Cette fonction prévoit la coupure de la haute tension lorsque le câble correspondant est débranché de la pompe.

La pompe peut être installée dans toutes les positions. Par facilité elle est généralement montée en position verticale avec bride d'entrée en partie haute, ou en position horizontale.

Les pompes peuvent également être suspendues dans toutes les positions à l'aide de leur bride d'entrée.

Pour plus de détails sur l'installation de la pompe, consulter l'appendice "Technical Information".

Utilisation

Toutes les instructions pour le fonctionnement correct de la pompe sont fournies dans la notice de l'unité de contrôle.

Il est conseillé de lire attentivement cette notice avant d'utiliser la pompe.

Il est recommandé de porter la pompe à une pression de 1×10^{-4} mbars de façon à obtenir un démarrage plus rapide. Pour ce faire il est préférable de ne pas utiliser une pompe mécanique scellée à l'huile mais au cas où il serait nécessaire d'utiliser une telle pompe, il est recommandé d'introduire un dispositif de retenue le long de la ligne de vide afin de réduire la pression due aux vapeurs d'eau et d'huile engendrées par la pompe mécanique. Essayer de réduire au minimum le temps où la pompe mécanique est ouverte vers le système et vers la pompe ionique car ses vapeurs se répandent dans le système à des pressions inférieures à 1×10^{-1} mbars causant une contamination de celui-ci. Dans les systèmes où l'huile doit être complètement absente, il est préférable d'effectuer ladite opération à l'aide d'une pompe Turbo.

Des dépôts hygroscopiques et l'absorption d'hydrogène dans le composé de titane peuvent provoquer, par effet du vieillissement de la pompe, l'allongement du temps de démarrage. Pendant l'exposition à l'air, le dépôt du composé de titane absorbe de la vapeur d'eau et cette action a pour conséquence qu'au démarrage suivant le chauffage de la pompe entraînera la dispersion de la vapeur et d'une partie de l'hydrogène pompée précédemment ce qui aura pour effet d'allonger le temps de démarrage.

Procédure d'utilisation

Observer la procédure suivante pour l'utilisation de la pompe:

1. A l'aide d'une pompe de pré-vide, porter le système à une pression de 1×10^{-4} mbars ou inférieure.
2. Brancher l'unité de contrôle à une source d'alimentation appropriée et l'allumer.
3. Laisser la soupape de pré-vide ouverte pendant le démarrage de la pompe ionique tant qu'une pression de démarrage appropriée n'a pas été atteinte. Si la tension de la pompe ionique descend après la fermeture de la soupape, la rouvrir pour un pré-pompage supplémentaire. Dès que la pression diminue, la tension remonte et la soupape de prévide doit être refermée.
4. Lorsque l'on porte la pompe à la pression atmosphérique, utiliser de l'azote sec de façon à éviter l'absorption de vapeur aqueuse par les parois de la pompe.

AVERTISSEMENT!



Lorsque la pompe est utilisée pour le pompage de gaz toxiques, inflammables ou radioactifs, suivre les procédures appropriées à chaque gaz. Ne pas utiliser la pompe en présence de gaz explosifs.

ATTENTION!

Ne pas approcher de dispositifs électroniques de la pompe. Le champ magnétique environnant cette dernière peut entraîner des dysfonctionnements desdits dispositifs.

Maintenance

Les pompes de la série 2 l/s ne demandent aucun entretien. Toute intervention doit être exécutée par un personnel agréé.

AVERTISSEMENT! Avant toute intervention sur la pompe, la débrancher de la haute tension.



En cas de mise au rebut d'une pompe, procéder à son élimination dans le respect des normes nationales en vigueur.

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.

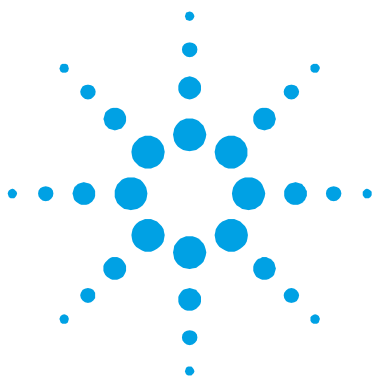


Pour plus d'informations, rendez-vous à l'adresse:

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

3 Mode d'emploi

Mise au rebut



4

Instructions for Use

General Information	38
Preparation for Installation	39
Installation	40
Use	42
Operating Procedure	43
Maintenance	44
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.

The 2 l/s series pumps are ion pumps commonly used to create ultra-high vacuum, due to their cleanliness, ability to pump different gases, and maintenance- and vibration-free operation.

