

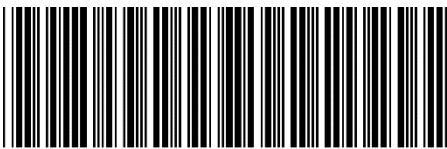


**Agilent Technologies**

*Vacuum Products Division*

***IDP 7 / IDP 10  
VPI Valve Minor  
Replacement  
Kit***

*INSTRUCTION MANUAL*



\*X3807-90015\*

Manual No. X3807-90015  
Rev. A  
February 2017

# 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 7 / IDP 10 VPI Valve Minor Replacement Kit

---

### IDP 7 / IDP 10 VPI Valve Minor Replacement Kit

Part Name	Qty
Rubber thermal protector	1
Orifice	1
Diaphragm retainer plate	1
Piston	1
O-ring 2-151	1
O-ring 2-121	1
O-ring 2-008	5
Screw	8

Table 1-1

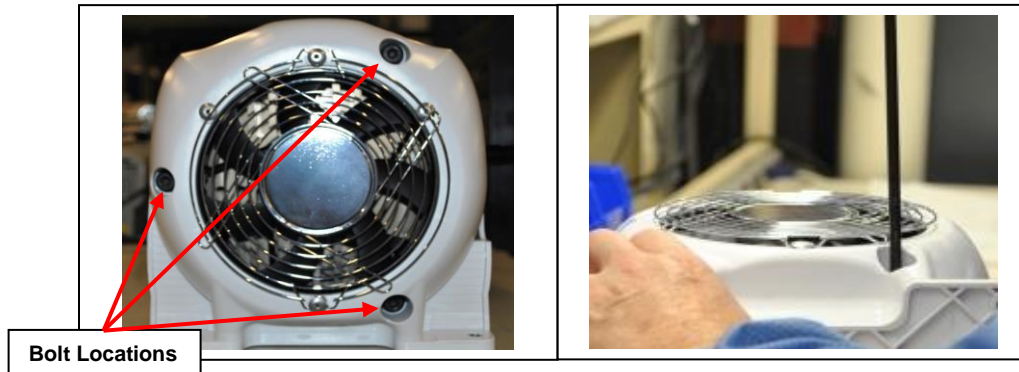
### Installation :



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

## IDP 7 / IDP 10 VPI Valve Minor Replacement Kit

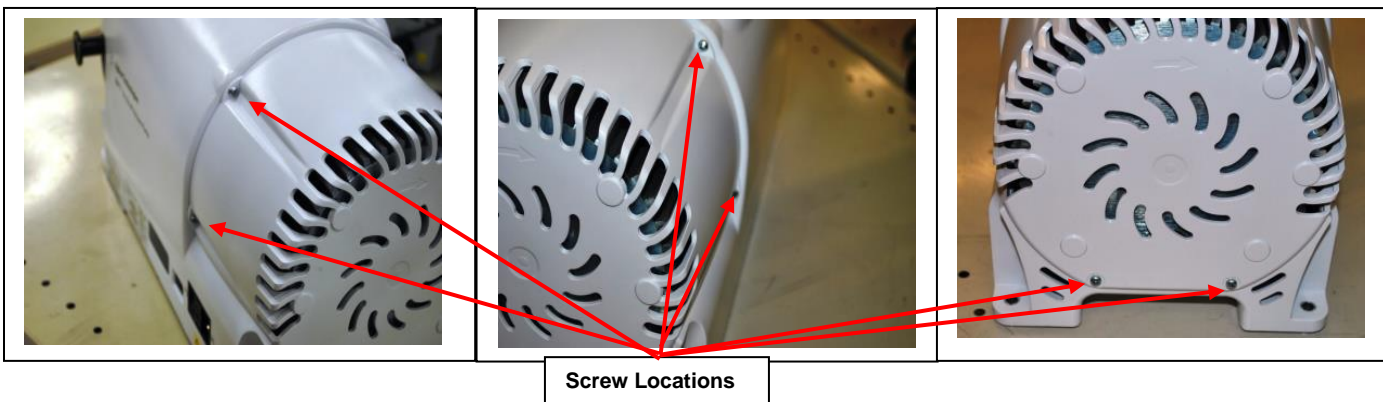
1. Detach the front cowling from the front of the pump by removing three M8x20mm flange bolts using a 6mm T handled hex key.



