



Agilent Technologies

Vacuum Products Division

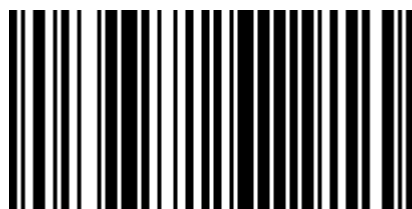
VS PD03 and G8610 Scroll Pump and Components Replacement

*FIELD INSTALLATION
INSTRUCTIONS*

Part Number 699910017

Rev. B

July 2017



699910017

This page intentionally left blank.




Agilent Field Instruction Sheet

VS PD03 and G8610 Scroll Pump and Components Replacement

Preface










Documentation Standards

This manual uses the following documentation standards:

NOTE	Notes contain important information.
	
CAUTION	Cautions appear before instructions, which if not followed, could cause damage to the equipment or data loss.
	
WARNING	Warnings appear for a particular procedure or practice which, if not followed correctly, could lead to serious injury or death.
	

Hazard and Safety Information

The common international symbols used in this manual and on the equipment are defined below.

	OFF Supply (Power)		Earth (Ground) Terminal
	ON Supply (Power)		Caution, Hot Surface
	AC – Alternating Current		Caution, Risk of Electrical Shock
	Warning, Risk of danger		Protective Conductor Terminal
	Frame or chassis Terminal		



Operators and service personnel must be aware of all hazards associated with this equipment. They must know how to recognize hazardous and potentially hazardous conditions, and know how to avoid them. The consequences of unskilled, improper, or careless operation of the equipment can be serious. Every operator or service person must read and thoroughly understand operation/maintenance manuals and any additional information provided by Agilent. All warning and cautions must be read carefully and strictly observed. Consult local, state, and national agencies regarding specific requirements and regulations. Address any safety, operation, and/or maintenance questions to your nearest Agilent office.

Solvents

WARNING



The mechanical components of leak detectors may be cleaned with one of the recommended solvents. When heated, sprayed, or exposed to high-temperature equipment, these solvents become flammable and explosive, causing serious injury or death. Do not use these solvents near a high-temperature source. Ventilate the working area with a blower and work in a large, well-ventilated room.

Solvents are irritants, narcotics, depressants and/or carcinogens. Their inhalation and/or ingestion may produce serious side effects. Prolonged or continued contact with the skin results in absorption through the skin and moderate toxicity. Always ensure that cleaning operations are carried out in large, well-ventilated rooms, and wear eye shields, gloves, and protective clothing.

Due to the effective cleaning nature of VacuSolv solvent and its residue-free properties, Agilent' Component and Spectrometer Cleaning Kit (Part Number 670029096), used in accordance with the kit instructions, is recommended for cleaning spectrometer components. The kit can also be used for fine cleaning of other parts in the leak detector's vacuum system such as valves and fittings. No rinsing steps or high-temperature drying is required following cleaning with VacuSolv. Although appropriate precautions are advised, VacuSolv is compatible with most materials and does not contain toxic chemicals or CFCs (chlorofluorocarbons). Other acceptable solvents are isopropyl alcohol (IPA) or Dow Corning[®] OS-20.

To clean the leak detector plastic enclosure, the LCD display and Front Panel buttons, use only a soft cloth slightly dampened with water or a mild soap.

Do NOT use excess water or cleaning solvents of any kind.

Avoid splashing any cleaning solvents into the unit through the ventilation openings or Front Panel buttons. Wipe the surface with a dry lint-free cloth.

Vacuum Equipment and Cleanliness

Cleanliness is vital when servicing the leak detector or any vacuum equipment. There are some techniques that are more important in leak detector servicing than in general vacuum work:

CAUTION



Wear non-powdered, ESD-safe Nitride or equivalent gloves to prevent skin oils from getting on spectrometer internal components.



O-ring Care

When removing, checking or replacing O-rings, keep in mind the following:

NOTE



Agilent recommends replacing all O-rings during routine maintenance or during any maintenance procedure requiring that O-rings be removed.

CAUTION



Remove O-rings carefully with your fingers. Do not use metal tools for this task; this prevents scratching of any sealing surfaces.

- Wipe all O-rings clean with a lint-free cloth before installation to ensure that no foreign matter is present to impair the seal.
- Do not use grease or any other substance on O-rings that will come in contact with the vacuum surfaces.
- Do not use alcohol, methanol or other solvents on O-rings. Doing so causes deterioration and reduces their ability to hold a vacuum.
- Agilent does not recommend the use of vacuum grease. If applicable, apply a small amount of Apiezon[®] L grease and wipe the O-rings shiny dry.

Metal Seal Care

CAUTION



Metal Seals must be replaced any time a spectrometer is opened. All fasteners must be installed and torqued per assembly procedure specifications. Remove Metal Seals carefully with your fingers or a soft tool. Metal tools scratch sealing surfaces.

- Metal Seals are supplied in pre-cleaned condition. No cleaning is required. If necessary, Metal Seals can be cleaned using the recommended solvents. Wipe Metal Seals clean with a lint-free cloth before installation to ensure that no foreign matter impairs the seal.
- Do not use grease or any other substance on Metal Seals that will come in contact with the spectrometer.

Spectrometer

CAUTION



Store the Ion Source/Preamplifier sub-assembly in a cool, dry area in a tightly sealed, ESD protected container. Wear non-powdered, ESD-safe Nitride or equivalent gloves when handling the spectrometer. Wash hands thoroughly after handling the spectrometer filaments and especially before smoking or eating.

The spectrometer and PCB's are static sensitive devices. Wear a grounding strap when performing any maintenance on these units and especially when performing maintenance of static sensitive parts.

CAUTION



The spectrometer operates at a very high vacuum produced by the high vacuum turbomolecular pump. Service of the spectrometer requires that this vacuum be vented to the atmosphere.

This manual contains instructions for VS (Part One) and G8610 (Part Two)

Part One: VS Model - Scroll Pump Replacement

Equipment Required

- Slotted screw driver
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. If a VS unit (top), using an extended length M5 Allen wrench.

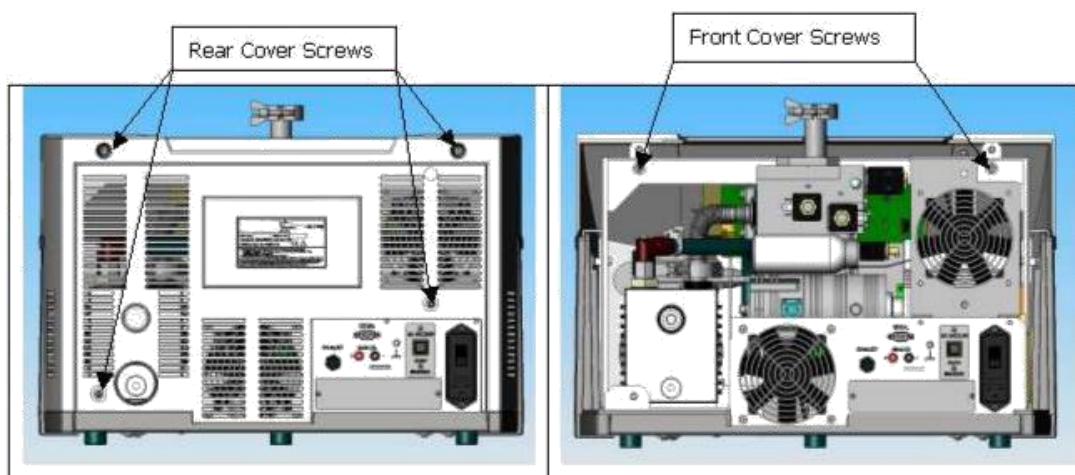


Figure 1: Covers

4. Disconnect the NW 16 connection (Figure 2: NW 16 Connection).



Figure 2: NW 16 Connection

5. Tilt the unit and unscrew the four connections underneath that hold the pump (Figure 3: Scroll Pump Mounting Connections). Place the unit flat again and remove the whole pump assembly.



