

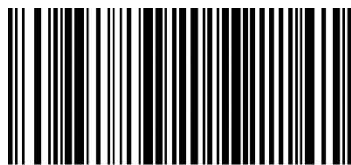


**Agilent Technologies**

*Vacuum Products Division*

# ***IDP-3 Tip Seal Replacement***

*INSTRUCTION MANUAL*



X3803-90002

Part No. X3803-90002

Rev. B

April 2017

# ***IDP-3 Tip Seal Replacement***

# Warranty

Products manufactured by Seller are warranted against defects in materials and workmanship for twelve (12) months from date of shipment thereof to Customer, and Seller's liability under valid warranty claims is limited, at the option of Seller, to repair, replacement, or refund an equitable portion of the purchase price of the Product. Items expendable in normal use are not covered by this warranty. All warranty replacement or repair of parts shall be limited to equipment malfunctions which, in the sole opinion of Seller, are due or traceable to defects in original materials or workmanship. All obligations of Seller under this warranty shall cease in the event of abuse, accident, alteration, misuse, or neglect of the equipment. In-warranty repaired or replaced parts are warranted only for the remaining unexpired portion of the original warranty period applicable to the repaired or replaced parts. After expiration of the applicable warranty period, Customer shall be charged at the then current prices for parts, labor, and transportation.

When products are used with toxic chemicals, or in an atmosphere that is dangerous to the health of humans, or is environmentally unsafe, it will be the responsibility of the Customer to have the product cleaned by an independent agency skilled and approved in handling and cleaning contaminated materials before the product will be accepted by Agilent, Inc. for repair and/or replacement.

Reasonable care must be used to avoid hazards. Seller expressly disclaims responsibility for loss or damage caused by use of its Products other than in accordance with proper operating procedures.

Except as stated herein, Seller makes no warranty, express or implied (either in fact or by operation of law), statutory or otherwise; and, except as stated herein, Seller shall have no liability under any warranty, express or implied (either in fact or by operation of law), statutory or otherwise. Statements made by any person, including representatives of Seller, which are inconsistent or in conflict with the terms of this warranty shall not be binding upon Seller unless reduced to writing and approved by an officer of Seller.

## Warranty Replacement and Adjustment

All claims under warranty must be made promptly after occurrence of circumstances giving rise thereto, and must be received within the applicable warranty period by Seller or its authorized representative. Such claims should include the Product serial number, the date of shipment, and a full description of the circumstances giving rise to the claim. Before any Products are returned for repair and/or adjustment, written authorization from Seller or its authorized representative for the return and instructions as to how and where these Products should be returned must be obtained. Any Product returned to Seller for examination shall be prepaid via the means of transportation indicated as acceptable by Seller. Seller reserves the right to reject any warranty claim not promptly reported and any warranty claim on any item that has been altered or has been returned by non-acceptable means of transportation. When any Product is returned for examination and inspection, or for any other reason, Customer shall be responsible for all damage resulting from improper packing or handling, and for loss in transit, notwithstanding any defect or non-conformity in the Product. In all cases, Seller has the sole responsibility for determining the cause and nature of failure, and Seller's determination with regard thereto shall be final.

If it is found that Seller's Product has been returned without cause and is still serviceable, Customer will be notified and the Product returned at its expense; in addition, a charge for testing and examination may be made on Products so returned.

## IDP-3 Tip Seal Replacement

---

*This page intentionally left blank.*

# Preface

## Documentation Conventions

This manual uses the following documentation conventions:

**WARNING**



*Warnings indicate a particular procedure or practice, which if not followed correctly, could lead to serious injury.*

**CAUTION**



*Cautions indicate a particular procedure or practice, which if not followed, could cause damage to the equipment.*

**NOTE**



*Notes contain important information.*

Before operating or servicing equipment, read and thoroughly understand all operation/maintenance manuals provided by Agilent. Be aware of the hazards associated with this equipment, know how to recognize potentially hazardous conditions, and how to avoid them. Read carefully and strictly observe all cautions and warnings. The consequences of unskilled, improper, or careless operation of the equipment can be serious.

In addition, consult local, state, and national agencies regarding specific requirements and regulations. Address any safety, operation, and/or maintenance questions to your nearest Agilent office.