2. Detach the fan cable from the front cowling and remove the cowling from the pump.



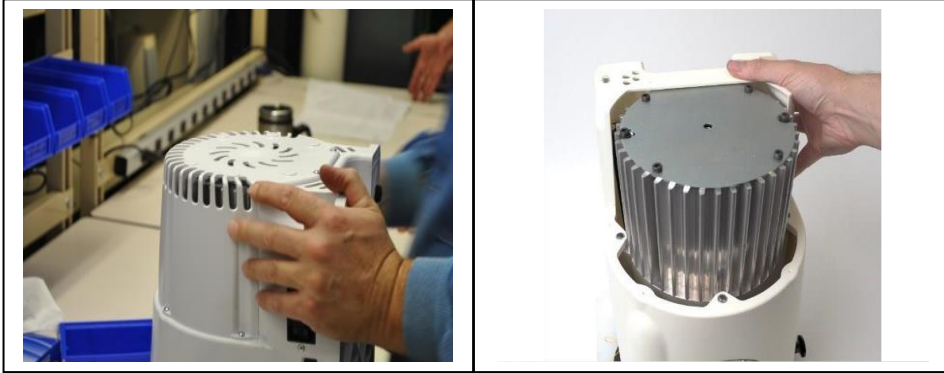
3. Disassemble the rear cowling by remove the six self-tapping screws and two washers that hold the rear cowling to the chassis and front cowling using a Philips head screwdriver.



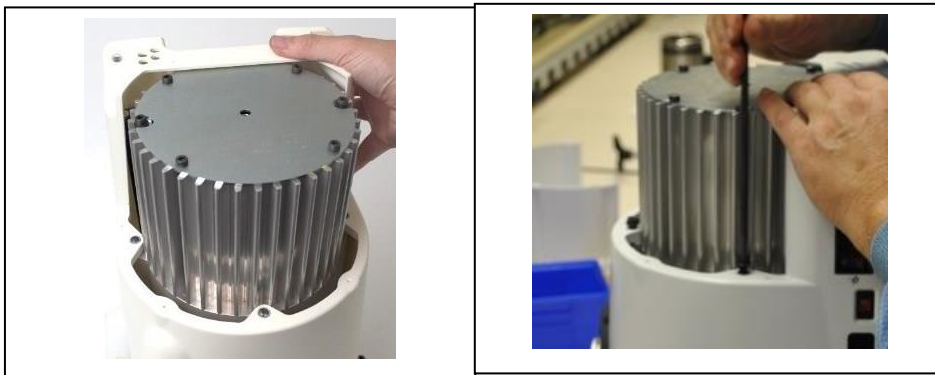
## ***IDP 7 / IDP 10 VPI Valve Minor Replacement Kit***

