



# Rough Pump Communications Guide (RS-232/RS-485/USB)

## Addendum

# Notices

## Manual Part Number

D0144169

Edition A.00 February 2026

## Copyright

© Agilent Technologies, Inc. 2026

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](http://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

Validity	4
Window Protocol Description	4
Communication Format	4
Communication Protocol	4
Byte Type Format:	4
Examples	8

## Validity

This addendum lists a minimum set of information necessary to understand the operations of the Rough Pump Communications.

## Window Protocol Description

### Communication Format

- 8 data bits
- no parity
- 1 stop bit
- baud rate: 9600 default, programmable from 600 to 38400 depending on model.

### Communication Protocol

The communication protocol is a MASTER/MINION type where:

- Host PC/PLC/MCU = MASTER
- Inverter, Motor, or Pump = MINION.

The communication is performed in the following way:

- 1 The MASTER sends a COMMAND to the MINION.
- 2 The MINION returns a RESPONSE to the MASTER.

### Byte Type Format:

- 1 ASCII character bytes are shown between quotes. Example: '3'  
ASCII control character bytes will be shown underlined in italics. Example: *STX*
- 2 Hex bytes are shown with the prefix 0x \_\_ . Example: 0x03

Decimal bytes are shown as numerical values with no additional formatting.  
Example: 3

**NOTE**

The Agilent Pump Serial Module is a Python-based library built for RS-232 and RS-485 serial communication and is available at Agilent.com, ready for use. This library is complete with documentation that can aid the programmer in understanding and implementing this protocol in their application.

## Sending

The COMMAND is a byte string with the following format:

<STX>+<ADDR>+<WIN>+<COM>+<DATA>+<ETX>+<CRC>

Where:

- <STX> (Start of transmission) = STX = 0x02 = 2
- <ADDR> (Unit address) = 0x80 + device number (0 to 31)  
(For RS 232 the device number is always 0, so the <ADDR> is 0x80)
- <WIN> (Window) Similar to a register or memory location.  
The three digit window number (from '000' to '999').  
Each digit encoded as an ASCII character.  
For the meaning of each window number see the specific pump manual.
- <COM> (Command type)  
To read from a window use '0' = 0x30 = 48  
To write into the window use '1' = 0x31 = 49
- <DATA>  
An ASCII string of the data to be written into the window.  
This field is not present when using a read command.

The <DATA> field length and contents vary according to the data type as shown in the following table:

**Table 1**

Data Type	Field Length (Bytes)	Valid Characters
Logic (L)	1	'0' = FALSE '1' = TRUE
Numeric (N)	6	',' ','0' ... '9' right justified with '0'
Alphanumeric (A)	10 or greater	from blank to '_' (ASCII) left justified with space " "

- <ETX> (End of transmission) = ETX = 0x03 = 3
- <CRC> (Cyclic Redundancy Check) Is an XOR of all bytes from <ADDR> to <ETX>. The XOR result is first encoded to hexadecimal, then each of the two hex digits are sent as sperate ASCII characters.

Example: XOR = 0xB3 -> CRC = "B" & "3" = 0x42 & 0x33 = 66 & 51.

(**Note:** if Hex digits A-F are present, they **must be uppercase**).

## Receiving

The addressed MINION will reply with a RESPONSE whose structure depends on the MASTER COMMAND command type (<COM>).

When the MASTER COMMAND <COM> is a read type ('0'), the MINION will reply by transmitting a RESPONSE byte string similar in structure to the COMMAND.

<STX>+<ADDR>+<WIN>+<COM>+<DATA>+<ETX>+<CRC>

Where the content and format of the <DATA> bytes depend on the type of data requested. See the table below.

When the MASTER COMMAND <COM> is a write type ('1'), or if an error is encountered, the MINION will reply with a simplified RESPONSE byte string. The <WIN> & <COM> fields are removed, and the <REPLY> field is a single byte.

<STX>+<ADDR>+<REPLY>+<ETX>+<CRC>

This simplified RESPONSE format can be recognized by its short 6 byte length.

After a successful write command, the <REPLY> is an acknowledgment = ACK = 0x06 = 6.

If the MASTER COMMAND was invalid, the MINION will provide a simplified RESPONSE, with the <REPLY> byte corresponding to the encountered error as seen in the table below.

