



Agilent Technologies

Vacuum Products Division



Vacuum Gauge Controller

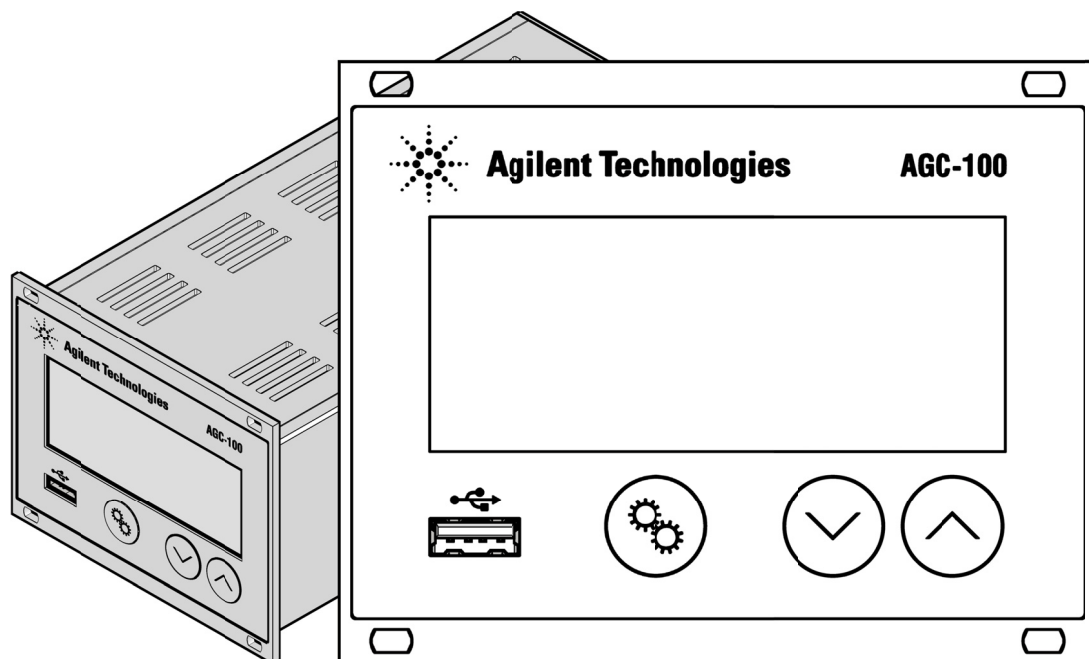
AGC-100

Operating Instructions

Manual No. tqna96e1
Revision -
February 2016

Vacuum Gauge Controller



AGC-100



Contents

Product Identification	6
Validity	6
Intended Use	7
Scope of Delivery	7
1 Safety	8
1.1 Symbols Used	8
1.2 Personnel Qualifications	8
1.3 General Safety Instructions	9
1.4 Liability and Warranty	9
2 Technical Data	10
3 Installation	13
3.1 Installation, Setup	13
3.1.1 Rack Installation	13
3.1.2 Installation in a control panel	16
3.1.3 Use as Desk-Top Unit	17
3.2 Mains Power Connector	17
3.3 Gauge Connector <i>CH 1</i>	18
3.4 <i>CONTROL</i> Connector	20
3.5 Interface Connector USB Type B	21
3.6 Interface Connector USB Type A	21
3.7 Interface Connector Ethernet	21
4 Operation	22
4.1 Front panel	22
4.2 Turning the AGC-100 On and Off	23
4.3 Operating Modes	23
4.4 Measurement Mode	25
4.5 Parameter Mode	27
4.5.1 Switching Function Parameters	28
4.5.2 Gauge parameters	32
4.5.3 Gauge Control	40
4.5.4 General Parameters	42
4.5.5 Test Parameters	47
4.5.6 Data Logger Mode	50
4.5.7 Parameter Transfer Mode	52
5 Communication Protocol (Serial Interface)	54
5.1 Data Transmission	54
5.2 Communication Protocol	55
5.3 Mnemonics	56
5.4 Measurement Mode	58
5.5 Switching Function Parameters	60
5.6 Gauge Parameters	61
5.7 Gauge Control	66
5.8 General Parameters	67
5.9 Data Logger Parameters	72
5.10 Parameter Transfer	73
5.11 Test Parameters	73
5.12 Further	77
5.13 Example	78
6 Maintenance	79
7 Troubleshooting	80
8 Repair	81
9 Accessories	81
10 Storage	81
11 Disposal	81

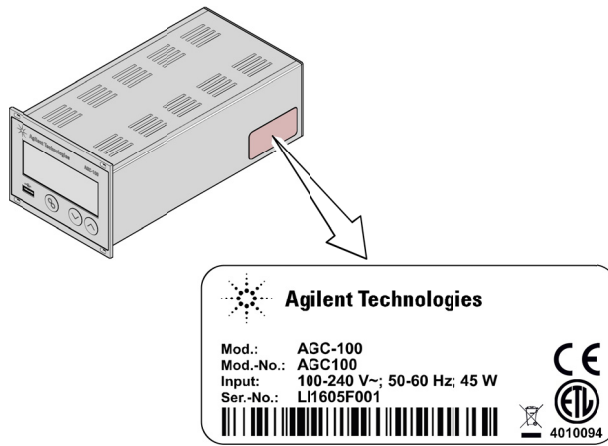
Appendix	82
A: ConversionTables	82
B: Firmware Update	83
C: Ethernet Configuration	86
C 1: Connect the AGC-100 to a Network	86
C 2: Connect the AGC-100 to a Computer	87
C 3: Ethernet Configuration Tool	87
D: Literature	90
ETL Certification	91

For cross-references within this document, the symbol (→  XY) is used, for cross-references to further documents and data sources, the symbol (→  [Z]).

Vacuum Gauge Controller AGC-100

Product Identification

In all communications with Agilent Technologies, please specify the information on the product nameplate:



Specimen nameplate

Validity

This document applies to products with part number AGC100.
The part number (Mod.-No.) can be taken from the product nameplate.

This manual is based on firmware version V1.02.
If your unit does not work as described in this document, please check that it is equipped with the above firmware version (→ 47).

We reserve the right to make technical changes without prior notice.

All dimensions are indicated in mm.

Vacuum Gauge Controller AGC-100

Intended Use

The Vacuum Gauge Controller AGC-100 is used together with Agilent gauges for total pressure measurement. All products must be operated in accordance with their respective operating manuals.

Scope of Delivery

The scope of delivery consists of the following parts:

- 1× Vacuum Gauge Controller
- 1× Power cord
- 1× Rubber bar
- 2× Rubber feet
- 4× Collar screws
- 4× Plastic sleeves
- 1× CD-ROM (manuals, tools, ...)
- 1× EU Declaration of Conformity
- 1× Installation Manual

1 Safety

1.1 Symbols Used


Symbols for residual risks

 **DANGER**

Information on preventing any kind of physical injury.






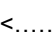
 **WARNING**

Information on preventing extensive equipment and environmental damage.


 **Caution**

Information on correct handling or use. Disregard can lead to malfunctions or minor equipment damage.

Further symbols

-  The lamp / display is lit.
-  The lamp / display flashes.
-  The lamp / display is off.
-  Press the key (example: parameter key).
-  Do not press any key.
-  Labeling


1.2 Personnel Qualifications


 **Skilled personnel**

All work described in this document may only be carried out by persons who have suitable technical training and the necessary experience or who have been instructed by the end-user of the product.

1.3 General Safety Instructions

Adhere to the applicable regulations and take the necessary precautions for all work you are going to do and consider the safety instructions in this document.



**DANGER**

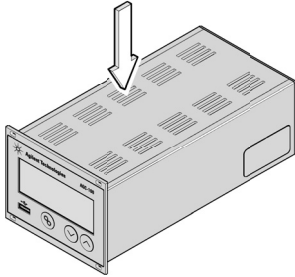


DANGER: mains voltage

Contact with live parts is extremely hazardous when any objects are introduced or any liquids penetrate into the unit.

Make sure no objects enter through the louvers and no liquids penetrate into the equipment.

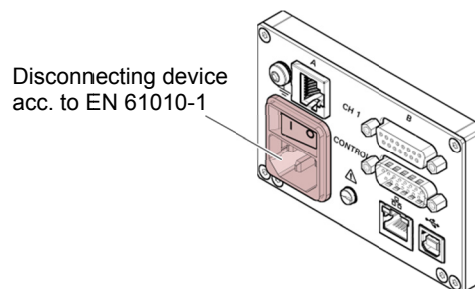




Disconnecting device

The disconnecting device must be readily identifiable by and easily reached by the user.

To disconnect the unit from the mains supply, you must unplug the mains cable.



Communicate the safety instructions to all other users.

1.4 Liability and Warranty






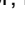
AgilentTechnologies assumes no liability and the warranty is rendered null and void if the end-user or third parties

- disregard the information in this document
- use the product in a non-conforming manner
- make any kind of interventions (modifications, alterations etc.) on the product
- use the product with accessories not listed in the corresponding product documentation.

