

## Vacuum Products Division

# IDP-3 Dry Scroll Vacuum Pump

INSTRUCTION MANUAL

Manual No. 699904350 Revision G October 2017



## 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, to replace, or refund of 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.

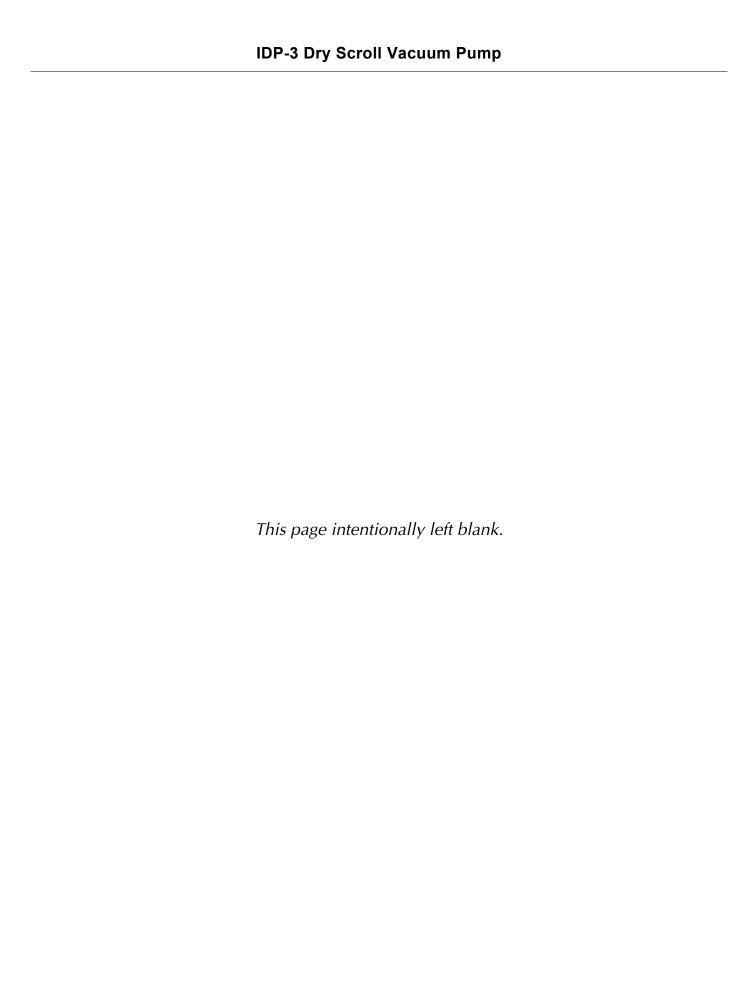
Reasonable care must be used to avoid hazards. Seller expressly disclaims responsibility for loss or damage caused by use of its Products other than in accordance with proper operating procedures.

Except as stated herein, Seller makes no warranty, expressed 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, expressed 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

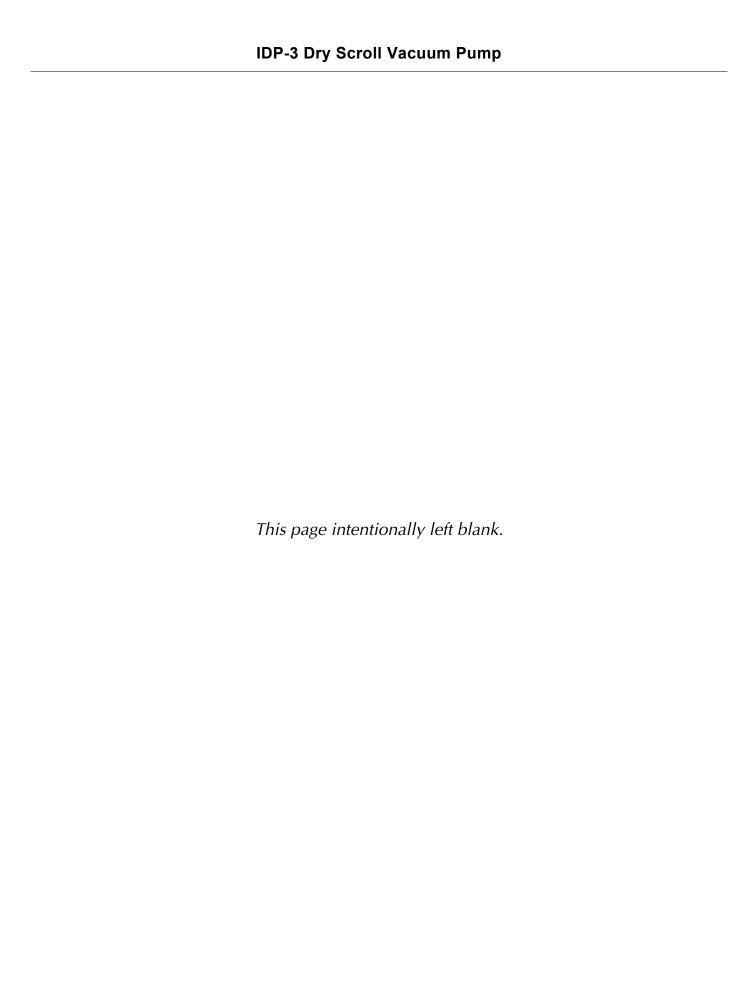
All claims under warranty must be made promptly after occurrence of circumstances giving rise thereto, and must be received within the applicable warranty period by Seller or its authorized representative. Such claims should include the Product serial number, the date of shipment, and a full description 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 the Customer's expense; in addition, a charge for testing and examination may be made on Products so returned.



## **Contents**

Warranty iii	List	of Figures
Warranty Replacement and Adjustment iii  Instructions for Use	Figure	Caption Page
General Information	1	Interface Drawing with Dimensions (AC Version
Storage		
Installation2	2	Interface Drawing with Dimensions (DC Version
Use		
Maintenance	3	Outline Drawing and Principal Items 6
Tb-:	4	24 VDC Power Cable 8
Technical Information	5	Shield Cable 8
Unpacking and Inspection6	6	RPM. vs. 0-10 V Speed Control Signal 9
Installation	7	Grounding Plug and Outlet 10
Safety	8	ON/OFF Power Switch
Startup	9	IDP-3 with Inlet Valve
Electrical Connections	10	Mount Inlet Valve with IDP-3 Pump 14
Run Currents	11	24 VDC Rev Inlet Valve Kit 15
Start Current	12	Tip Seal Seating
Mechanical Connections	13	Exploded View of Pump Body 19
Operation	14	Module Replacement - Exploded View 20
Cleaning the Pump	Table	Title Page
Startup Procedure	Table	3
Shutdown Procedure	1	Specifications
IDP-3 Inlet Valve Kit	2	IDP-3 Power Cord Selection
Model Selection	3	IDP-3 Typical Run Currents (A) 11
Installation	4	Typical IDP-3 Maximum Starting Current and
Troubleshooting		Duration
Maintenance	5	Inlet Valve Kit Selection
Kits and Service Options	6	IDP-3 Model Numbers with Inlet Valve 13
Cleaning	7	Technical Specifications
Tip Seal Replacement	8	Troubleshooting Chart
Pump Module Replacement19	9	Tip Seal Replacement Kit 17
Request for Return Health and Safety Certification		



## Instructions for Use

## **General Information**

This equipment is designed 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 that occur due to non-compliance 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 IDP-3 is a hermetic, dry scroll vacuum pump. This pump is suitable for pumping air or inert gases. The pump is not intended to pump corrosive, explosive, or particulate-forming gases.