The following paragraphs contain all the information necessary to guarantee the safety of the operator when using the equipment. Detailed information is supplied in the appendix "Technical Information".

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.

Preparation for Installation

The pump is supplied in a special protective packing. If this shows signs of damage which may have occurred during transport, contact your local sales office.

When unpacking the pump, be sure not to drop it and avoid any kind of sudden impact or shock vibration to it.

Do not dispose of the packing materials in an unauthorized manner. The material is 100 % recyclable and complies with EEC Directive 85/399.

CAUTION!

In order to prevent outgassing problems, do not use bare hands to handle components which will be exposed to vacuum. Always use gloves or other appropriate protection.

NOTE

Normal exposure to the environment cannot damage the pump. Nevertheless, it is advisable to keep it closed until it is installed in the system, thus preventing any form of pollution by dust.

4 Instructions for Use Installation

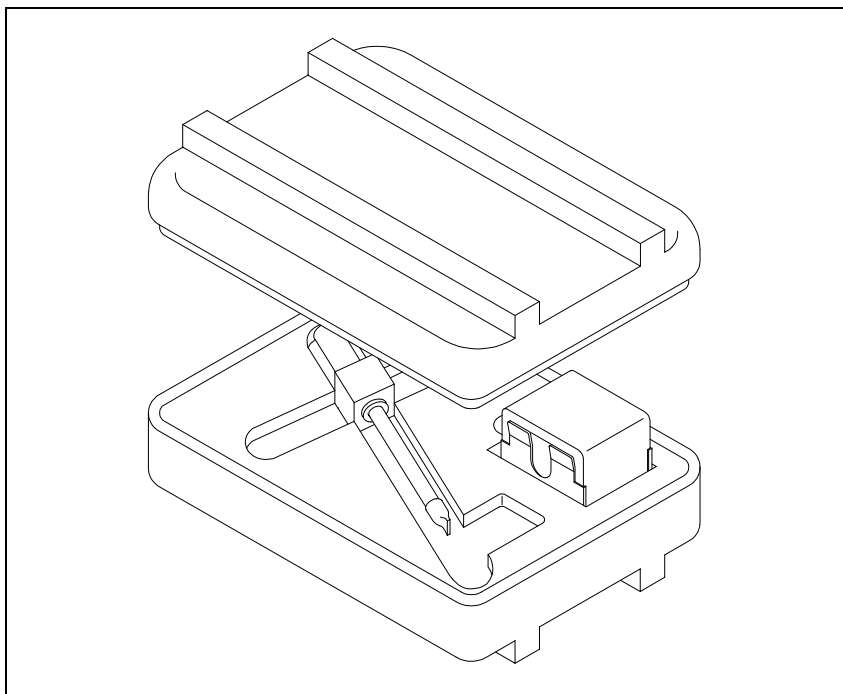


Figure 1

Installation

Do not install or use the pump in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk. During operation, to obtain the declared functioning specification, the ambient temperature must be between 0 °C and +85 °C.

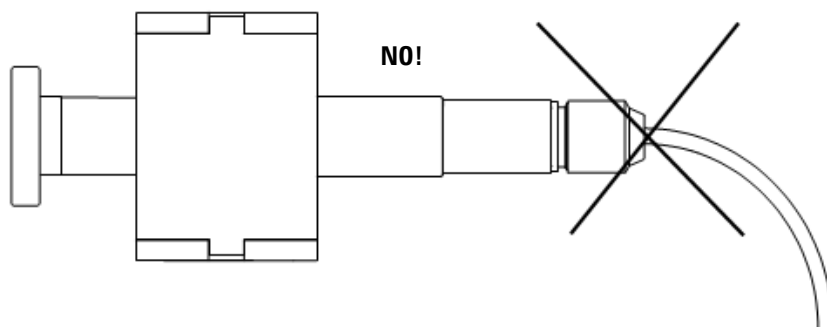
CAUTION!

The pump should be kept sealed until it is ready for attachment to the vacuum system.

WARNING!



To avoid injury, never connect the high voltage to the pump before it is installed into the system and all the inlet flanges are properly connected or blanked off.



CAUTION!

The safety specifications agreement using the pump is guaranteed using the Agilent controller and cable only.

CAUTION!

The High Voltage cable feedthrough is designed for use with cable Agilent (P/N 9290706).
Only in this case the “H.V. Cable Safety Interlock” function is available. This function switch off the high voltage when the relevant cable is disconnected from the pump.

4 Instructions for Use

Use

The pump can be installed in any position. For convenience, a pump is usually mounted vertically with the inlet up, or placed horizontally.

Pumps can be supported by the mounting flange in any position.

For detailed information about the pump installation, see the appendix “Technical Information”.