---

4. Remove the Rear Cowling from the pump assembly.



5. Loosen the three M6x20mm that secure the top part of the base to the Pump Head.

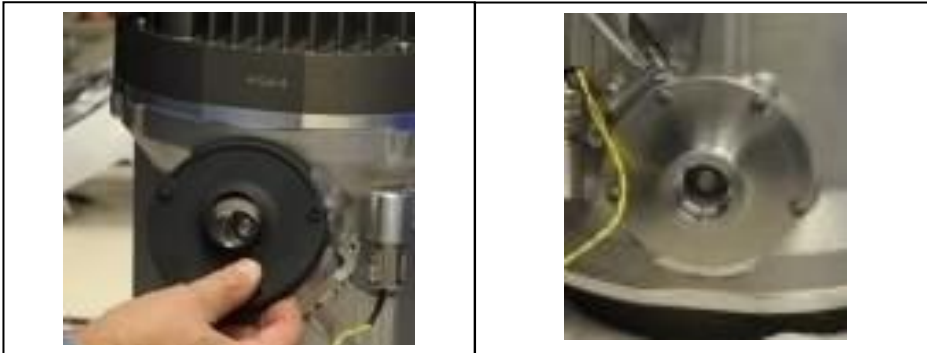


6. Loosen and remove the two M6x16 mm bolts on the bottom of the base to the pump module using a 5mm hex key. Move the base assembly up enough to facilitate removal of the Rubber Thermal Protector.



## IDP 7 / IDP 10 VPI Valve Minor Replacement Kit

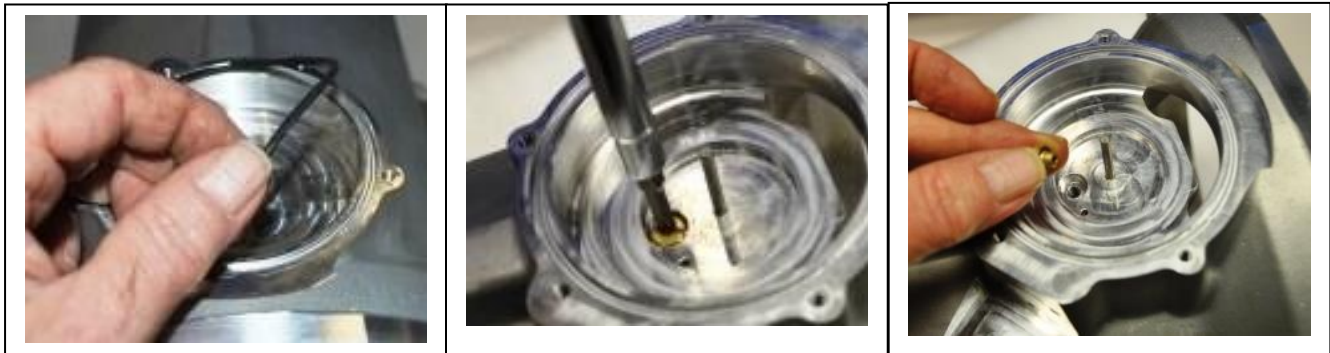
7. Remove the thermal rubber protector from the inlet assembly and discard.



8. Loosen the four M4x10mm screws holding the inlet flange on to the frame. Remove the inlet flange from the top of the frame assembly. Remove the VPI spring and diaphragm/piston from the frame assembly. Discard the piston assembly.



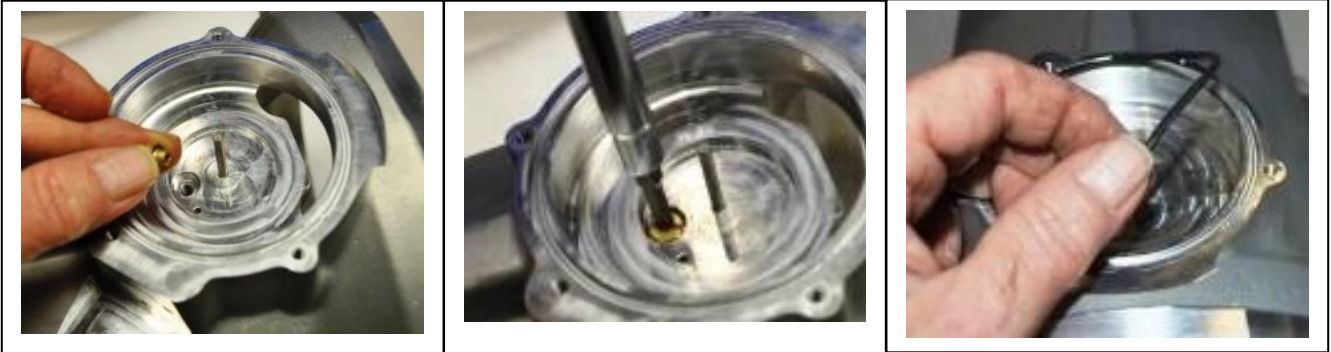
9. Remove the O-ring from the assembly and discard. Remove the 10-32 threaded orifice and discard.



