



Vacuum Products Division



Field Installation Instructions

Hot Filament Replacement

for Agilent Technologies Helium
Leak Detection (G8610, G8611,
or G8612)

or for VS Series Leak Detector

Part Number G8600-90005

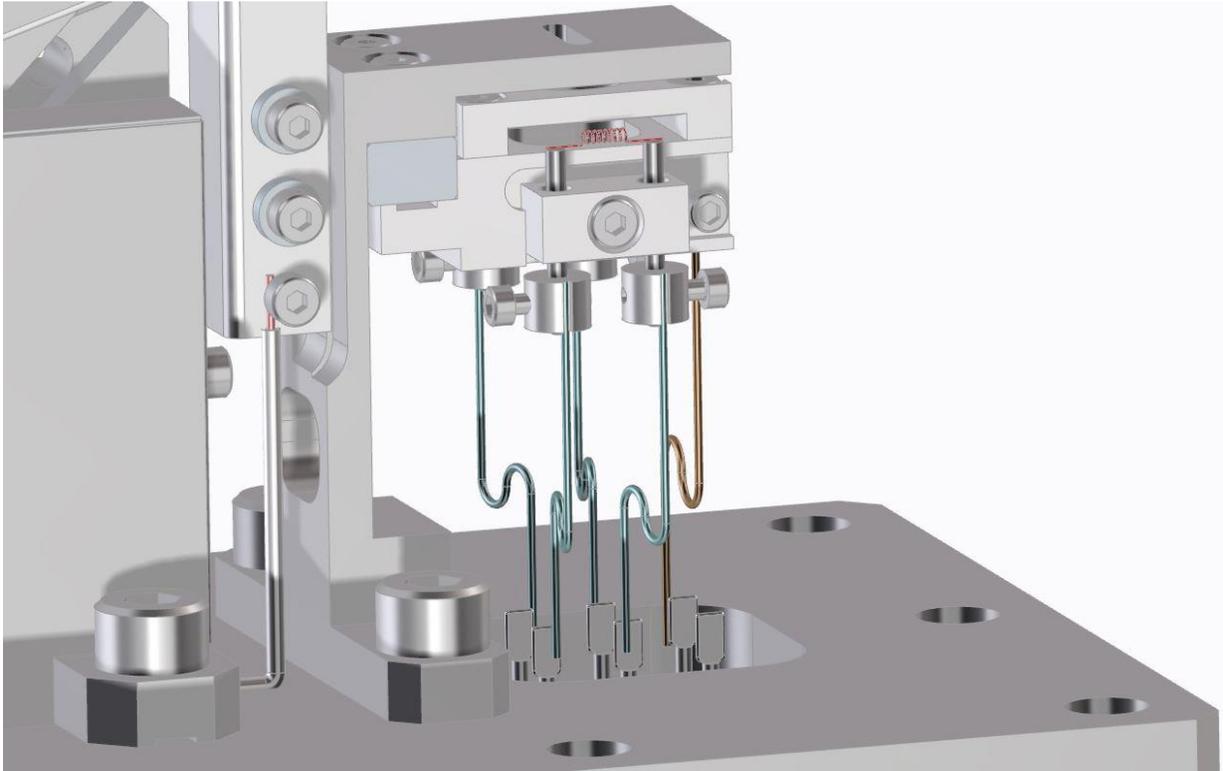
Rev. A

May 2018



G8600-90005

Filament Replacement for Agilent Helium Leak Detector



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

Warranty Replacement and Adjustment

All claims under warranty must be made promptly after occurrence of circumstances giving rise thereto, and must be received within the applicable warranty period by Seller or its authorized representative. Such claims should include the Product serial number, the date of shipment, and a full description of the circumstances giving rise to the claim. Before any Products are returned for repair and/or adjustment, written authorization from Seller or its authorized representative for the return and instructions as to how and where these Products should be returned must be obtained. Any Product returned to Seller for examination shall be prepaid via the means of transportation indicated as acceptable by Seller. Seller reserves the right to reject any warranty claim not promptly reported and any warranty claim on any item that has been altered or has been returned by non-acceptable means of transportation. When any Product is returned for examination and inspection, or for any other reason, Customer shall be responsible for all damage resulting from improper packing or handling, and for loss in transit, notwithstanding any defect or non-conformity in the Product. In all cases, Seller has the sole responsibility for determining the cause and nature of failure, and Seller's determination with regard thereto shall be final. If it is found that Seller's Product has been returned without cause and is still serviceable, Customer will be notified and the Product returned at Customer's expense; in addition, a charge for testing and examination may be made on Products so returned.

This page intentionally left blank

Table of Contents

Warranty	iii
Warranty Replacement and Adjustment	iii
1 Preface.....	1
1.1 Documentation Standards	1
1.1.1 Symbols.....	1
1.2 Hazard and Safety Information.....	2
1.2.1 Solvents	3
1.2.2 Vacuum Equipment and Cleanliness	3
1.2.3 O-ring Care.....	4
1.2.4 Metal Seal Care	4
1.2.5 Spectrometer.....	5
2 Filament Installation: Agilent Helium Leak Detector Model G8610, G8611, or G8612	6
2.1 Equipment Required for G8610, G8611, or G8612	6
2.2 Installation Procedure for G8610, G8611, or G8612.....	6
3 Filament Installation: Agilent Legacy VS Series Helium Leak Detector	16
3.1 Equipment Required for Agilent Legacy VS Series Helium Leak Detector	16
3.2 Installation Procedure for Agilent Legacy VS Series Helium Leak Detector.....	17

Table of Figures

Figure 2-1. Covers.....	6
Figure 2-2. Spectrometer access.....	7
Figure 2-3. Spectrometer exploded view	9
Figure 2-4. Filament removal and replacement	9
Figure 2-5. Filament positioning	11
Figure 2-6. Ion support assembly	12
Figure 2-7. Filament alignment	13
Figure 2-8. Spectrometer alignment.....	14
Figure 2-9. Tightening spectrometer header screws.....	15
Figure 2-10. Ion source header pin schematic.....	15
Figure 2-11. Ion source header pin locations.....	16
Figure 3-1. Rear screws	17
Figure 3-2. Fan assembly	17
Figure 3-3. Cable locations.....	18
Figure 3-4. Spectrometer access.....	19
Figure 3-5. Spectrometer exploded view	21
Figure 3-6. Filament removal and replacement	21
Figure 3-7. Filament positioning	23
Figure 3-8. Filament alignment	23
Figure 3-9. Spectrometer alignment.....	25
Figure 3-10. Tightening spectrometer header screws.....	25
Figure 3-11. Ion source header pin schematic.....	26
Figure 3-12. Ion source header pin locations.....	26

List of Tables

Table 2-1. Equipment required for G8610, G8611, or G8612 models.....	6
Table 3-1. Equipment required for VS models	16

1 Preface

1.1 Documentation Standards

1.1.1 Symbols

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.

1.2 Hazard and Safety Information

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.

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

	OFF Supply (Power)		Earth (Ground) Terminal
	ON Supply (Power)		Protective Conductor Terminal
	Warning, Risk of danger Refer to documentation		Frame or Chassis Terminal
	Alternating Current		Caution, Hot Surface
	Direct Current		Caution, Risk of Electrical Shock
	Do Not Place in Trash		

1.2.1 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.

1.2.2 Vacuum Equipment and Cleanliness

Caution



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

1.2.3 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.*

1.2.4 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.*

1.2.5 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.

2 Filament Installation: Agilent Helium Leak Detector Model G8610, G8611, or G8612

The same filament replacement kit is compatible with any model G8610, G8611, or G8612, as well as legacy Agilent model VS series leak detectors.

2.1 Equipment Required for G8610, G8611, or G8612

The following tools are suggested for assembly.

Table 2-1. Equipment required for G8610, G8611, or G8612 models

Item	Note
Filament Service Kit (PN: VSFLDHFR)	--
Slotted screw driver	Use to remove lead detector covers
Metric Allen wrench set	--
Needle nose pliers or tweezers	--
M3 Phillips head screw driver	--
Torque wrench (must be adjustable to 40 in-lbs (4.5 N-m) and 90 in-lbs (10.2 N-m))	--
Digital multimeter (Fluke 187 or equivalent)	For resistance and continuity measurements

2.2 Installation Procedure for G8610, G8611, or G8612

For clarity, some items have been omitted from views.