Use

All the instructions for the correct use of the pump are contained in the control unit manual.

Read the manual carefully before using the pump.

Rough pumping down to 1×10^{-4} mbar is recommended for the most rapid starting. Roughing with an oil-sealed mechanical pump is not desirable, but when used, a trap in the roughing line is recommended to reduce pressure due to water vapor and oils from the mechanical pump. Be careful to minimize the time that this pump is open to the system and ion pump, since mechanical pump vapors will start diffusing into the system at pressures below 1×10^{-1} mbar and cause contamination. In systems where oils must be completely eliminated, turbopump roughing pumps should be used.

Hygroscopic deposits and hydrogen absorption into titanium may cause starting times to increase with age. During exposure to air, the deposits of titanium compound absorb water vapor. In subsequent start ups, pump heating causes release of the water vapor and some previously pumped hydrogen; thus, the starting time may be lengthened.

Operating Procedure

Follow the procedure below when operating the pump:

1. With a clean roughing pump, establish a roughing pressure of 1×10^{-4} mbar or lower in the vacuum system.
2. Plug the control unit into a suitable power source and switch the power ON.
3. Allow the roughing valve to remain open after turning on the ion pump until an adequate starting pressure is reached. If the ion pump voltage drops after closing the roughing valves, reopen the valve for additional rough pumping. As the pressure decreases, the voltage again will rise, and the roughing valve should be closed.
4. When venting the pump, use dry nitrogen. This will avoid water vapor absorption on the pump walls.

WARNING!



When employing the pump for pumping toxic, flammable, or radioactive gases, please follow the required procedures for each gas disposal. Do not use the pump in presence of explosive gases.

CAUTION!

Do not put any electronic device near the pump otherwise the magnetic field around it may cause a device malfunctioning.

Maintenance

The 2 l/s series pumps do not require any maintenance. Any work performed on the pump must be carried out by authorized personnel.

WARNING!



Before carrying out any work on the pump, disconnect it from the High Voltage supply.

If a pump is to be scrapped, it must be disposed of in accordance with the specific national standards.

Disposal

Meaning of the "WEEE" logo found in labels

The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive.

This symbol (**valid only in countries of the European Community**) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system.

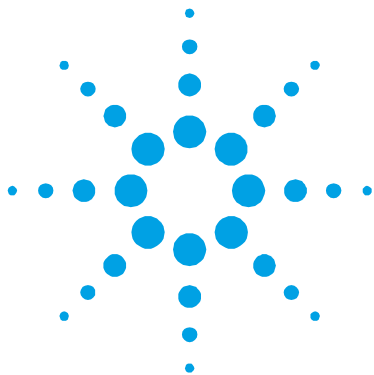
The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.



For more information refer to:

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

4 Instructions for Use Disposal



5

Technical Information

Description of the Pump	48
Operating Notes	50
Technical Specification	51
Outline Drawing	53
Installation	56
Pump Assembling	56
Pump Installation	57
Control Unit Connection	58
Bakeout	59
Unusual Operating Conditions	59
Pump Troubleshooting	61
2 l/s Pump Replacement Parts and Accessories	62

Original Instructions



Description of the Pump

The Agilent VacIon 2 l/s pumps are ion pumps and are available in the following versions:

Tab. 1

PART NUMBER	MODEL
9190520	2 l/s pump mini CFF
9190521	2 l/s pump ¾" O.D. 180 ° SST tube
9190522	2 l/s pump 3/8" O.D. 180 ° copper tube
9190523	2 l/s pump ¾" O.D. 180 ° SST tube vacuum proc.
9190524	2 l/s pump ¾" O.D. 90 ° SST tube tee style

NOTE

The part numbers in the table are relevant to the pump without the magnet. The magnet part numbers is 9190038.

The 2 l/s pump is a pump which operates on the principle of ion gettering in the pressure range from 10^{-4} to less than 10^{-9} mbar. A roughing pump is needed to lower the system pressure from atmosphere to below 10^{-4} mbar.