2 Technical Data

Mains specifications	Voltage	100 ... 240 V (ac) $\pm 10\%$
	Frequency	50 ... 60 Hz
	Power consumption	≤ 45 W
	Overvoltage category	II
	Protection class	1
	Connection	European appliance connector IEC 320 C14
Ambience	Temperature	
	Storage	$-20 \dots +60$ °C
	Operation	$+ 5 \dots +50$ °C
	Relative humidity	$\leq 80\%$ up to $+31$ °C, decreasing to 50% at $+40$ °C
	Use	indoors only max. altitude 2000 m NN
	Pollution degree	II
Gauge connections	Degree of protection	IP30
	Number	1
	Gauge connections per channel	RJ45 (FCC68), 8-pin (\rightarrow 19) D-Sub, 15-pin, female (\rightarrow 19) (connected in parallel)
	Compatible gauges	
	Pirani	PVG-500, PVG-502, PVG-550, PVG-552
	Pirani / Capacitance	PCG-750, PCG 752
	Cold cathode	IMG-500
	Cold cathode / Pirani	FRG-700, FRG-702
	Hot ionization / Pirani	FRG-720, FRG-730
	Capacitance	CDG-500
Gauge supply	Voltage	$+24$ V (dc) $\pm 5\%$
	Ripple	$< \pm 1\%$
	Current	0 ... 1 A (per channel)
	Power	25 W (per channel)
	Fuse protection	1.5 A (per channel) with PTC element, self-resetting after turning the unit off or disconnecting the gauge. The supply conforms to the grounded protective extra low voltage requirements.
Operation	Front panel	via 3 keys
	Remote control	via USB type B interface via Ethernet interface

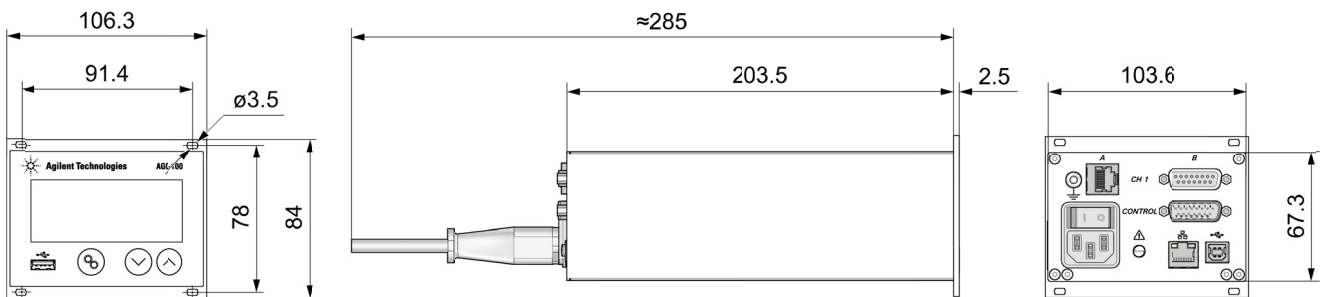
Vacuum Gauge Controller AGC-100

Measurement values	Measurement ranges	depending on gauges (→  [1] ... [8])
	Measurement error analog	
	Gain error	≤0.01% F.S. (typical) ≤0.10% F.S. (over temperature range, time)
	Offset error	≤0.01% F.S. (typical) ≤0.10% F.S. (over temperature range, time)
	Measurement rate analog	≥100 / s
	Display rate	≥10 / s
	Filter time constant	
	Slow	8 s ($f_g = 0.02$ Hz)
	Normal	800 ms ($f_g = 0.2$ Hz)
	Fast	160 ms ($f_g = 1$ Hz)
	Measurement units	mBar, hPa, Torr, Pa, Micron, V
	Offset correction	for linear gauges
	Calibration factor	0.10 ... 10.00
	A/D converter	resolution 0.001% F.S. (the measurement values of FRG-720, FRG-730, CDG-500 are transmitted digitally)
Switching functions	Number	2
	Reaction delay	≤10 ms, if switching threshold close to measurement value (for larger differences consider filter time constant)
	Adjustment range	depending on gauge (→  30, 31)
	Hysteresis	≥1% F.S. for linear gauges, ≥10% of measurement value for logarithmic gauges
Switching function relays	Contact type	floating changeover contact
	Load max.	60 V(dc), 30 W (ohmic) 30 V(ac), 1 A (ohmic)
	Service life	
	Mechanical	1×10 ⁸ cycles
	Electrical	1×10 ⁵ cycles (at max. load)
	Contact positions	→  20)
Error signal	Connector (<i>CONTROL</i>)	D-Sub appliance connector, male, 15-pin (pin assignment →  20)
Error signal	Number	1
	Reaction time	≤10 ms
Error signal relay	Contact type	floating normally open contact
	Load max.	60 V(dc), 0.5 A, 30 W (ohmic) 30 V(ac), 1 A (ohmic)
	Service life	
	Mechanical	1×10 ⁸ cycles
	Electrical	1×10 ⁵ cycles (at max. load)
	Contact positions	→  20
Error signal relay	Connector (<i>CONTROL</i>)	D-Sub appliance connector, male, 15-pin (pin assignment →  20)

Vacuum Gauge Controller AGC-100

Analog outputs	Number	1
	Voltage range	−5 ... +14.5 V (dc) If no gauge is connected, +14.5 V (dc) is output
	Deviation from display value	±20 mV
	Output resistance	<50 Ω
	Measuring signal vs. pressure	depending on gauge (→ [1] ... [8])
	CONTROL connector	D-Sub appliance connector, male, 15-pin (pin assignment → [20])
USB Type A interface	Protocol	FAT file system file handling in ASCII format
USB Type B interface	Protocol	ACK/NAK, ASCII with 3-character mnemonics
	Data format	bi-directional data flow, 1 start bit, 8 data bits, 1 stop bit, no parity bit, no handshake
	Transmission rate	9600, 19200, 38400, 57600, 115200
Ethernet interface	Protocol	ACK/NAK, ASCII with 3-character mnemonics
	Data format	bi-directional, 1 start bit, 8 data bits, 1 stop bit, no parity bit, no handshake
	Transmission rate	9600, 19200, 38400, 57600, 115200
	IP Address	DHCP (default) or manual setting (→ [86])
	MAC Address	readable via "MAC" parameter


Dimensions [mm]



Use For incorporation into a rack or control panel or as a desk-top unit

Weight AGC-100 0.85 kg


3 Installation



Skilled personnel

The unit may only be installed by persons who have suitable technical training and the necessary experience or who have been instructed by the end-user of the product.

The unit is suited for incorporation into a 19" rack or a control panel or for use as a desk-top unit.




STOP DANGER

Putting a product which is visibly damaged into operation can be extremely hazardous. If the product is visibly damaged do not put it into operation and make sure it is not inadvertently put into operation.

3.1 Installation, Setup

3.1.1 Rack Installation

The unit is designed for installation into a 19" rack chassis adapter according to DIN 41 494. For this purpose, four collar screws and plastic sleeves are supplied with it.



STOP DANGER

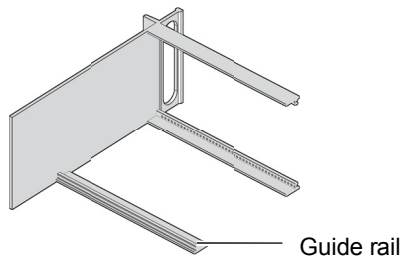
DANGER: protection class of the rack

If the product is installed in a rack, it is likely to lower the protection class of the rack (protection against foreign bodies and water) e.g. according to the EN 60204-1 regulations for switching cabinets.

Take appropriate measures for the rack to meet the specifications of the protection class.

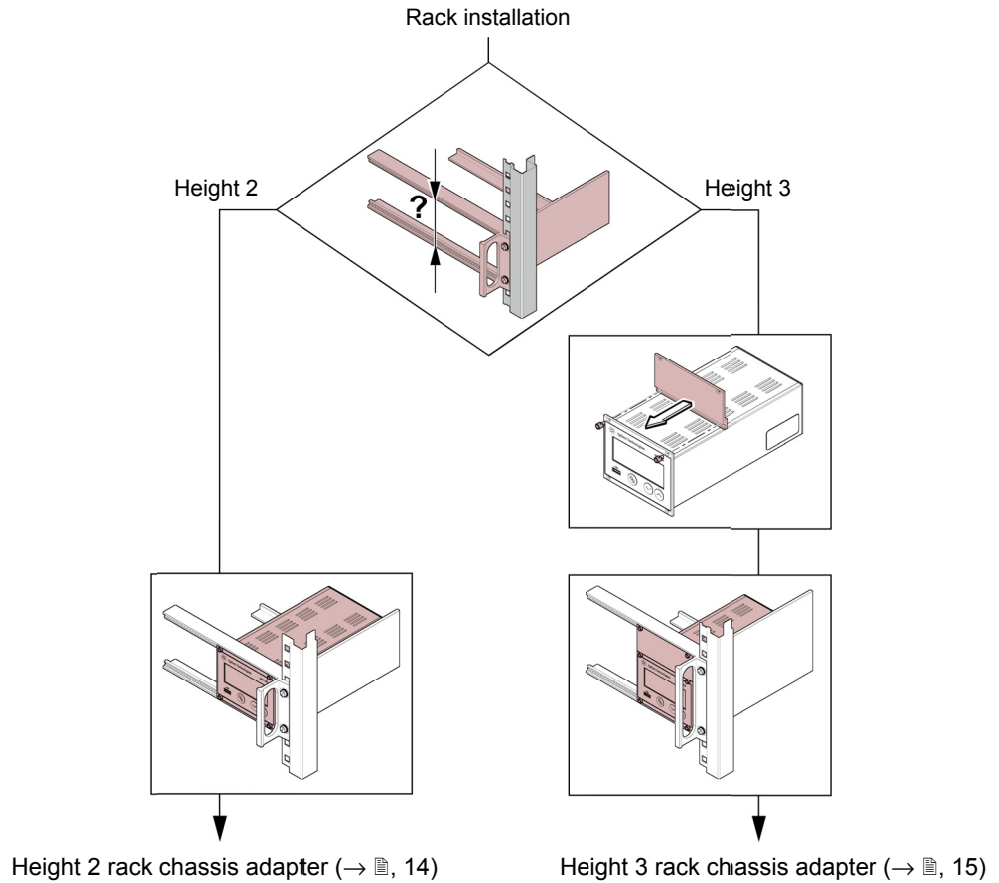
Guide rail

In order to reduce the mechanical strain on the front panel of the AGC-100, preferably equip the rack chassis adapter with a guide rail.