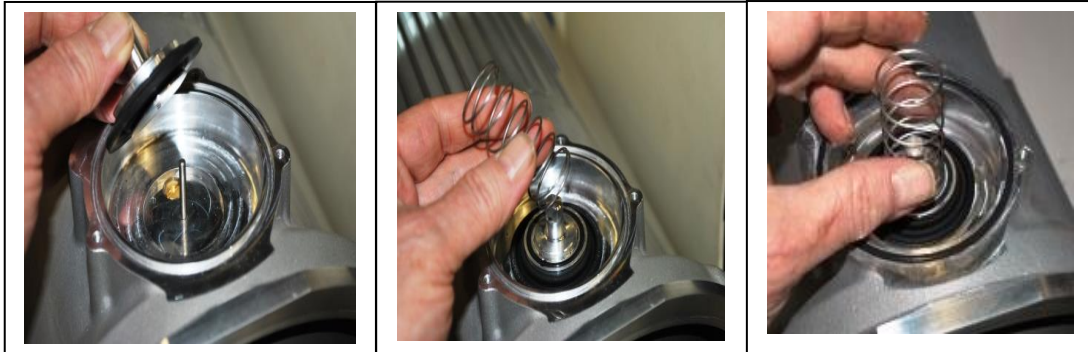
## *IDP 7 / IDP 10 VPI Valve Minor Replacement Kit*

---

10. Install a new 10-32 threaded orifice. Apply LVP grease to a new O-ring and install it into the O-ring groove.



11. Install a new O-ring from the kit into the diaphragm/piston assembly. Install the new diaphragm/piston assembly from the kit into the frame. Install the spring onto the diaphragm/piston assembly.



12. Install the Inlet Flange into the frame assembly and secure with four M4x10 mm screws. Torque to 1.6 N.m if a torque wrench is available.



## ***IDP 7 / IDP 10 VPI Valve Minor Replacement Kit***

---

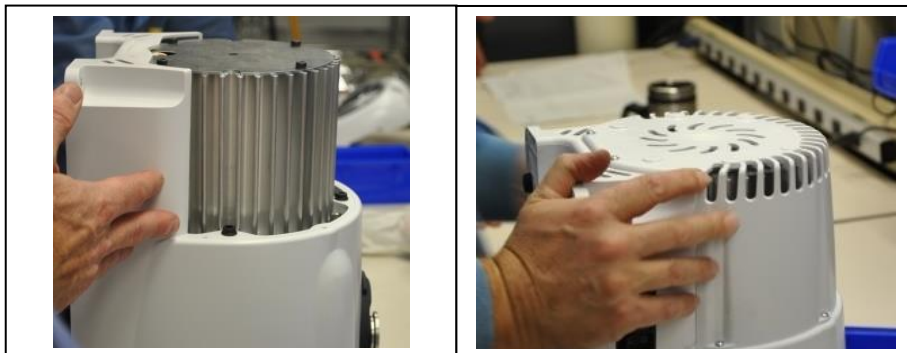
13. Install a new Thermal Rubber Protector onto the inlet assembly.



14. Reseat the base assembly. Verify that all connections are in place and that there are no pinched wires. Tighten the five screws that hold the base assembly in place.