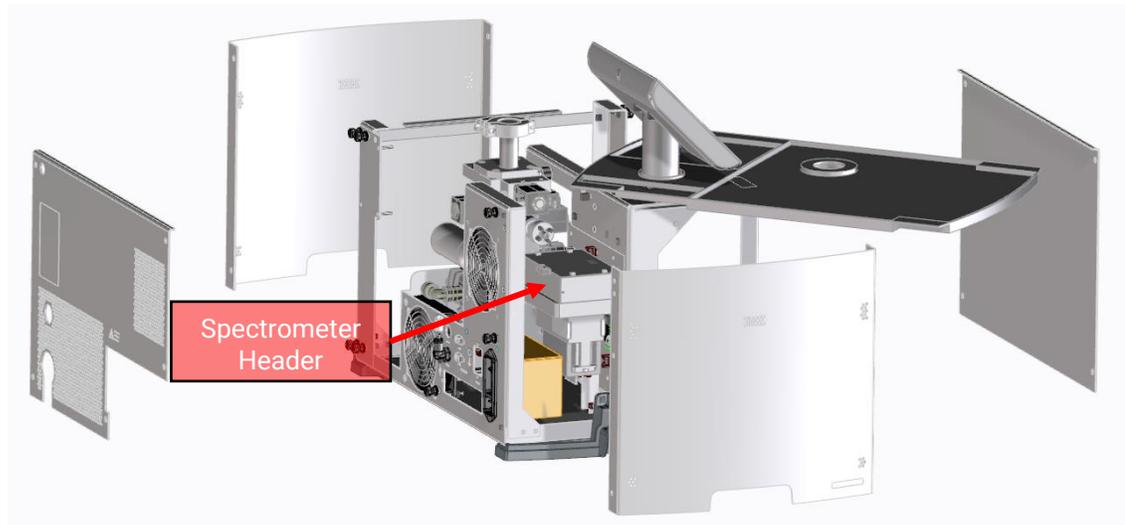


Figure 2-1. Covers

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, unfasten the captive screws holding each of the four covers. If the unit is equipped with a handle, remove the handle.

4. Detach all four covers and rotate the top lid about the display, as to give access to the spectrometer.

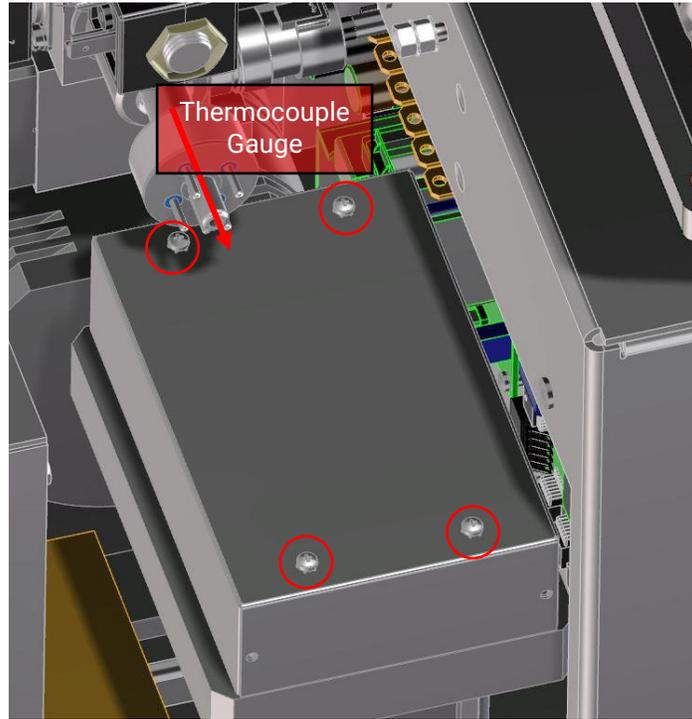


Figure 2-2. Spectrometer access

5. To remove the spectrometer top cover, it may be necessary to first pull straight out the thermocouple gauge (shown in Figure 2-2 with the red arrow. It is shown without the outside black plastic cover). Leave the thermocouple gauge connected to its wiring. Use an M3 Phillips head screw driver to remove the four screws securing the aluminum spectrometer cover. The screw shown in the top right of Figure 2-2 will be securing a grounding strap. Leave the ground strap hanging from the leak detector frame. Disconnect the two electrical cables plugged into the spectrometer board, the pre-amp RJ45 style connector on the turbo side of the spectrometer, and the ion source multi-pin connector (Cables not shown for clarity: refer to Figure 2-3). Remove the top cover, set aside with the screws.

CAUTION



Static sensitive device, ensure that personnel are properly grounded before proceeding.

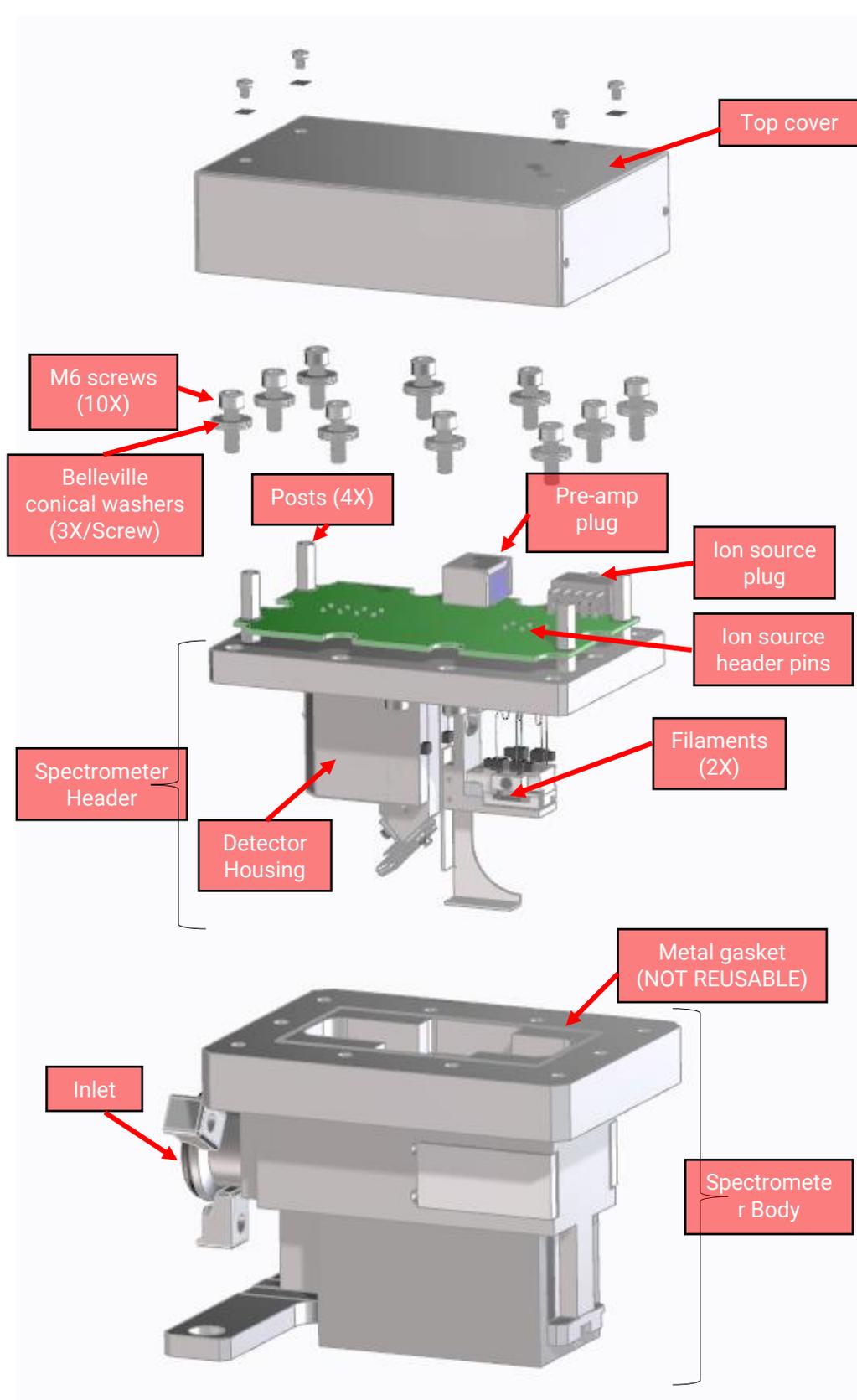


Figure 2-3. Spectrometer exploded view

CAUTION



Wear non-powdered, ESD-safe Nitrile or equivalent gloves (not included in kit) to prevent skin oils from getting on vacuum surfaces.

