



Agilent Wireless Remote Base Unit for HLD series G8610, G8611 or G8612 Leak Detectors (G8610-63000) and VS Series (G8600-68001)

Instructions for Installation and Technical Specification



Notices

Manual Part Number

699910002

Edition E.00, September 2024

Copyright

© Agilent Technologies, Inc. 2024

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies Italia S.p.A.
Vacuum Products Division
Via F.lli Varian, 54
10040 Leini (TO)
ITALY

www.agilent.com

Instrument Manufacturing

Agilent Technologies LDA Malaysia
Sdn Bhd Bayan Lepas
Free Industrial Zone 11900 Bayan
Lepas, Penang Malaysia

Printed in Italy

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Contents

1	Instructions for Installation and Technical Specification	4
	About this manual	5
	Validity	5
	Definitions and terminology	6
	Definition of Caution, Warning and Note	6
	Warning Symbols	7
	Safety	8
	Installation on HLD series G8610, G8611 or G8612 Leak Detector (G8610-63000)	9
	Installation Procedure	9
	Installation on VS Series Leak Detector (G8600-68001)	18
	Installation Procedure	18
	Disposal	26
	Service	27



1

Instructions for Installation and Technical Specification

About this manual	5
Validity	5
Definitions and terminology	6
Definition of Caution, Warning and Note	6
Warning Symbols	7
Safety	8
Installation on HLD series G8610, G8611 or G8612 Leak Detector (G8610-63000)	9
Installation Procedure	9
Installation on VS Series Leak Detector (G8600-68001)	18
Installation Procedure	18
Disposal	26
Service	27

About this manual

Validity

This manual lists the instructions for the users of the Wireless Remote Base Unit, with particular reference to the notions relating to safety, operation and first level maintenance, limited to maintenance operations for which the user is responsible.

The maintenance operations, illustrated in the specific sections, with specific provisions relating to the higher level of maintenance (personnel specifically trained for maintenance operations) must not be carried out by the user.

NOTE

- 1 This manual contains useful information so that all personnel using the Wireless Remote Base Unit can operate it safely and guarantee perfect efficiency, for its entire life span.
 - 2 Keep this manual, together with all the related publications, in an accessible place known to all operators/maintenance personnel.
-

Definitions and terminology

Definition of Caution, Warning and Note

Some important references of this manual are highlighted and framed in contrasting color.

CAUTION

Caution messages are displayed before procedures which, if not observed, could cause damage to the equipment.

WARNING

Warning messages draw the operator's attention to a specific procedure or practice which, if not performed correctly, could result in serious personal injury.








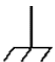


NOTE

Notes are intended to call attention to important information and provide more detail regarding specific steps.

Warning Symbols

The following is a list of symbols that appear in conjunction with warnings on the Wireless Remote Base Unit. The hazard they describe is also shown. A triangular symbol indicates a warning. The meanings of the symbols that may appear alongside warnings in the documentation are as follows.

The following symbol may be used on warning labels attached to the instrument. When you see this symbol, refer to the relevant operation or service manual for the correct procedure referred to by that warning label.

	Generic hazard
	Dangerous voltages
	Hot surfaces
	Alternate current (AC)
	Protective Ground Terminal
	Power ON
	Power OFF
	Frame or chassis Terminal
	CE certification
	Waste Electrical and Electronic Equipment

Safety

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.

NOTE

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 liquids into the unit through the ventilation openings or Front Panel buttons. Wipe the surface with a dry lint-free cloth.

CAUTION

The 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.

This manual includes two sections:

- the first gives instructions for installation with systems G8610, G8611 or G8612 HLD Series Leak Detectors (G8610-63000);
- the second section gives instructions for installation with the VS Series Leak Detector (G8600-68001).

Installation on HLD series G8610, G8611 or G8612 Leak Detector (G8610-63000)

Equipment Required	Kit Includes
Slotted Screw Driver	Wireless Base PCB
Metric Allen Wrench Set	Wireless Remote Label
M3 Philips Head Screw Driver	RS232 Cable
Adjustable Wrench	Antenna
	M3 Screw (Qty. 4)

Installation Procedure

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.