Vacuum Gauge Controller AGC-100

Mounting height

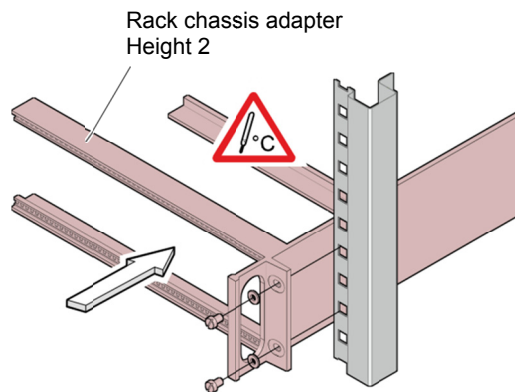


Height 2 rack chassis adapter

- 1 Secure the rack chassis adapter in the rack frame.

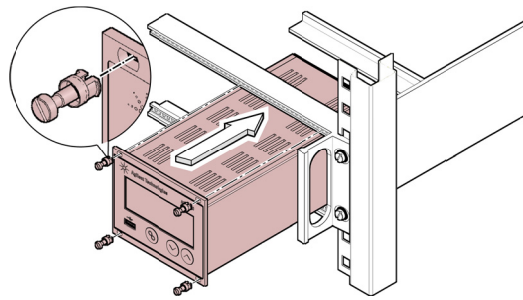


The maximum admissible ambient temperature (→ 10) must not be exceeded and the air circulation must not be obstructed.



Vacuum Gauge Controller AGC-100

- 2 Slide the AGC-100 into the adapter ...



... and fasten the AGC-100 to the rack chassis adapter using the screws supplied with it.

Height 3 rack chassis adapter

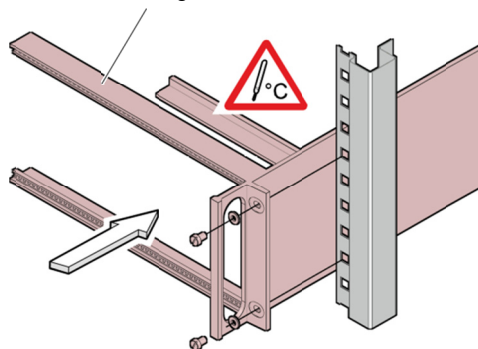
For incorporation into a 19" rack chassis adapter, height 3, an adapter panel (incl. two collar screws and plastic sleeves) is available (Accessories → 81).

- 1 Secure the rack adapter in the rack frame.



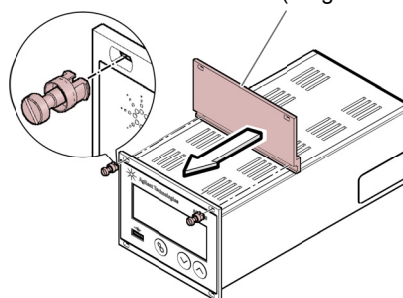
The maximum admissible ambient temperature (→ 10) must not be exceeded and the air circulation must not be obstructed.

Rack chassis adapter
Height 3



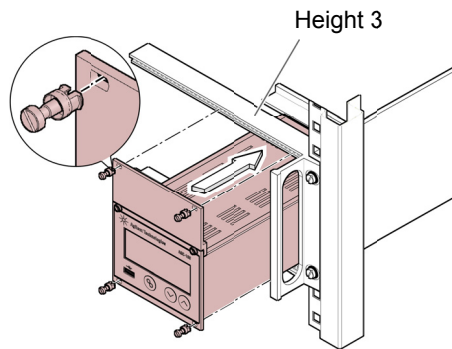
- 2 Mount the adapter panel as upper extension to the front panel of the AGC-100 using the screws supplied with the adapter panel.

Adapter panel
(Height 2 to height 3)



Vacuum Gauge Controller AGC-100

- Slide the AGC-100 into the rack chassis adapter ...



... and fasten the adapter panel to the rack chassis adapter using the screws supplied with the AGC-100.

3.1.2 Installation in a control panel



DANGER

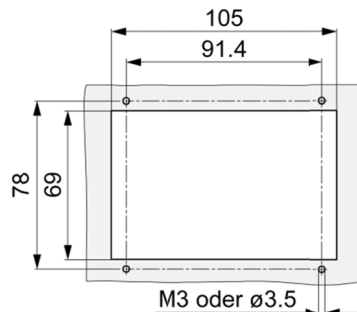


DANGER: protection class of the rack

If the product is installed in a rack, it is likely to lower the protection class of the rack (protection against foreign bodies and water) e.g. according to the EN 60204-1 regulations for switching cabinets.

Take appropriate measures for the rack to meet the specifications of the protection class.

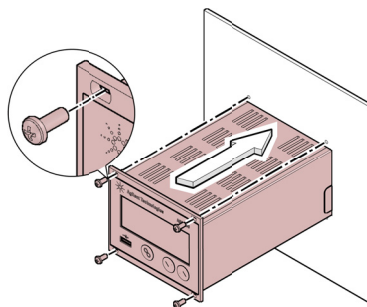
For mounting the AGC-100 into a control panel, the following cut-out is required:



The maximum admissible ambient temperature (→ 10) must not be exceeded and the air circulation must not be obstructed.

For reducing the mechanical strain on the front panel of the AGC-100, preferably support the unit.

- Slide the AGC-100 into the cut-out of the control panel ...



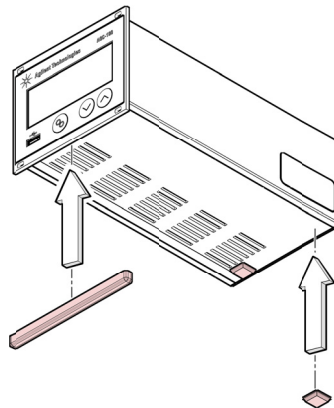
... and secure it with four M3 or equivalent screws.

Vacuum Gauge Controller AGC-100

3.1.3 Use as Desk-Top Unit

The AGC-100 may also be used as a desk-top unit. For this purpose, two self-adhesive rubber feet and a slip-on rubber bar are supplied with it.

- 1 Stick the two supplied rubber feet to the rear part of the bottom plate ...



Select a location where the admissible maximum ambient temperature is not exceeded (e.g. due to sun irradiation) (→ 10).

... and slip the supplied rubber bar onto the bottom edge of the front panel.

3.2 Mains Power Connector



DANGER



DANGER: line voltage

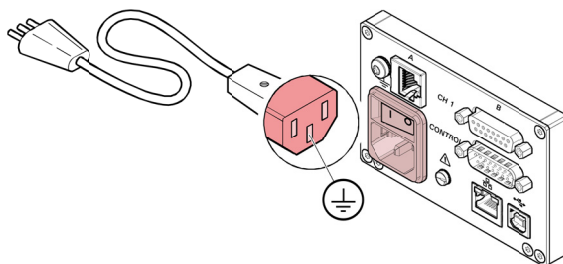
Incorrectly grounded products can be extremely hazardous in the event of a fault.

Use only a 3-conductor power cable with protective ground. The mains power connector may only be plugged into a socket with a protective ground. The protection must not be nullified by an extension cable without protective ground.

The unit is supplied with a power cord. If the mains connector is not compatible with your system, use your own, suitable cable with protective ground ($3 \times 1.5 \text{ mm}^3$).



The socket must be fuse-protected with $10 \text{ A}_{\text{max}}$



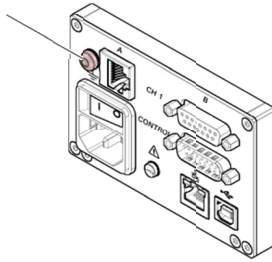
If the unit is installed in a switching cabinet, the mains voltage should be supplied and turned on via a central distributor.

Vacuum Gauge Controller AGC-100

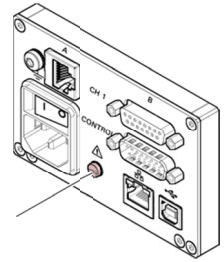
Ground Connection

On the rear of the unit is a screw enabling the AGC-100 where necessary to be connected via a ground conductor, e.g. with the protective ground of the pump stand.

Protective
grounding




Do not unfasten
this screw
(internal ground
protection)



3.3 Gauge Connector *CH 1*

For each channel there are two connections available which are connected in parallel:

- one RJ45 appliance connector, female, 8-pin (CH A)
- one D-Sub appliance connector, female, 15-pin (CH B)



Connect the gauge to the *CH 1* connector via a sensor cable set available from us (→ sales literature) or your own, screened (electromagnetic compatibility) sensor cable. Use compatible gauges (→ 10).

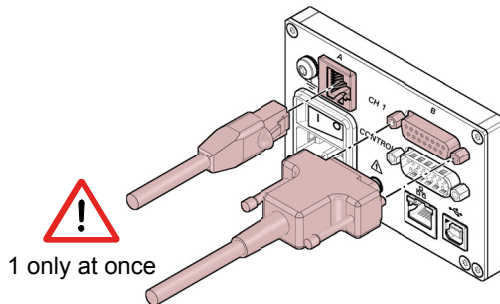


Caution



Caution: Multiple connection

Only one sensor may be connected to each of the channels (connection CH A or CH B). Otherwise the connected sensors may be damaged.



DANGER



DANGER: Hazardous voltage

According to EN 61010, voltages exceeding 30 V (ac) or 60 V (dc) are hazardous.

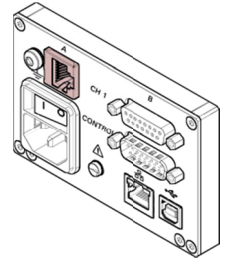
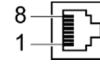
Only connect a protective low voltage (PELV).