Refer to Figure 2-3 for disassembly of the spectrometer header.

6. Loosen screws (10X) and remove spectrometer header from spectrometer body. The vacuum system vents to atmosphere as the screws are loosened. Retain the socket head cap screws and Belleville washers. Maintain correct orientation of Belleville (conical) washers when removing.
7. Remove the spectrometer header. Rest upside-down on the four posts sticking up from the printed circuit board.
8. Remove metal gasket and discard. Do not scratch the mating surface when removing the gasket. To prevent scratching of any sealing surface, do not use metal tools for this task. Do not attempt to reuse the gasket.
9. Examine the spectrometer cavity for discolored areas. If the inside of the spectrometer body appears dirty, clean it with Vac-u-solv spectrometer cleaning kit (Agilent part number 670029096). Ensure no lint is present after cleaning.
10. On a clean work surface, remove the defective filament(s) from the spectrometer by loosening the two coupler screws. Then loosen the M2 screw and washer and remove the defective filament(s). See Figure 2-4.

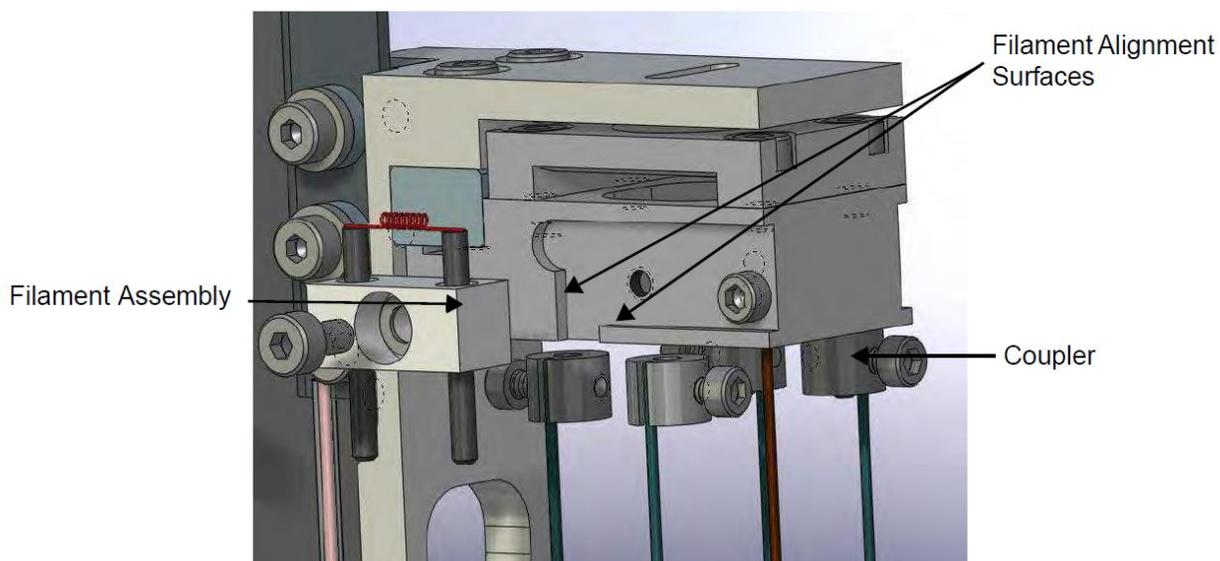


Figure 2-4. Filament removal and replacement

11. Inspect the new filament assembly for deformation, damage, and flaking of the coating prior to installation. Do not install if damaged.

12. Insert new filament assembly into the coupler.

13. Guide the ceramic against the bottom and side surfaces of the ion chamber (see Figure 2-5).

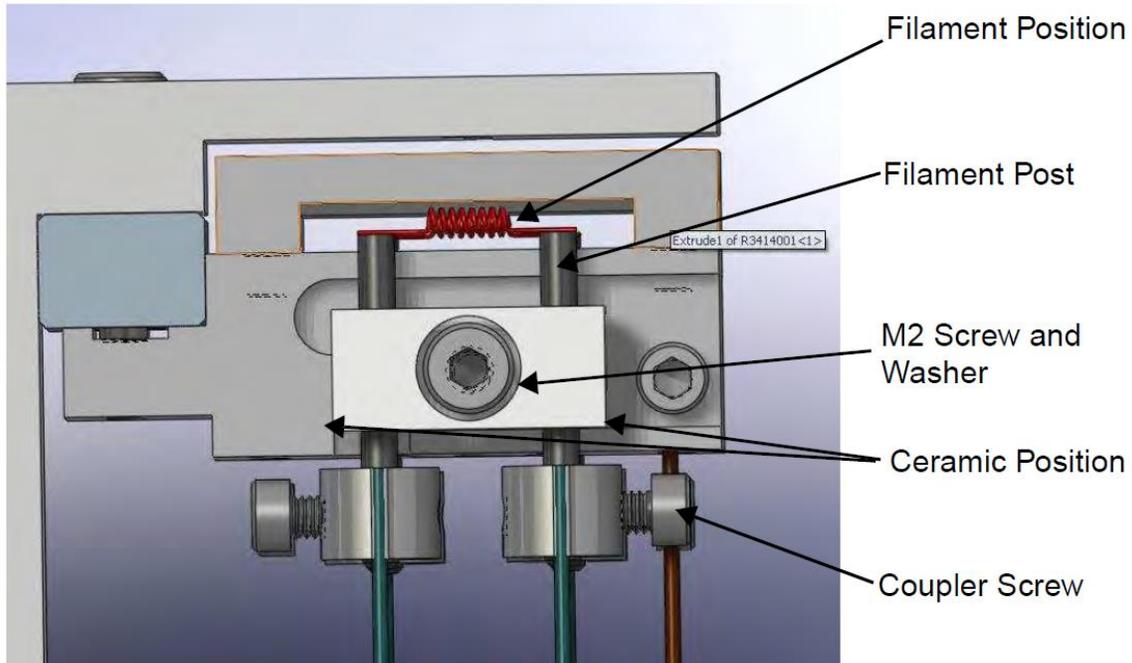


Figure 2-5. Filament positioning

14. While holding the filament in position, insert the M2 screw and washer and tighten using an Allen key.
15. Tighten the coupler screws onto the filament posts. Using small pliers, pull lightly on the extractor and chamber wires to ensure that the coupler screws are secured tightly.