Instructions for Installation and Technical Specification

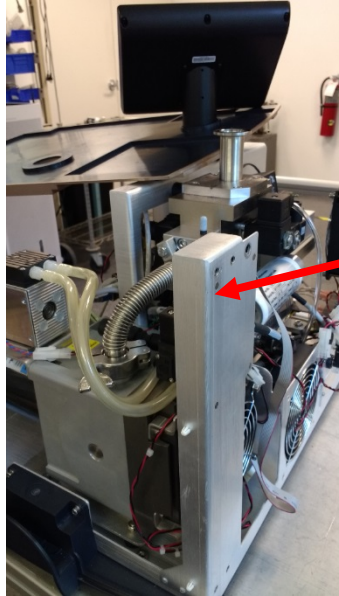
- 1 Turn off the power switch located on the back of the unit and unplug.
- 2 Wait 30 seconds for the high voltage to dissipate.



Figure 1 Orientation

- 3 Using slotted screw driver, unfasten the four screws holding the front, rear and right side covers (Figure 1: Covers) and detach. The left side cover is secured with hex-head screws.

Instructions for Installation and Technical Specification



PCB to mount to opposite side of this aluminum support

Figure 2 Covers

- 4 Rotate the lid 90° to give better access to the area where the antenna will be installed.

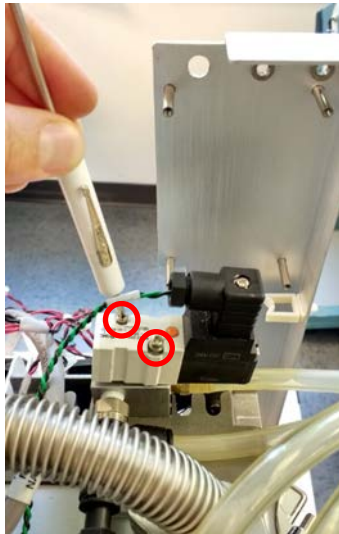


Figure 3 Removing Valve

Instructions for Installation and Technical Specification

- 5 If the leak detector configuration has a valve in front of the bottom left post, move that valve aside by unscrewing the two screws circled (Figure 3: Removing valve). Keeping the wiring connected, set valve aside. This valve is only used on certain models of leak detectors. If the valve is not present, ignore this step.
- 6 Mount the printed circuit board to the four posts, noting the correct orientation. Fasten with screws (Figure 4: Base Board Assembly with Posts).



Figure 4 Base Board Assembly with Posts

Instructions for Installation and Technical Specification

- Using the provided washer and nut, hand tighten the antenna to the outside of the leak detector wall, and then apply a $\frac{1}{2}$ turn using an adjustable wrench (Figure 5: Wireless Antenna).