Vacuum Gauge Controller AGC-100

Pin assignment *CH 1*

Appliance socket RJ45

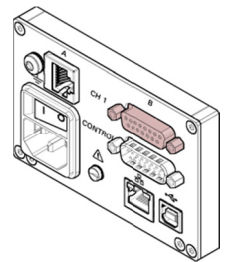
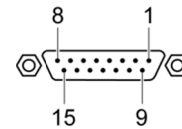
Pin assignment of the female 8-pin RJ45 appliance connectors:



Pin	Signal
1	Supply +24 V (dc)
2	Supply common GND
3	Signal input (measuring signal 0 ... +10 V (dc))
4	Identification
5	Signal common
6	Status
7	HV_L
8	HV_H / HV_EMI

Appliance socket D-Sub

Pin assignment of the female 15-pin D-Sub appliance connectors:



Pin	Signal
1	EMI status
2	Signal input (measuring signal 0 ... +10 V (dc))
3	Status
4	HV_H / HV_EMI
5	Supply common GND
6	n.c.
7	Degas
8	Supply +24 V (dc)
9	n.c.
10	Identification
11	Supply +24 V (dc)
12	Signal common
13	RxD
14	TxD
15	Chassis

3.4 CONTROL Connector

This connector allows the user to read the measuring signal, evaluate the state of the floating contacts of the error relay.



Connect the peripheral components to the **CONTROL** connector on the rear of the unit using your own, screened (electromagnetic compatibility) cable.



DANGER



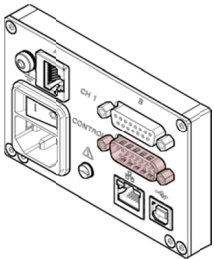
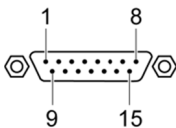
DANGER: Hazardous voltage

According to EN 61010, voltages exceeding 30 V (ac) or 60 V (dc) are hazardous.

Only connect a protective low voltage (PELV).

Pin assignment

Pin assignment of the male 15-pin D-Sub appliance connector:



Pin	Signal
1	Analog output $-5 \dots +13 \text{ V (dc)}$
2	Analog output GND
Switching function 1	
3	Pressure above threshold or power supply turned off
4	
5	
6	HV_H on +24 V off 0 V
7	+24 V (dc), 200 mA
8	Chassis = GND
Fuse-protected at 300 mA with PTC element, self-resetting after power off or pulling the CONTROL connector. Meets the requirements of a grounded protective extra low voltage.	
Error signal	
9	No error
10	
11	
Switching function 2	
12	Pressure above threshold or power supply turned off
13	
14	
15	Chassis = GND

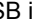


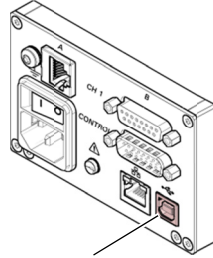
The analog output (pin 1) differs from the displayed value by no more than $\pm 20 \text{ mV}$.

3.5 Interface Connector USB Type B

The USB Type B interface connector facilitates direct communication with the AGC-100 via a computer (e.g. firmware update, parameter saving (read/write)).



Connect the USB interface connector to the  connector on the rear of the unit using a screened (electromagnetic compatibility) cable.




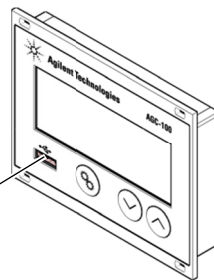
USB Type B

3.6 Interface Connector USB Type A

The USB Type A interface connector with master functionality is situated on the front of the unit and is used for the connection of a USB memory stick (e.g. firmware update, parameter saving (read/write), data logger).



Connect the USB memory stick to the connector  on the front of the unit.

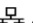


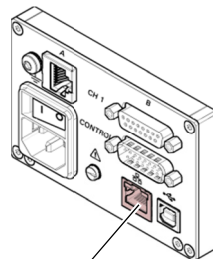
USB Type A

3.7 Interface Connector Ethernet

The Ethernet interface allows direct communication with the AGC-100 via a computer.



Connect the Ethernet cable to the connector  on the rear of the unit.



Ethernet

Green LED

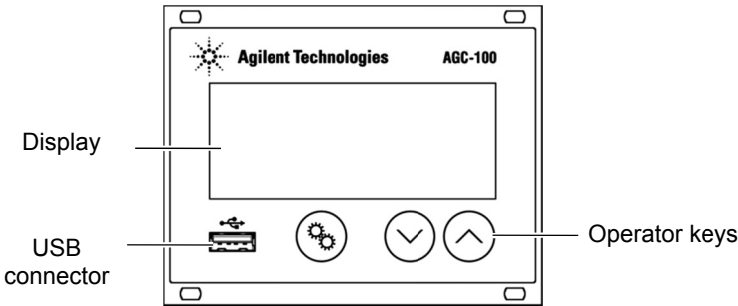
Link or transmit LED. Indicates that a hardware connection has been established.

Yellow LED

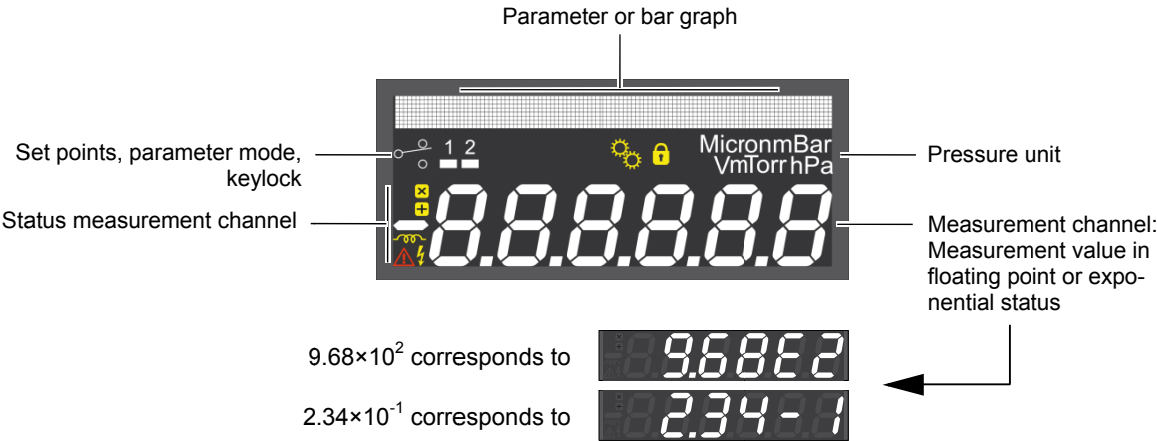
Status or packet detect LED. Indicates the status of the transmission. When this LED flashes or flickers, data are being transmitted.

4 Operation

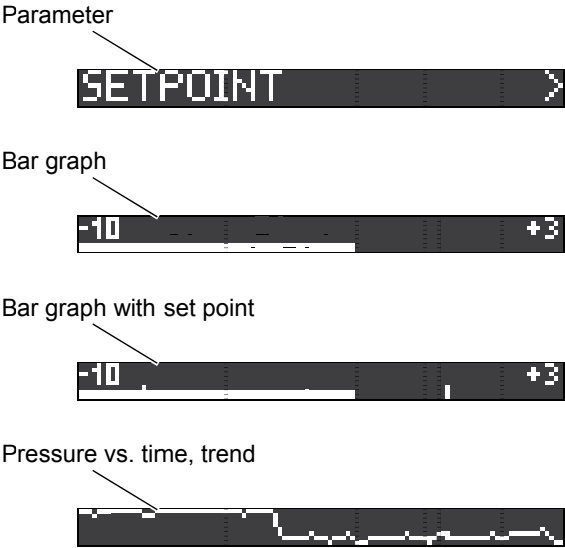
4.1 Front panel



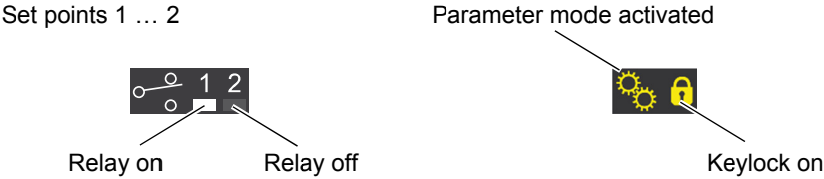
Display



Parameter, bar graph

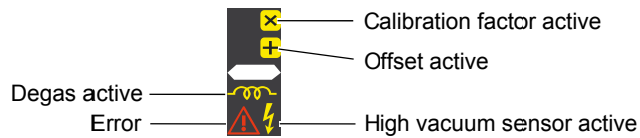


Set points, parameter mode, keylock



Vacuum Gauge Controller AGC-100

Measurement channel
specifically

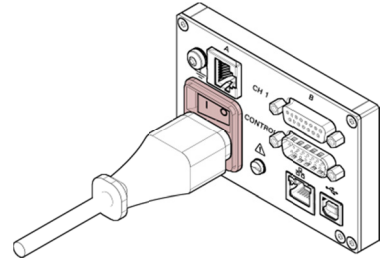


4.2 Turning the AGC-100 On and Off

Turning the AGC-100 on

The power switch is on the rear of the unit.

Turn the AGC-100 on with the power switch (or centrally, via a switched power distributor, if the unit is incorporated in a rack).



After power on, the AGC-100...

- automatically performs a self-test
- identifies the connected gauges
- activates the parameters that were in effect before the last power off
- switches to the Measurement mode
- adapts the parameters if required (if a different gauge was previously connected).

Turning the AGC-100 off

Turn the AGC-100 off with the power switch (or centrally, via a switched power distributor, if the unit is incorporated in a rack).