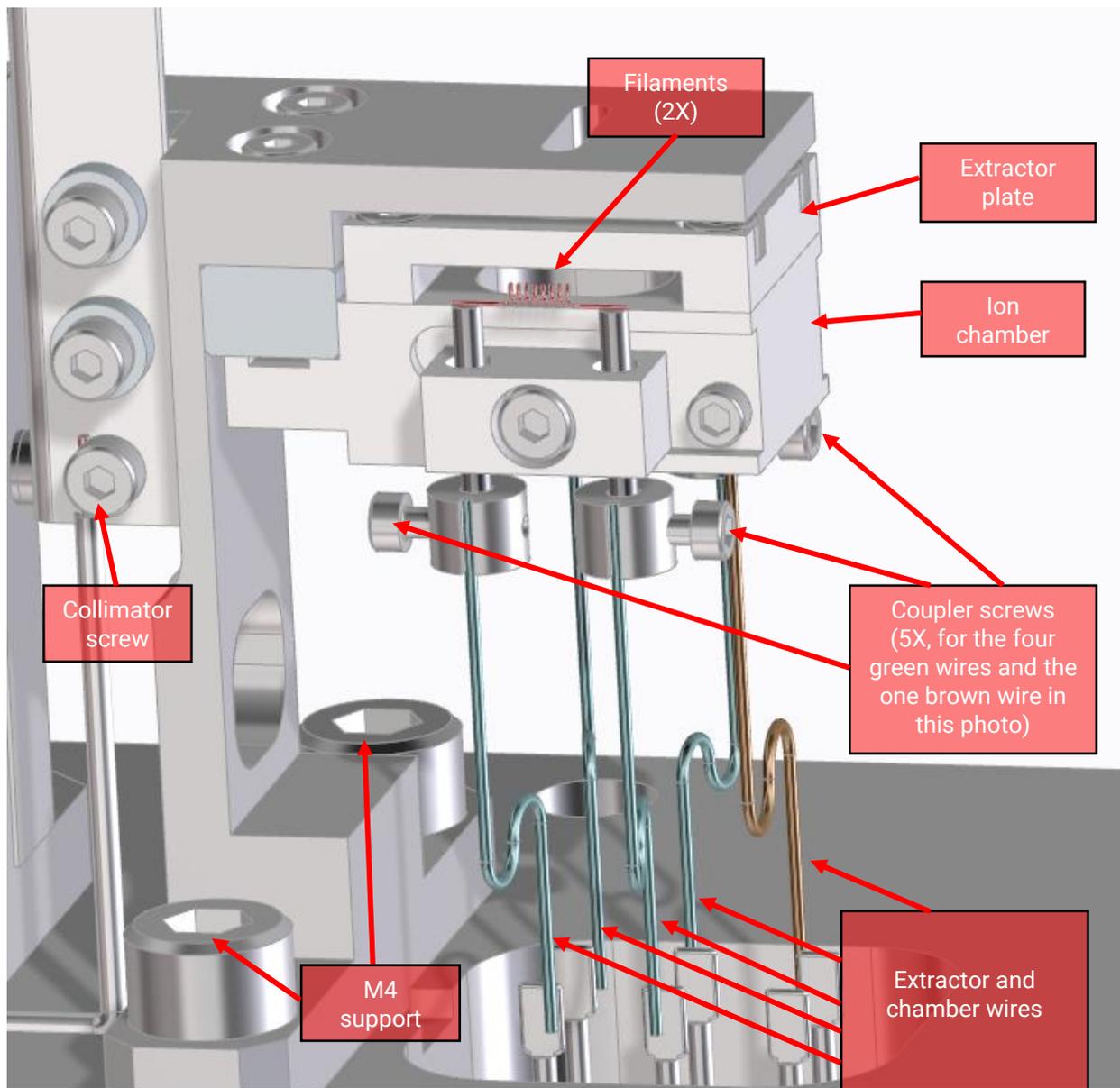


Figure 2-6. Ion support assembly

16. Visually verify ceramic is tight against the bottom and left-side ribs. Ensure that the filament is positioned in the slot of the ion source as shown in Figure 2-7.

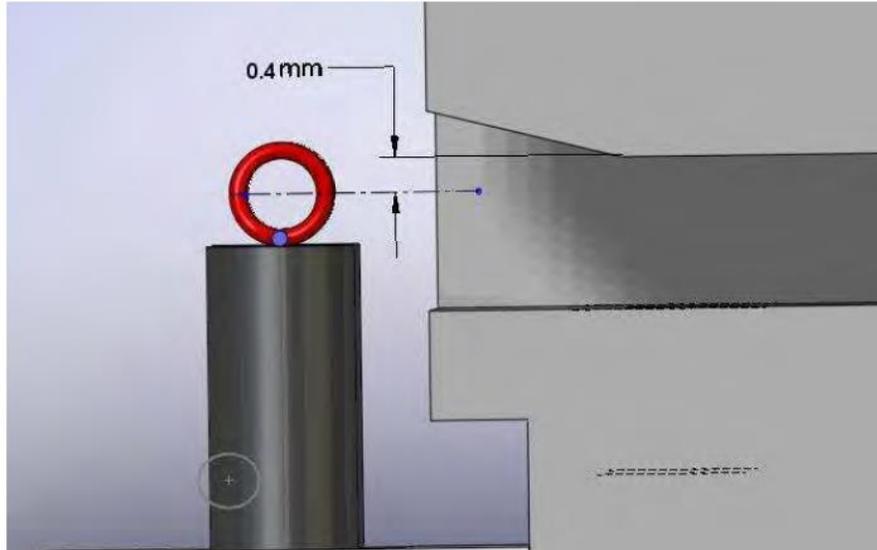


Figure 2-7. Filament alignment

17. Clean the mating surfaces of the spectrometer body and header with isopropyl alcohol and a clean lint free wipe. Agilent recommends use of the Vac-u-solv spectrometer cleaning kit (Agilent part number 670029096). Be careful to ensure no debris or lint remains in the spectrometer body after cleaning
18. Center the replacement metal gasket inside the bolt pattern and outside of the body cavity. To prevent scratching of any sealing surface, do not use metal tools for this task.
19. Guide the spectrometer header into the pocket with the detector housing closest to the inlet.

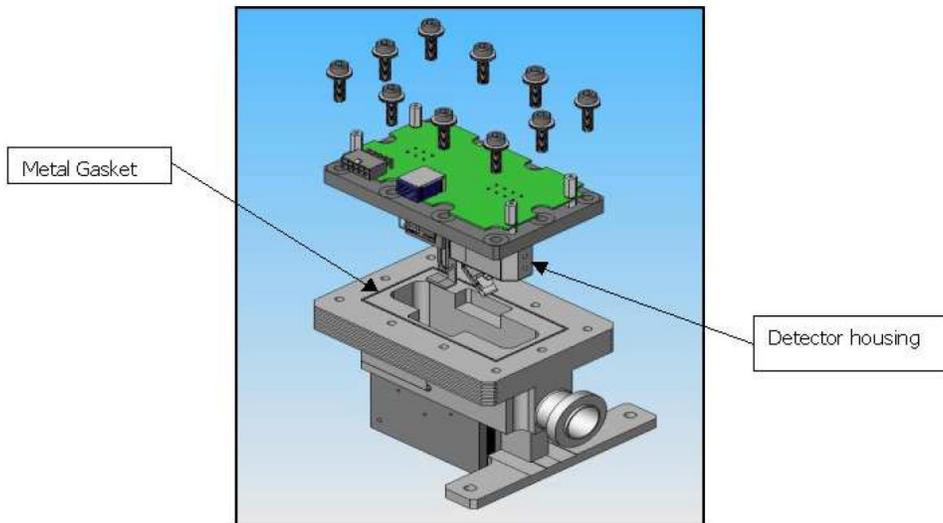


Figure 2-8. Spectrometer alignment

20. Insert a screw with three Belleville washers into each hole and finger tighten (see Figure 2-9).
21. Follow the torquing pattern in Figure 2-9. Torque screws to 40 in-lbs (4.5 N-m).
22. Re-torque screws to 90 in-lbs (10.2 N-m) following the same pattern. Go through the entire torque pattern twice to ensure the metal gasket is firmly compressed.
23. Wait a minimum of five minutes, then torque the screws again in the same pattern (Figure 2-9) to a value of 90 in-lbs (10.2 N-m). This wait period allows the washers to decompress.