**Table 2**

Response Type	Response Length	Response Value	Description
Logic	1 byte	-	After a read command of a logic window
Numeric	6 bytes	-	After a read command of a numeric window
Alphanumeric	10 or greater bytes	-	After a read command of an alphanumeric window
ACK	1 byte	(0x6)	The command execution has been successfully completed
NACK	1 byte	(0x15)	The command execution has been failed
Unknown Window	1 byte	(0x32)	The specified window in the command is not a valid window
Data Type Error	1 byte	(0x33)	The data type specified in the command (Logic, Numeric or Alphanumeric) is not accorded with the specified Window
Out of Range	1 byte	(0x34)	The value expressed during a write command is out of range value of the specified window
Win Disabled	1 byte	(0x35)	The specified window is Read Only or temporarily disabled (for example) you can't write the Soft Start when the Pump is running)

## Examples

### Example 1: Start the pump

Pump address: 0

Window: "000" (Start/Stop)

Command Type: Write

Data: '1' to start (Logic data type)

MESSAGE Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<WIN>1	'0'	0x30	48	00110000	10110000	
<WIN>2	'0'	0x30	48	00110000	10000000	
<WIN>3	'0'	0x30	48	00110000	10110000	
<COM>	'1'	0x31	49	00110001	10000001	
<DATA>	'1'	0x31	49	00110001	10110000	
<ETX>	<u>ETX</u>	0x03	3	00000011	10110011	0xB3
<CRC> 1	'B'	0x38	56	00111000		
<CRC> 2	'3'	0x35	53	00110101		

### Step-by-step COMMAND construction

<STX> The COMMAND begins with an ASCII "start of transmission" byte.

<STX> = STX = 0x02 = 2.

<ADDR> In this example the pump address is 0, so the <ADDR>= is 0x80+ 0.

<ADDR> = 0x80= 128

<WIN> In this example the Start/Stop window is 0. This would be found in the specific pump manual. The <WIN> field is three ASCII digits, right justified with leading zeros. In this case, <WIN> = '0', '0', '0' =0x30, 0x30, 0x30 = 48, 48, 48.

<COM> This example is writing a value to a window, so the write command

<COM> = 0x31 = "1" = 49, is used.

<DATA> The start/stop window will accept a logical '1' or '0' to start or stop the pump. Logic <DATA> is formatted as a single ASCII character byte. In this case a '1' to start the pump. <DATA> = '1' = 0x31 = 49.

<ETX> The COMMAND is ended with an ASCII "end of transmission" byte.  
 <ETX> = ETX = 0x03 = 3

<CRC> The two byte CRC is generated from the XOR of all bytes from the <ADDR> to <ETX>. The XOR is shown in the above table to the right of the divider. The XOR result is encoded to Hexadecimal, in this case 0xB3. The CRC is the ASCII character bytes for each Hexadecimal digit, in this case "B" and "3"

The resulting <CRC> is two bytes = 'B', '3' = 0x42, 0x33 = 66, 51.

### **Step by step RESPONSE deconstruction**

The pump replies with an acknowledgement in the following format.

RESPONSE Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<REPLY>	<u>ACK</u>	0x06	6	00000110	10000110	
<ETX>	<u>ETX</u>	0x03	3	00000011	10000101	0x85
<CRC> 1	'8'	0x38	56	00111000		
<CRC> 2	'5'	0x35	53	00110101		

### **Reading the serial line**

When reading the serial line, listen for the <STX> byte to begin recording the byte string. Continue recording bytes while listening for the <ETX> byte, signaling the end of the COMMAND. Then record the following two CRC bytes for comparison.

Note: If you're using a multi-address RS485 network, the <ADDR> address byte should be verified.

### Verifying the CRC

When receiving a RESPONSE It is good practice to recompute the CRC and verify it against the received CRC. In this case the XOR from <ADDR> to <EXT> is 0x85, resulting in a CRC of the two ASCII characters '8', '5'. The recomputed CRC matches the received CRC, so it is safe to interpret the reply.

### Extracting the Reply

Since the RESPONSE length is 6 bytes long, it is the simplified reply format. Reading the 3<sup>rd</sup> byte or removing the first two bytes (<STX>, <ADDR>), and the last three bytes (<ETX>, <CRC>1, <CRC>2) will isolate the reply. In this case the <REPLY> is an acknowledgement = ACK = 0x06 = 6.