Figure 2 Model 9190520

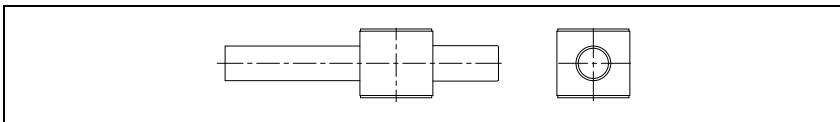


Figure 3 Model 9190521

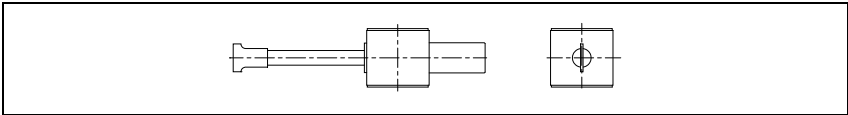


Figure 4 Model 9190522

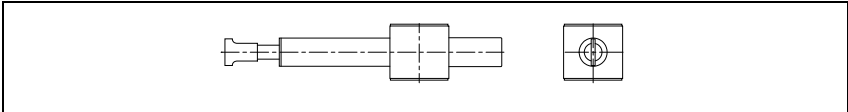


Figure 5 Model 9190523

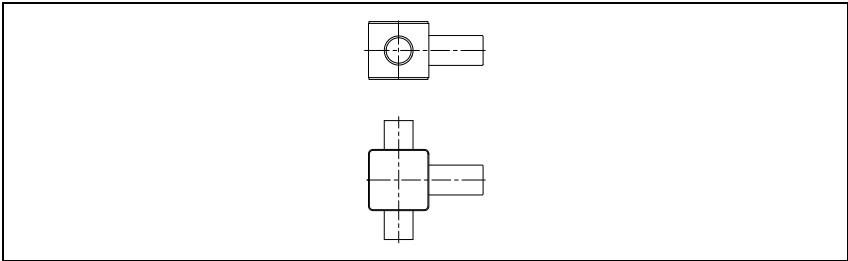


Figure 6 Model 9190524

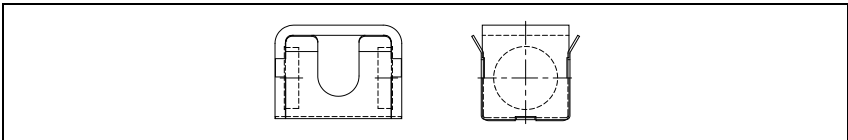


Figure 7 Magnet model 9190038

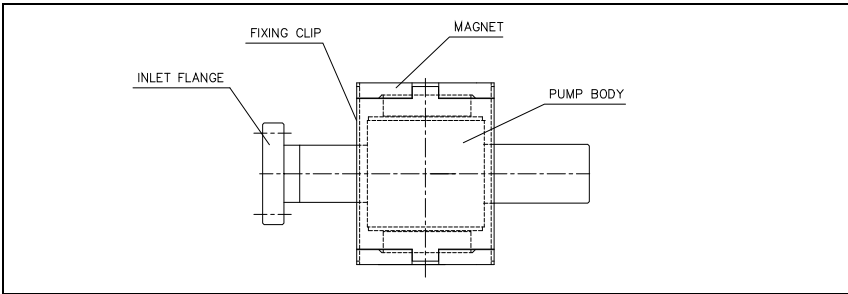


Figure 8 Pump main assemblies

Operating Notes

VacIon pumps operate as a single-cell Penning discharge. A potential of about 3 kV is applied to the anode cell; the cathodes are at ground potential. Electrons are constrained from going directly to the anode by the presence of an axial magnetic field of about 1200 gauss. A self-shielding insulator isolates the anode from ground potential. When the VacIon pump is operating properly (with magnet installed), the current drawn is proportional to pressure (see the figure in the following page).

After the system is rough-pumped to below 10^{-4} mbar, the pump can be started by simply switching on the VacIon pump control unit. A current draw on the control unit will indicate that the pump discharge has started.

The pump can operate in any position and can be supported on the mounting flange.

NOTE

Always wait at least 30 seconds after turning off the high voltage switch before disconnecting the power supply leads. This should allow the power supply output capacitor to discharge adequately.

Technical Specification

The following table details the main technical specifications of the 2 l/s Ion Pump.

Tab. 2

Nominal pumping speed for Nitrogen (l/s)	2
Operating life at 5x10 ⁻⁶ mbar (hours)	8000
Operating voltage	+5000 Vdc
Maximum starting pressure (mbar)	1x10 ⁻⁴
Ultimate pressure (mbar)	10 ⁻⁹
Internal volume (litres)	0.064
Maximum baking temperature (°C)	400 (without magnet) 150 (with magnet)
Material: Body	AISI 304
Cathode	Titanium
Anode	AISI 304
Magnet	Sm-Co
Weight, (kg)	0.3 net 0.6 with magnet