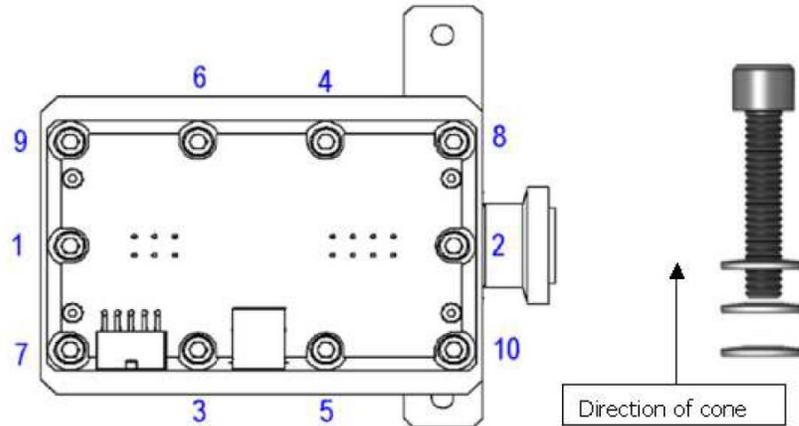


Figure 2-9. Tightening spectrometer header screws

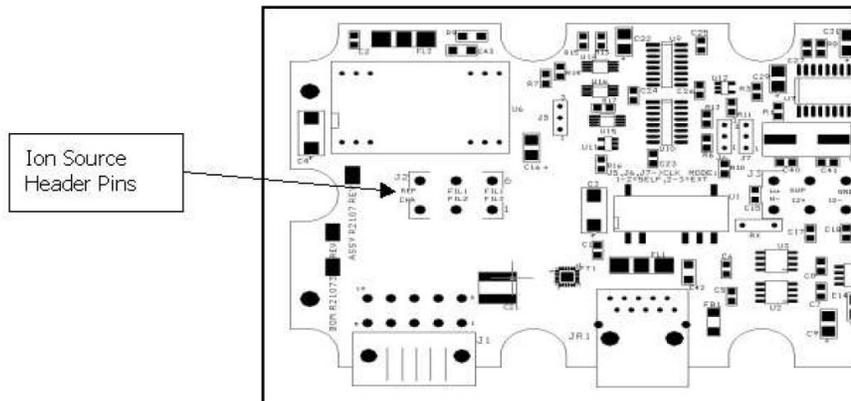


Figure 2-10. Ion source header pin schematic

24. On the PCB, use the resistance meter to verify an open circuit between any two of the six ion source header pins (except FIL-1 to FIL-1 and FIL-2 to FIL-2 which should be 0.3 Ohms or less). Also verify an open circuit between the body of the spectrometer and any of the ion source header pins (Figure 2-10). If there is continuity (short circuit) at any of the points, remove the header and inspect for shorting.

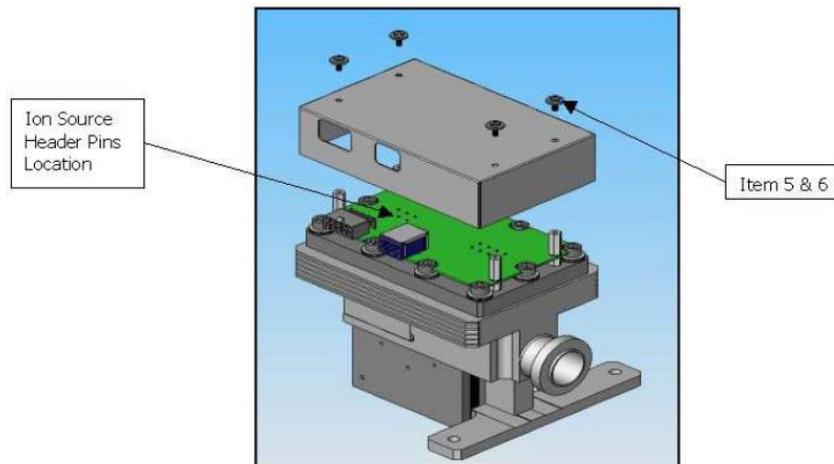


Figure 2-11. Ion source header pin locations

25. Place the spectrometer cover over the header and align the holes (Figure 2-11).
26. Install screws and washers (items 5 & 6), then tighten using a M3 Phillips screw driver (Figure 2-11). Tighten one screw over the ground strap connector.
27. Re-plug in both the pre-amp and ion source cable.
28. Reinstall the leak detector covers and fasten hardware.
29. Power up the leak detector.
30. Let the unit run for 20-30 minutes, then perform a calibration (via the I/O, front panel display, or RS232), per the user manual, to validate a successful installation. Leak check the spectrometer to ensure a leak free seal by spraying helium around the spectrometer metal seal while the leak detector is running in Test, blanked off, after calibration.

3 Filament Installation: Agilent Legacy VS Series Helium Leak Detector

The same filament replacement kit is compatible with any model G8610, G8611, or G8612, as well as legacy Agilent model VS series leak detectors.

3.1 Equipment Required for Agilent Legacy VS Series Helium Leak Detector

The following tools are suggested for assembly.

Table 3-1. Equipment required for VS models

Item	Note
Filament Service Kit (PN: VSFLDHFR)	--
Extended Length M5 Allen Wrench	--
Metric Allen wrench set	--

Needle nose pliers or tweezer	--
M3 Phillips head screw driver	--
Torque wrench (must be adjustable to 40 in-lbs (4.5 N-m) and 90 in-lbs (10.2 N-m))	--
Digital multimeter (Fluke 187 or equivalent)	For resistance and continuity measurements

3.2 Installation Procedure for Agilent Legacy VS Series Helium Leak Detector

For clarity, some items have been omitted from views.

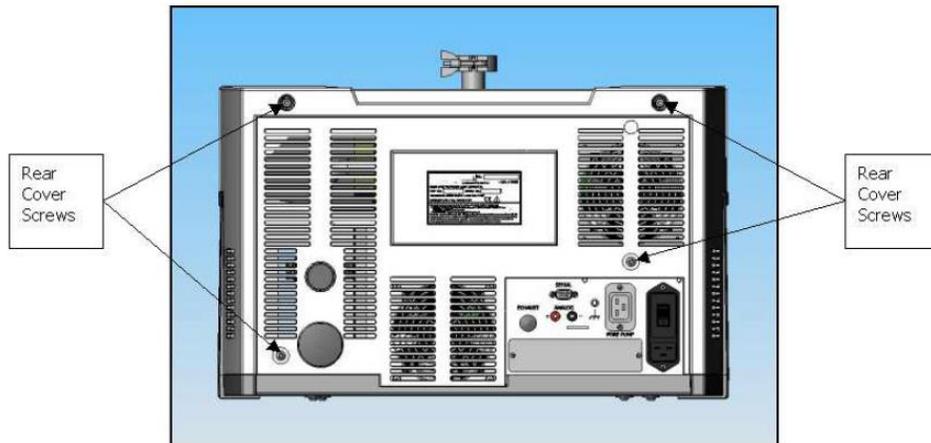


Figure 3-1. Rear screws

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 an extended length M5 Allen wrench, remove the four screws holding the rear plastic cover (Figure 3-1).
4. Detach the rear plastic cover from the unit.

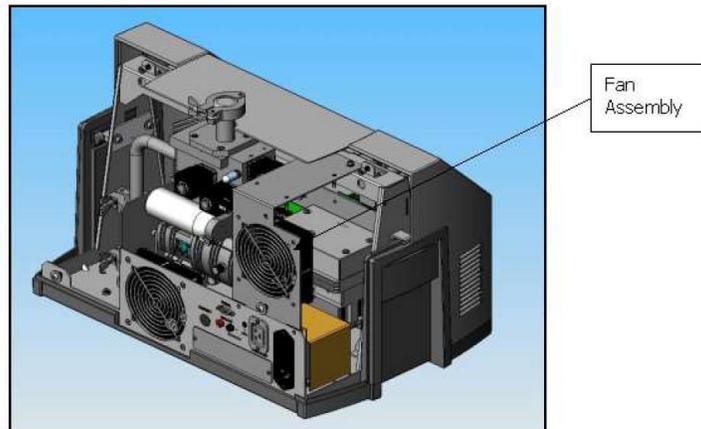


Figure 3-2. Fan assembly

5. Disconnect the fan cable and remove the fan assembly by unfastening two wing nuts and one M4 socket head cap screw (Figure 3-2).