### **Example 2: Set Pump Speed**

Pump address: 0  
 Window: "120" (Pump Speed (Hz))  
 Command Type: Write  
 Data: "60" (Numerical (Hz))

COMMAND Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<WIN>1	'1'	0x31	49	00110001	10110001	
<WIN>2	'2'	0x32	50	00110010	10000011	
<WIN>3	'0'	0x30	48	00110000	10110011	
<COM>	'1'	0x31	49	00110001	10000010	
<DATA>1	'0'	0x30	48	00110000	10110010	
<DATA>2	'0'	0x30	48	00110000	10000010	
<DATA>3	'0'	0x30	48	00110000	10110010	
<DATA>4	'0'	0x30	48	00110000	10000010	
<DATA>5	'6'	0x36	54	00110110	10110100	
<DATA>6	'0'	0x30	48	00110000	10000100	
<ETX>	<u>ETX</u>	0x03	3	00000011	10000111	0x87
<CRC> 1	'8'	0x38	56	00111000		
<CRC> 2	'7'	0x37	55	00110111		

### Step by step COMMAND construction

<STX> The COMMAND begins with the ASCII “start of transmission” byte.

<STX> = STX = 0x02 = 2.

<ADDR> In this example the pump address is 0, so the <ADDR> is 0x80+ 0 = 0x80.

<WIN> In this example the “target speed” window is 120. This would be found in the specific pump manual. The <WIN> field is three ASCII digits, right justified with leading zeros. In this case <WIN> = ‘1’, ‘2’, ‘0’ =0x31, 0x32, 0x30 = 49, 50, 48.

<COM> This example is writing a value to the window, so the write command

<COM> = 0x31 = “1” = 49 is used.

<DATA> Since numerical data is being sent, the field needs to be formatted to 6 bytes long, right justified, padded with ‘0’s. In this case we are sending a 60Hz command so our <DATA> byte string would be ‘000060’.

<ETX> The COMMAND is ended with the ASCII “end of transmission” byte.

<ETX> = ETX = 0x03 = 3

<CRC> The two byte CRC is generated from the XOR of all bytes from <ADDR> to <ETX>. The XOR is shown in the above table, to the right of the divider. The XOR result is encoded to Hexadecimal, in this case 0x87. The CRC is the ASCII character bytes for each Hexadecimal digit, in this case “8” and “7”

The resulting <CRC> is two bytes = ‘8’, ‘7’ = 0x38, 0x37 = 56, 55.

### Pump RESPONSE

The pump replies with an acknowledgement in the following format.

RESPONSE Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<REPLY>	<u>ACK</u>	0x06	6	00000110	10000110	
<ETX>	<u>ETX</u>	0x03	3	00000011	10000101	0x85
<CRC> 1	'8'	0x38	56	00111000		
<CRC> 2	'5'	0x35	53	00110101		

This is the same response as Example 1, see that example for the step-by-step deconstruction.

### Example 3: Read the pump's onboard pressure gauge

Pump address: 0

Window: "224" (Gauge Pressure)

Command Type: Read

Expected Data: Alphanumerical Scientific Notation ('X.XXE±XX')

COMMAND Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<WIN>1	'2'	0x32	50	00110010	10110010	
<WIN>2	'2'	0x32	50	00110010	10000000	
<WIN>3	'4'	0x34	52	00110100	10110100	
<COM>	'0'	0x30	48	00110000	10000100	
<ETX>	<u>ETX</u>	0x03	3	00000011	10000111	0x87
<CRC> 1	'8'	0x38	56	00111000		
<CRC> 2	'7'	0x35	53	00110101		

#### **Step by step COMMAND construction**

<STX> The COMMAND begins with the ASCII "start of transmission" byte.

<STX> = STX = 0x02 = 2.

<ADDR> In this example the pump address is 0, so the <ADDR> is 0x80+ 0 = 0x80

<WIN> In this example the pump pressure gauge window is 224. This would be found in the specific pump manual. The <WIN> field is three ASCII digits, right justified with leading zeros. In this case <WIN> = '2', '2', '4' =0x32, 0x32, 0x34 = 50, 50, 52.

<COM> This example is reading the data from a window, so the read command, <COM> = "0"=0x30 = 49 is used.

<DATA> Since this is a read command no <DATA> bytes are sent.

<ETX> The COMMAND is ended with the ASCII "end of transmission" byte.