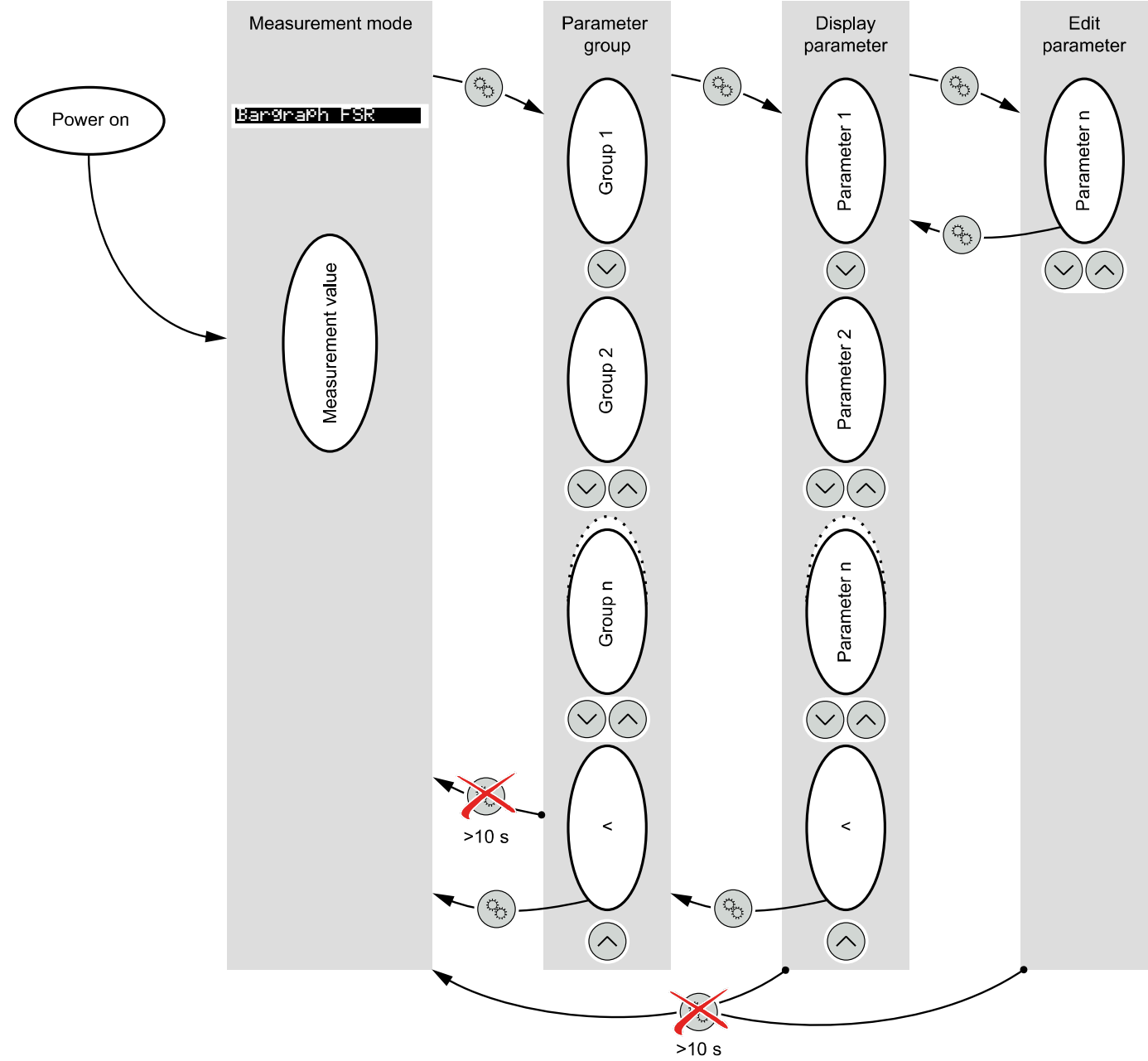
Wait at least 10 s before turning the AGC-100 on again in order for it to correctly initialize itself.

4.3 Operating Modes

The AGC-100 works in the following operating modes:

- Measurement mode
for displaying measurement values or statuses (→ 25)
- Parameter mode
for displaying and editing parameters (→ 27)
 - Switching function parameter group **SETPOINT** >
for entering and displaying thresholds (→ 28)
 - Gauge parameter group **SENSOR** >
for entering and displaying gauge parameters (→ 32)
 - Gauge control group **SENSOR-CONTROL** >
for entering and displaying gauge control parameters (→ 40)
 - General parameter group **GENERAL** >
for entering and displaying general parameters (→ 42)
 - Test program group **TEST** >
for running internal test programs (→ 47)
 - Data logger mode **DATA LOGGER** >
for logging measurement data (→ 50)
 - Parameter transfer mode **SETUP** >
for saving (read/write) parameters (→ 52)

Vacuum Gauge Controller AGC-100



4.4 Measurement Mode

Measurement mode is the standard operating mode of the AGC-100 with display of

- a bar graph (if required)
- a measurement value for each measurement channel
- status messages for each measurement channel

Adjusting bar graph

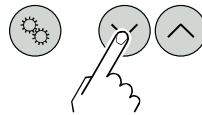
If required a bar graph may be displayed (→ 44).

Turning the gauge on/off

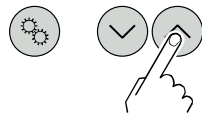
Certain gauges can be turned on and off manually, provided the gauge control is set to **S-ON HAND** (→ 41).

Available for the following gauges:

- | | |
|--|--------------------|
| <input type="checkbox"/> Pirani | (PVG) |
| <input type="checkbox"/> Pirani / Capacitance | (PCG) |
| <input checked="" type="checkbox"/> Cold cathode | (IMG) |
| <input type="checkbox"/> Cold cathode / Pirani | (FRG-70x) |
| <input type="checkbox"/> Hot ionization / Pirani | (FRG-720, FRG-730) |
| <input type="checkbox"/> Capacitance | (CDG) |



⇒ Press key for >1 s:
Gauge switches off. Instead of a measurement value the word OFF is displayed.



⇒ Press key for >1 s:
Gauges switches on. Instead of the measurement value a status message may be displayed.

Switching the emission on / off

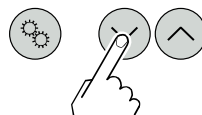
For certain gauges the emission can be switched on and off manually, provided the sensor parameter is set to **EMISSION HAND** (→ 39).



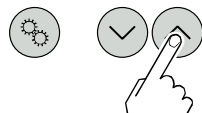
Switching on the emission is only possible if the pressure is below 2.4×10^{-2} mbar.


Available for the following gauges:

- | | |
|---|----------------|
| <input type="checkbox"/> Pirani | (PVG) |
| <input type="checkbox"/> Pirani / Capacitance | (PCG) |
| <input type="checkbox"/> Cold cathode | (IMG) |
| <input type="checkbox"/> Cold cathode / Pirani | (FRG-70x) |
| <input checked="" type="checkbox"/> Hot ionization / Pirani | (FRG-730 only) |
| <input type="checkbox"/> Capacitance | (CDG) |



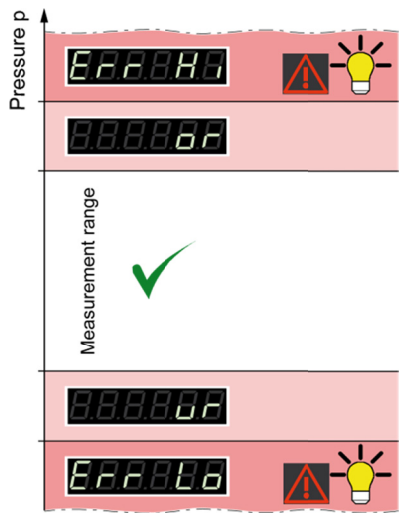
⇒ Press key for >1 s:
The emission is switched off. The measurement value of the Pirani sensor is displayed instead of the measurement value of the hot cathode ionization sensor.



⇒ Press key for >1 s:
The emission is switched on. The measurement value of the hot cathode ionization sensor is displayed and  lid solid.

Vacuum Gauge Controller AGC-100

Measurement range

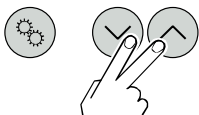


If the unit is operated with linear gauges (CDG), negative pressures may be indicated.

Possible causes:

- negative drift
- activated offset correction.

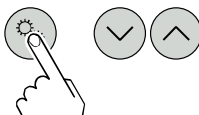
Displaying the gauge identification



⇒ Press keys for >0.5 ... 1 s:
For the measurement channel in question the type of the connected gauge is automatically identified and displayed for 6 s:

Pirani gauge (PVG-500, PVG502, PVG-550, PVG552)	PVG5xx	
Pirani / Capacitance gauge (PCG-750, PCG-752)	PCG75x	
Cold cathode gauge (IMG-500)	IMG500	
Cold cathode / Pirani gauge (FRG-700, FRG-702)	FRG70x	
Hot ionization / Pirani gauge (FRG-720)	FRG720	
(FRG-730)	FRG730	
Linear gauge (capacitance, digital) (CDG-500)	CDG500 Ux.xx	Version during 3 s, then
	CDG500 1000MBAR	FSR during 3 s
No gauge connected	noSENSOR	
Gauge connected, but not identifiable	noIDENT.	

Changing to the Parameter mode



→ 27

4.5 Parameter Mode

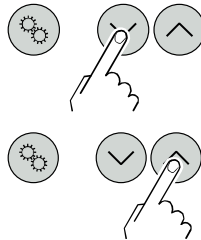
The Parameter mode is used for displaying, editing and entering parameter values as well as for testing the AGC-100 and for saving measurement data. For ease of operation the individual parameters are divided into groups.



Unit switches from measurement mode to parameter mode. The respective parameter group is displayed in place of the bar graph.



Selecting a parameter group



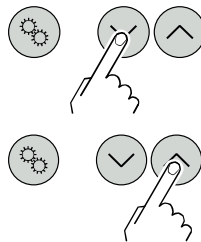
Select group

- ⇒ Switching function parameters → 28
- Gauge parameters → 32
- Gauge control → 40
- General parameters → 42
- Test parameters → 47
- Data logger → 50
- Parameter transfer → 52

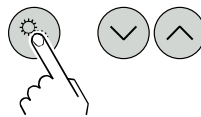


Confirm group

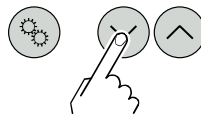
Reading a parameter in a parameter group



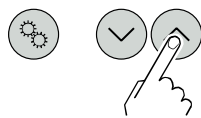
Editing and saving a parameter in a parameter group



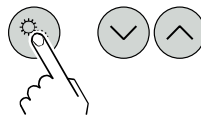
Confirm the parameter. The value flashes and can now be edited.