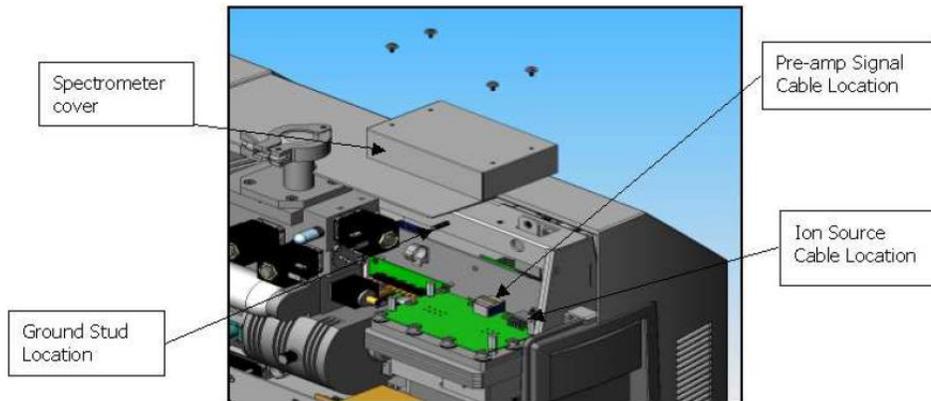


Figure 3-3. Cable locations

6. Disconnect the spectrometer ground cable, the preamp signal cable and the ion source cable from the spectrometer (Figure 3-3).

NOTE *Cables not shown for clarity.*



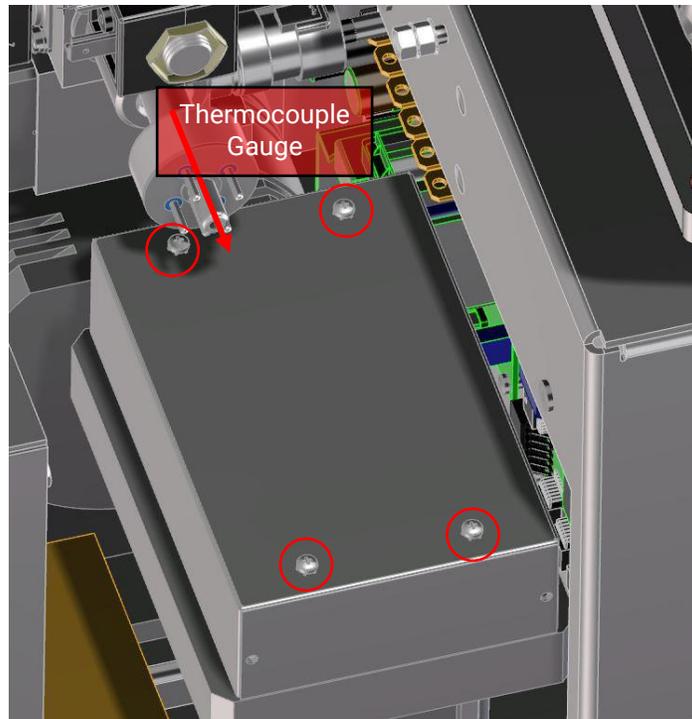


Figure 3-4. Spectrometer access

7. To remove the spectrometer top cover, it may be necessary to first pull straight out the thermocouple gauge (shown in Figure 3-4 with the red arrow. It is shown without the outside black plastic cover). Leave the thermocouple gauge connected to its wiring. Use an M3 Phillips head screw driver to remove the four screws securing the aluminum spectrometer cover. The screw shown in the top right of Figure 3-4 will be securing a grounding strap. Leave the ground strap hanging from the leak detector frame. Disconnect the two electrical cables plugged into the spectrometer board, the pre-amp RJ45 style connector on the turbo side of the spectrometer, and the ion source multi-pin connector (Cables not shown for clarity; refer to Figure 3-5). Remove the top cover, set aside with the screws.

CAUTION



Static sensitive device, ensure that personnel are properly grounded before proceeding.

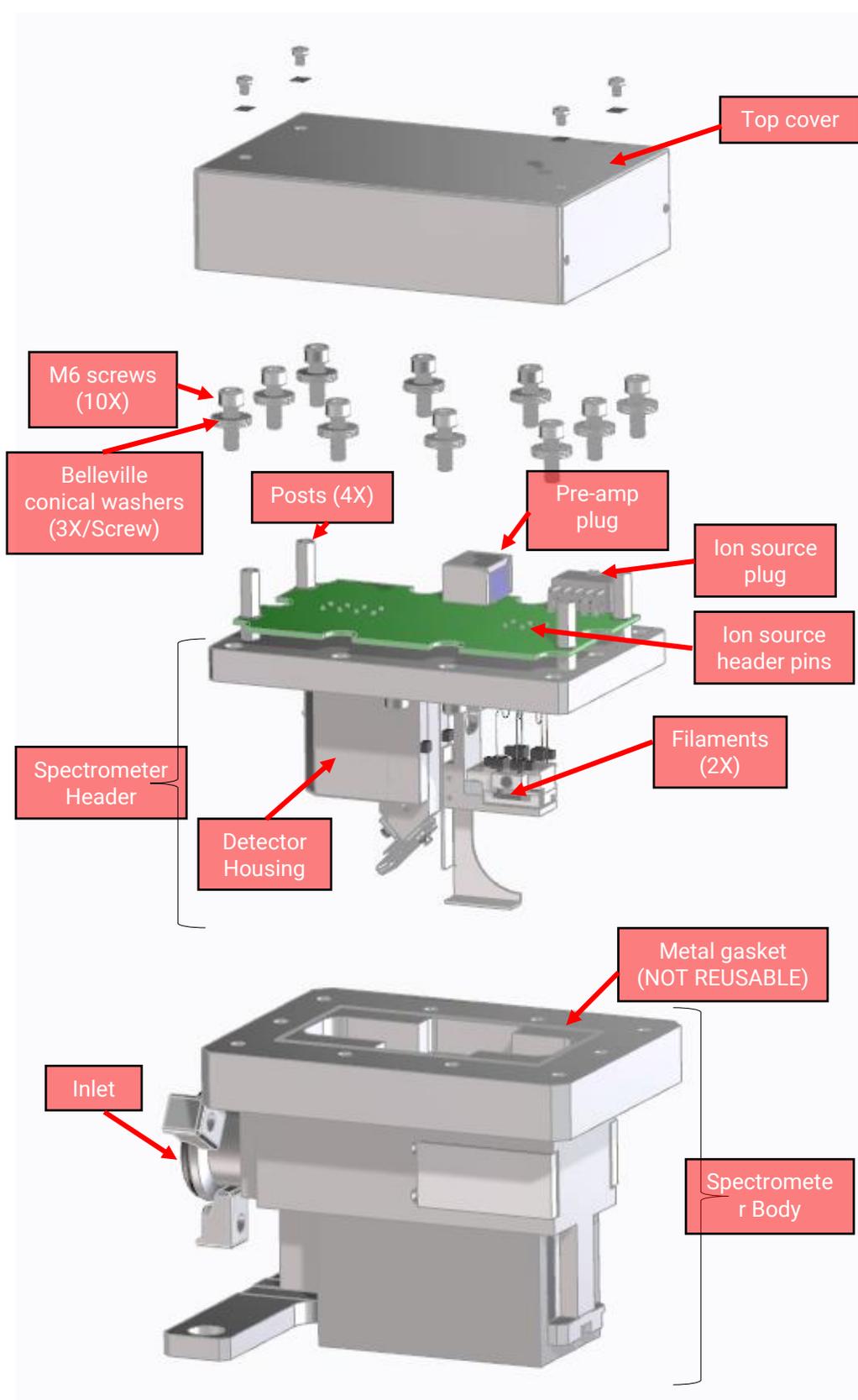


Figure 3-5. Spectrometer exploded view

CAUTION



Wear non-powdered, ESD-safe Nitrile or equivalent gloves (not included in kit) to prevent skin oils from getting on vacuum surfaces.