5 Technical Information
Technical Specification

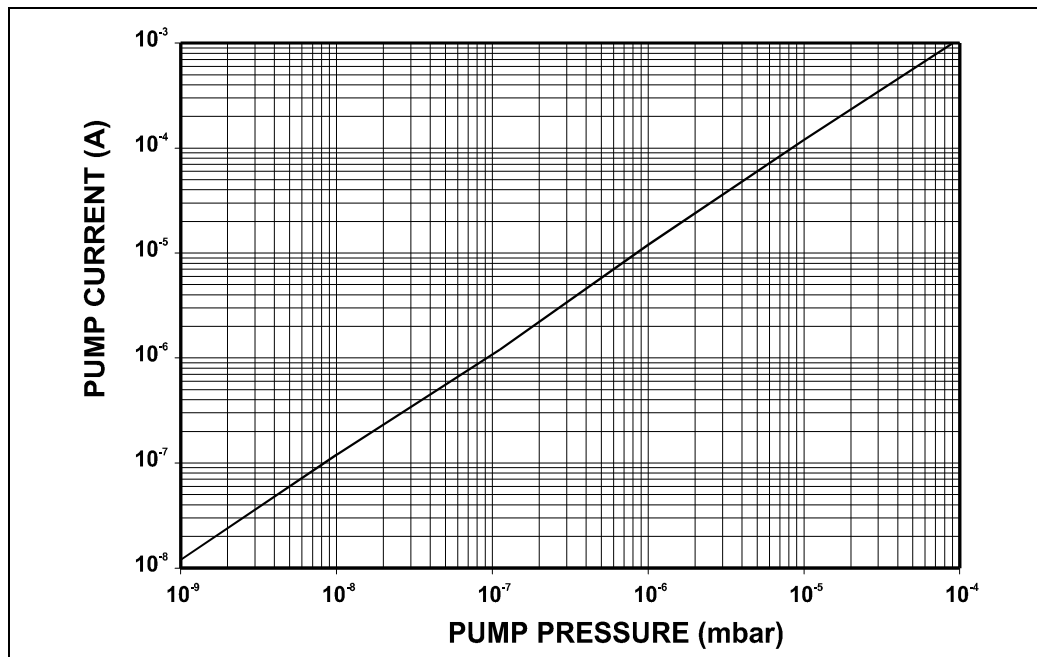


Figure 9 Pressure vs Current curve

Outline Drawing

The following figure shows the outline drawing for the 2 l/s pump and the magnet (all dimensions are in inches [mm]).