<ETX> = ETX = 0x03 = 3

<CRC> The two byte CRC is generated from the XOR of all bytes from <ADDR> to <ETX>. The XOR is shown in the above table, to the right of the divider. The XOR result is encoded to Hexadecimal, in this case 0x87. The CRC is the ASCII character bytes for each Hexadecimal digit, in this case "8" and "7"

The resulting <CRC> is two bytes = '8', '7' = 0x38, 0x37 = 56, 55.

### **Step by step RESPONSE deconstruction**

The pump responds with the following RESPONSE.

RESPONSE Byte String					CRC XOR	
Byte Field	ASCII	HEX	DEC	BIN	BIN XOR	HEX XOR
<STX>	<u>STX</u>	0x02	2	00000010	Not Used	
<ADDR>	N/A	0x80	128	10000000	10000000	
<WIN>1	'2'	0x32	50	00110010	10110010	
<WIN>2	'2'	0x32	50	00110010	10000000	
<WIN>3	'4'	0x34	52	00110100	10110100	
<COM>	'0'	0x30	48	00110000	10000100	
<DATA>1	'3'	0x33	51	00110011	10110111	
<DATA>2	':'	0x2E	46	00101110	10011001	
<DATA>3	'6'	0x36	54	00110110	10101111	
<DATA>4	'5'	0x35	53	00110101	10011010	
<DATA>5	'E'	0x45	69	01000101	11011111	
<DATA>6	':'	0x2D	45	00101101	11110010	
<DATA>7	'0'	0x30	48	00110000	11000010	
<DATA>8	'3'	0x33	51	00110011	11110001	
<DATA>9	''	0x20	32	00100000	11010001	
<DATA>10	''	0x20	32	00100000	11110001	
<DATA>11	''	0x20	32	00100000	11010001	
<ETX>	<u>ETX</u>	0x03	3	00000011	11010010	0xD2
<CRC> 1	'D'	0x44	68	01000100		
<CRC> 2	'2'	0x32	50	00110010		

### **Reading the serial and CRC validation**

See Example #1.

Since this RESPONSE is greater than 6 bytes, it is the full-length standard format.

### **Confirming the Window (optional)**

In a full length RESPONSE, bytes 3, 4, & 5 should contain the same window number as the requested read window. In this case the received window '224' matches the requested window 224.

### **Extracting the Reply**

Removing the first five bytes (<STX>, <ADDR>, <WIN>1, <WIN>2, <WIN>3, <COM>), and the last three bytes (<ETX>, <CRC>1, <CRC>2) will isolate the <DATA>. Since the remaining <DATA> is 11 bytes long, it is an alphanumeric data type (>=10 bytes). The expected format is ASCII characters, left justified with spaces.

The isolated <DATA> ASCII byte string reads '3.65E-03 '

In this case it is gauge pressure in scientific notation. A separate window typically controls the reported pressure units. See the specific pump manual for more details.

### **Common Pitfalls**

- 1** No response – Verify Serial/Baud rate settings
- 2** Python's default Hex literal output is lower case. The CRC must use capital ASCII Characters for the HEX digits.  
If the CRC uses Lowercase ASCII letters, you will get a simplified RESPONSE with <REPLY> = NACK = 0x15 = 21.
- 3** Since the RESPONSE lengths vary significantly by data type, and there is no line terminator, typical serial read line functions will timeout. Seeking the *ETX* character is a more responsive way to read the serial RESPONSE.

#### **NOTE**

Python Library for WIN Protocol exists on Agilent.com.

---

## 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](mailto: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](mailto: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](mailto: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](mailto:vacuum_cnmarketing@agilent.com)  
[vpc-customerservice@agilent.com](mailto: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](mailto: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](mailto: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.Ili 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](mailto: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](mailto: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](mailto: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](mailto: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](mailto: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](mailto: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.Ili 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](mailto:vpt-customer-care@agilent.com)

### Customer Support & Service

#### **NORTH AMERICA:**

Toll Free: 800 882 7426  
[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

#### **EUROPE:**

Toll Free: 00 800 234 234 00  
[vpt-customer-care@agilent.com](mailto: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](http://www.agilent.com/chem/vacuum)  
Representatives in most countries

# In This Book

The Addendum describes the following:

- Rough Pump Communications Guide

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026

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



D0144169

Edition 02/26  
Issue A.00