8. Refer to Figure 3-5 for reference in disassembly of the spectrometer header. Loosen screws (10X) and remove spectrometer header from spectrometer body. The vacuum system vents to atmosphere as the screws are loosened. Retain the socket head cap screws and Belleville washers. Maintain correct orientation of Belleville (conical) washers when removing.
9. Remove the spectrometer header. Rest upside-down on the four posts sticking up from the printed circuit board.
10. Remove metal gasket and discard. Do not scratch the mating surface when removing the gasket. To prevent scratching of any sealing surface, do not use metal tools for this task. Do not attempt to reuse the gasket.
11. Examine the spectrometer cavity for discolored areas. If the inside of the spectrometer body appears dirty, clean it with Vac-u-solv spectrometer cleaning kit (Agilent part number 670029096).
12. On a clean work surface, remove the defective filament(s) from the spectrometer by loosening the two coupler screws. Then loosen the M2 screw and washer and remove the defective filament(s). See Figure 3-6.

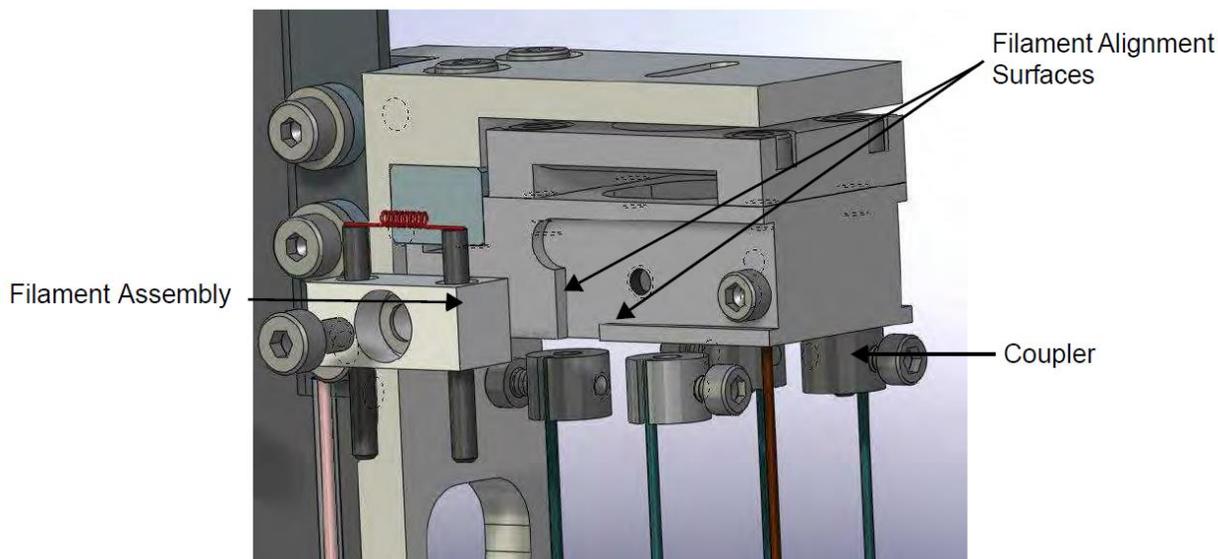


Figure 3-6. Filament removal and replacement

13. Inspect the new filament assembly for deformation, damage, and flaking of the coating prior to installation. Do not install if damaged.
14. Insert new filament assembly into the coupler.

15. Guide the ceramic against the bottom and side surfaces of the ion chamber (see Figure 3-7).

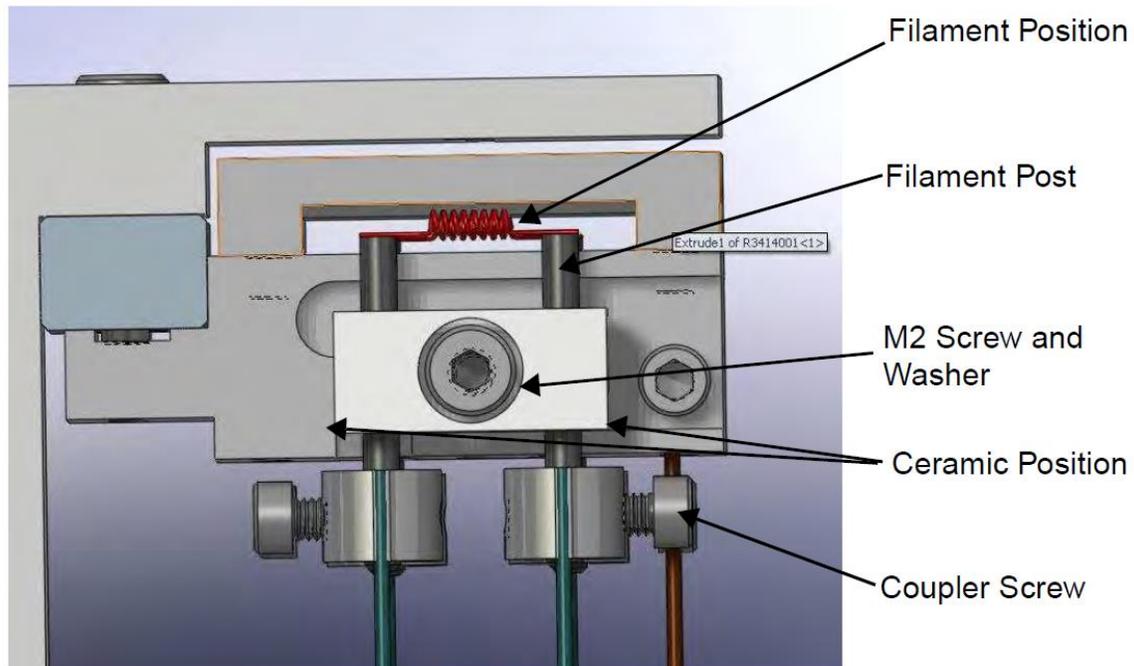


Figure 3-7. Filament positioning

16. While holding the filament in position, insert the M2 screw and washer and tighten using an Allen key.
17. Tighten the coupler screws onto the filament posts. Using small pliers, pull lightly on the extractor and chamber wires to ensure that the coupler screws are secured tightly.
18. Visually verify ceramic is tight against the bottom and left-side ribs. Ensure that the filament is positioned in the slot of the ion source as shown in Figure 3-7.

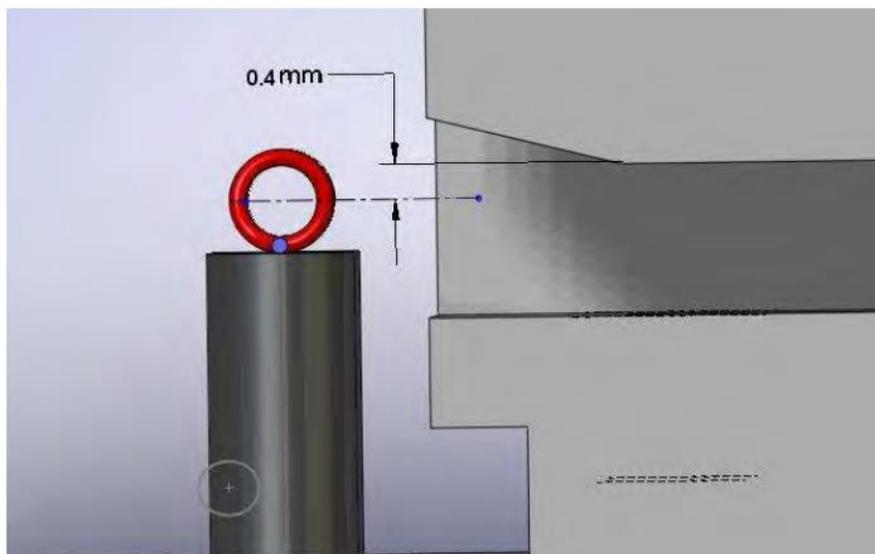


Figure 3-8. Filament alignment

19. Clean the mating surfaces of the spectrometer body and header with isopropyl alcohol and a clean lint free wipe. Agilent recommends use of the Vac-u-solv spectrometer cleaning kit (Agilent part number 670029096). Be careful to ensure no debris or lint remains in the spectrometer body after cleaning
20. Center the replacement metal gasket inside the bolt pattern and outside of the body cavity. To prevent scratching of any sealing surface, do not use metal tools for this task.
21. Guide the spectrometer header into the pocket with the detector housing closest to the inlet.