Figure 3: Scroll Pump Mounting Connections

6. Disconnect the exhaust hose (Figure 4: Exhaust Hose).



Figure 4: Exhaust Hose

- 7. Disconnect the diaphragm, fan and motor power connectors.
- 8. Remove the fan assembly bracket and lay it aside (Figure 5: Fan Assembly Mounts).

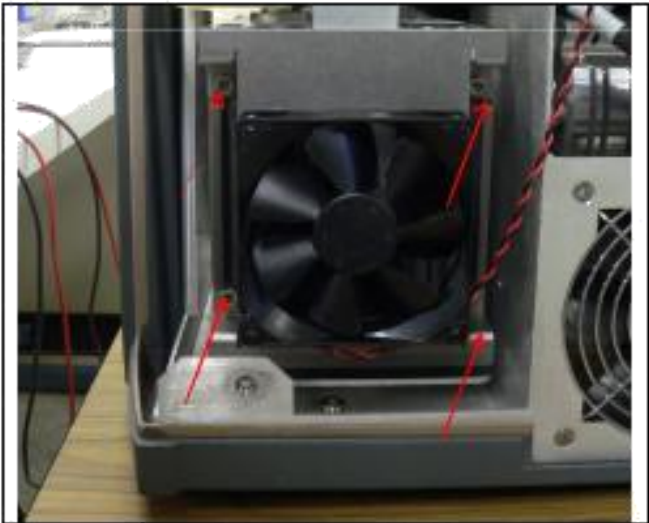


Figure 5: Fan Assembly Mounts

9. Unscrew the four M2 screws and remove the manifold block (Figure 6: Manifold Block Connections).

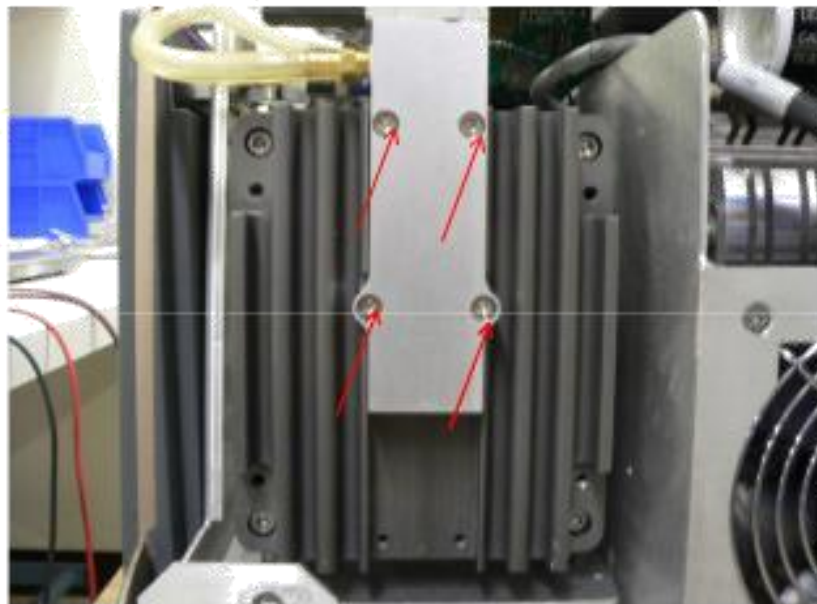


Figure 6: Manifold Block Connections

10. Remove the exhaust valve unit. See “Exhaust Valve Replacement” on page 10.
11. Remove the diaphragm pump. See “Diaphragm Pump Replacement” on page 12.
12. Mount the new scroll pump using the four connections seen in “Figure 3: Scroll Pump Mounting Connections” on page 5.
13. Reattach the diaphragm pump. See “Diaphragm Pump Replacement” on page 12.
14. Reattach the exhaust valve. See “Exhaust Valve Replacement” on page 10.
15. Reattach the four M2 screws to remount the manifold block.
16. Reattach the fan assembly.
17. Reconnect the diaphragm, fan and motor power connectors.
18. Reconnect the exhaust hose.
19. Reattach covers using existing hardware and fasten.
20. Connect the power cord and power up the unit.
21. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
22. Refer to the operator's manual if the system fails to reach the *System Ready* mode. Agilent recommends a full calibration of the unit prior to leak test operations.

Part One: VS Model - Clippard (Vent) Valve Replacement

Equipment Required

- Extended Length M5 Allen Wrench
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. If a VS unit (top), using an extended length M5 Allen wrench (VS)

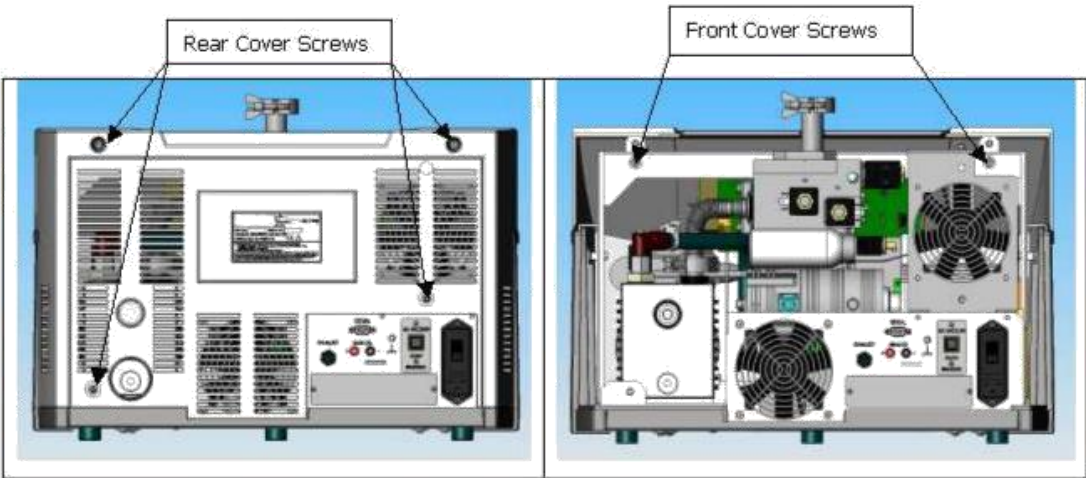


Figure 7: Covers

4. Disconnect the two spade connectors (red arrow in Figure 8: Spade Connectors):

- Red/black
- Black/black

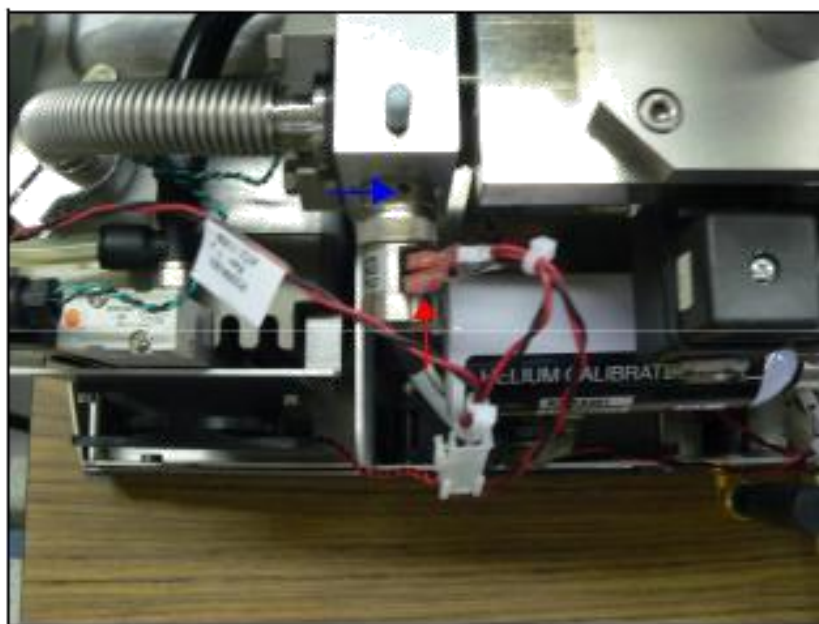


Figure 8: Spade Connectors

4. Use an Allen wrench to loosen valve from housing (blue arrow in Figure 8: Spade Connectors).

5. Mount the new valve and tighten.

6. Reconnect the two spade connections.

7. Attach the all covers using existing hardware.

8. Fasten covers using Allen wrench

9. Connect the power cord and power up the unit.

10. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.