Edit the value.



Save the change and return to read mode.




4.5.1 Switching Function Parameters

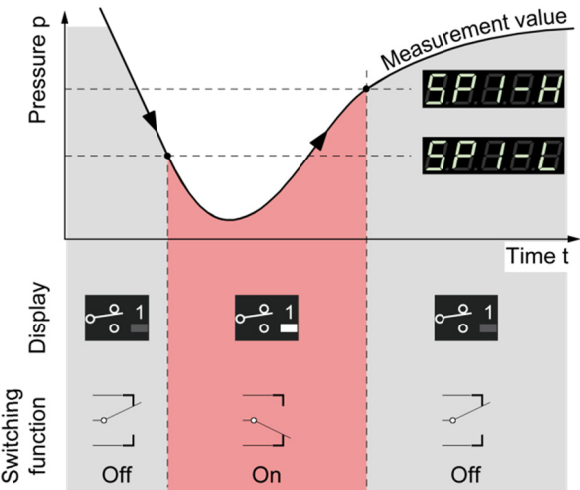
SETPOINT 

The switching function parameter group is used for displaying, editing and entering threshold values and assigning the two switching functions to a measurement channel.

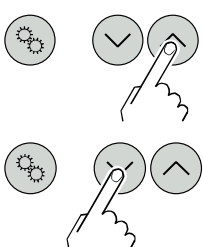
Parameters in this group

SP1-CH	Configuration of switching function 1
SP1-L	Switching function 1 lower threshold
SP1-H	Switching function 1 upper threshold
SP2-CH	Configuration of switching function 2
SP2-L	Switching function 2 lower threshold
SP2-H	Switching function 2 upper threshold
<	One level back

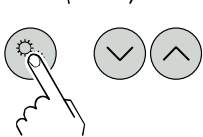
The AGC-100 has two switching functions with two adjustable thresholds each. The status of the switching functions is displayed on the front panel and can be evaluated via the floating contacts at the *CONTROL*, respectively *RELAY* connector (→  20).



Selecting a parameter



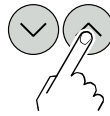
⇒ The name of the parameter and the currently valid parameter value are displayed.
e. g.: **SP1-CH DISABLED**
Switching function 1 turned off



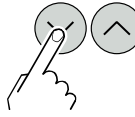
⇒ Select parameter. The value flashes and can now be edited.

Vacuum Gauge Controller AGC-100

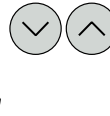
Editing and saving the parameter



⇒ Press key for <1 s:
The value is increased/decreased by 1 increment.



Press key for >1 s:
The value is increased/decreased continuously.



⇒ Save the change and return to read mode.



We recommend setting the upper threshold $\frac{1}{2}$ decade above the lower, or the lower threshold $\frac{1}{2}$ decade below the upper, threshold limit.

Vacuum Gauge Controller AGC-100

Configuring a switching function

	Value
SP1-CH	Configuring a switching function.
SP1-CH DISABLED	⇒ Switching function 1 is factory-deactivated
SP1-CH ENABLED	⇒ Switching function 1 is always turned on
SP1-CH 1	⇒ Switching function 1 is assigned to channel 1



The lower and the upper threshold of a switching function are always assigned to the same channel. The last assignment is valid for both thresholds.

Limits of the lower switching thresholds

	Value																
SP1-L	The lower threshold (Setpoint low) defines the pressure at which the switching function is activated when the pressure is dropping.																
e.g.: SP1-L 5.00-4	⇒ gauge dependent. If another gauge type is connected, the AGC-100 automatically adjusts the switching threshold if required.																
	<table><tr><th>SPx-L min.</th><th>SPx-L max.</th></tr><tr><td>PVG5xx</td><td>2×10^{-3} *)</td></tr><tr><td>PCG75x</td><td>2×10^{-3} *)</td></tr><tr><td>IMG500</td><td>1×10^{-9}</td></tr><tr><td>FRG70x</td><td>1×10^{-9}</td></tr><tr><td>FRG720</td><td>1×10^{-8}</td></tr><tr><td>FRG730</td><td>1×10^{-8}</td></tr><tr><td>CDG500</td><td>F.S. / 1000</td></tr></table>	SPx-L min.	SPx-L max.	PVG5xx	2×10^{-3} *)	PCG75x	2×10^{-3} *)	IMG500	1×10^{-9}	FRG70x	1×10^{-9}	FRG720	1×10^{-8}	FRG730	1×10^{-8}	CDG500	F.S. / 1000
SPx-L min.	SPx-L max.																
PVG5xx	2×10^{-3} *)																
PCG75x	2×10^{-3} *)																
IMG500	1×10^{-9}																
FRG70x	1×10^{-9}																
FRG720	1×10^{-8}																
FRG730	1×10^{-8}																
CDG500	F.S. / 1000																

all values in mbar, GAS=nitrogen

*) 2×10^{-4} mbar if RNG-EXT (Pirani range extension) is activated (→ 43)



The minimum hysteresis between the upper and lower switching threshold amounts to at least 10% of the lower threshold (logarithmic gauges) or 1% of the full scale value (linear gauges). The upper threshold is if necessary automatically adjusted to a minimum hysteresis. This prevents unstable states.

Vacuum Gauge Controller AGC-100

Limits of the upper switching thresholds

	Value										
SP1-H	<p>The upper switching threshold (Setpoint high) defines the pressure at which the switching function is deactivated when the pressure is rising.</p> <p>⇒ gauge dependent.</p> <p>If another gauge type is connected, the AGC-100 automatically adjusts the threshold if required.</p>										
e.g.: SP1-H 1500											
	<table> <tr> <th>SPx-H min.</th><th>SPx-H max.</th></tr> <tr> <td rowspan="7">= SPx-L min.</td><td>1×10^3</td></tr> <tr> <td>1.5×10^3</td></tr> <tr> <td>1×10^{-2}</td></tr> <tr> <td>1×10^3</td></tr> <tr> <td>1×10^3</td></tr> <tr> <td>1×10^3</td></tr> <tr> <td>F.S.</td></tr> </table>	SPx-H min.	SPx-H max.	= SPx-L min.	1×10^3	1.5×10^3	1×10^{-2}	1×10^3	1×10^3	1×10^3	F.S.
SPx-H min.	SPx-H max.										
= SPx-L min.	1×10^3										
	1.5×10^3										
	1×10^{-2}										
	1×10^3										
	1×10^3										
	1×10^3										
	F.S.										
PUG5xx											
PCG75x											
IMG500											
FRG70x											
FRG720											
FRG730											
CDG500											

all values in mbar, GAS=nitrogen



The minimum hysteresis between the upper and lower switching threshold amounts to at least 10% of the lower threshold (logarithmic gauges) or 1% of the full scale value (linear gauges). This prevents unstable states.

4.5.2 Gauge parameters

SENSOR

The sensor parameter group is used for displaying, entering and editing parameters of the connected gauges.

Parameters in this group

DEGAS

Cleaning the electrode system.

FSR

Measurement range linear gauges.

FILTER

Measurement value filter.

OFFSET

Offset correction.

GAS

Correction factor for other gases.

COR

Offset correction.

HV-CTRL

Activating / deactivating high vacuum measurement circuit.

EMISSION

Emission.

FILAMENT

Filament selection.

DIGITS

Display resolution.

<

One level back.

Some parameters are not available for all gauges and thus not always displayed.

→ 33 34 35 36 38 38 39 39 39 40

	DEGAS	FSR	FILTER	OFFSET	GAS	COR	HV-CTRL	EMISSION	FILAMENT	DIGITS
PUG5xx	-	-	✓	-	✓	✓	-	-	-	✓
PCG75x	-	-	✓	-	✓	✓	-	-	-	✓
IMG500	-	-	✓	-	✓	✓	✓	-	-	✓
FRG70x	-	-	✓	-	✓	✓	-	-	-	✓
FRG720	✓	-	✓	-	✓	✓	-	-	-	✓
FRG730	✓	-	✓	-	✓	✓	-	✓	✓	✓
CDG500	-	✓	✓	✓	-	✓	-	-	-	✓

Degas

Contamination deposits on the electrode system of hot cathode gauges may cause instabilities of the measurement values. The degas function facilitates cleaning of the electrode system.



The degas process works only at pressures below 7.2×10^{-6} mbar.



FRG730 gauges: The Degas function acts only upon the active filament.

Available for the following gauges:

- | | |
|---|------------------|
| <input type="checkbox"/> Pirani | (PVG) |
| <input type="checkbox"/> Pirani / Capacitance | (PCG) |
| <input type="checkbox"/> Cold cathode | (IMG) |
| <input type="checkbox"/> Cold cathode / Pirani | (FRG-70x) |
| <input checked="" type="checkbox"/> Hot ionization / Pirani | (FRG-720 / -730) |
| <input type="checkbox"/> Capacitance | (CDG) |

	Value	
DEGAS		
DEGAS OFF	⇒ Normal operation (Degas blocked)	
DEGAS ON	⇒ Degas: The electron collection grid is heated to ≈700 °C by electron bombardment and the electrode system is thus cleaned. Duration = 180 s.	

Editing and saving a parameter



⇒ Start Degas. Duration of the Degas function 180 seconds (may also be aborted).



Abort Degas.



⇒ Save change and return to read mode.

Vacuum Gauge Controller AGC-100

Measuring range (F.S.) of linear gauges

The full scale (F.S.) is automatically recognized.