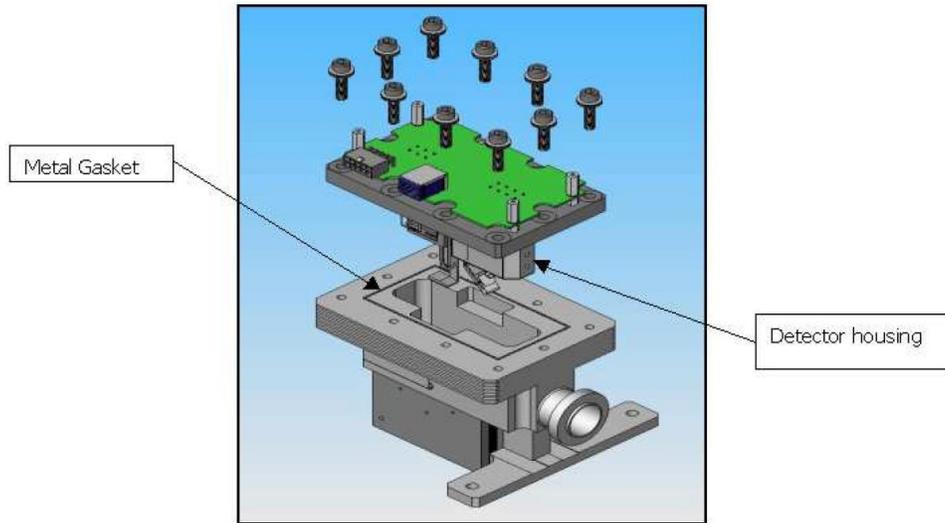


Figure 3-9. Spectrometer alignment

22. Insert a screw with three Belleville washers into each hole and finger tighten (see Figure 3-10).
23. Follow the torquing pattern in Figure 2-9. Torque screws to 40 in-lbs (4.5 N-m).
24. Re-torque screws to 90 in-lbs (10.2 N-m) following the same pattern. Go through the entire torque pattern twice to ensure the metal gasket is firmly compressed.
25. Wait a minimum of five minutes, then torque the screws again in the same pattern (Figure 3-10) to a value of 90 in-lbs (10.2 N-m). This wait period allows the washers to decompress.

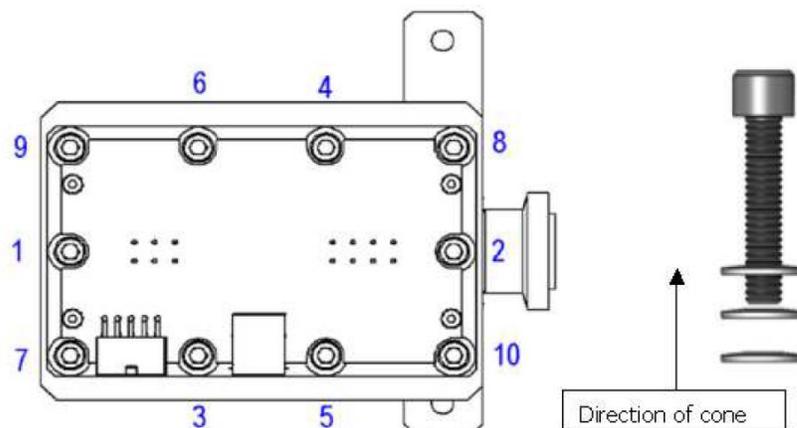


Figure 3-10. Tightening spectrometer header screws

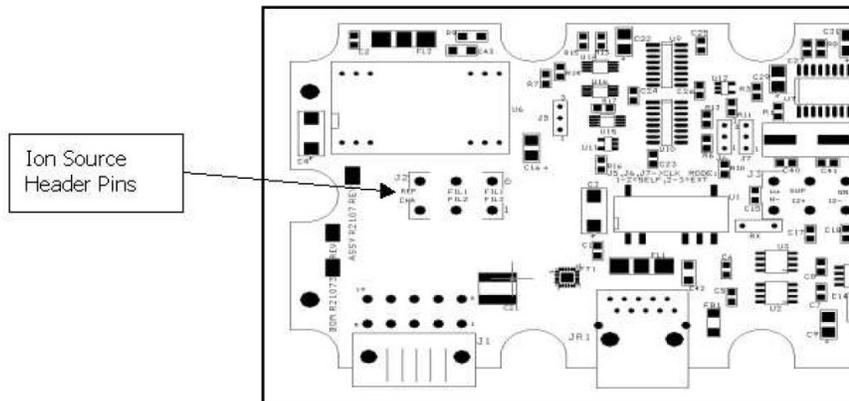


Figure 3-11. Ion source header pin schematic

26. On the PCB, use the resistance meter to verify an open circuit between any two of the six ion source header pins (except FIL-1 to FIL-1 and FIL-2 to FIL-2 which should be 0.3 Ohms or less). Also verify an open circuit between the body of the spectrometer and any of the ion source header pins (Figure 3-11). If there is continuity (short circuit) at any of the points, remove the header and inspect for shorting.

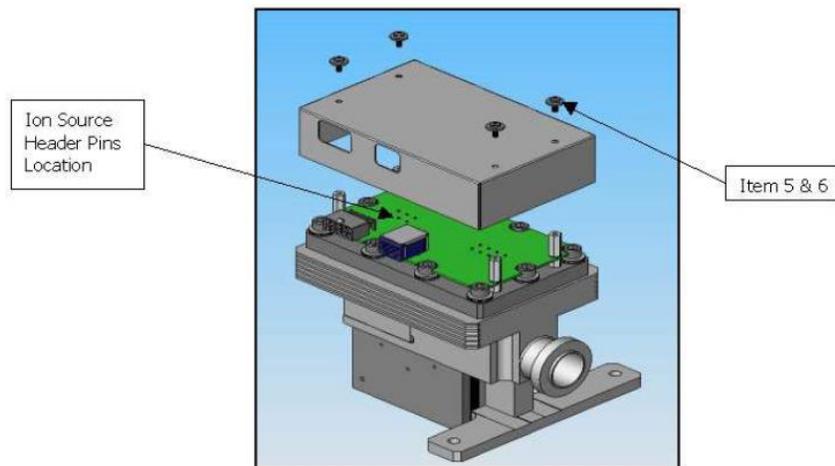


Figure 3-12. Ion source header pin locations

27. Place the spectrometer cover over the header and align the holes (Figure 3-12).
28. Install screws and washers (items 5 & 6), then tighten using a M3 Phillips screw driver (Figure 3-12). Tighten one screw over the ground strap connector.
29. Re-plug in both the pre-amp and ion source cable.
30. Reinstall the leak detector covers and fasten hardware.
31. Power up the leak detector.
32. Let the unit run for 20-30 minutes, then perform a calibration (via the I/O, front panel display, or RS232), per the user manual, to validate a successful installation. Leak check the spectrometer to ensure a leak free seal by spraying helium around the spectrometer metal seal while the leak detector is running in Test, blanked off, after calibration.

This page intentionally left blank.

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, etc.).

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 nonreturned/ 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.



Agilent Vacuum Products Division

Instructions for returning products

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

Equipment			
Product Description:	Agilent Part No.:	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				
Sensitive	<input type="checkbox"/> YES <input type="checkbox"/> NO				
Other Dangerous Substance	<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? APPLICABLE	<input type="checkbox"/> YES <input type="checkbox"/> NO <input type="checkbox"/> NOT
I confirm to place this declaration on the outside of the shipping box.	

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, Sao Paulo CEP: 06460-040
Toll-Free: +0800 728 1405

China

Agilent Technologies (China) Co.
Ltd
No.3, Wang Jing Bei Lu, Chao Yang
District
Beijing, 100102, China
Tel. +86 (0)10 64397888
Fax. +86 (0)10 64392765
Toll free: 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

© Agilent Technologies, Inc. 2018