Refer to the operator's manual if the system fails to reach the *System Ready* mode.

Part One: VS Model - Exhaust Valve Replacement

Equipment Required

- Extended Length M5 Allen Wrench
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING *Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.*



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. If a VS unit (top), using an extended length M5 Allen wrench

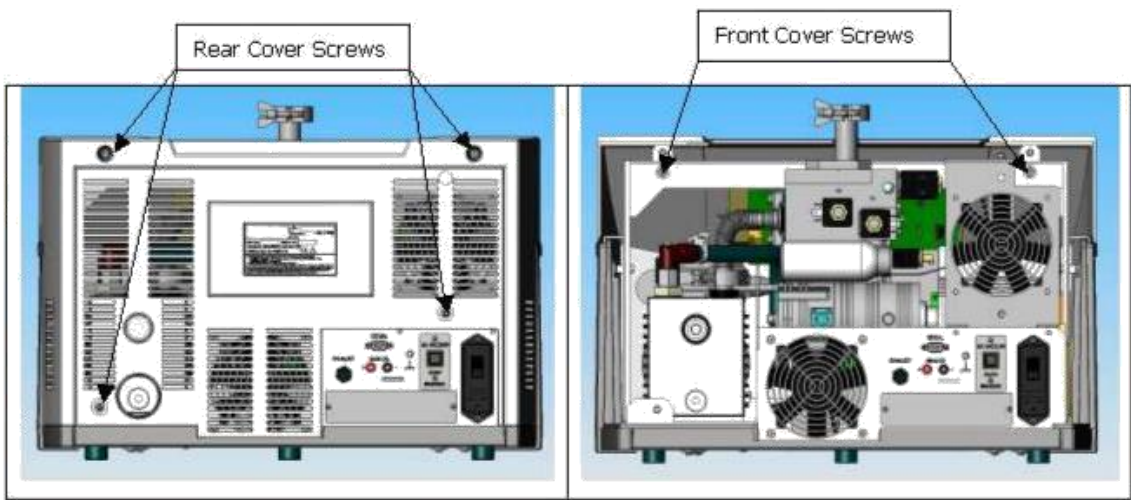


Figure 9: Covers

4. Unscrew the black cover using the Philips head screw, pull up the cover and disconnect (red arrow in Figure 10: Unit Screws)

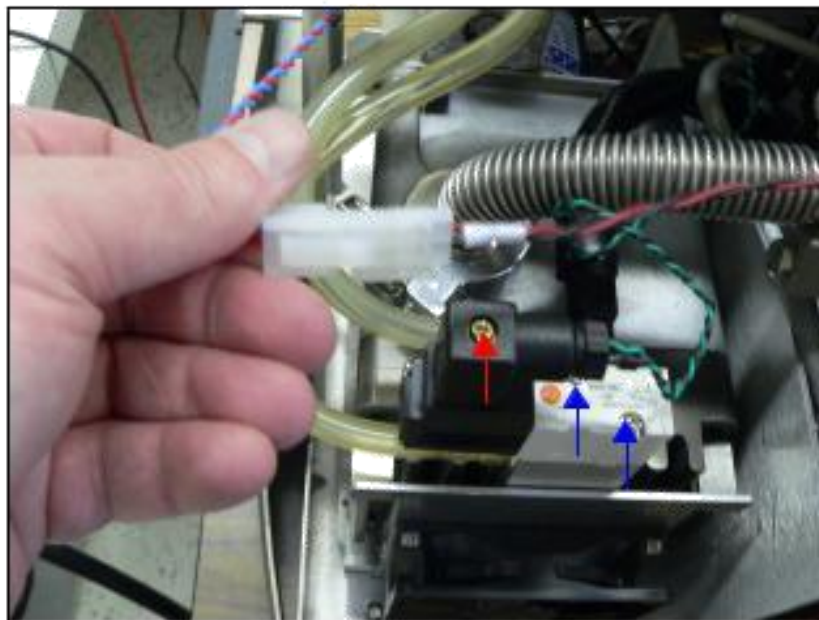


Figure 10: Unit Screws

5. Unscrew the two Philips head screws (blue arrow in “Figure 10: Unit Screws”) and pull the complete black/grey body up and out.
6. Mount new unit and replace the three Philips screws.
7. Attach the all covers using existing hardware.
8. Fasten covers using Allen wrench.
9. Connect the power cord and power up the unit.
10. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
Refer to the operator's manual if the system fails to reach the *System Ready* mode.
11. Agilent recommends a full calibration of the unit prior to leak test operations.

Part One: VS Model - Diaphragm Pump Replacement

Equipment Required

- Extended Length M5 Allen Wrench
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver
- Adjustable Wrench

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. If a VS unit (top), using an extended length M5 Allen wrench (Figure 11: Covers).