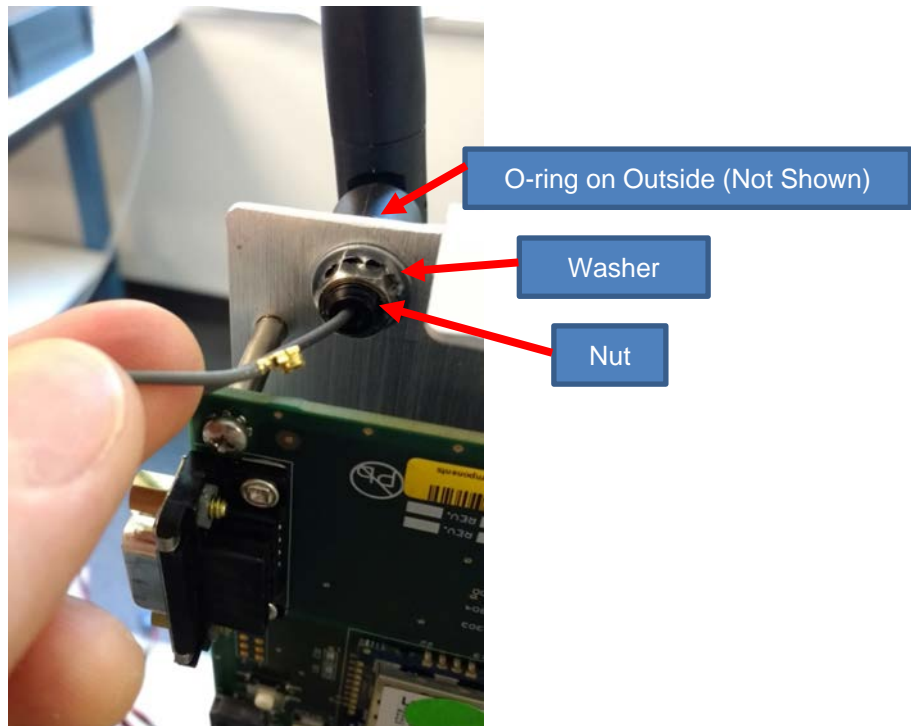


Figure 5 Wireless Antenna

Instructions for Installation and Technical Specification

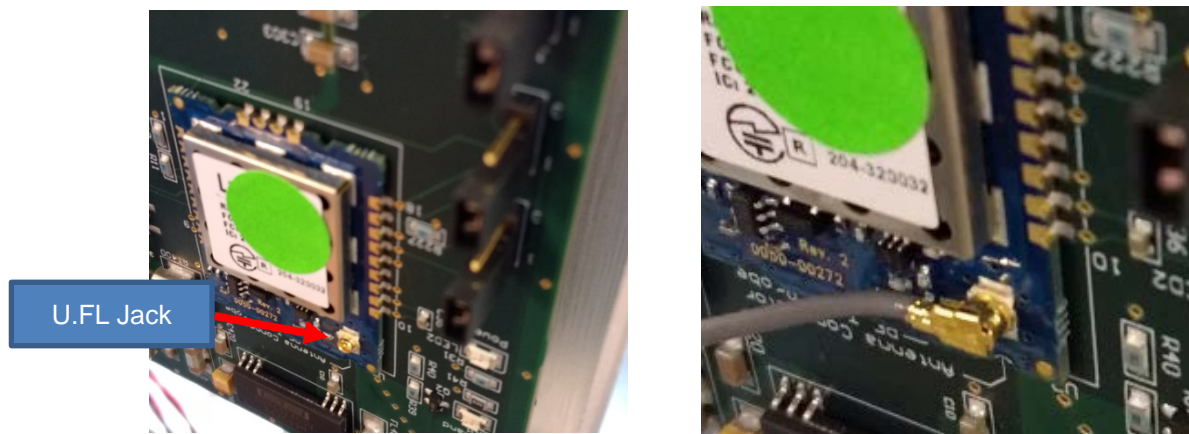


Figure 6 Antenna Connector

CAUTION

Use proper ESD safety techniques when performing a PCB card install.

- 8 Connect the antenna cable to the U.FL on the wireless base board (Figure 6: Antenna Connector). Carefully center the connector on the jack and press the connector onto the jack until it snaps into place. Rotate the antenna cable to ensure the connector is firmly seated in the jack.

Instructions for Installation and Technical Specification

- 9 Connect the RS-232 on the wireless board using its two slotted screws (Figure 7: RS232 Connection, left side). Plug in the other side of the ribbon cable to the green port on the CPU board, located on the front side of the leak detector, away from the antenna's printed circuit board (Figure 7: RS232 Connection, right side).

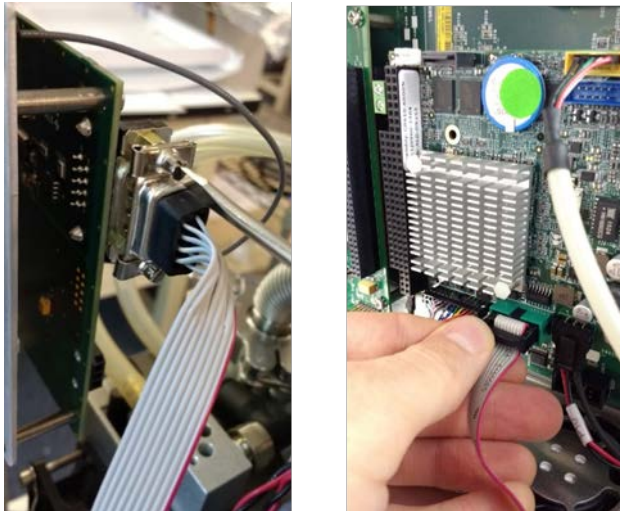


Figure 7 RS232 Connection

Instructions for Installation and Technical Specification

10 Unfurl the black/red wire (P1-RF) with the 24 V three-pin connector and the ground lug (Figure 8: 24 V Power Cable), which is tie rapped at the base of the leak detector, and:

- Plug the 24 V connector labeled P1-RF into the three-pin connector on the wireless board.
- Attach the 24 V ground lug to one of the bracket screws.