The following paragraphs contain all the information necessary to guarantee the safety of the operator when using the equipment. Detailed information is supplied in "Technical Information" on page 3.

This manual uses the following standard safety 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.

## Storage

When transporting and storing the pump, the following environmental requirements should not be exceeded:

Temperature:  $-20 \,^{\circ}\text{C}$  to  $+60 \,^{\circ}\text{C}$  ( $-4 \,^{\circ}\text{F}$  to  $140 \,^{\circ}\text{F}$ ) Relative humidity: 0 to 95% (non-condensing)

## **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.

Total weight of the packing, IDP-3 pump included, is approximately 10.5 kg (23 lbs).

## **WARNING**



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

## **NOTE**



Normal exposure to the environment cannot damage the pump. Nevertheless, it is advisable to keep the pump inlet closed until the pump is installed in the system.

## 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.

If placing the IDP-3 pump inside an enclosure, provide ample room to supply ambient air to both the front and rear air intakes of the pump.

During operation, the following environmental conditions must be respected:

Temperature: +5 °C to +40 °C (41 °F to 104 °F)

Relative humidity: 0 to 95% (non-condensing)

There are four versions of the IDP-3:

□ 100 VAC, 50/60 Hz

□ 115 VAC, 60 Hz

□ 220-230 VAC, 50/60 Hz

□ 24 VDC

## CAUTION



Be certain that your electrical mains power voltage corresponds to that indicated on the rear of the pump.

#### For the:

- □ AC version, connect the pump to the power supply using an IEC-320 style power cord of at least 10 A capacity.
- □ DC version, connect to the power supply using the power cord supplied with the pump.

## **CAUTION**



Never disturb the two hex head bolts on either side of the pump. Disturbing these bolts will cause loss of performance and/or pump damage.

## Use

In order to reach maximum vacuum, the pump must be left running for about an hour with the inlet sealed.

There are no special instructions for starting the pump; it need only be switched on using the On/Off switch.

#### WARNING



The pump is designed for operation with neutral or noncorrosive fluids. It is absolutely forbidden to use it with potentially explosive or inflammable substances.

There are no special instructions for stopping the pump; it need only be disconnected from the electric power source by the On/Off switch.

## **Maintenance**

Personnel responsible for pump operation and maintenance must be well-trained and aware of the accident prevention rules.

## **WARNING**



- □ Death may result from contact with high voltages. Always take extreme care and observe the accident prevention regulations in force.
- □ When machine is powered up, be careful of moving parts and high voltages.
- ☐ If you have to perform maintenance on the pump after a considerable time in operation, allow it to cool as the temperature of the outer surface may be in excess of 60 °C.
- ☐ Always disconnect your power supply to the pump before beginning maintenance work.

## NOTE



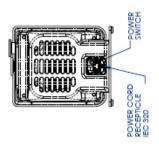
Before returning the pump to the factory for repair, the "Health and Safety" sheet attached to this instruction manual must be completed and sent to the local sales office. A copy of the sheet must be inserted in the pump package before shipping.

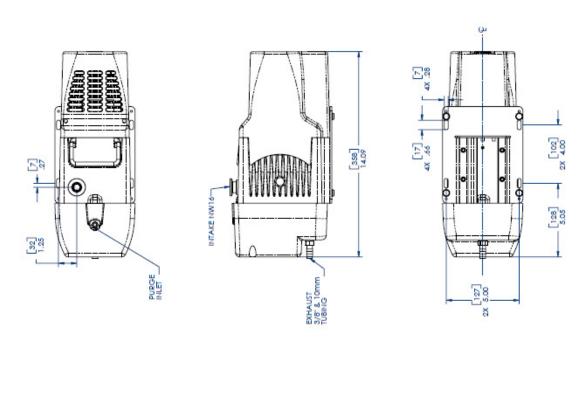
If a pump is to be discarded, it must be disposed of in accordance with specific national standards.

## **Technical Information**

Table 1 Specifications

Model	IDP-3 Dry Scroll Single Hermetic Vacuum Pump	
Interface dimensions	See Figure 1 on page 4	
Peak pumping speed	□ 50 Hz: 50 L/m, 3.0 m <sup>3</sup> /hr (1.8 cfm) □ 60 Hz: 60 L/m, 3.6 m <sup>3</sup> /hr (2.1 cfm) □ 24 VDC: 60 L/m, 3.6 m <sup>3</sup> / hr (2.1 cfm) at full speed setting	
Media	No corrosive, explosive or particulate forming gases	
Ultimate pressure	$2.5 \times 10^{-1}$ Torr (3.3 x $10^{-1}$ mbar)	
Maximum inlet pressure	1.0 atmosphere (0 psig)	
Maximum outlet pressure	6.5 psig	
Inlet connection	NW16	
Exhaust connection	Female 1/4" National Pipe Thread (10 mm hose barb provided)	
Gas ballast	Female 1/8" National Pipe Thread (adapter provided)	
Ambient operating temperature	5 °C to 40 °C (41 °F to 104 °F)	
Storage temperature	-20 °C to 60 °C (-4 °F to 140 °F)	
Motor rating	☐ AC: 0.16 HP (0.12 kW) ☐ DC: 0.16 HP (0.12 KW) Peak rating: 0.27 HP (0.20 KW)	
Operating voltages	□ 1 phase/ 100 VAC, ±10%, 50/60 Hz, 2.31/2.39 FLA □ 1 phase/ 115 VAC, ±10%, 60 Hz, 2.41 FLA □ 1 phase/ 220-230 VAC, ±10%, 50/60 Hz, 1.17/1.26 FLA □ 24 VDC, ±10%, 7.5 FLA	
Run current	See Table 3 on page 11	
Motor thermal protection	Automatic thermal protection	
Operating speed	□ 60 Hz: 3200 RPM, 50 Hz: 2600 RPM □ 24V DC: variable speed, factory setting@ 3200 RPM	
Cooling system	Air-cooled	
Weight	Pump only: 9.5 kg (21 lbs) Shipping weight: 10.5 kg (23 lbs)	
Leak rate (with exhaust and gas ballast sealed)	<1 x 10 <sup>-6</sup> sccs helium	
Vibration level at inlet, per ISO 10816-1	≤ 1.5 mm/sec	
Noise level, per ISO 11201	≤ 55 dBA	
Installation	<ul> <li>□ Pollution Degree: 2</li> <li>□ Installation Category: II</li> <li>□ Altitude 2000 m, indoor usage and storage</li> </ul>	