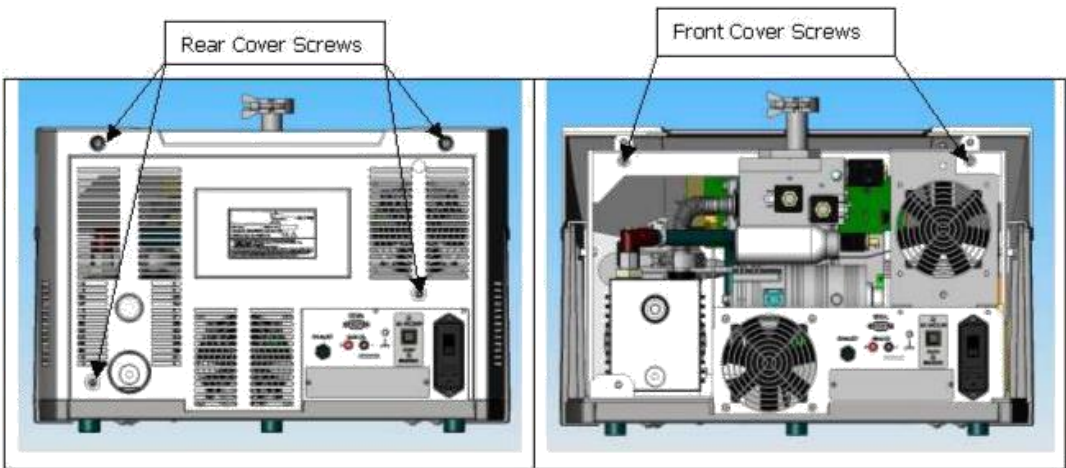


Figure 11: Covers

4. Disconnect the R3359-301 Molex connector to the Clippard valve (Figure 12: Molex Connector).

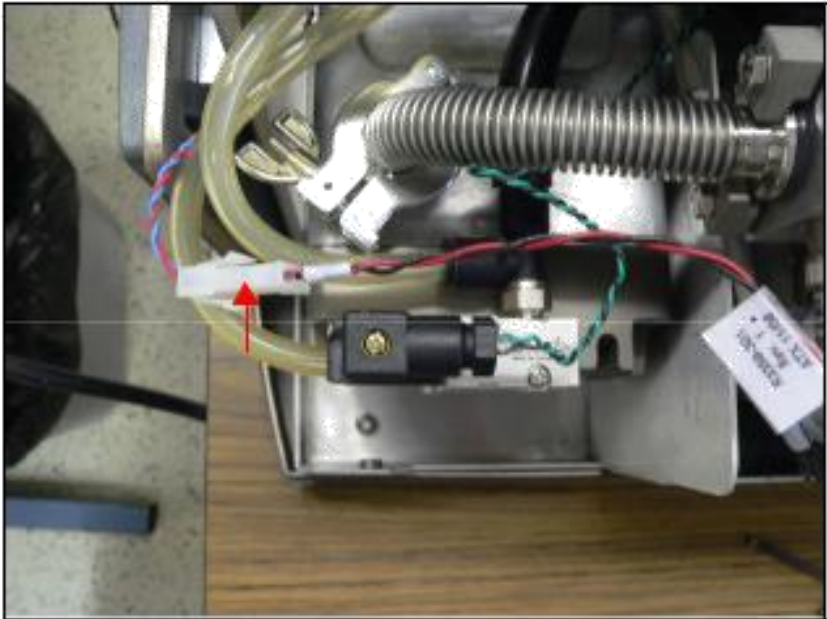


Figure 12: Molex Connector

5. Remove the ground connection to the system body (red arrow in Figure 13: Ground Connection).

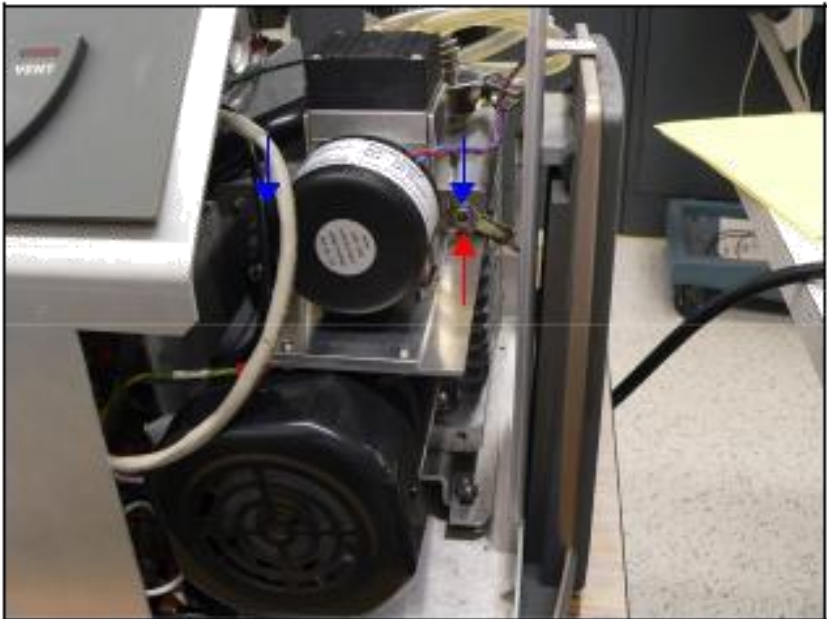


Figure 13: Ground Connection

6. Disconnect the two M4 screws mounting the plate to the scroll pump assembly (blue arrow in Figure 13: Ground Connection).

7. Remove the diaphragm pump assembly, turn it over and unscrew the four M2 screw holding the pump to the plate (Figure 14: Pump to Plate Connections - two shown).

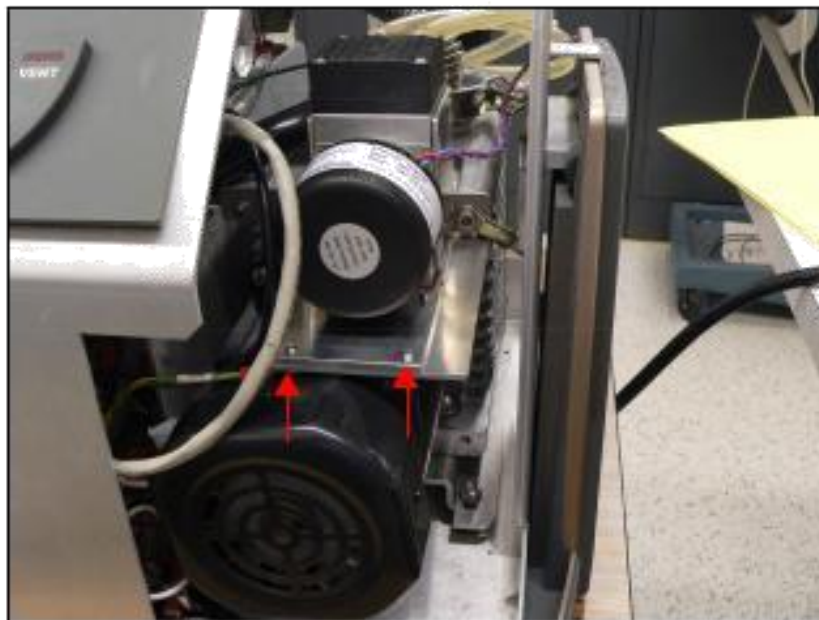


Figure 14: Pump to Plate Connections

8. Reattach the new pump to the plate.
9. Reattach the plate to the scroll pump.
10. Reattach the ground and Molex connections.
11. Attach the front cover and secure to the frame using existing hardware.
12. Attach the rear cover and secure to the unit using existing hardware.
13. Connect the power cord and power up the unit.
14. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
Refer to the operator's manual if the system fails to reach the *System Ready* mode.
15. Agilent recommends a full calibration of the unit prior to leak test operations.

Part One: VS Model - Diaphragm and Seal Replacement

Equipment Required

- Extended Length M5 Allen Wrench
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver
- Phillips Head Screw Driver, #1
- Felt Tip Pen

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. If a VS unit (top), using an extended length M5 Allen wrench (Figure 15: Covers).

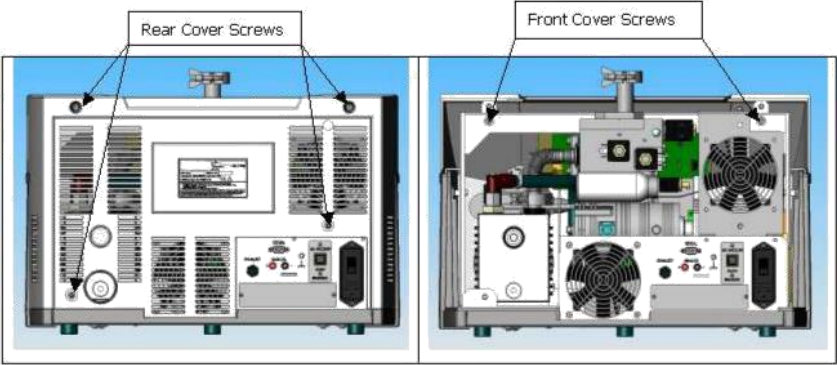


Figure 15: Covers

NOTE



Always change the valve plates, diaphragm and sealing rings at the same time.