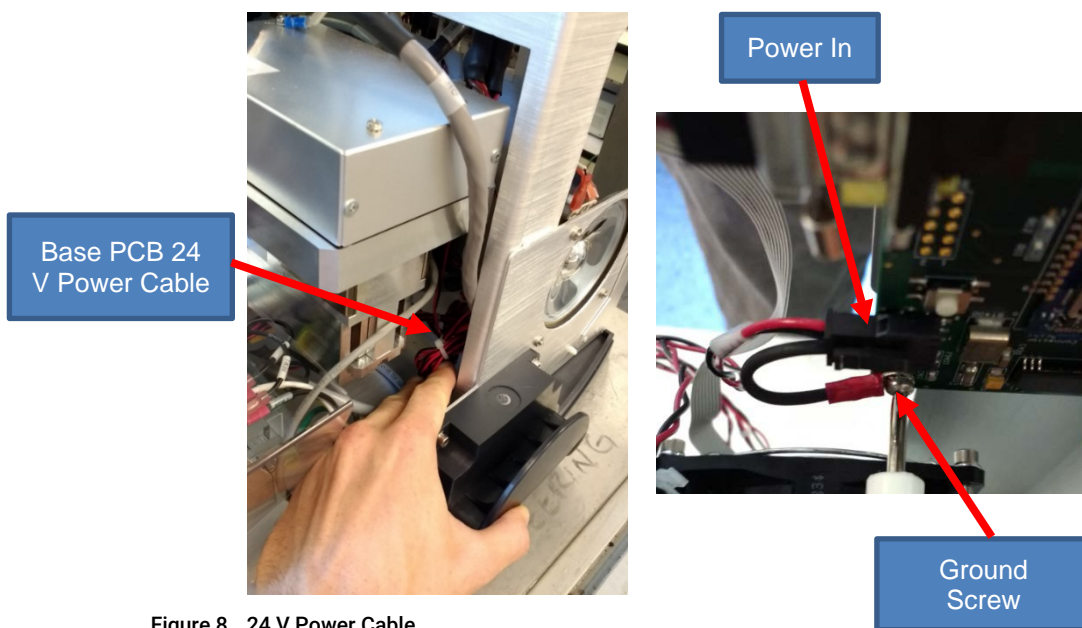


Figure 8 24 V Power Cable

- 11** Reinstall the valve and dress internal wires inside of leak detector (Figure 3: Removing Valve).
- 12** Reattach front and side (Figure 2: Covers).
- 13** Rotate the antenna to its horizontal locked position.

Instructions for Installation and Technical Specification

- 14 Guide the antenna through its location hole in the rear cover. Then secure the rear cover to the frame using existing hardware.



Figure 9 Label

- 15 Attach the label to the rear cover in location shown (Figure 9: Label). This new label will help service identify the equipment used. Do not cover the existing label on the leak detector.
- 16 Connect the power cord and power up the unit.
- 17 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.
- 18 After installation refer to the Leak Detector manual section for leak detector activation of the wireless remote.
- 19 Refer to the wireless remote user's manual (P/N # 699909945) for operational instructions.
- 20 Agilent recommends a full calibration of the unit prior to leak test operations.

Installation on VS Series Leak Detector (G8600-68001)

Equipment Required	Kit Includes
Extended Length M5 Allen Wrench	Wireless Base PCB
Metric Allen Wrench Set	Wireless Remote Label
M3 Philips Head Screw Driver	RS232 Cable
Adjustable Wrench	Antenna
	M3 Screw (Qty. 4)

Installation Procedure

For clarity, some items have been omitted from views.

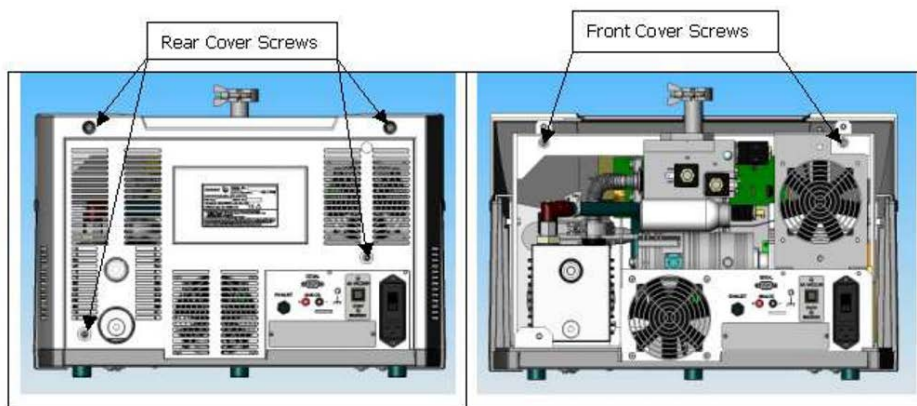


Figure 10 Rear and Front Cover Screws

WARNING



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

NOTE

Prior to wireless PCB installation, ensure that the software is revision 4.02 or higher by navigating from the Home screen menu to System Information. If not, contact Agilent Customer Service. See the back cover of this manual for a listing of our sales and service offices.