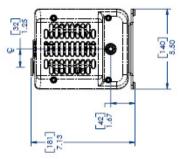


Figure 1 Interface Drawing with Dimensions (AC Version)



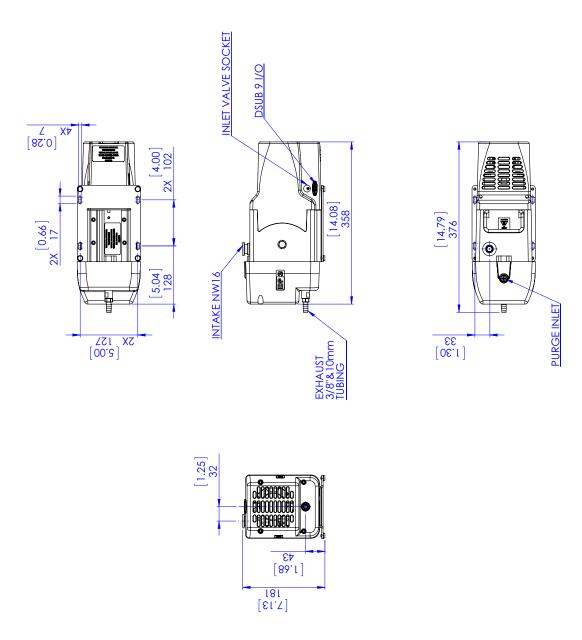
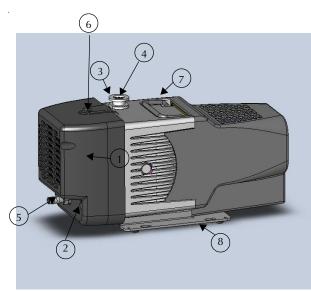
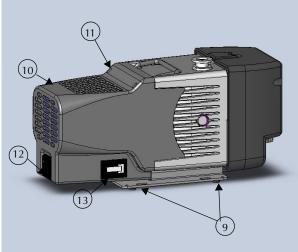


Figure 2 Interface Drawing with Dimensions (DC Version)





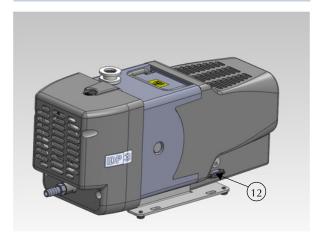


Figure 3 Outline Drawing and Principal Items

- 1. Front Cowling Screws; M5 (4)
- 2. Front Cowling
- 3. Inlet (NW16)
- 4. Inlet Screen
- 5. 10 mm Hose Barb
- 6. Gas Ballast Port (1/8" NPT adapter provided)
- 7. Frame
- 8. Base
- 9. Mounting Holes; (4) for 1/4" or M6 Hardware
- 10. Rear Cowling
- 11. Rear Cowling Screws: M5 (4)
- 12. On/Off Switch/Power Connection (IEC-320) or power cord with free leads for DC version
- 13. Hour Meter

## **Unpacking and Inspection**

- 1. Orient the shipping container with *This End Up* on top.
- 2. Open the box and carefully lift the IDP-3 and foam blocks out of the box. Remove the foam blocks.
- 3. Save the carton and all packing materials.
- 4. Inspect the pump for damage. If there is shipping damage, contact the freight carrier and your local Agilent sales office immediately.

## Installation

## Safety

Do not remove or modify any safety or insulating equipment from the pump. To do so may create a serious safety hazard and may void the warranty.

## **WARNING**



- ☐ This pump is capable of pumping and exhausting air and inert gases only. It is not designed to pump explosive, flammable, corrosive or particulate forming gases. They can cause bodily injury, explosion, or fire.
- ☐ Install in an area that is not exposed to rain, steam, or excessive humidity. They can cause electric shock, short circuits, and severe bodily injury.
- ☐ Before inspecting or servicing the pump, be sure the electrical supply is disconnected.
- ☐ The gas ballast must be sealed whenever pumping any gas not intended to be vented to the atmosphere.
- ☐ An exhaust silencer option is available for those applications in which it is desirable to diminish the sound level produced by the pump.
- ☐ To prevent bodily injury, avoid exposing any part of the human body to vacuum.

#### CAUTION



Although the pump can pump trace particulates normally found in the atmosphere, it is not designed to process solids, chemicals, powders, solvents, condensates, or other particulates. They can damage the equipment, degrade its performance, or shorten its useful life.

There are four versions of the IDP-3:

- □ 100 VAC, 50/60 Hz
- □ 115 VAC, 60 Hz
- □ 220-230 VAC, 50/60 Hz
- □ 24 VDC

## **CAUTION**

Ensure that the pump version conforms to the supply voltage before using.



## Startup

1. Check that the inlet screen is installed before beginning operation.

## **WARNING**



Do not insert a finger or any foreign object in the path of the fan; serious personal injury may result or the pump may be damaged.

2. Operate the pump at an ambient temperature of 5 °C to 40 °C (41 °F to 104 °F), otherwise damage to the pump or shortened operating life may result.

## **CAUTION**



Do not block the fan ducts. Blocking these ducts can cause pump overheating. A pump surface temperature in excess of 65 °C (150 °F) is potentially damaging. If such conditions are observed, turn the pump off and allow it to cool. Disassemble, inspect for damage, and repair, if necessary.

## **Electrical Connections**

## **Power Cord**

Several power cord options are available from your Agilent dealer. Descriptions of the available power cords and their ordering numbers are given in Table 2.

## **NOTE**



For high voltage operation, the pump must be connected to the power supply using a high voltage IEC-320 type power cord of at least 10 A capacity.

## **CAUTION**



Be certain that your electrical mains power voltage corresponds to that indicated on the rear of the pump.