## IDP-3 Tip Seal Replacement

---

*This page intentionally left blank.*

# IDP-3 Tip Seal Replacement

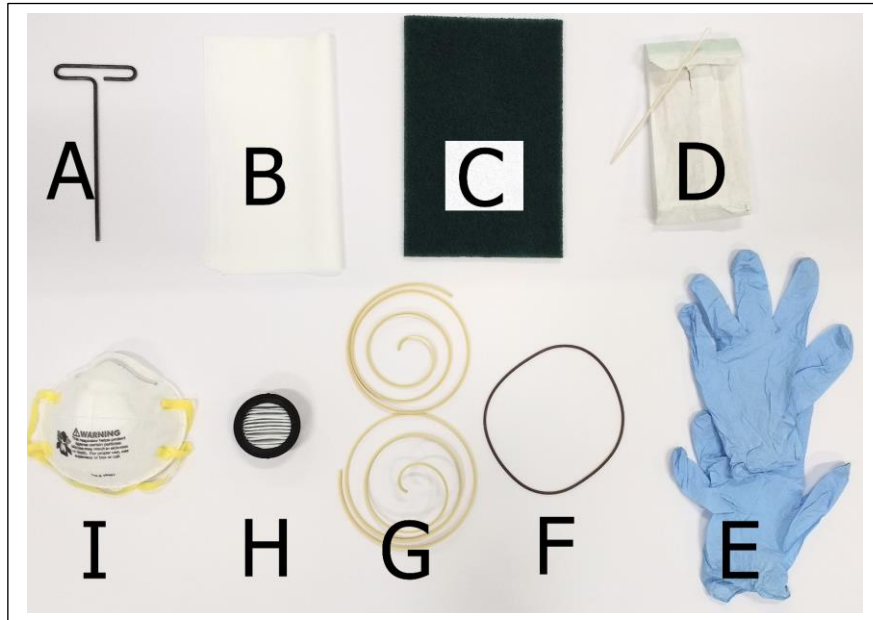


Figure 1 IDP-3 Tip Seal Replacement Kit

Table 1 IDP-3 Tip Seal Replacement Kit Contents

Part Label	Part Name	Qty.
A	4mm T-handle Allen wrench	1
B	Low lint cleaning wipes	1 pkg.
C	Scotch Brite™ 9" by 6" piece	1
D	Low Lint Cleaning Swabs	1 pkg.
E	Nitrile gloves (color may vary)	1 pr.
F	O-ring	1
G	Tip seal	1 set (2 pc.)
H	Silencer Filter	1
I	Dust mask	1

To prevent any dust inhalation, Agilent recommends a dust mask to be used during IDP-3 maintenance. Additional tools and supplies which may be used (razor, needle nose pliers, flat-head screw driver, adjustable wrench, and isopropyl alcohol) are not included.

## IDP-3 Tip Seal Replacement

### Installation

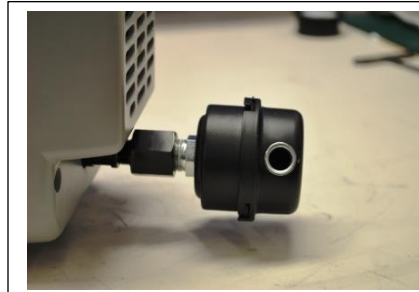
**NOTE**

*To ensure proper operation and prevent damage, the kit must be installed and operated only as specified.*

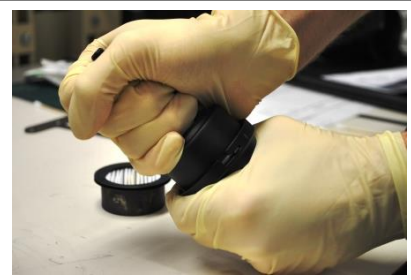
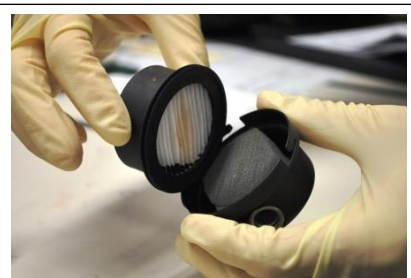


To install:

1. Remove the muffler assembly from the front of the IDP-3 if muffler assembly option is equipped. If an optional barb fitting is equipped instead, remove it and set aside.