- 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 10: Rear and Front Cover Screws) and detach the rear plastic cover from the unit.
- 4 Remove the four screws holding the front plastic cover and detach the cover from the unit. Two screws are situated at the front of the unit (not shown) and two screws are positioned inside the unit (Figure 10: Rear and Front Cover Screws).

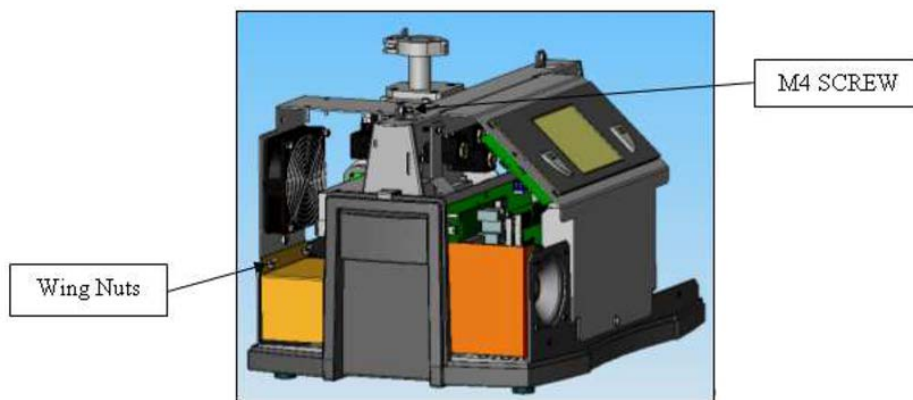


Figure 11 Fan Assembly and Socket Head Cap Screw

- 5 Disconnect the fan connector from the 24 V cable harness.
- 6 Detach the fan assembly by removing the M4 socket head cap screw and loosening the two wing nuts (Figure 11: Fan Assembly and Socket Head Cap Screw). Place the fan assembly on a flat surface on its back.

Instructions for Installation and Technical Specification

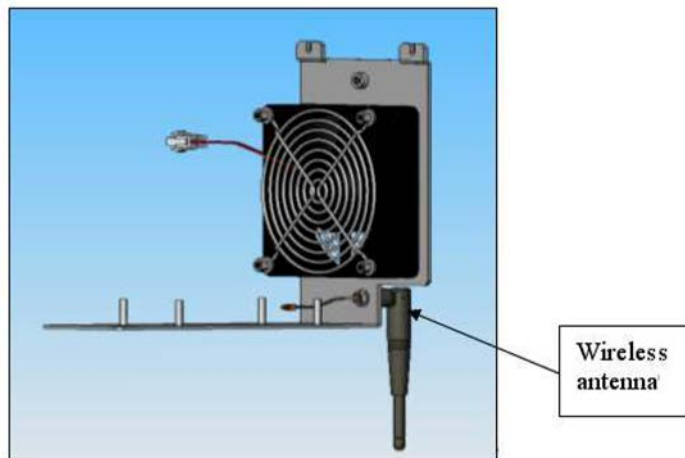


Figure 12 Wireless Antenna

- 7 Hand tighten the antenna to the fan bracket, and then apply a $\frac{1}{2}$ turn using an adjustable wrench (Figure 12: Wireless Antenna).