Table 2 IDP-3 Power Cord Selection

Country	Power Cord Specification	Order
Europe	10 A / 220-230 VAC, 2.5 m	656494220
Denmark	10 A / 220-230 VAC, 2.5 m	656494225
Switzerland	10 A / 230 VAC, 2.5 m	656494235
UK/Ireland	13 A / 230 VAC, 2.5 m	656494250

Table 2 IDP-3 Power Cord Selection (Continued)

Country	Power Cord Specification	Order
India	10 A / 220-250 VAC, 2.5 m	656494245
Israel	10 A / 230 VAC, 2.5 m	656494230
Japan	12 A / 100 VAC, 2.3 m	656494240
North	15 A / 125 VAC, 2.0 m	656458203
America	10 A / 230 VAC, 2.5 m	656494255

## Power Cord (24 VDC)

The interface for the 24VDC power entry is a 9-pin D-sub connector (Figure 4). On this connector pins 4, 5 and 9 are all connected 24V+. Pins 2, 7 and 8 are all connected to 24V-. It is recommended that the Power Cord (Agilent Part Number X3803-60033) included with the pump is used when connecting to the power supply. To correctly connect the power cord to a power supply, attach the 16Awg red wire to 24V+ and attach the Black wire to 24V-.

Pin 1	Control voltage +
Pin 2	24V -
Pin 3	N/A
Pin 4	24V +
Pin 5	24V +
Pin 6	Control voltage -
Pin 7	24V -
Pin 8	24V -
Pin 9	24V +

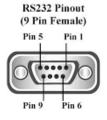


Figure 4 PIN Information

The IDP3 24 VDC model also offers variable speed control. Apply 0-10 VDC through Pin1 and Pin 6 of the 9-pin D-sub connector, or white (+) and blue (-) 20AWG wires on the Agilent power cord supplied. The motor runs at full speed if there is no control signal supplied.

If a power supply cable other than Agilent Part Number X3803-60033 is used to supply power to the IDP3 24VDC Dry Scroll Pump, the user becomes responsible for ensuring that the pump and cable combination complies with any EMC requirements that might exist in that application. All tests performed to ensure compliance to the standards listed in the user manual were done using Agilent Part Number X3803-60033 power cord. If a customer supplied power cable is used, Agilent recommends using a shielded cable with length no longer than 1.5 meters.

## **NOTE**



If a customer supplied cable is being used, ensure that the construction of the cable has a wire/conductor connecting each of the 24V+ and 24V- pins to the power supply. This is required to ensure proper handling of the in-rush and running currents.

#### **NOTE**



There must be good contact between the metallic case and the external shield of the wire. Moreover, this contact must be made at least on the controller side. In this way, it is possible to minimize the influence of any external noise that may impact the ability to satisfy the EMC compliance requirements. Figure 5 shows a correctly connected cable assembly.

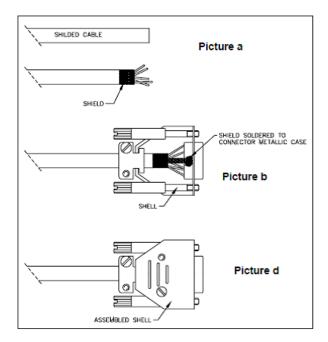


Figure 5 Shield Cable

The minimum rotation speed of the motor is 2200 rpm  $\pm 10\%$ \* when the control is between 0-2.6 V. The motor reaches the full speed when the signal is 8.2 to 10 V. Speed increases proportionally between 2.6 and 8.2 V (Figure 6).

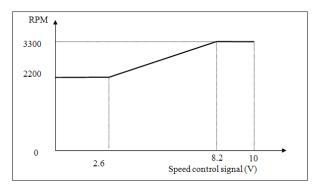


Figure 6 RPM. vs. 0-10 V Speed Control Signal

\*RPM varies with power supply's voltage. Generally, the higher the power voltage is, the higher rotating speed.

Base pressure decreases and power consumption increases as the RPM increases.

## **Grounding Instructions**

This product should be grounded. In the event of an electrical short circuit, grounding reduces the risk of electric shock by providing an escape wire for the electric current. This pump is equipped with a power cord that has a grounding wire with an appropriate grounding plug. The plug must be inserted into an outlet that is properly installed and grounded in accordance with all local codes and ordinances.

**DANGER** 



Improper installation of the grounding plug can result in a risk of electrical shock.

For United States and Canadian installations:

- □ When this product is configured for use on a nominal 120 V circuit, it must be used with a grounding plug that looks like the plug illustrated in Figure 7.
- □ If repair or replacement of the cord or plug is necessary, connect the grounding wire to the grounding terminal only.
- ☐ The grounding wire is insulated and its outer surface is green. It may or may not have yellow stripes.
- □ When this product is configured for use on a nominal 220 V circuit, it must be used with a factory supplied cord and plug that permits connection to the proper electric circuit. See "Electrical Connections" on page 7 for proper rating and type of cord set.

WARNING



Check with a qualified electrician or serviceman if the grounding instructions are not completely understood, or if you are in doubt as to whether the product is properly grounded.

Do not modify the plug provided; if it does not fit the outlet, have the proper outlet installed by a qualified electrician.

Connect the product only to an outlet that has the same configuration as the plug.

Do not use an adapter with this product.

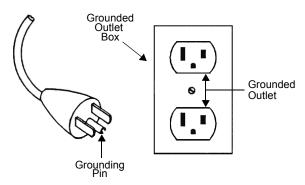


Figure 7 Grounding Plug and Outlet

**WARNING** 



If the product must be reconnected for use on a different type of electric circuit, the connector should be replaced by qualified service personnel.

## **Grounding Instructions (DC version)**

**NOTE** 



Ensure the 24 VDC power supply and 0-10 V power supply are grounded according to the correct practice.

## **Extension Cords**

If you must use an extension cord with this product:

- □ For this product, Agilent recommends using only extension cords with a minimum of 16-gauge wire and a maximum length of 25 feet (7.6 m).
- ☐ Use only a 3-wire extension cord that will accept the plug.
- ☐ Make sure your extension cord is in good condition.
- ☐ Be sure the extension cord is rated high enough to carry the current your product will draw. An undersized cord will cause a drop in line voltage resulting in loss of power and overheating.

## **Run Currents**

Typical run currents listed in Table 3 are approximately constant from minimum to maximum intake pressure.

Table 3 IDP-3 Typical Run Currents (A)

Frequency	100 V Nominal		11	5 V Nomi	nal	2	220 - 230	V Nomina	I	
	90 V	100 V	110 V	103.5 V	115 V	126.5 V	198 V	220 V	230 V	253 V
60 Hz	2.3	2.2	2.2	2.1	2.0	2.0	1.1	1.1	1.1	1.1
50 Hz	1.7	1.7	1.9	N/A	N/A	N/A	0.8	0.8	0.8	1.0
DC	24	V Nomin	ial							
Version	21.6 V	24 V	26.4 V							
	6.8	7.0	8.2							

## **Start Current**

Table 4 lists the typical IDP-3 maximum starting current and duration.