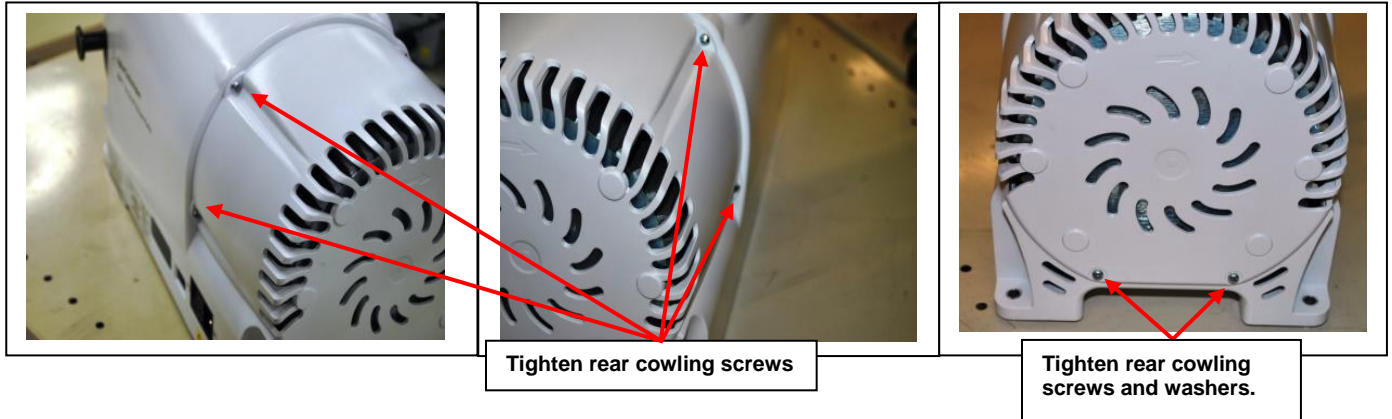
15. Reinstall the rear cowling on to the pump assembly.



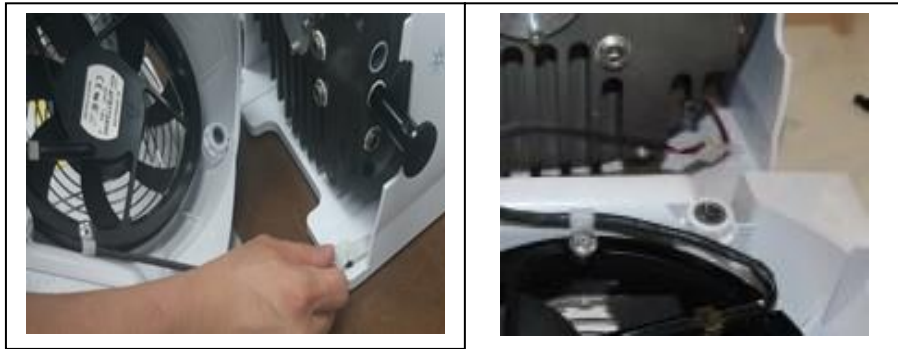


## IDP 7 / IDP 10 VPI Valve Minor Replacement Kit

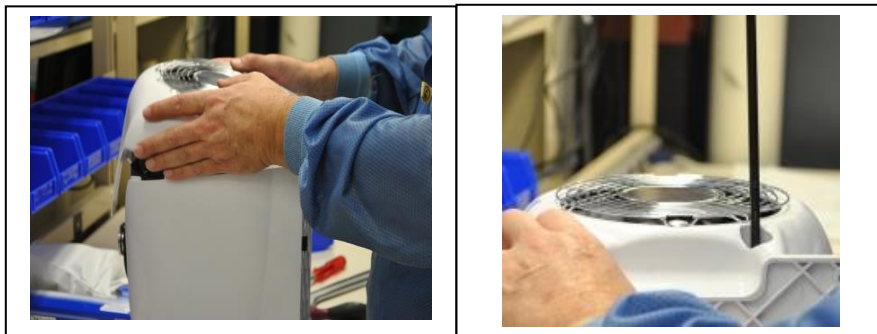
16. Install the six self-tapping screws and two washers that hold the rear cowling to the chassis and front cowling.