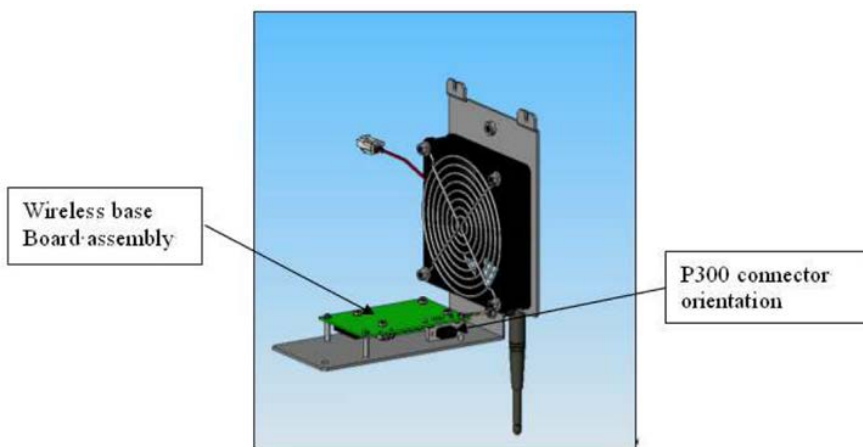


Figure 13 Base Board Assembly with Posts

Instructions for Installation and Technical Specification

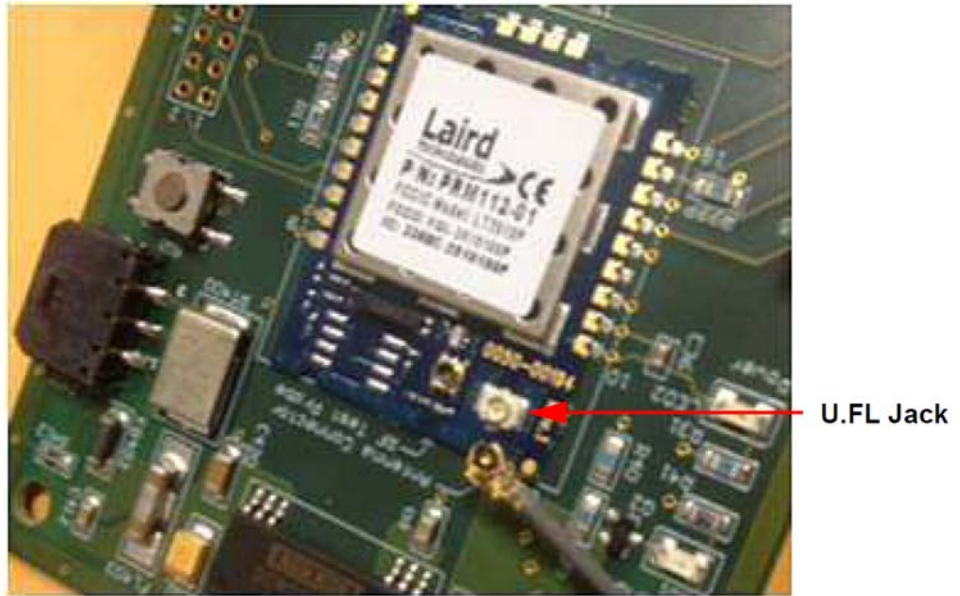


Figure 14 Antenna Connector

- 8 Place the wireless board assembly on the four fan bracket standoffs. Note the correct orientation of the wireless base board P300 connector (Figure 4: Base Board Assembly with Posts).

CAUTION

Use proper ESD safety techniques when performing a PCB card install.

Instructions for Installation and Technical Specification

- 9 Connect the antenna cable to the U.FL (Color: silver) on the wireless base board (Figure 14: Antenna Connector). Carefully center the connector on the jack and press the connector onto the jack until it snaps into place. Rotate the antenna cable to ensure the connector is firmly seated in the jack.
- 10 Secure the wireless base board to the fan bracket using the four M3 Phillips screws provided with the kit.
- 11 Unfurl the black/red wire (P1-RF) with the 24 V three-pin connector and the ground lug (Figure 15: 24 V Power Cable), which is tie rapped at the base of the leak detector, and:
 - Plug the 24 V connector labeled P1-RF into the three-pin connector on the wireless board.
 - Attach the 24 V ground lug to one of the bracket screws.
 - Reinstall the fan assembly by tightening the M4 socket head cap screw and the two wing nuts (Figure 11: Fan Assembly and Socket Head Cap Screw)