Table 4 Typical IDP-3 Maximum Starting Current and Duration

	50 Hz		60 Hz		DC		
Mains Voltage	Current (A)	Time (ms)	Current (A)	Time (ms)	Current (A)	Time (ms)	
100 V	2.8	300	2.8	500	N/A	N/A	
115 V	N/A	N/A	3.3	500	N/A	N/A	
220 V	1.3	300	1.3	500	N/A	N/A	
230 V	1.4	300	1.4	300	N/A	N/A	
24 V	N/A	N/A	N/A	N/A	15	300 ms	

## **Mechanical Connections**

## Pump Inlet

Use NW16, or larger, clean vacuum hardware with as short a length as practical between the pump and vacuum chamber.

Insert a bellows between the pump and vacuum chamber to provide both vibration isolation and strain relief.

## **Pump Exhaust**

A female 1/4" National Pipe Thread exhaust fitting is located on the front of the pump. Additionally, a 10 mm (3/8") hose barb fitting is provided. To avoid overheating the pump, do not restrict the exhaust flow with long lengths of small diameter tubing. Use as short as practical lengths of 10 mm ID, or larger, diameter hardware.

#### **Gas Ballast**

The pump incorporates an automatic gas ballast to prevent water and other condensates from accumulating within the pump. The standard configuration has a seal screw installed in the 1/4-20 thread in the top of the exhaust manifold (item 6 on Figure 3 on page 6). This configuration can be used for relatively dry applications. When appreciable amounts of water, solvents, etc., are being pumped, remove the seal screw and replace it with the provided adapter and 1/8" NPT sintered filter plug. Do not exceed 25 in-lb of torque when installing the adapter into the manifold.

For applications where the ingress of atmospheric air is undesirable, dry nitrogen at a flow rate of approximately 5 l/min can be provided to the gas ballast. Remove the 1/8" NPT sintered filter plug from the adapter and replace it with an appropriate fitting (supplied by end user).

#### **WARNING**



The gas ballast must be sealed whenever pumping any gas not intended to be vented to atmosphere.

## **Operation**

## **Cleaning the Pump**

Unlike conventional oil-sealed pumps, Agilent dry scroll pumps do not contain fluid for the cleansing of accumulated dust and debris. Run the pump periodically at atmosphere for a minute or two to flush it out. Until experience is gained on your specific process, flush the pump regularly and adjust this schedule according to your specific conditions.

## **Startup Procedure**

1. Make sure that the pump is configured for the mains voltage to which the pump is connected.

## NOTE



The pump ON/OFF switch is a rocker type switch that has symbols in accordance with IEC Publication 417 to represent the ON and OFF positions. Figure 8 shows a switch in the ON position.

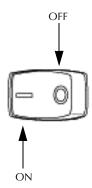


Figure 8 ON/OFF Power Switch

2. Switch the pump ON using the pump ON/OFF switch for AC versions. The 24 VDC pump switches on once the 24 VDC power is connected to the pump and is switched on.

## Shutdown Procedure

To shutdown the pump:

□ Switch the pump OFF for AC versions. For a DC version, switch off the 24 VDC power or disconnect the power cord.

## **IDP-3 Inlet Valve Kit**

The IDP-3 scroll pump is provided with an exhaust valve to prevent rapid leak-up to atmosphere when turned off. However, any vacuum pump generates a small increase in inlet line pressure when turned off. A small amount of pump debris can be carried back into a vacuum line or chamber when the pump is turned off. For applications where the process is sensitive to pump debris, to prevent this, install a fast acting, automatic, normally closed valve. For your convenience, Agilent offers an optional inlet valve and operator kit for the IDP-3 (Figure 9).



Figure 9 IDP-3 with Inlet Valve

This kit includes upper and lower brackets, delay on made timer, solenoid valve with wire harness to the pump and a cooling fan, centering O-ring, quick clamp, nuts and screws for installation, a cord clamp, and a ½" flexible screw driver.

- ☐ The opening of this valve occurs a minimum of 10 s after pump startup.
- □ Valve closing occurs  $\leq$  25 ms after pump shut off.

## **Model Selection**

The inlet valve kit can be obtained separately for existing standard IDP-3 pumps. Refer to Table 5 for model selection, and "Installation" on page 14 for detailed procedures.

Table 5 Inlet Valve Kit Selection

Part Number	Model	Input AC
VPI16IDP220	IDP3A01	220-230 V 50/60 Hz
VPI16IDP115	IDP3B01	115 V 60 Hz
VPI16IDP100	IDP3C01	100 V 50/60 Hz
VPI16IDP24DC	IDP3D01	24 VDC

It is strongly recommended to order the IDP-3 pump with the inlet valve factory-installed to prevent malfunctions due to incorrect installation (see Table 6).

Table 6 IDP-3 Model Numbers with Inlet Valve

Part Number	Input AC
IDP3A21	220-230V 50/60Hz
IDP3B21	115V 60Hz
IDP3C21	100V 50/60Hz
IDP3D21	24 VDC

The technical specification is shown in Table 7.

**Table 7 Technical Specifications** 

Leak Rate:	
□ Body	< 1X10 <sup>-9</sup> atm cc/sec He
□ Seal	<1X10 <sup>-9</sup> atm cc/sec He
Closing time (after solenoid valve power loss)	25 ms
Closing burst	0.01 torr-liter
Power	7.7 W (AC), 9.5 W (DC)
Materials:	
□ Body	Aluminum (6061-T6)
□ Seal	Viton
Temperature Range:	
☐ Valve Body	0~100 °C
□ Solenoid	0~50 °C

## Installation

## AC

Mount the valve body on the pump inlet with centering ring and quick clamp provided in the pump package of IDP3A21 B21 or C21. Observe valve body orientation, otherwise the valve will not function properly (Figure 10).

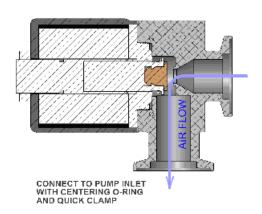




Figure 10 Mount Inlet Valve with IDP-3 Pump

**NOTE** 

If VPI16IDP100, VPI16IDP115, or VPI16IDP220 is purchased separately from the IDP3 AC rev pump, follow the installation procedure in the IDP3 inlet valve kit instruction manual (PN# 699904370).



## DC

24VDC inlet valve kit for IDP series (VPI16IDP24DC) is shown in Figure 11. It is included in the package if a pump with part number IDP3D21 is purchased. To install:



- 1. Plug the connector at the end of solenoid valve to the DC jacket on the rear cowling.
- 2. Follow the same instructions to mount the valve body to the pump inlet as AC rev.



Figure 11 24 VDC Rev Inlet Valve Kit

## **Troubleshooting**

Use the Troubleshooting chart in Table 8 to assist in defining a problem, determining a possible cause, and defining action steps to remedy the situation.