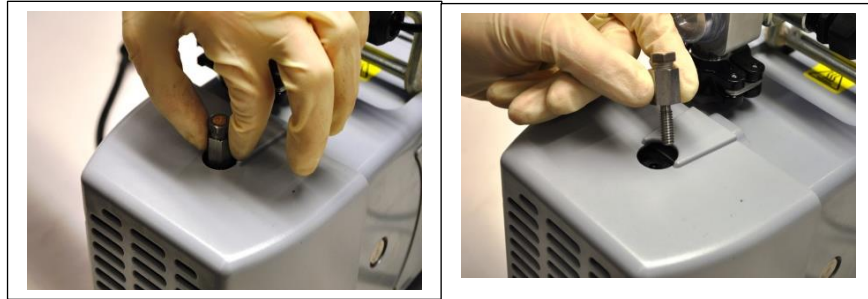
2. Remove the Filter from the muffler assembly (if equipped) and replace it with the new filter.





### IDP-3 Tip Seal Replacement

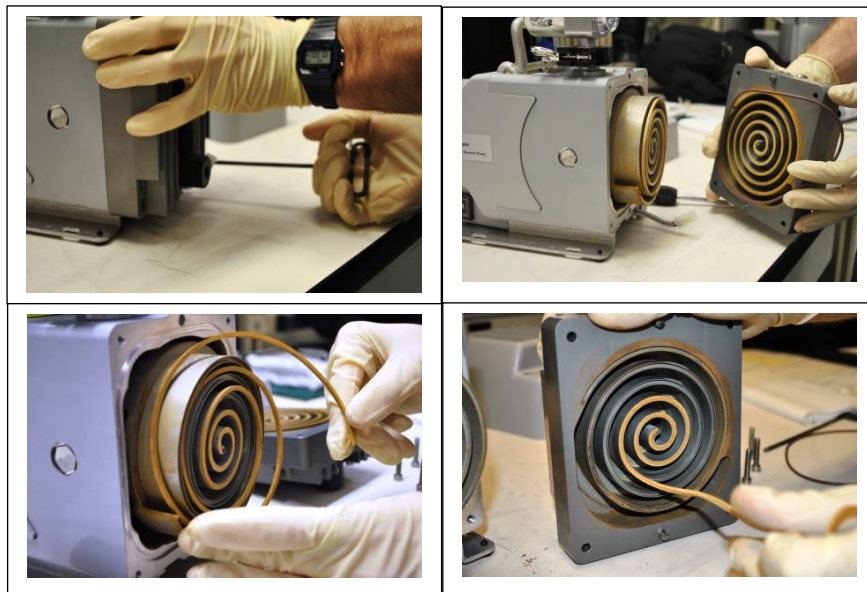
3. Unscrew and remove the sintered plug from the top of the IDP-3. An adjustable wrench may be used if necessary.



4. Using the 4mm T handled hex key, remove the four M5 screws that hold the front cowling on to the IDP-3 and disconnect the cable that powers the fan.