17. Attach the fan cable from the front cowling to the motor.



18. Attach the front cowling to the pump using three M8x20mm flange bolts. Tighten with 6mm Hex Key. Torque screws to 11.3 Nm if a torque wrench is available.



## ***IDP 7 / IDP 10 VPI Valve Minor Replacement Kit***

---

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





## **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:
  - 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 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: 00390119979  
330  
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: 800 882 7426, Option 3  
[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)

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



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>		USA/Canada only: <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. Checkboxes 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):

<b>I understand and agree to the terms of Section 6, Page 3/3.</b>		
<b>Print Name:</b>	<b>Authorized Signature:</b> .....	<b>Date:</b>



**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: ..... OPERATING TIME:

**ON PUMPS/CONTROLLERS**

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

**VALVES/COMPONENTS**

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

**LEAK DETECTORS**

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

**INSTRUMENTS**

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

**SCROLL AND ROTARY VANE PUMPS**

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

**DIFFUSION PUMPS**

- Heater failure	- Electrical problem
- Doesn't reach vacuum	- Cooling coil damage
- Vacuum leak	- 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.

- 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.

# Agilent Technologies

## Vacuum Product Division

### United States & Canada

*Agilent Technologies*  
*Vacuum Products Division*  
121 Hartwell Avenue  
Lexington, MA 02421 USA  
Tel: +1 781 861 7200  
Toll-Free: +1 800 882 7426  
Fax: +1 871 860 5437

### Benelux

*Agilent Technologies*  
*Vacuum Products Division*  
Herculesweg 8  
4338 PL Middelburg  
THE NETHERLANDS  
Tel: +31 118 671570  
Fax: +31 118 671569

### China

*Agilent Technologies*  
*Vacuum Products Division*  
Room 1648  
Central Tower South Wing  
Beijing Junefield Plaza  
No. 10 XuanWuMenWai Street  
Beijing 100052 P.R. CHINA  
Tel.: +86 (10) 6310 8550  
Toll-Free: 800 820 6556  
Fax: +86 (10) 6310 0141

### France

*Agilent Technologies*  
*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

### Germany & 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

### India

*Agilent Technologies*  
*Vacuum Product Division*  
205-A, "A" wing of Galleria,  
2nd floor, Hiranandani Gardens,  
Powai, Mumbai-400 076, India  
Tel.: +91 22-2570 8595 / 8597  
Fax: +91 22- 2570 8599

### Italy

*Agilent Technologies*  
*Vacuum Products Division*  
via F.lli Varian 54  
10040 Leini, (Torino) ITALY  
Tel.: +39 011 997 9111  
Toll-Free: 00 800 234 234 00  
Fax: +39 011 997 9350

### Japan

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

### Korea

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

### 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

### Southeast Asia

*Agilent Technologies*  
*Vacuum Products Division*  
South East Asia (SEA) - Alex Ho  
H/P: +601 2213 1253  
Fax: +603 6733 8121

### Singapore

*Agilent Technologies*  
*Vacuum Products Division*  
Singapore  
Unit 10-04 Helios  
Biopolis @ one-north  
11 Biopolis Way, 138667  
Singapore  
H/P.: +65 92364988  
Fax: +65 64789603

### Taiwan

*Agilent Technologies*  
*Vacuum Products Division*  
14F-6, No. 77, Hsin Tai Wu Road,  
Sec. 1  
Hsi chih, Taipei Hsien, Taiwan,  
R.O.C.  
Tel.: +886 2 2698 9555  
Toll Free: 0800 051 342  
Fax: +886 2 2698 9678

### UK & Ireland

*Agilent Technologies*  
*Vacuum Products Division*  
6 Mead Road  
Oxford Industrial Park Tel.:  
+44 (0) 1865 291570  
Fax: +44 (0) 1865 291



**Agilent Technologies**