**Table 8 Troubleshooting Chart** 

Problem	Possible Causes	Corrective Actions
Will not start	Motor thermal protector open	Allow motor to cool. Identify cause of overload.
	Excessive voltage drop	Check size and length of cable.
	Defective motor	Inspect. Contact Agilent .
Poor ultimate	System leak	Locate and repair leak.
pressure	Water in pump	Flush pump with air or dry nitrogen.
	Solvent in pump	Flush pump with air or dry nitrogen. Install trap or filter.
	Seals worn out	Replace tip seals.
	Poor conductance to pump	Replumb with shorter and/or larger diameter tubing.
Hammering noise	Pump overheated	Check ventilation to pump. Check ambient temperature.
	Debris in pump	Check intake screen. Flush pump. Disassemble pump and inspect.
Pump runs intermittently	Motor thermal protector is cycling open and closed.*	Allow the motor to cool.  Identify the cause of the overload.

<sup>\*</sup> The IDP-3 is equipped with an auto-reset thermal motor protector. This protector automatically shuts down the pump when it detects an overload condition, and automatically restarts the pump when the motor has cooled to within an acceptable temperature range.

## **Maintenance**

## Kits and Service Options

Agilent pumps will provide many years of trouble-free service if the maintenance procedures and intervals are observed. Cleaning and tip seal replacement are recommended when pump base pressure has risen to an unacceptably high level for your application. If your pump exhibits humming or grinding noises from the bearings, a major overhaul should be done by Agilent personnel or an authorized rebuild center. Advance exchange pumps are available to minimize downtime.

The parts needed for tip seal replacement on the IDP-3 are available in the kit described in Table 9. This kit contains seals and O-rings, and can be obtained from your Agilent dealer.

Table 9 Tip Seal Replacement Kit

Part Number	Description	Contents
IDP3TS	Replacement Tip Seal Set	Replacement Tip Seals and O-rings for IDP-3 pumps

## Cleaning

## **Exterior**

The exterior surfaces of the IDP-3 may be cleaned with alcohol or mild detergents only.

#### Interior

Run the pump periodically at atmosphere for a minute or two to flush it out. For more information, see "Cleaning the Pump" on page 12.

## Tip Seal Replacement

The parts and tools required to replace tip seals are listed below:

- ☐ Tip Seal Replacement Kit P/N IDP3TS
- 4 mm Allen wrench
- □ Razor blade or side-cutting pliers
- ☐ Compressed air (optional)

**WARNING** 



If dangerous gases were being pumped, flush the pump with air or inert gas for at least 10 minutes prior to disassembly. Figure 13 on page 19 shows the various components involved in a tip seal replacement procedure. Refer to Figure 13 as you follow the procedure.

To remove the worn tip seals:

- 1. Disconnect the pump from electrical power.
- 2. Remove (4) M5 socket head bolts (item 1).
- 3. Remove the front cowling and disconnect the electrical connector and set the cowling aside.
- 4. Remove (4) M5 bolts (item 4).
- 5. Remove the outboard housing axially off the frame (item 5).
- 6. Remove and discard the worn tip seals (item 6) and the main O-ring (item 7).
- 7. If compressed air is available, blow any remaining seal debris off the scroll parts. If seal debris is attached to the sides, use a razor or Exacto knife to scrape the debris off.

To install the new seals and O-rings:

- Unpack the tip seals.
   Two tip seals are provided in the kit; one fits the orbiting scroll and the other fits the outboard housing scroll.
- 2. Install the correct tip seal into the groove of the orbiting scroll (item 8) by:
  - a. Facing the cutting burrs on the edges of the seal upwards, inserting the tightly curled end into the center of the orbiting scroll, until the seal is well seated in the groove and the anchor inside the groove protrudes into the seal's side (Figure 12). The seal may need to be pushed firmly to sit well into the groove.

The seal is correctly installed when it sits in its entirety into the groove, sticking out just slightly over the top.



Figure 12 Tip Seal Seating

If the seal's curl seems to have lost some of its curl radius and does not fit into the center of the scroll, re-form the seal end using a pair of round nose pliers to make it fit into the groove. Tip-Seal kit instructions indicate which face should be up (valid for Service Kits only).

- b. Cutting the tip seal off about 1/8" (3 mm) from the outer end of the groove.
- 3. Install the correct tip seal into the groove of the outboard housing scroll (item 4) as in step 3.
- 4. Place the new main O-ring into the groove in the frame (item 10). Ensure that the groove is clean.
- 5. Ensure the sealing face of the outboard housing is clean. Carefully replace the outboard housing by lining up the locating pins. Ensure the tip seal has not fallen out of its groove.
- 6. Reinstall (4) M5 bolts (item 4) and torque the (4) M5 bolts to 5.6 N-m (50 in-lb).
- 7. Reconnect the electrical connector at the front cowling.
- 8. Place the front cowling in place and replace the M5 bolts.
- 9. Reconnect the pump to the electrical power mains.

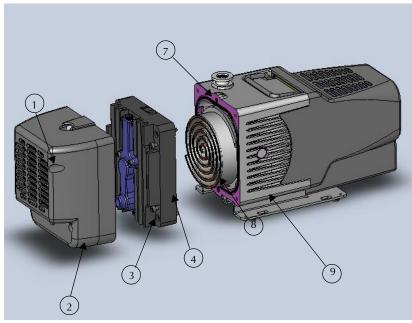


Figure 13 Exploded View of Pump Body

- 1. Front Cowling Bolts; M5 (4)
- 3. Frame Bolts; M5 (4)
- 5. Tip Seals (Not shown)
- 7. Locating Pins (2)
- 9. Frame

## To test the pump:

1. Run the pump for about 5 seconds. Verify that the front fan is running.

If you hear loud noises or observe labored operation, this indicates that the tip seal or main O-ring are possibly out of place.

2. Disassemble and repair as necessary.

The pump is now ready to return to service.

#### NOTE



Newly installed tip seals may require several hours of run time to seat properly and enable the pump to meet speed and base pressure specifications.

- 2. Front Cowling
- 4. Outboard Housing
- 6. Main O-ring Parker No. 2-160 (Not shown)
- 8. Orbiting Scroll

## **Pump Module Replacement**

The parts and tools required to replace the pump module include:

☐ Replacement Module Kit P/N IDP3

☐ 4 mm Allen wrench

Figure 14 on page 20 shows the various components involved in pump module replacement.

To replace the pump module:

- 1. Disconnect the pump from electrical power and disconnect the power cord from the power cord socket.
- 2. Stand the pump on the fan cowling (cover the work surface to protect the pump finish).
- 3. Loosen the four M5 captive screws (item 1) that hold the motor cowling in place (item 2), but do not remove the motor cowling.
- 4. Stand the pump on the motor cowling (cover the work surface to protect the pump finish).
- 5. Remove the four M5 socket head screws (item 3) that secure the fan cowling (item 4).