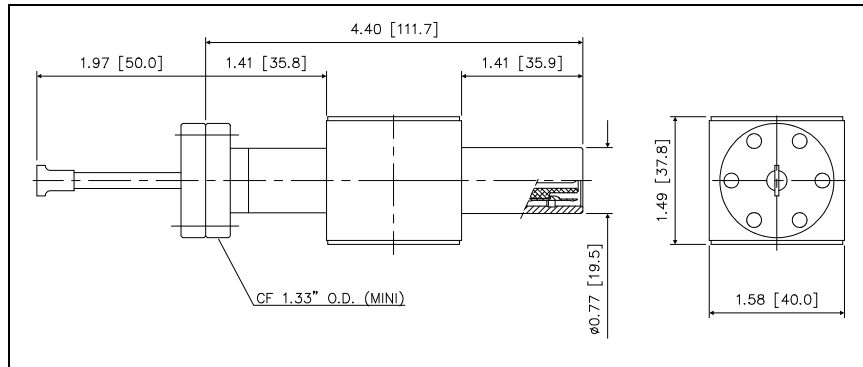


Figure 10 Model 9190520

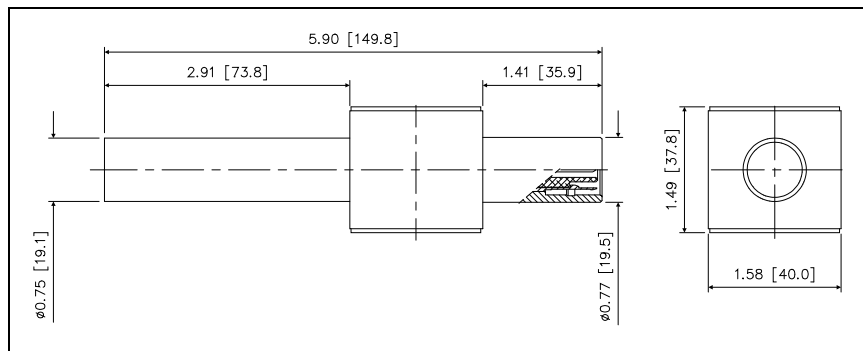


Figure 11 Model 9190521

5 Technical Information

Outline Drawing

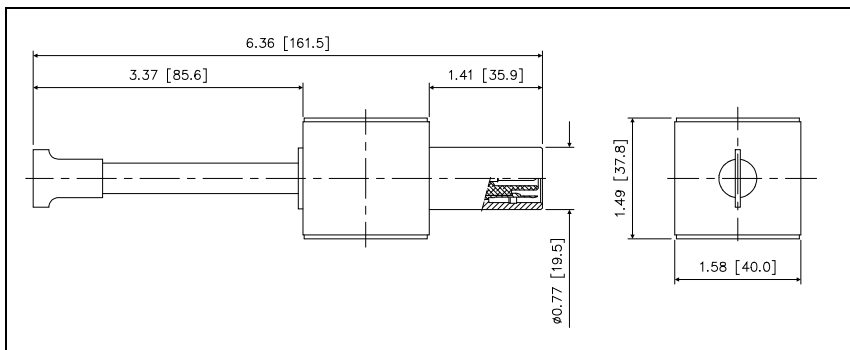


Figure 12 Model 9190522

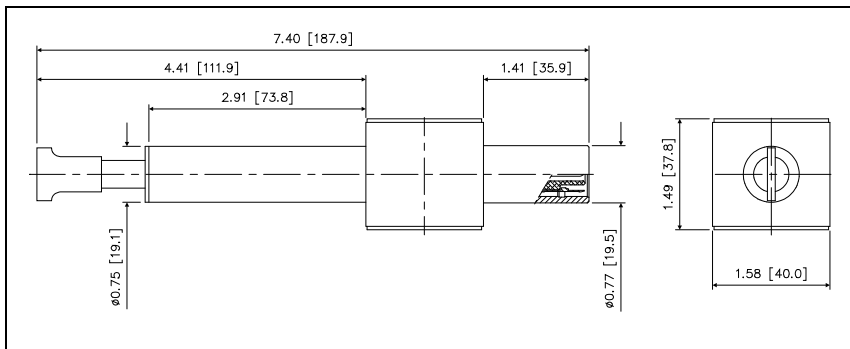


Figure 13 Model 9190523

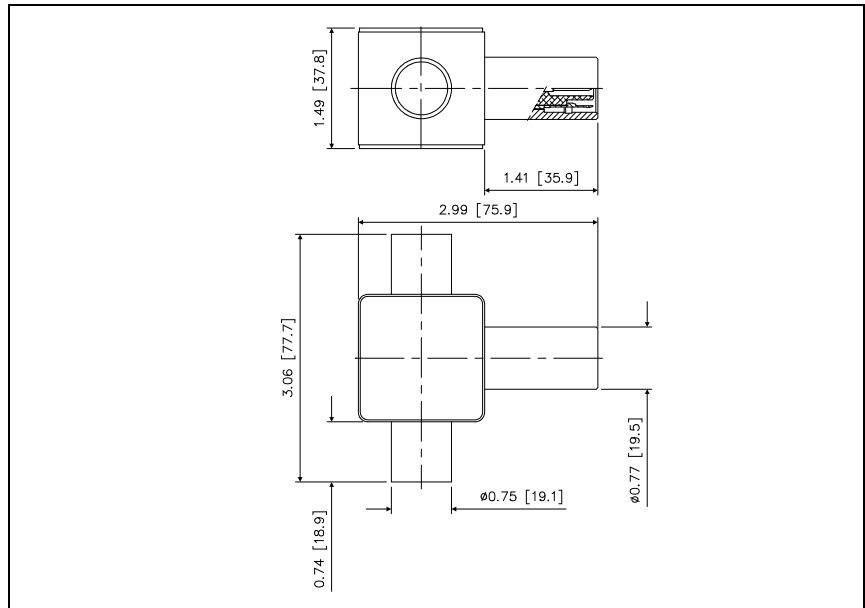


Figure 14 Model 9190524

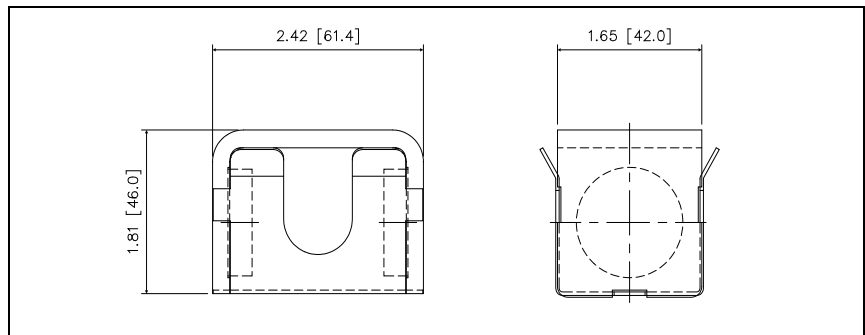


Figure 15 Magnet model 9190038

Installation

Pump Assembling

Before installing the pump into the system, the pump must be assembled.