4. Remove the pump head by (Figure 16: Pump Head Removal):

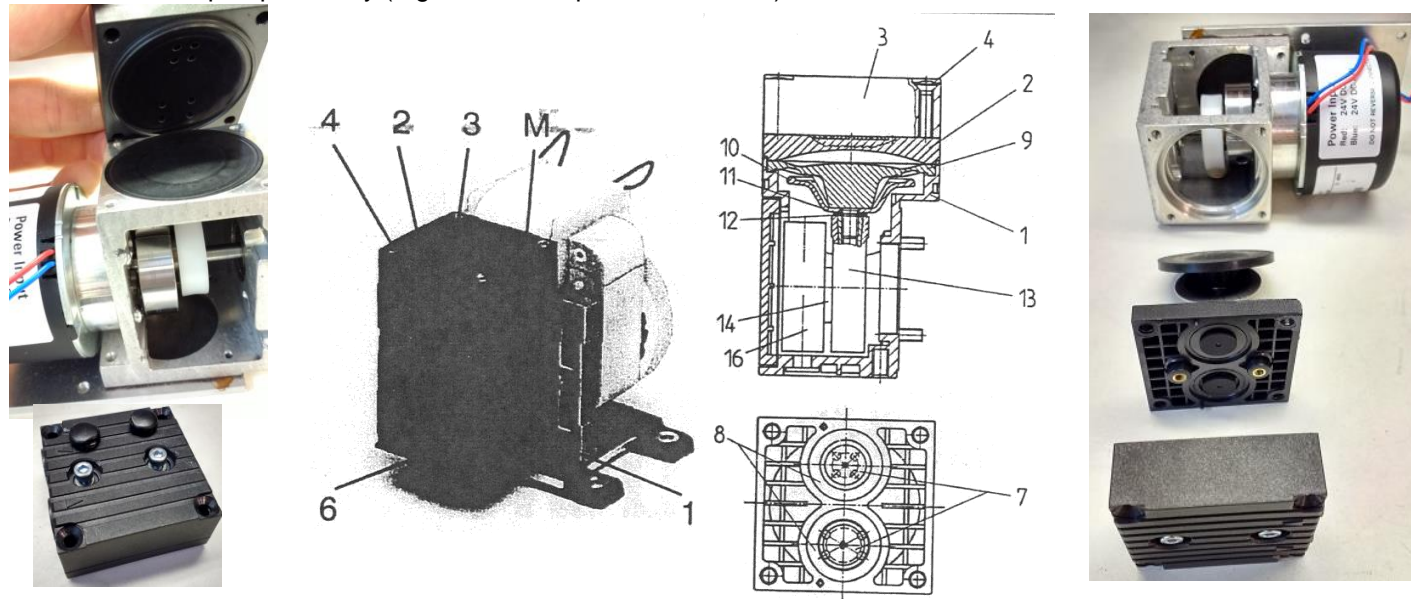


Figure 16: Pump Head Removal

- a. Making a mark (M) on the head plate (3), intermediate plate (2), and housing (1) with a felt-tip pen.
- b. Undoing the 4 screws (4) in the head plate and lifting the head plate, with the intermediate plate, off the pump housing.
- c. *For models with:*
 - *DC motors (no cooling fan):* Removing the cover (6) from the pump housing (1) by inserting a screwdriver into the slots on the sides, and prying the cover off (with care the adhesive gasket between cover and housing may be reused).
 - *No fan:* Reference is made to turning or holding the cooling fan, the necessary operations must be carried out by turning or holding the counterweight (16).

5. Change the diaphragm by (see Figure 16: Pump Head Removal):

- a. Turning the fan to bring the structured diaphragm (9) to top dead center.
- b. Lifting the edge of the diaphragm and, gripping it on opposite sides, unscrewing it by turning counterclockwise.

CAUTION



Ensure that the disc spring (12) and diaphragm spacers (11) on the threaded portion of the diaphragm do not fall into the housing.

- c. Taking the disc spring (12), diaphragm spacer(s) (11), and diaphragm support (10) off the threaded portion of the diaphragm and retaining them.
- d. Ensuring that all parts are free from dirt and clean them, if necessary.
- e. Placing the diaphragm support, diaphragm spacer(s), and disc spring, in that order, on the thread part of the new diaphragm. The concave side of the disc spring must be towards the diaphragm.

- f. Turning the fan until the connecting rod (13) is at top dead center.
 - g. Screwing the diaphragm, complete with diaphragm support, diaphragm spacer(s), and disc spring, into the connecting rod (clockwise) and tightening it by hand.
 - 6. Change the valve plates by (see Figure 16: Pump Head Removal):
 - a. Separating the head plate (3) from the intermediate plate (2).
 - b. Removing the valve plates (7) and sealing rings (8) from the intermediate plate.
 - c. Checking that the valve seats in the head plate and the intermediate plate are clean. If scratches, distortion, or corrosion are evident, replace that part.
 - d. Laying the new valve plates in the intermediate plate recesses. The valve plates for suction and pressure sides are identical, as are upper and lower sides of the plates.
 - e. Ensuring the valve plates are not deformed by moving them gently sideways in their recesses.
 - f. Laying the sealing rings on the intermediate plate.
 - 7. Refit the pump heads by (see Figure 16: Pump Head Removal):
 - g. Turning the fan to bring the diaphragm to top dead center.
 - h. Placing the intermediate plate (2), with valve plates (7), sealing rings (8), and head plate (3) on the housing, in the position indicated by the (M).
 - i. Checking that the head plate is centered by moving it gently sideways.
 - j. Gently tightening the screws (4), evenly and diagonally.
 - k. Turning the fan to check that the pump rotates freely.
 - l. Turning the fan again to bring the diaphragm to top dead center.
 - m. Tightening the screws (4).
 - n. *For DC versions (no cooling fan):* Reattaching the cover (6) to housing (1).
 - 8. Attach the front cover and secure to the frame using existing hardware.
 - 9. Attach the rear cover and secure to the unit using existing hardware.
 - 10. Connect the power cord and power up the unit.
 - 11. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
- Refer to the operator's manual if the system fails to reach the *System Ready* mode.

Part Two: G8610 Model - Scroll Pump Replacement

Equipment Required

- Slotted screw driver
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING



Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.

1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. Using a slotted screw driver, remove all four covers and rotate the top about the screen, which does not need to be removed (Figure 1: Covers).

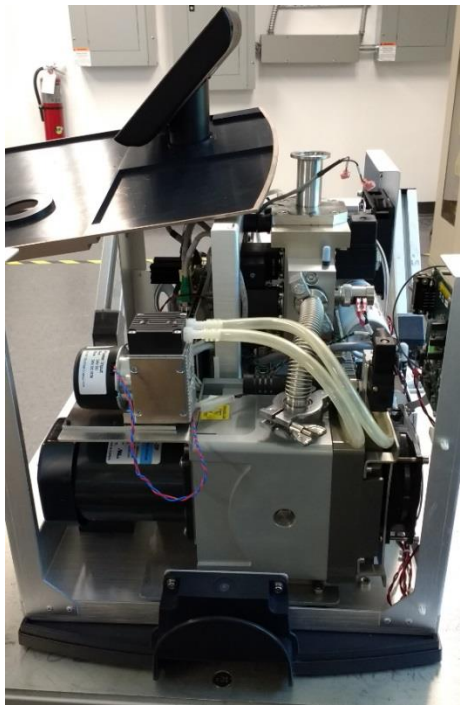


Figure 1: Covers

4. Disconnect the NW 16 connection (Figure 2: NW 16 Connection).



Figure 2: NW 16 Connection

5. Tilt the unit and unscrew the four connections underneath the leak detector that hold the scroll pump (Figure 3: Scroll Pump Mounting Connections). Place the unit flat again and remove the whole pump assembly.

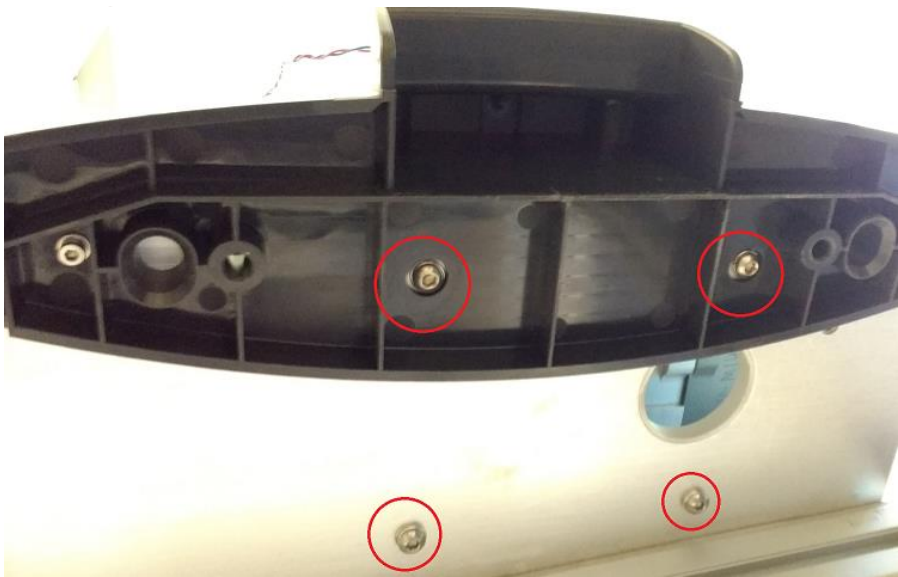


Figure 3: Scroll Pump Mounting Connections

6. Disconnect the exhaust hose (Figure 4: Exhaust Hose).



Figure 4: Exhaust Hose

- 7. Disconnect the diaphragm, fan and motor power connectors.
- 8. Remove the fan assembly bracket and lay it aside (Figure 5: Fan Assembly Mounts).