- 6. Remove the fan cowling, disconnect the electrical connector, and set the cowling aside.
- 7. Remove the four M5 socket head screws (item 5) that secure the base plate (item 6); remove the base plate, and set aside.
- 8. Lay the pump on its side and slide the motor cowling off the pump. Disconnect the electrical connector and set the cowling aside.
- 9. Remove the four M6 button head screws (item 7) that secure the motor to the pump module and set the motor aside.
- 10. Install the new coupling spider (item 8) to the coupling (motor side or pump module side) and install the motor to the replacement pump module with the label facing up (in the same direction as the inlet flange of the pump module). Secure the motor with four M6 button head screws.
- 11. Reconnect the electrical connector on the motor cowling to the motor.

- 12. Tuck the electrical connector behind the aluminum plate mounted to the motor cowling (item 9) and slide the motor cowling into place.
- 13. Holding the motor cowling in place, stand the pump on the end of the motor cowling with the exhaust fitting pointing up.
- 14. Reinstall the base plate with four M5 socket head screws.
- 15. Connect the electrical connector on the fan cowling to the wiring harness and reinstall the fan cowling with four M5 socket head screws.
- 16. Holding the motor cowling in place, stand the pump on the fan cowling.
- 17. Tighten the four M5 captive screws to secure the motor cowling.
- 18. Reattach the power cord and confirm proper operation.

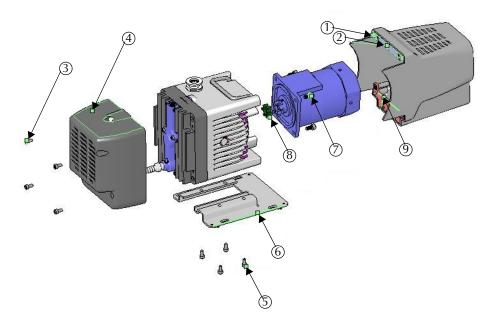


Figure 14 Module Replacement - Exploded View



# 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, eq).

- 3) Important steps for the shipment of returning product:
  - Remove all accessories from the core product (e.g. inlet screens, vent valves).
  - Prior to shipment, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
  - If ordering an Advance Exchange product, <u>please use the packaging from the Advance Exchange to return the defective</u> product.
  - Seal the product in a plastic bag, and package product carefully to avoid damage in transit. You are responsible for loss or damage in transit.
  - Agilent Technologies is not responsible for returning customer provided packaging or containers.
  - Clearly label package with RA number. Using the shipping label provided will ensure the proper address and RA number
    are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will
    be returned.
- 4) Return only products for which the RA was issued.
- 5) Product being returned under a RA must be received within 15 business days.
- 6) Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information. Customer is responsible for freight charges on returning product.
- 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:

 Fax:
 00 39 011 9979 330

 Fax Free:
 00 800 345 345 00
 Fax:
 1 781 860 9252
 please visit our website for individual office information

 Toll Free:
 00 800 234 234 00
 Toll Free: 800 882 7426, Option 3
 office information

 vpt-customercare@agilent.com
 vpl-ra@agilent.com
 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.

) CUSTOMER INFORMATION			
Company Name:		Contact Name:	
Tel:	Email:	Fax:	
Customer Ship To:		Customer Bill To:	
Europe only: VAT reg. Numb	er:	USA/Canada only: Tax	kable Non-taxable
) PRODUCT IDENTIFICATION			
Product Description	Agilent P/N	Agilent S/N	Original Purchasing Reference
<u> </u>			
_			
TYPE OF RETURN (Choose on	e from each row and supply	/ Purchase Order if requesting a billa	ble service)
		f (hard copy must be submitted with	
			Evaluation Return for Credit
<b>3B</b> . Exchange Repair	UpgradeConsign	iment/ DemoCalibration!	Evaluationneturn for Credit
) HEALTH and SAFETY CERTIFIC	CATION		
AGILENT TECHNOLOGIES CAN	NOT ACCEPT ANY PRODUC	CTS CONTAMINATED WITH BIOLOG	ICAL OR EXPLOSIVE HAZARDS,
RADIOACTIVE MATERIAL, OR			
Call Agilent Technologies to d	iscuss alternatives if this re	equirement presents a problem.	
The equipment listed above (c			
		toxic or hazardous materials. OR wing toxic or hazardous materials. I	f this hav is chacked the following
	-	boxes for all materials to which pro	<u> </u>
Toxic Corrosive		lammable Explosive	Biological Radioactive
		_ ,	•
LIST all toxic/ nazardous matei	als. Include product name	, chemical name, and chemical syn	iboi or tormuia:
			ed, <b>the customer will be held responsible</b> for all
costs incurred to ensure the safe hand exposure to toxic or hazardous materia	= -	r any harm or injury to Agilent employees as	well as to any third party occurring as a result of
Print Name:	•	nature:	Date:
) FAILURE INFORMATION:			
Failure Mode (REQUIRED FIELD	See next name for summes	tions of failure terms):	
Detailed Description of Malfun		·	
·	, ,	ioi ilicosaye)	
Application (system and model	):		
I understand and agree to the t	terms of Section 6 Dags 2/	3	
Print Name:		o. nature:	Date:



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

## Please use these Failure Mode to describe the concern about the product on Page 2.

#### TURBO PUMPS and TURBO CONTROLLERS

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

## ION PUMPS/CONTROLLERS

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

## LEAK DETECTORS

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

#### **SCROLL AND ROTARY VANE PUMPS**

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

## VALVES/COMPONENTS

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

#### **INSTRUMENTS**

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

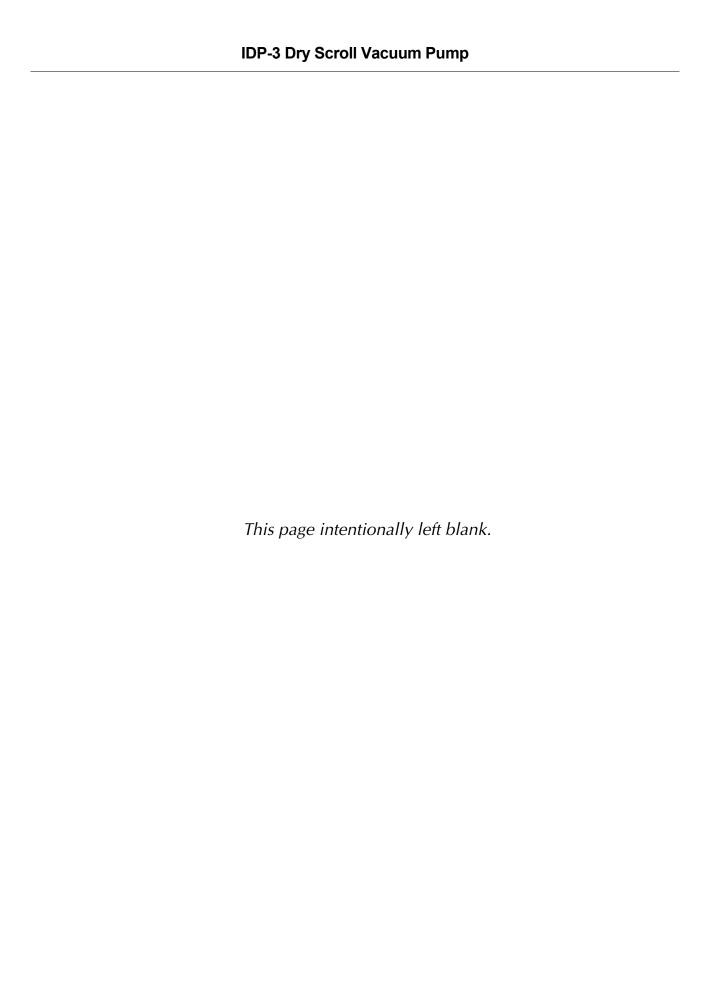
#### **DIFFUSION PUMPS**

- 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.

- Customer is responsible for the freight charges for the returning product. Return shipments must comply with all
  applicable Shipping Regulations (IATA, DOT, etc.) and carrier requirements.
- Customers receiving an Advance Exchange product agree to return the defective, rebuildable part to Agilent Technologies
  within 15 business days. Failure to do so, or returning a non-rebuildable part (crashed), will result in an invoice for the
  non-returned/non-rebuildable part.
- Returns for credit toward the purchase of new or refurbished Products are subject to prior Agilent approval and may incur
  a restocking fee. Please reference the original purchase order number.
- Units returned for evaluation will be evaluated, and a quote for repair will be issued. If you choose to have the unit
  repaired, the cost of the evaluation will be deducted from the final repair pricing. A Purchase Order for the final repair price
  should be issued within 3 weeks of quotation date. Units without a Purchase Order for repair will be returned to the
  customer, and the evaluation fee will be invoiced.
- A Special Cleaning fee will apply to all exposed products per Section 4 of this document.
- If requesting a calibration service, units must be functionally capable of being calibrated.



## **Service & Support**

#### **North America**

Agilent Technologies Vacuum Products Division 121 Hartwell Avenue Lexington, MA 02421 USA Tel.: +1 781 861 7200

Toll-Free: +1 800 882 7426 Fax: +1 781 860 5437

#### **Benelux**

Agilent Technologies Netherlands B.V. Vacuum Products Division Herculesweg 8 4338 PL Middelburg The Netherlands Tel: +31 118 671570 Fax: +31 118 671569

Toll free: 00 800 234 234 00

#### Canada

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

#### China

Agilent Technologies (China) Co. Ltd Vacuum Products Division No.3, Wang Jing Bei Lu, Chao Yang District

Beijing, 100102 China Tel.: +86 (0)10 64397888 Fax: +86 (0)10 64391318 Toll-Free: 800 820 3278

#### France

Agilent Technologies France Vacuum Products Division 3 avenue du Canada, Parc Technopolis Z.A. Courtaboeuf - CS 90263 91978 Les Ulis cedex France

Tel.: +33 (0)1 64 53 61 15 Fax: +33 (0)1 64 53 50 01 Toll free: 00 800 234 234 00 e-Mail: vpf.sales@agilent.com

## Germany & Austria

Agilent Technologies Sales & Services GmbH & Co. KG Vacuum Products Division Lyoner Str. 20 60 528 Frankfurt am Main Germany

Tel.: +49 69 6773 43 2230 Fax: +49 69 6773 43 2250

This information is subject to change without notice.

© Agilent Technologies, Inc., 2017

#### India

Agilent Technologies India Pvt. Ltd. Vacuum Products Division G01. Prime corporate Park, 230/231, Sahar Road, Opp. Blue Dart Centre, Andheri (East), Mumbai – 400 099. India

Tel: +91 22 30648287/8200 Fax: +91 22 30648250 Toll Free: 1800 113037

#### Italy

Agilent Technologies Italia S.p.A. Vacuum Products Division Via F.Ili Varian 54 10040 Leini, (Torino) ITALY Tel.: +39 011 997 9111 Fax: +39 011 997 9350 Toll-Free: 00 800 234 234 00

#### Japan

Agilent Technologies Japan, Ltd.
Vacuum Products Division
8th Floor, Sumitomo Shibaura Building
4-16-36 Shibaura Minato-ku
Tokyo 108-0023 JAPAN
Tel.: +81 3 5232 1253
Toll-Free: 0120 655 040

#### Korea

Agilent Technologies Vacuum Products Division Shinsa 2nd Bldg. 2F 966-5 Daechi-dong Kangnam-gu, Seoul KOREA 135-280 Tel.: +82 (0)2 3452 2455

Toll-Free: 080 222 2452 Fax: +82 (0)2 3452 3947

Fax: +81 3 5232 1710

## Singapore

Agilent Technologies Singapore Pte. Ltd Vacuum Products Division Agilent Technologies Building, No.1 Yishun Avenue 7 Singapore 768923

Tel: +65 6215 8045 Fax: +65 6754 0574

#### Southeast Asia

Agilent Technologies Sales Sdn Bhd Vacuum Products Division Unit 201, Level 2 uptown 2, 2 Jalan SS21/37, Damansara Uptown 47400 Petaling Jaya, Selangor, Malaysia

Tel: +603 7712 6106 Fax: +603 6733 8121

#### Taiwan

Agilent Technologies Taiwan Limited Vacuum Products Division (3F) 20 Kao-Shuang Rd., Pin-Chen City, 32450 Taoyuan Hsien, Taiwan, R.O.C. Tel. +886 34959281 Toll Free: 0800 051 342

#### **UK & Ireland**

Agilent Technologies Vacuum Products Division 6 Mead Road Oxford Industrial Park Yarnton, Oxford OX5 1QU UK Tel.: +44 (0) 1865 291570 Fax: +44 (0) 1865 291571

Toll free: 00 800 234 234 00

#### **Other Countries**

Agilent Technologies Italia S.p.A. Vacuum Products Division Via F.Ili Varian 54 10040 Leini, (Torino) ITALY Tel.: +39 011 997 9111

Fax: +39 011 997 9111 Fax: +39 011 997 9350 Toll-Free: 00 800 234 234 00

## **Customer Service and Support**

North America

Toll Free: 800 882 7426, Option 3

vpl-ra@agilent.com

#### Europe

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

#### Pacific Rim:

Please visit our website for individual office information http://www.agilent.com

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

www.agilent.com Representatives in most countries