To assemble the pump, proceed as follows (see the following figure):

1. Remove the pump parts from the packaging.
2. Remove the fixing clip (3) from the magnet (2).
3. Insert the pump (1) into the magnet.
4. Fix the pump to the magnet by means of the fixing clip.

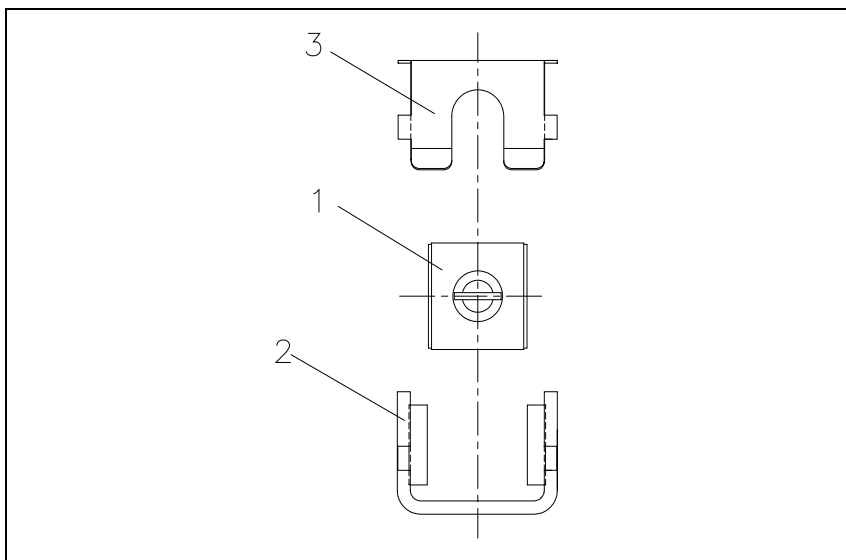


Figure 16 Pump assembly

Pump Installation

To install the pump into the system proceed as follows:

1. During installation, keep the pump and system clean. Avoid touching surfaces which will ultimately be under vacuum. When the VacIon pump is open to the atmosphere, take care to avoid introducing foreign particles, oil, solder flux, or any volatile substances. Before attaching the pump, clean the system thoroughly. Chemical cleaning is the most complete cleaning method for the vacuum system; however, do not attempt it when pockets and crevices exist that will prevent thorough rinsing. Degreasing with an acetone or methanol followed by water-rinsing is usually sufficient.
2. To achieve maximum pumping speed, connect the pump to the system with a tube that has the shortest length and the largest diameter possible.
3. For ultrahigh vacuum work, joints should be tungsten inert gas welded, hydrogen brazed, or vacuum brazed. Do not use flux in making joints for ultrahigh vacuum.
4. Gaskets in the system should be of soft metal such as annealed copper, aluminum, or gold.
5. After the pump is sealed in place, leak-check the entire system.
6. Assemble the magnet to the pump.

Control Unit Connection

WARNING!



The high voltage present in the high voltage cable which connects the control unit to the ion pump, can cause severe injury or death. Before mounting the high voltage connector of the cable on the pump high voltage feedthrough, or before removing it, be sure the main power is removed from the control unit.

WARNING!



To avoid injury, never connect the high voltage to the pump before it is installed into the system and all the inlet flanges are properly connected or blanked off. Make sure that the pump is well connected to the grounded vacuum system.

Connect the control unit to the ion pump with the coaxial high voltage cable assembly as follows:

1. Push the female end of the cable connector over the high voltage feedthrough.
 2. Push the male end of the cable connector into the socket on the control unit rear panel. (Refer to the control unit instruction manual).
-

WARNING!



Before removing the high voltage connector of the cable from the control unit, be sure the main power is removed from the control unit. Wait at least 10 seconds after removing the main power from the control unit, to allow capacitors to discharge completely.

To disconnect the coaxial high voltage cable from the controller, slide the safety locking sleeve (very little sleeve travel is required) from the control unit and at the same time pull on the male end of the cable connector to remove it from the socket on the control unit.

Bakeout

When a 2 l/s ion pump does not reach the desired base pressure, and there are no leaks, it is necessary to bake the system. This can be done by heating the pump and all the components in the system, and is generally required to achieve less than 10^{-8} mbar base pressures.

- The pump body can be baked up to 400 °C if the magnet and the H.V. cable connector have been removed.
- If the bakeout is done with magnets and H.V. cable connected, the maximum allowable temperature is 150 °C and the pump control unit can be left on in order to be able to monitor the pressure variations. During the bakeout cycle the pressure must not increase above 5×10^{-5} mbar, if this value is reached, turn the bakeout off and then on again when low pressure is restored.
- Bake the 2 l/s pump for 24 hours. Longer bakeout periods are recommended when the pump has been used with heavy gas load applications.
- Wait until the pump cools down to room temperature and recovers its initial pressure before using it in the application.