Figure 5: Fan Assembly Mounts

9. Unscrew the four M2 screws and remove the manifold block (Figure 6: Manifold Block Connections).



Figure 6: Manifold Block Connections

10. Remove the exhaust valve unit. See “Exhaust Valve Replacement” on page 24.
11. Remove the diaphragm pump. See “Diaphragm Pump Replacement” on page 26.
12. Mount the new scroll pump using the four connections seen in “Figure 3: Scroll Pump Mounting Connections” on page 19.
13. Reattach the diaphragm pump. See “Diaphragm Pump Replacement” on page 26.
14. Reattach the exhaust valve. See “Exhaust Valve Replacement” on page 24.
15. Reattach the four M2 screws to remount the manifold block.
16. Reattach the fan assembly.
17. Reconnect the diaphragm, fan and motor power connectors.
18. Reconnect the exhaust hose.
19. Reattach covers using existing hardware and fasten.
20. Connect the power cord and power up the unit.
21. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
22. Refer to the operator's manual if the system fails to reach the *System Ready* mode. Agilent recommends a full calibration of the unit prior to leak test operations.

Part Two: G8610 Clippard (Vent) Valve Replacement

Equipment Required

- Slotted screw driver
- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING *Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.*



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. Using a slotted screw driver, remove the only back cover (Figure 7: Covers).



Figure 7: Covers

4. Disconnect the two spade connectors (red arrow in Figure 8: Spade Connectors):

- Red/black
- Black/black

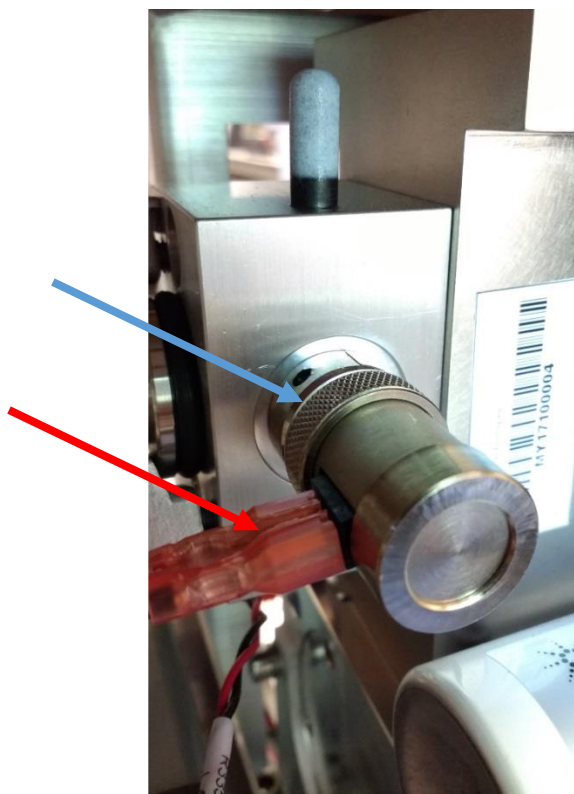


Figure 8: Spade Connectors

5. Use an Allen wrench to loosen valve from housing (blue arrow in Figure 8: Spade Connectors).

6. Mount the new valve and tighten.

7. Reconnect the two spade connections.

8. Attach the all covers using existing hardware.

9. Fasten covers using a slotted screw driver.

10. Connect the power cord and power up the unit.

11. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.

Refer to the operator's manual if the system fails to reach the *System Ready* mode.

Part Two: G8610 Exhaust Valve Replacement

Equipment Required

- Metric Allen Wrench Set
- Slotted Screw Driver
- M3 Phillips Head Screw Driver

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. Using a slotted screw driver, remove all four covers and rotate the lid about the screen, without removing the screen itself (Figure 9: Covers).

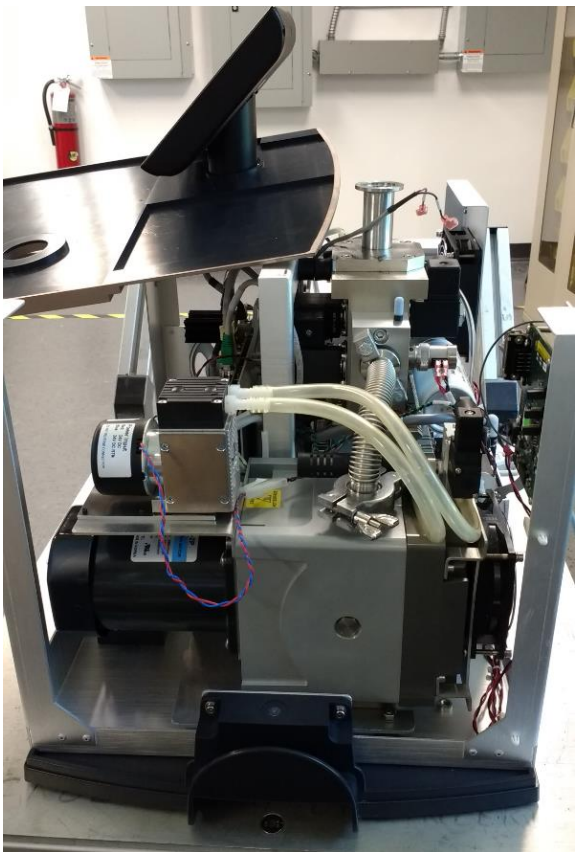


Figure 9: Covers

4. Unscrew the black cover using the Philips head screw, pull up the cover and disconnect (red arrow in Figure 10: Unit Screws)

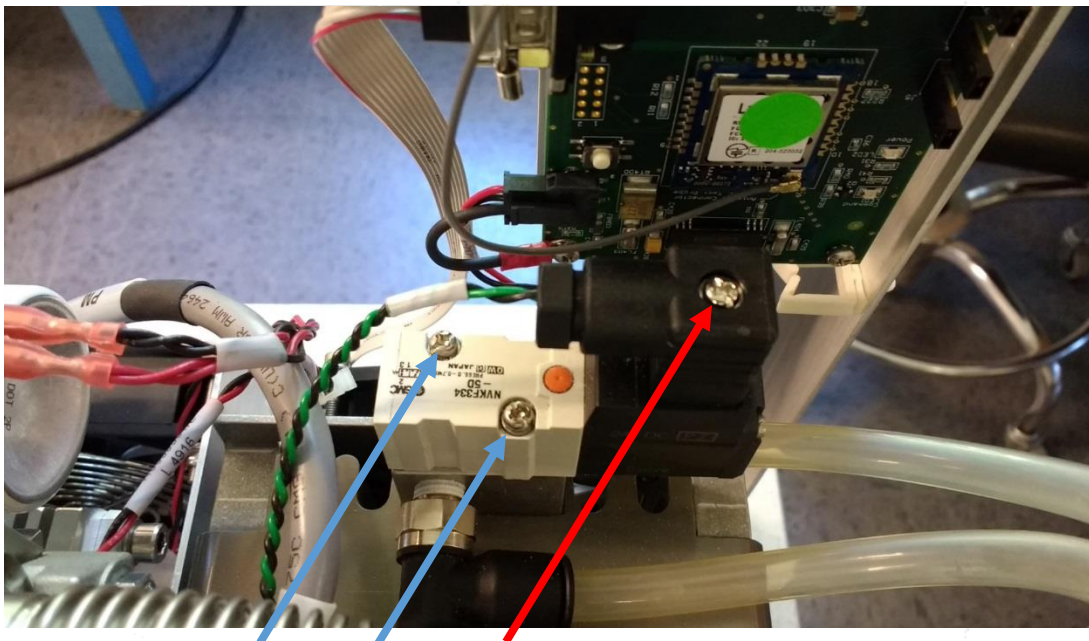


Figure 10: Unit Screws

5. Unscrew the two Philips head screws (blue arrow in “Figure 10: Unit Screws”) and pull the complete black/grey body up and out.
 6. Mount new unit and replace the three Philips screws.
 7. Attach the all covers using existing hardware.
 8. Fasten covers using a slotted screw driver.
 9. Connect the power cord and power up the unit.
 10. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
- Refer to the operator's manual if the system fails to reach the *System Ready* mode.
11. Agilent recommends a full calibration of the unit prior to leak test operations.