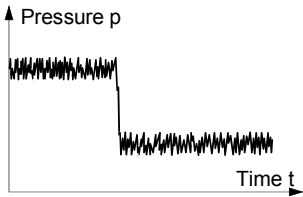
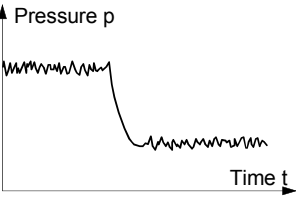
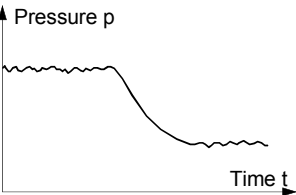
Available for the following gauges:

- | | |
|--|------------------|
| <input type="checkbox"/> Pirani | (PVG) |
| <input type="checkbox"/> Pirani / Capacitance | (PCG) |
| <input type="checkbox"/> Cold cathode | (IMG) |
| <input type="checkbox"/> Cold cathode / Pirani | (FRG-70x) |
| <input type="checkbox"/> Hot ionization / Pirani | (FRG-720 / -730) |
| <input checked="" type="checkbox"/> Capacitance | (CDG) |

	Value
FSR	
e.g. FSR 1000 MBAR	⇒ 0.01 mbar, 0.02 mbar, 0.05 mbar 0.01 Torr, 0.02 Torr, 0.05 Torr 0.10 mbar, 0.25 mbar, 0.50 mbar 0.10 Torr, 0.25 Torr, 0.50 Torr 1 mbar, 2 mbar, 5 mbar 1 Torr, 2 Torr, 5 Torr 10 mbar, 20 mbar, 50 mbar 10 Torr, 20 Torr, 50 Torr 100 mbar, 200 mbar, 500 mbar 100 Torr, 200 Torr, 500 Torr 1000 mbar, 1100 mbar 1000 Torr 2 bar, 5 bar, 10 bar, 50 bar A conversion table can be found in the Appendix (→ 82).

Measurement value filter

The measurement value filter permits a better evaluation of unstable or disturbed measuring signals.

	Value
<div>FILTER</div> <div>FILTER OFF</div> <div>FILTER FAST</div>	<div>⇒ No measurement value filter</div> <div>⇒ Fast: The AGC-100 responds quickly to fluctuations in the measurement value. As a result, it will respond faster to interference in the measured values.</div> <div></div>
<div>FILTER NORMAL</div>	<div>⇒ Normal (factory setting): Good relationship between response and sensitivity of the display and the switching function to changes in the measured values.</div> <div></div>
<div>FILTER SLOW</div>	<div>⇒ Slow: The AGC-100 does not respond to small changes in measured values. As a result, it will respond more slowly to changes in the measured values.</div> <div></div>

Offset correction of the controller




The offset value is displayed and readjusted according to the actual measurement value.

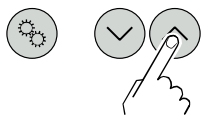
Available for the following gauges:

- ☐ Pirani (PVG)
- ☐ Pirani / Capacitance (PCG)
- ☐ Cold cathode (IMG)
- ☐ Cold cathode / Pirani (FRG-70x)
- ☐ Hot ionization / Pirani (FRG-720 / -730)
- ☒ Capacitance (CDG)

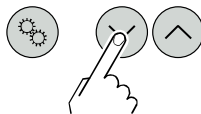
The offset correction affects:

- ☒ the displayed measurement value
- ☐ the displayed threshold value of the switching functions
- ☐ the analog outputs at the *CONTROL* connector (→ 20)

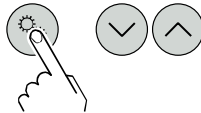
	Wert	
OFFSET		
OFFSET OFF	⇒ Offset correction factory-deactivated	
e.g. OFFSET 9.53	⇒ Offset correction activated (display in the relevant units of measurement)	



- ⇒ Press key for >1.5 s:
The offset value is readjusted (the actual measurement value is accepted as new offset value).



Reset the offset value.



- ⇒ Save change and return to read mode.

When offset correction is activated, the saved offset value is subtracted from the actual measurement value. This allows measuring relative to a reference pressure.

Vacuum Gauge Controller AGC-100

Zero adjustment of a digital CDG



First adjust the gauge and then the controller.

Available for the following gauges:

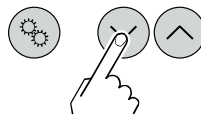
- | | |
|--|------------------|
| <input type="checkbox"/> Pirani | (PVG) |
| <input type="checkbox"/> Pirani / Capacitance | (PCG) |
| <input type="checkbox"/> Cold cathode | (IMG) |
| <input type="checkbox"/> Cold cathode / Pirani | (FRG-70x) |
| <input type="checkbox"/> Hot ionization / Pirani | (FRG-720 / -730) |
| <input checked="" type="checkbox"/> Capacitance | (CDG) |



When the zero of the gauge is readjusted, the offset correction must be deactivated.

	Value	
OFFSET		
e.g. OFFSET OFF	⇒ Zero adjustment deactivated	

Lit solid after >1.5 s and as long as key remains pressed



⇒ Press >1.5 s:
Zero adjustment of the digital CDG.



After adjusting the zero point, a zero value is displayed. Due to the measuring resolution of the CDG (noise, drift), a zero with plus/minus several digits are displayed.

Vacuum Gauge Controller AGC-100

Correction factor GAS

The correction factor GAS allows

- the measured value to be calibrated for the preset gases N₂, Ar, H₂, He, Ne, Kr and Xe, or
- manual input of the correction factor for other gases (COR).

→ Characteristic curves in  [1] ... [8].




This parameter is not available for the unit of measurement: Volt.

Available for the following gauges:

<input checked="" type="checkbox"/> Pirani	(PVG)	Only for pressures
<input checked="" type="checkbox"/> Pirani / Capacitance	(PCG)	<1 mbar
<input checked="" type="checkbox"/> Cold cathode	(IMG)	<1 mbar
<input checked="" type="checkbox"/> Cold cathode / Pirani	(FRG-70x)	<1×10 ⁻³ mbar
<input checked="" type="checkbox"/> Hot ionization / Pirani	(FRG-720 / -730)	<1×10 ⁻³ mbar
<input type="checkbox"/> Capacitance	(CDG)	

	Value
GAS	
GAS N2	⇒ Gas: nitrogen / air (factory setting)
GAS AR	⇒ Gas: argon
GAS H2	⇒ Gas: hydrogen
GAS HE	⇒ Gas: helium
GAS NE	⇒ Gas: neon
GAS KR	⇒ Gas: krypton
GAS XE	⇒ Gas: xenon
GAS COR	⇒ Calibration factor for other gases by manually entering parameter COR

Calibration factor COR

The calibration factor COR allows the measured value to be calibrated for other gases (→ characteristic curve in  [1] ... [8]). This parameter is effective in the entire measurement range of the gauge.



Precondition: Parameter "GAS COR" is set (except capacitance gauges).



This parameter is not available with the measurement unit: Volt.


Available for the following gauges:

<input checked="" type="checkbox"/> Pirani	(PVG)
<input checked="" type="checkbox"/> Pirani / Capacitance	(PCG)
<input checked="" type="checkbox"/> Cold cathode	(IMG)
<input checked="" type="checkbox"/> Cold cathode / Pirani	(FRG-70x)
<input checked="" type="checkbox"/> Hot ionization / Pirani	(FRG-720 / -730)
<input checked="" type="checkbox"/> Capacitance	(CDG)

	Value	
COR		
e.g. COR 1.00	⇒ No correction	
e.g. COR 1.53	⇒ Measurement value corrected by a factor of 0.10 ... 10.00	



Vacuum Gauge Controller AGC-100

Turning the gauge on / off

Activating / deactivating the high vacuum measurement circuit (→ also  [25]).

Available for the following gauges:

- ☐ Pirani (PVG)
- ☐ Pirani / Capacitance (PCG)
- ☒ Cold cathode (IMG)
- ☐ Cold cathode / Pirani (FRG-70x)
- ☐ Hot ionization / Pirani (FRG-720, FRG-730)
- ☐ Capacitance (CDG)

	Value	
HV-CTRL		
HV-CTRL ON	⇒ High vacuum measurement circuit activated	
HV-CTRL OFF	⇒ High vacuum measurement circuit deactivated	

Emission


Switching the emission on and off.

Available for the following gauges:

- ☐ Pirani (PVG)
- ☐ Pirani / Capacitance (PCG)
- ☐ Cold cathode (IMG)
- ☐ Cold cathode / Pirani (FRG-70x)
- ☒ Hot ionization / Pirani (FRG-730 only)
- ☐ Capacitance (CDG)

	Value	
EMISSION		
EMISSION AUTO	⇒ The emission is switched on and off automatically by the gauge	
EMISSION HAND	⇒ The emission is switched on and off by the user	



The symbol  lid solid, if the emission is switched on.

Filament

Means of selection.

Available for the following gauges:

- ☐ Pirani (PVG)
- ☐ Pirani / Capacitance (PCG)
- ☐ Cold cathode (IMG)
- ☐ Cold cathode / Pirani (FRG-70x)
- ☒ Hot ionization / Pirani (FRG-730 only)
- ☐ Capacitance (CDG)

	Value	
FILAMENT		
FILAMENT AUTO	⇒ The gauge automatically alternates between the filaments	
FILAMENT FIL 1	⇒ Filament 1 active	
FILAMENT FIL 2	⇒ Filament 2 active	

Display resolution

Display resolution of measured values.