Unusual Operating Conditions

If the pump current reads lower than normal for the pressure, the discharge may have been extinguished.

To check for the pump being "out of strike" (discharge extinguished), remove the magnet and observe the residual current. If the current is the same with and without the magnet in place, no discharge exists in the pump.

5 Technical Information

Bakeout

1. To restart the discharge, tap the pump body with a penny held between the thumb and forefinger, or heat the pump body locally with a small soldering iron. This should liberate surface gas within the pump and restart the discharge. NEVER strike the pump with hand tools.
2. If the magnet strength is below its rated level, the pump discharge and speed will be reduced and the discharge may extinguish at a pressure below 1×10^{-7} mbar.

Pump Troubleshooting

In the event of a power failure (momentary or long period) the controller is switched off. When power is restored, the controller will automatically restart.

Tab. 3

SYMPTOM	POSSIBLE CAUSE	CORRECTION PROCEDURE
1. Slow starting (more than 30 minutes).	Air leaks which limit pressure to above 10^{-6} mbar and cause longer starting time.	Leak check the vacuum system with a helium leak detector.
2. Slow pump-down due to absorption of vapours on pump and system walls.	Vapours and gases admitted to a system are absorbed on the walls of the system and pump. Subsequent reduction in pressure depends on the rate of depletion of this vapour. Heavy hydrocarbons are most troublesome because of their relative low vapour pressure and are very difficult to remove, even by baking.	Heat the system walls, thereby accelerating the desorption process. Baking mobilizes the vapours so they can be cracked and pumped by discharge.
3. Slow starting or slow pump-down.	High voltage feedthrough is leaking.	Replace the feedthrough.
4. Current higher than expected at any given pressure.	Ion pump leakage current causing higher pressure reading.	Highpot the pump.

2 l/s Pump Replacement Parts and Accessories

Tab. 4

PART NUMBER					
Basic pump	9190520	9190521	9190522	9190523	9190524
Magnet Cobalt-Samarium			9190038		
CABLES					
HV Bakeable cable, radiation resistant, 13' (4 m) long with safety interlock	9290706				

For a complete overview of Agilent's extensive vacuum product line,
please refer to the Agilent Vacuum Catalogue.



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.Illi 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.

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

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

Product preparation

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

Shipping

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

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

EUROPE:

Fax: 00 39 011 9979 330

Fax Free: 00 800 345 345 00

Toll Free: 00 800 234 234 00

vpt-customerccare@agilent.com

NORTH AMERICA:

Fax: 1 781 860 9252

Toll Free: 800 882 7426

vpl-ra@agilent.com

PACIFIC RIM:

please visit our website for individual
office information

<http://www.agilent.com>



TERMS AND CONDITIONS

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

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

Vacuum Products Division Request for Return Form

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

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

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

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

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

I declare that the above information is true and complete to the best of my knowledge and belief. I understand and agree to the terms and conditions on page 2 of this document.	
Name:	Authorized Signature:
Position:	
Date:	
NOTE: If a product is received at Agilent which is contaminated with a toxic or hazardous material that was not disclosed, the customer will be held responsible for all costs incurred to ensure the safe handling of the product, and is liable for any harm or injury to Agilent employees as well as to any third party occurring as a result of exposure to toxic or hazardous materials present in the product.	

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies

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

Netherlands

Agilent Technologies Netherlands B.V.

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

Belgium

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

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

Brazil

Agilent Technologies Brasil

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

China

Agilent Technologies (China) Co. Ltd

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

France

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

Southeast Asia

Agilent Technologies Sales Sdn Bhd

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

India (Sales)

Agilent Technologies India Pvt. Ltd.

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

Italy

Agilent Technologies Italia S.p.A.

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

Japan

Agilent Technologies Japan, Ltd.

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

Singapore

Agilent Technologies Singapore Pte. Ltd,

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

Korea

Agilent Technologies Korea, Ltd.

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

UK and Ireland

Agilent Technologies LDA UK Limited

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

© Agilent Technologies, Inc. 2019
06/2019

India (Service)

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

Taiwan

Agilent Technologies Taiwan Limited

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

Germany and Austria

Agilent Technologies

Sales & Services GmbH & Co. KG

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

Mexico

Agilent Technologies

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

Other Countries

Agilent Technologies Italia S.p.A.

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

Customer Support & Service

NORTH AMERICA:

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

EUROPE:

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

PACIFIC RIM:

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

**Worldwide Web Site, Catalog and Order On-
line:**

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