Part Two: G8610 Diaphragm Pump Replacement

Equipment Required

- Metric Allen Wrench Set
- M3 Phillips Head Screw Driver
- Adjustable Wrench

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. Using a slotted screw driver, remove the covers (Figure 11: Covers).

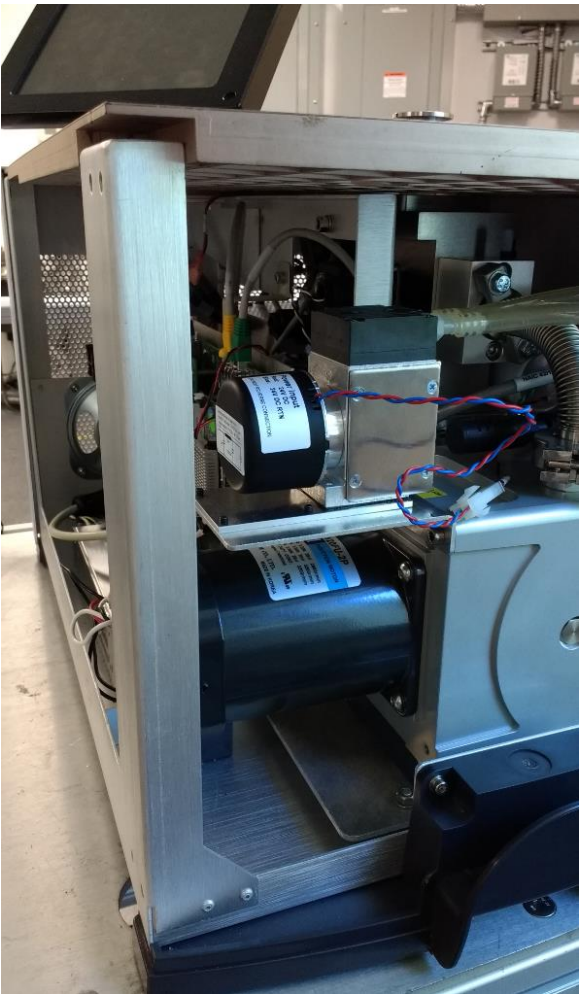


Figure 11: Covers

4. Disconnect the R3359-301 Molex connector to the Clippard valve (Figure 12: Molex Connector).

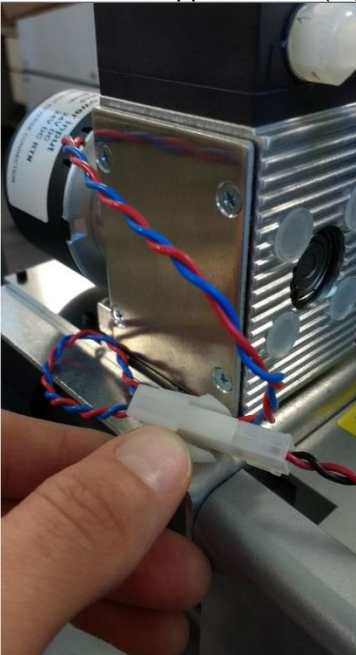


Figure 12: Molex Connector

5. Remove the ground connection to the system body (red arrow in Figure 13: Ground Connection).

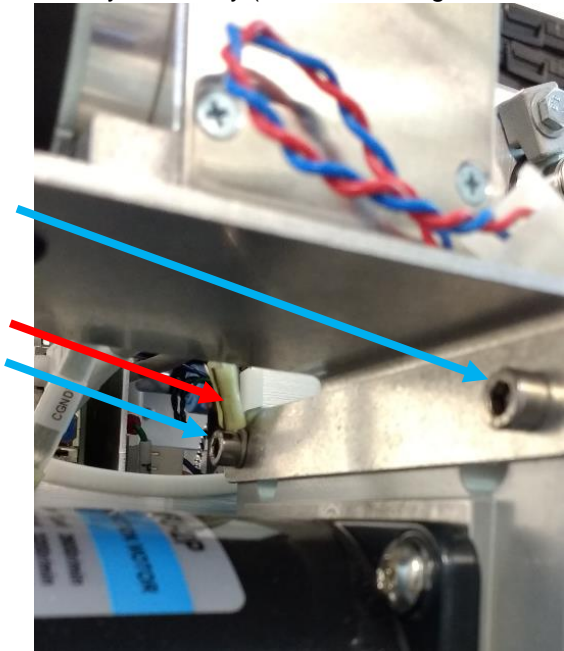


Figure 13: Ground Connection

6. Disconnect the two M4 screws mounting the plate to the scroll pump assembly (blue arrow in Figure 13: Ground Connection).

7. Remove the diaphragm pump assembly, turn it over and unscrew the four M2 screw holding the pump to the plate (Figure 14: Pump to Plate Connections).

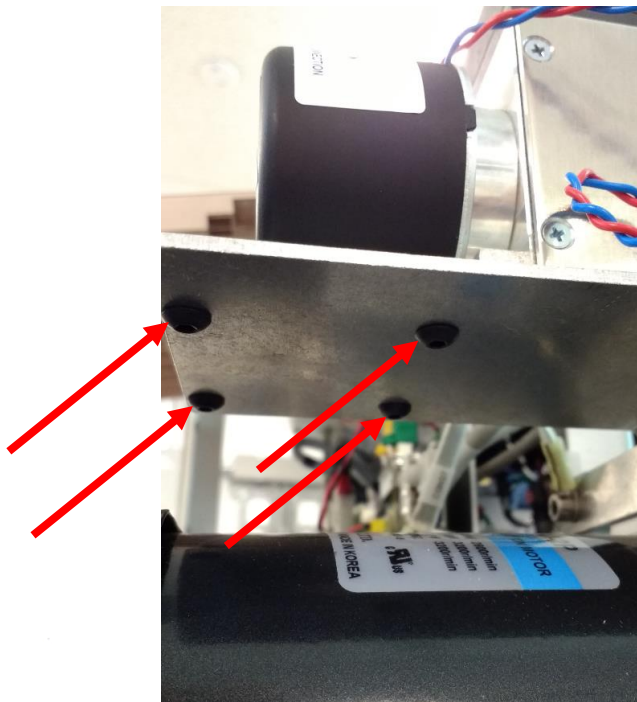


Figure 14: Pump to Plate Connections

- 8. Reattach the new pump to the plate.
- 9. Reattach the plate to the scroll pump.
- 10.Reattach the ground and Molex connections.
- 11.Attach the front cover and secure to the frame using existing hardware.
- 12.Attach the rear cover and secure to the unit using existing hardware.
- 13.Connect the power cord and power up the unit.
- 14.Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
Refer to the operator's manual if the system fails to reach the *System Ready* mode.
- 15.Agilent recommends a full calibration of the unit prior to leak test operations.

Part Two: G8610 Diaphragm and Seal Replacement

Equipment Required

- Phillips Head Screw Driver, #1
- M3 Phillips Head Screw Driver
- Metric Allen Wrench Set
- Slotted Screw Driver
- Felt Tip Pen

For clarity, some items have been omitted from views.

WARNING

Disconnect power from the unit before performing any maintenance procedure that requires physically disconnecting any part of the system.



1. Turn off the power switch located on the back of the unit and unplug.
2. Wait 30 seconds for the high voltage to dissipate.
3. Using a slotted screw driver, remove the covers (Figure 15: Covers).

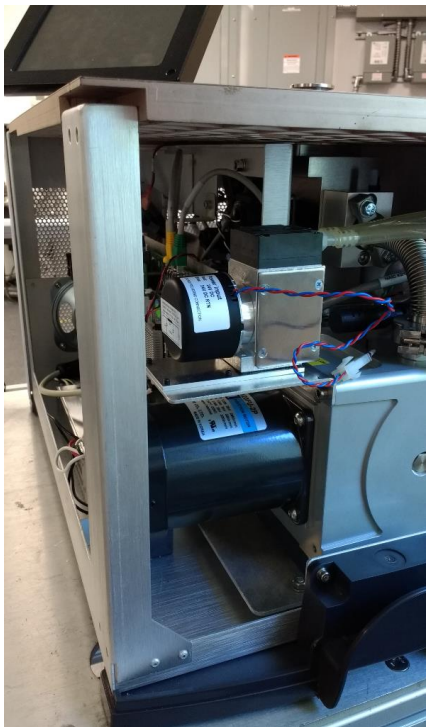


Figure 15: Covers

Always change the valve plates, diaphragm and sealing rings at the same time.