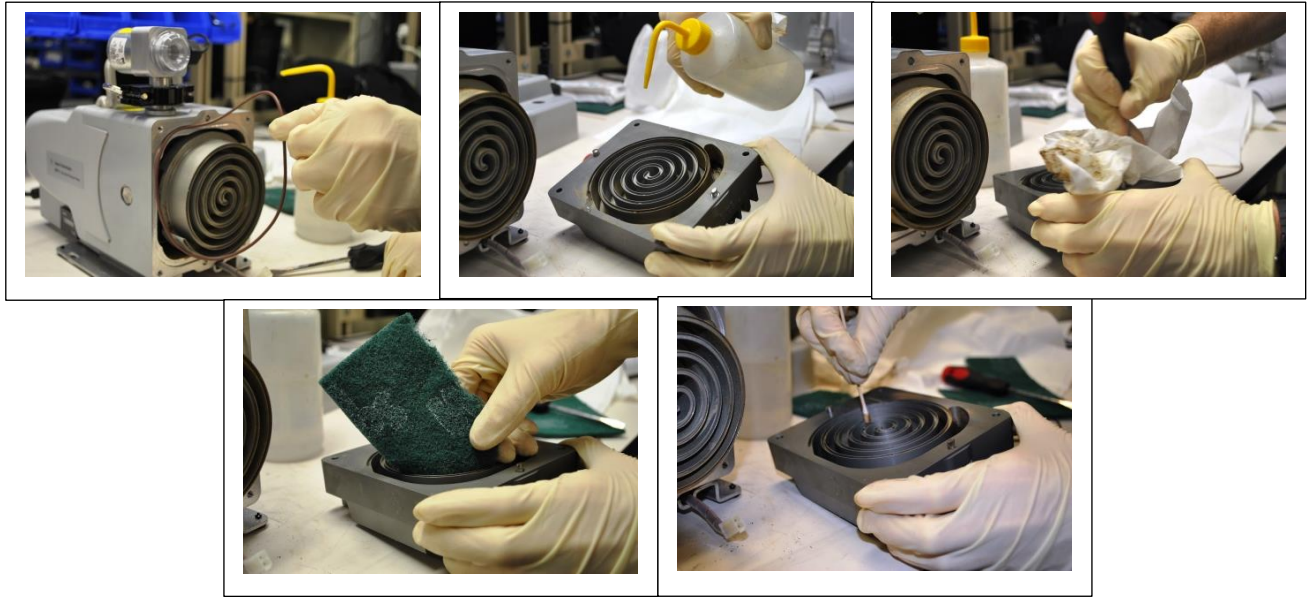


5. Using a 4mm hex key, remove the four M5 screws that hold on the Outboard Housing and remove the housing from the IDP-3. Remove the tip seal from both the outboard housing and the orbiting plate.

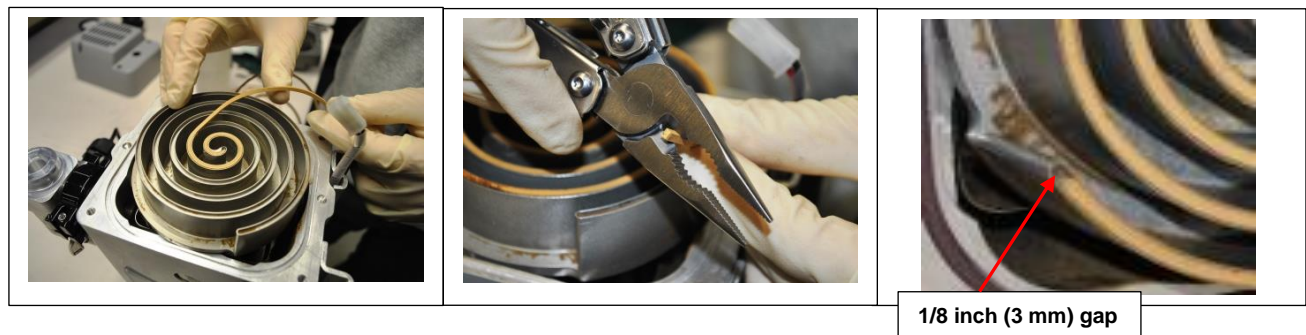


## IDP-3 Tip Seal Replacement

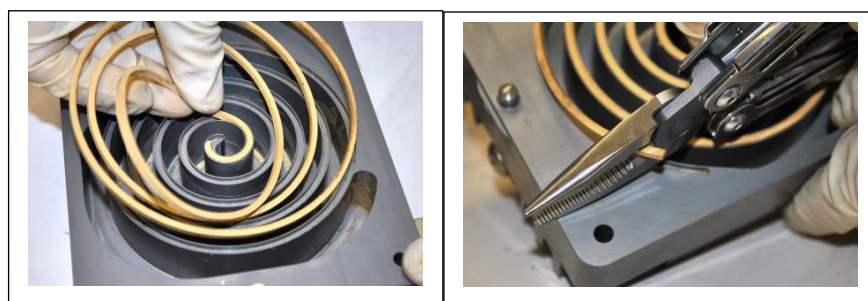
- Remove the O-ring that provides the seal between the orbiting plate and the outboard housing from the pump frame. Clean the grooves and interior of both the Outboard Housing and the Orbital Plate with a Kim Wipe or equivalent, lint free cotton swab or equivalent, Scotch-Brite as required and isopropyl alcohol (use if necessary or desired: not included) to remove tip seal debris. If in a cleanroom environment, do not use compressed air or canned air to remove tip seal dust from the interior of the pump. Do not use water, detergent, or home/industrial cleaning products.



- Position the pump with the orbiting plate facing up. If possible, using a pair of needle nose pliers, form fit the tip seal to match the tip seal groove at the center of the orbiting plate. Install the tip seal from the center of the Orbiting plate and work it outward. When the tip seal groove is filled, cut off the excess, leaving 1/8<sup>th</sup> of an inch (3 mm) of open space in the groove to allow for expansion of the tip seal.



- Repeat the installation process with the outboard housing.