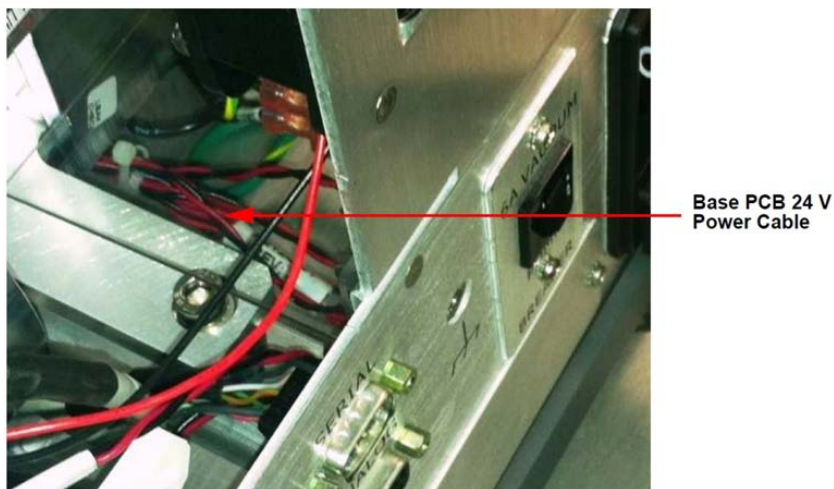


Figure 15 24 V Power Cable

Instructions for Installation and Technical Specification

- 12 Connect the RS-232 cable end labeled J300-1 to P300 on the wireless board using its two slotted screws (Figure 14: Antenna Connector).

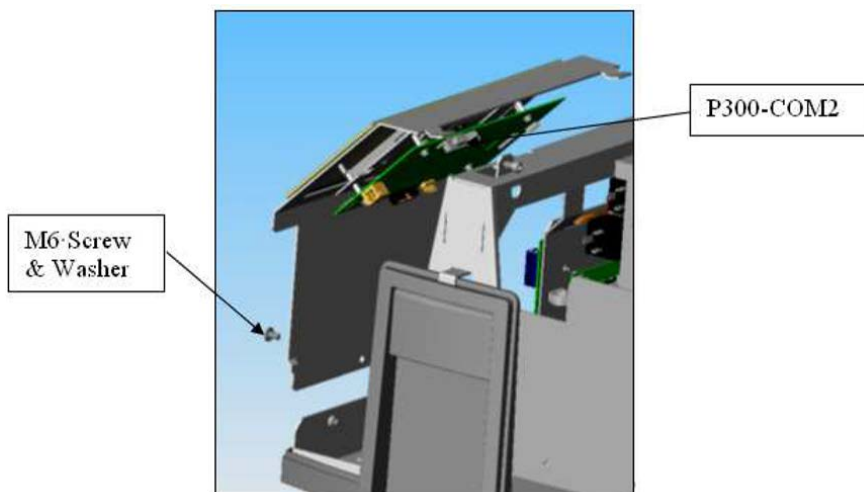


Figure 16 Front Panel Display

- 13 Feed the RS-232 the RS-232 cable through the leak detector to the front panel display.
- 14 Remove the M6 screw and washer from the front panel display and detach the display (Figure 16: Front Panel Display).

Instructions for Installation and Technical Specification

- 15 Connect the RS-232 cable end labeled J300-2 to P300-COM2 on the front panel display board using its two slotted screws (Figure 17: RS232 Cable).

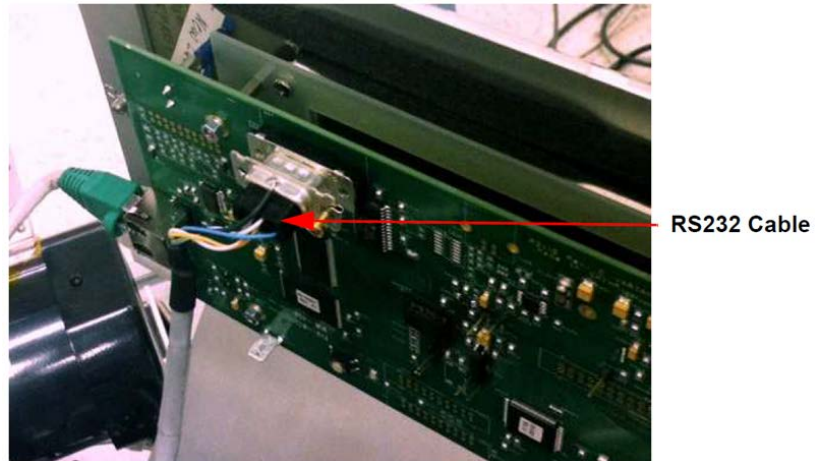


Figure 17 RS232 Cable

- 16 Reattach the display by inserting the tabs into their slotted positions and secure using the M6 screw and washer.
- 17 Attach the front cover and secure to the frame using existing hardware.
- 18 Rotate the antenna to its horizontal locked position.
- 19 Guide the antenna through its location hole in the rear cover. Then secure the rear cover to the frame using existing hardware.