NOTE



4. Remove the pump head by (Figure 16: Pump Head Removal):

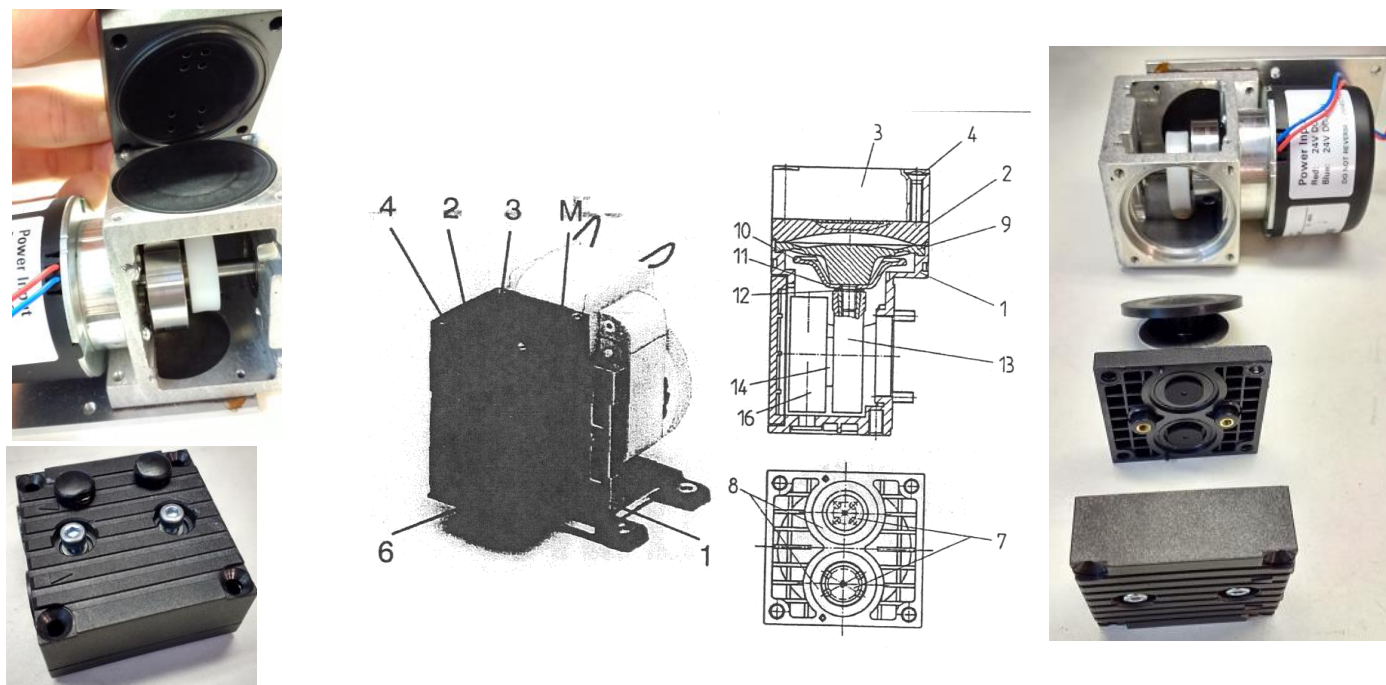


Figure 16: Pump Head Removal

- d. Making a mark (M) on the head plate (3), intermediate plate (2), and housing (1) with a felt-tip pen.
 - e. Undoing the 4 screws (4) in the head plate and lifting the head plate, with the intermediate plate, lift off the pump housing.
 - f. *For models with:*
 - *DC motors (no cooling fan):* Removing the cover (6) from the pump housing (1) by inserting a screwdriver into the slots on the sides, and prying the cover off (with care the adhesive gasket between cover and housing may be reused).
 - *No fan:* Reference is made to turning or holding the cooling fan, the necessary operations must be carried out by turning or holding the counterweight (16).
5. Change the diaphragm by (see Figure 16: Pump Head Removal):
- a. Turning the fan to bring the structured diaphragm (9) to top dead center.
 - b. Lifting the edge of the diaphragm and, gripping it on opposite sides, unscrewing it by turning counterclockwise.

CAUTION



Ensure that the disc spring (12) and diaphragm spacers (11) on the threaded portion of the diaphragm do not fall into the housing.

- f. Taking the disc spring (12), diaphragm spacer(s) (11), and diaphragm support (10) off the threaded portion of the diaphragm and retaining them.
- g. Ensuring that all parts are free from dirt and clean them, if necessary.
- h. Placing the diaphragm support, diaphragm spacer(s), and disc spring, in that order, on the thread part of the new diaphragm. The concave side of the disc spring must be towards the diaphragm.

- h. Turning the fan until the connecting rod (13) is at top dead center.
 - i. Screwing the diaphragm, complete with diaphragm support, diaphragm spacer(s), and disc spring, into the connecting rod (clockwise) and tightening it by hand.
6. Change the valve plates by (see Figure 16: Pump Head Removal):
- a. Separating the head plate (3) from the intermediate plate (2).
 - b. Removing the valve plates (7) and sealing rings (8) from the intermediate plate.
 - c. Checking that the valve seats in the head plate and the intermediate plate are clean. If scratches, distortion, or corrosion are evident, replace that part.
 - d. Laying the new valve plates in the intermediate plate recesses. The valve plates for suction and pressure sides are identical, as are upper and lower sides of the plates.
 - e. Ensuring the valve plates are not deformed by moving them gently sideways in their recesses.
 - f. Laying the sealing rings on the intermediate plate.
7. Refit the pump heads by (see Figure 16: Pump Head Removal):
- a. Turning the fan to bring the diaphragm to top dead center.
 - b. Placing the intermediate plate (2), with valve plates (7), sealing rings (8), and head plate (3) on the housing, in the position indicated by the (M).
 - c. Checking that the head plate is centered by moving it gently sideways.
 - d. Gently tightening the screws (4), evenly and diagonally.
 - e. Turning the fan to check that the pump rotates freely.
 - f. Turning the fan again to bring the diaphragm to top dead center.
 - g. Tightening the screws (4).
 - h. *For DC versions (no cooling fan):* Reattaching the cover (6) to housing (1).
8. Attach the front cover and secure to the frame using existing hardware.
9. Attach the rear cover and secure to the unit using existing hardware.
10. Connect the power cord and power up the unit.
11. Watch the home screen to verify the *Spectube Pressure Wait* message progresses to *Stabilization Wait* and *System Ready* within ten minutes.
- Refer to the operator's manual if the system fails to reach the *System Ready* mode.



Vacuum Products Division Instructions for returning products

Dear Customer,

Please follow these instructions whenever one of our products needs to be returned.

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

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

Product preparation

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

Shipping

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

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

EUROPE:

Fax: 00 39 011 9979 330

Fax Free: 00 800 345 345 00

Toll Free: 00 800 234 234 00

vpt-customer@agilent.com

NORTH AMERICA:

Fax: 1 781 860 9252

Toll Free: 800 882 7426

vpl-ra@agilent.com

PACIFIC RIM:

please visit our website for individual
office information

<http://www.agilent.com>



TERMS AND CONDITIONS

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


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



Vacuum Products Division
Request for Return Form

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

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

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

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

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

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

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies
121 Hartwell Avenue
Lexington, MA 02421 - USA
Tel. +1 781 861 7200
Fax. +1 781 860 5437
Toll-Free: +1 800 882 7426
vpl-customer-care@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-care_belgium@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
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

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

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

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

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-care_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

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

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

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-care@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-care@agilent.com

Customer Support & Service

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

EUROPE:

Toll-Free: 00 800 234 234 00
vpt-customer-care@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
Representatives in most countries