## IDP-3 Tip Seal Replacement

9. Install the new O-ring from the kit into the O-ring groove on the IDP-3 pump frame. Re-mount the Outboard housing onto the IDP-3, using the guide pins on the housing. Make sure that the tip seal does not fall out of the groove and that the O-ring remains seated in the groove.



10. Using the 4mm hex key, re-install the four M5 screws that hold on the Outboard Housing to the pump frame. Torque to approximately 5 Nm if a torque wrench is available. Re-connect the fan connector that goes from the front cowling to the frame of the pump. Re-install the front cowling to the pump using the four M5 screws removed previously.



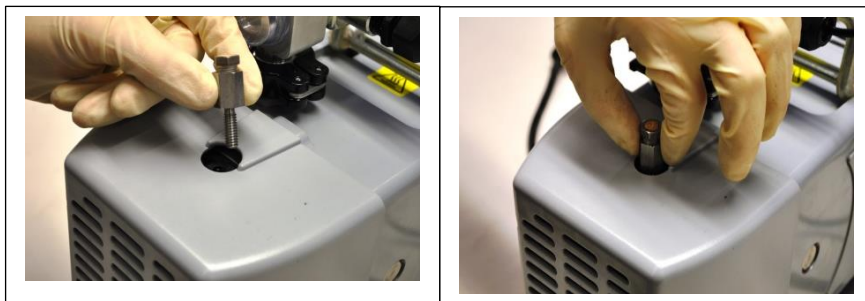
11. Re-install the muffler assembly into the port and hand tighten.



## IDP-3 Tip Seal Replacement

---

12. Re-install the sintered plug into the top of the pump.



13. Power the pump up to insure it is functioning correctly.



For more detailed instructions, search "**IDP3 TipSeal**" in YouTube and follow Agilent channels:





## Agilent Technologies

## 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 step for the shipment of returning product:
  - x Remove all accessories from the core product (e.g. inlet screens, vent valves).
  - x Prior to shipment, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
  - x If ordering an Advance Exchange product, please use the packaging from the Advance Exchange to return the defective product.
  - x 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.
  - x Agilent Technologies is not responsible for returning customer provided packaging or containers.
  - x 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: 00390119979330  
Fax Free: 00 800 345 345 00  
Toll Free: 0080023423400  
[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

**NORTH AMERICA:**  
Fax: 17818609252  
Toll Free: 8008827426, Option 3  
[vpl-ra@agilent.com](mailto: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**

<b>Company Name:</b>		<b>Contact Name:</b>	
<b>Tel:</b>	<b>Email:</b>	<b>Fax:</b>	
<b>Customer Ship To:</b>		<b>Customer Bill To:</b>	
<b>Europe only: VAT reg. Number:</b>		<b>USA/Canada only:</b> <input type="checkbox"/> <b>Taxable</b> <input type="checkbox"/> <b>Non-taxable</b>	

**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
<ul style="list-style-type: none"> <li>- Does not start</li> <li>- Does not spin freely</li> <li>- Does not reach full speed</li> <li>- Mechanical Contact</li> <li>- Cooling defective</li> </ul>	<ul style="list-style-type: none"> <li>- Noise</li> <li>- Vibrations</li> <li>-Leak</li> <li>-Over temperature (</li> <li>-Clogging</li> </ul>	<ul style="list-style-type: none"> <li>- Vertical</li> <li>-Horizontal</li> <li>-Upside-down</li> <li>-Other:</li> <li>.....</li> </ul>
		Power:                      Rotational Speed: Current:                   Inlet Pressure: Temp 1:                    Foreline Pressure: Temp 2:                    Purge flow: OPERATING TIME:

**ION PUMPS/CONTROLLERS**

<ul style="list-style-type: none"> <li>- Bad feedthrough</li> <li>- Vacuum leak</li> <li>- Error code on display</li> </ul>	<ul style="list-style-type: none"> <li>- Poor vacuum</li> <li>- High voltage problem</li> <li>- Other</li> </ul>
---	--

**VALVES/COMPONENTS**

<ul style="list-style-type: none"> <li>- Main seal leak</li> <li>- Solenoid failure</li> <li>- Damaged sealing area</li> </ul>	<ul style="list-style-type: none"> <li>- Bellows leak</li> <li>- Damaged flange</li> <li>-Other</li> </ul>
--	--