Instructions for Installation and Technical Specification

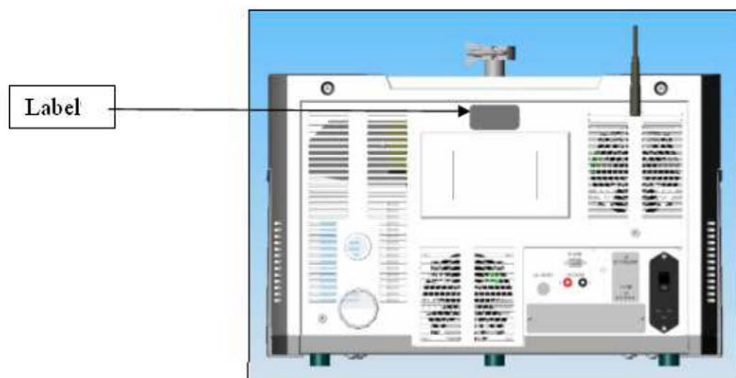


Figure 18 Label

- 20 Attach the label to the rear cover in location shown (Figure 9: Label).
- 21 Connect the power cord and power up the unit.
- 22 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.
- 23 After installation refer to the VS Series leak detector manual (P/N # 699909942) Section 3, for leak detector activation of the wireless remote.
- 24 Refer to the wireless remote user's manual (P/N # 699909945) for operational instructions.
- 25 Agilent recommends a full calibration of the unit prior to leak test operations.

Disposal

Meaning of the "WEEE" logo found in labels.

The following symbol is applied in accordance with the EC WEEE (Waste Electrical and Electronic Equipment) Directive.

This symbol (valid only in countries of the European Community) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system. The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.



Figure 19 Logo "WEEE"

For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

Service

Should a customer need an advanced exchange or repair service, please contact local distributor or directly e-mail to:

vpt-customer@agilent.com

vpl-customer@agilent.com

Completion of the "Request for Return" form is required to return your product to Agilent for service (provided at the end of this manual).

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
Toll Free: 00 800 234 234 00

vpt-customer@agilent.com

NORTH AMERICA:

Fax: 1 781 860 9252
Toll Free: 800 882 7426, Option 3

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 <input checked="" type="checkbox"/> 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"/> 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: <small>(by selecting 'YES' you MUST complete the table to the right)</small>	<small>* 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.</small>				
		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
Ph. +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426
vpl-customer-care@agilent.com

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

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 Penteadro de Ulhoa
Rodrigues, 939 - 6° andar
Castelo Branco Office Park
Torre Jacarandá - Tamboré
Barueri, Sao Paulo CEP: 06460-040
Toll free: 0800 728 1405

China

Agilent Technologies Co. Ltd (China)

No.3, Wang Jing Bei Lu, Chao Yang District
Beijing, 100102, China
Tel: +86 (0)10 64397888
Fax: +86 (0)10 64397666
Toll free: 400 8203278 (mobile)
Toll free: 800 8203278 (landline)
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

Southeast Asia

Agilent Technologies Sales Sdn Bhd

Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya, Selangor, Malaysia
Ph. +603 7712 6181
Fax: +603 7727 1239
Toll free: 1 800 880 805
vps-customerservice@agilent.com

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
Ph. +91 11 4623 7100
Fax: +91 4623 7105
Toll Free: 1 800 180 1517

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: 00 800 234 234 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

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

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
vpc-customerservice@agilent.com

UK and Ireland

Agilent Technologies LDA UK Limited

Lakeside Cheadle Royal Business Park
Cheadle, Cheshire SK8 3GR,
United Kingdom
Ph. +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

© Agilent Technologies, Inc. 2022
10/2022

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
Ph. +91 80 40614000
Fax: +91 80 40148991

Taiwan

Agilent Technologies Taiwan Limited

No. 20 Gao-shuang Road, Ping-zhen Dist
Tao-Yuan City
32450 Taiwan
Tel: +886 3 4959004
Toll free: 0800 018 768
vpw-customerservice@agilent.com

Germany and Austria

Agilent Technologies Sales & Services GmbH & Co. KG

Lyoner Str. 20
60 528 Frankfurt am Main
Germany
Tel: +49 69 6773 43 2230
Fax: +49 69 6773 43 2250

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

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 On- line:

www.agilent.com/chem/vacuum
Representatives in most countries

In This Book

The manual describes the following:

- Instructions for Installation and Technical Specifications

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024

Agilent Technologies Italia S.p.A.
Vacuum Products Division
Via F.lli Varian, 54
10040 Leini (TO)
ITALY



699910002

Edition 09/24
Issue E.00