**LEAK DETECTORS**

<ul style="list-style-type: none"> <li>- Cannot calibrate</li> <li>- Vacuum system unstable</li> <li>- Failed to start</li> </ul>	<ul style="list-style-type: none"> <li>-No zero/high background</li> <li>- Cannot reach test mode</li> <li>- Other</li> </ul>
---	---

**INSTRUMENTS**

<ul style="list-style-type: none"> <li>- Gauge tube not working</li> <li>- Communication failure</li> <li>- Error code on display</li> </ul>	<ul style="list-style-type: none"> <li>- Display problem</li> <li>- Degas not working</li> <li>- Other</li> </ul>
--	---

**SCROLL AND ROTARY VANE PUMPS**

<ul style="list-style-type: none"> <li>- Pump doesn't start</li> <li>- Doesn't reach vacuum</li> <li>- Pump seized</li> </ul>	<ul style="list-style-type: none"> <li>- Noisy pump (describe)</li> <li>- Over temperature</li> <li>- Other</li> </ul>
---	--

**DIFFUSION PUMPS**

<ul style="list-style-type: none"> <li>- Heater failure</li> <li>- Doesn't reach vacuum</li> <li>- Vacuum leak</li> </ul>	<ul style="list-style-type: none"> <li>- Electrical problem</li> <li>- Cooling coil damage</li> <li>- Other</li> </ul>
---	--

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.**

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

## IDP-3 Tip Seal Replacement

---

*This page intentionally left blank*



# Agilent Technologies Vacuum Product Division

## United States

Agilent Technologies  
121 Hartwell Avenue  
Lexington, MA 02421 - USA  
Tel. +1 781 861 7200  
Fax. +1 781 860 5437  
Toll-Free: +1 800 882 7426  
[vpl-customer@agilent.com](mailto:vpl-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](mailto:customer@agilent.com)

## Brazil

Agilent Technologies Brasil  
Avenida Marcos Penteado de Ulhoa  
Rodrigues, 939 - 6º andar  
Castelo Branco Office Park  
Torre Jacarandá - Tamboré  
Barueri, São Paulo CEP: 06460-040  
Toll-Free: +0800 728 1405

## China

Agilent Technologies (China) Co. Ltd  
No.3, Wang Jing Bei Lu, Chao Yang District  
Beijing, 100102, China  
Tel. +86 (0)10 64397888  
Fax. +86 (0)10 64392765  
Toll free: 800 820 6778 (Phone)  
Toll free: 400 820 6778 (Cell Phone)  
[vacuum.cnmarketing@agilent.com](mailto:vacuum.cnmarketing@agilent.com)  
[vpc-customerservice@agilent.com](mailto: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](mailto:vpf.sales@agilent.com)

## Germany and Austria

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

## 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  
Tel. +91 11 4623 7100  
Fax. +91 4623 7105

## 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  
Tel. +91 80 40614000  
Fax. +91 80 40148991

## 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: 00800 234 234 00  
[vpt-customerservice@agilent.com](mailto:vpt-customerservice@agilent.com)

## Japan

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

## 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

## 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@netherlands.agilent.com](mailto:customer@netherlands.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](mailto:vps-customerservice@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  
Tel. +603 7712 6181  
Fax. +603 7727 1239  
Toll-Free: +1 800 880 805  
[vps-customerservice@agilent.com](mailto:vps-customerservice@agilent.com)

## Taiwan

Agilent Technologies Taiwan Limited  
No. 20 Gao-shuang Road, Ping-zhen Dis  
Tao-Yuan City  
32450 Taiwan, R.O.C.  
Tel. +8863 3 4959004  
Toll-Free: 0800 018 768  
[vpw-customerservice@agilent.com](mailto:vpw-customerservice@agilent.com)

## UK and Ireland

Agilent Technologies LDA UK Limited

Lakeside Cheadle Royal Business Park  
Cheadle, Cheshire SK8 3GR,  
United Kingdom  
Tel. +44 01865291570  
Fax. +44 01865291571  
Toll-Free: 00 800 234 234 00  
Toll-Free fax.: 00 800 345 345 00  
[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

## 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](mailto:vpt-customer@agilent.com)

## Customer Support & Service

**NORTH AMERICA:**  
Toll Free: 800 882 7426  
[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

## **EUROPE:**

Toll-Free: 00 800 234 234 00  
[vpt-customer@agilent.com](mailto: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 Online:  
[www.agilent.com/chem/vacuum](http://www.agilent.com/chem/vacuum)  
Representatives in most countries