Available for the following gauges:

<input checked="" type="checkbox"/> Pirani	(PVG)
<input checked="" type="checkbox"/> Pirani / Capacitance	(PCG)
<input checked="" type="checkbox"/> Cold cathode	(IMG)
<input checked="" type="checkbox"/> Cold cathode / Pirani	(FRG-70x)
<input checked="" type="checkbox"/> Hot ionization / Pirani	(FRG-720 / -730)
<input checked="" type="checkbox"/> Capacitance	(CDG)

	Value
DIGITS	
DIGITS AUTO	⇒ Automatic ^{*)} (factory setting)
DIGITS 1	⇒ e.g. 2E-1 or 500
DIGITS 2	⇒ e.g. 2.5E-1 or 520
DIGITS 3	⇒ e.g. 2.47E-1 or 523
DIGITS 4	⇒ e.g. 2.473E-1 or 523.7

^{*)} The mantissa is dependent on the connected gauge and the currently valid pressure value.

With PVG and PCG gauges in the pressure range $p < 1.0E-4$ mbar and activated RNG-EXT (→ 43) the display is reduced by one decimal digit.

4.5.3 Gauge Control

SENSOR-CONTROL >

The sensor control group is used for displaying, entering and editing parameters which define how the connected gauges are activated/deactivated.




This group is available for the IMG gauges only.

Parameters in this group

S-ON	Gauge activation
S-OFF	Gauge deactivation
T-OFF	OFF threshold
<	One level back

Gauge activation


IMG gauges can be activated by different means.

	Value
S-ON	
S-ON HAND	⇒ Manual activation: The gauge is activated by pressing the  key.
S-ON HOTSTART	⇒ Hot start: The gauge is automatically activated when the AGC-100 is turned on. Measurement is thus automatically resumed after a power failure. Gauge deactivation → 41.

Vacuum Gauge Controller AGC-100

Gauge deactivation

IMG gauges can be deactivated by different means.

	Value
S-OFF	<div>⇒ Manual deactivation: The gauge is deactivated by pressing the  key.</div> <div>⇒ Self control: The subsequent parameter T-OFF is used to specify the switch-off threshold. The sensor is switched off when the pressure at the sensor exceeds the switch-off threshold.</div>
S-OFF HAND	
S-OFF SELF	

OFF threshold

Definition of the OFF threshold for the gauge to be deactivated by itself.
This parameter is only available if the sensor deactivation parameter is set to S-OFF SELF.

	Value
T-OFF	<div>⇒ The sensor is switched off when the pressure exceeds the switch-off threshold.</div>
e.g.: T-OFF 1.00-2	

4.5.4 General Parameters

GENERAL

The General parameters group is used for displaying, entering and editing generally applicable system parameters.

Parameters in this group

- UNIT
- Measurement unit
- BAUD USB
- Transmission rate USB interface
- RNG-EXT
- Pirani range extension
- AO-MODE
- Recorder output
- ERR-RELAY
- Error relay
- BARGRAPH
- Bar graph display
- BACKLIGHT
- Backlight
- SCREENSAVE
- Screensaver
- CONTRAST
- Contrast adjustment
- DEFAULT
- Factory settings
- LANGUAGE
- Language
- FORMAT
- Number format, measurement value
- END VAL
- Display of measurement range end value
- <
- One level back

Measurement unit

Unit of measured values, thresholds etc. (conversion table → 82).

	Value
UNIT	
UNIT MBAR	⇒ mBar (factory setting)
UNIT HPASCAL	⇒ hPa
UNIT TORR	⇒ Torr (only available if Torr lock is not activated → 48)
UNIT PASCAL	⇒ Pa
UNIT MICRON	⇒ Micron (= 0.001 Torr) (only available if Torr lock is not activated → 48)
UNIT VOLT	⇒ V

A change of the pressure unit influences also the pressure unit settings of the FRG-720 and FRG-730 gauges.

If the measurement unit micron is selected, automatic changeover to Torr occurs above 99000 micron. Below 90 Torr automatic changeover back to the measurement unit micron occurs.

Vacuum Gauge Controller AGC-100

Transmission rate

Transmission rate of the USB interface.

	Value
BAUD USB	
BAUD USB 9600	⇒ 9600 baud
BAUD USB 19200	⇒ 19200 baud
BAUD USB 38400	⇒ 38400 baud
BAUD USB 57600	⇒ 57600 baud
BAUD USB 115200	⇒ 115200 baud (factory setting)

Pirani range extension

The display and setpoint adjustment range can be extended (the setting only affects the control unit).



Use the parameter Pirani range extension only with Pirani and Pirani / Capacitance gauges with display / measurement range up to 5×10^{-5} mbar.

Available for the following gauges:

- ☒ Pirani (PVG)
- ☒ Pirani / Capacitance (PCG)
- ☐ Cold cathode (IMG)
- ☐ Cold cathode / Pirani (FRG-70x)
- ☐ Hot ionization / Pirani (FRG-720 / -730)
- ☐ Capacitance (CDG)

	Value
RNG-EXT	
RNG-EXT DISABLED	⇒ Deactivated (factory setting)
RNG-EXT ENABLED	⇒ Display extended to 5×10^{-5} mbar

Error relay

Switching behaviour of the error relay.

	Value
ERR-RELAY	
ERR-RELAY ALL	⇒ Switches for all errors (factory setting)
ERR-RELAY no SE	⇒ Only unit errors
ERR-RELAY CH 1	⇒ Error sensor 1 and unit error

Vacuum Gauge Controller AGC-100

Bar graph

In the dot matrix a bar graph or the measured pressure as a function of time ($p = f(t)$) may be shown.

During parameter setting the parameter and the parameter value may be displayed in place of this.

	Value
BARGRAPH	
BARGRAPH OFF	⇒ Factory setting.
BARGRAPH FSR	⇒ Bar graph covering full scale range.
BARGRAPH FSR h	⇒ Bar graph covering full scale range, high-level presentation.
BARGRAPH FSR+SP	⇒ Bar graph covering full scale range and setpoint threshold.
BARGRAPH DEC	⇒ Bar graph covering a decade according to current measurement value.
BARGRAPH DEC h	⇒ Bar graph covering a decade according to current measurement value, high-level presentation.
BARGRAPH DEC+SP	⇒ Bar graph covering a decade according to current measurement value and setpoint threshold.
BARGRAPH f(0.2s)	⇒ $p = f(t)$, autoscaled, 0.2 seconds / pixel For each measurement every 200 ms a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 20 seconds.
BARGRAPH f(1s)	⇒ $p = f(t)$, autoscaled, 1 second / pixel For each measurement every second a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 100 seconds.
BARGRAPH f(6s)	⇒ $p = f(t)$, autoscaled, 6 seconds / pixel For each measurement every 6 seconds a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 10 minutes.
BARGRAPH f(1min)	⇒ $p = f(t)$, autoscaled, 1 minute / pixel For each measurement every minute a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 100 minutes.

Backlight

	Value
BACKLIGHT	
e.g. BACKLIGHT 60%	⇒ Factory setting Adjustable from 0 ... 100% 100% = full brightness

Vacuum Gauge Controller AGC-100

Screensaver

The screensaver reduces the brightness of the backlight or switches it off completely (dark room).

	Value
SCREENSAVE	
SCREENSAVE OFF	⇒ factory setting
SCREENSAVE 10min	⇒ after 10 minutes
SCREENSAVE 30min	⇒ after 30 minutes
SCREENSAVE 1h	⇒ after 1 hour
SCREENSAVE 2h	⇒ after 2 hours
SCREENSAVE 8h	⇒ after 8 hours
SCREENSAVE DR	⇒ the backlight is switched off completely after 1 minute

Contrast



	Value
CONTRAST	
e.g. CONTRAST 40%	⇒ factory setting adjustable from 0 ... 100 % 100% = full contrast

Default parameter settings

All user parameter settings are replaced by the default values (factory settings).



Loading of the default parameter settings is irreversible.

	Value
DEFAULT	
DEFAULT ▼+▲ 2s	Press   keys at the same time for >2 s to start loading default values
DEFAULT SET	⇒ The default values are loaded

Language

Display language.

	Value
LANGUAGE	
LANGUAGE ENGLISH	⇒ English (factory setting)
LANGUAGE GERMAN	⇒ German
LANGUAGE FRENCH	⇒ French

Vacuum Gauge Controller AGC-100

Measurement value format

Measurement values in floating point or exponential format. If a measurement value cannot reasonably be expressed in the floating point format, it is automatically displayed in the exponential format.

	Value
FORMAT	
FORMAT X.X	⇒ Floating point format, if possible (factory setting)
FORMAT X.XESY	⇒ Exponential format

Display of measurement range end value

Display of underrange or overrange.

	Value
END VAL	
END VAL UR/OR	⇒ When an underrange or overrange occurs UR or OR is displayed (factory setting)
END VAL VALUE	⇒ When an underrange or overrange occurs the respective full scale value is displayed

4.5.5 Test Parameters

TEST

The Test parameter group is used for e.g. displaying the firmware version, entering and editing special parameter values, and for running test programs.



The group is only available if

- the key was pressed while the AGC-100 was turned on, or
- the key was pressed for 5 s while is displayed.

Parameters in this group

SOFTWARE	Firmware version
HARDWARE	Hardware version
MAC	MAC address
RUNHOURS	Operating hours
WATCHDOG	Watchdog control
TORR-LOCK	Torr lock
KEY-LOCK	Keylock
FLASH	FLASH test (program memory)
EEPROM	EEPROM test (parameter memory)
DISPLAY	Display test
I/O	I/O test
<	One level back

The parameters in this group are available for all gauges.

Firmware version

The firmware version (program version) is displayed.

	Version
e.g. SOFTWARE 1.00	This information is helpful when contacting Agilent Technologies

Hardware version

The hardware version is displayed.

	Hardware
e.g. HARDWARE 1.0	This information is helpful when contacting Agilent Technologies

MAC address

The MAC address is displayed.

	MAC address
e.g.: MAC 00A0410A0008	The address is displayed without any separators (e.g. 00-A0-41-0A-00-08)

Operating hours

The operating hours are displayed.

	Hours
e.g. RUNHOURS 24 h	⇒ Operating hours

Vacuum Gauge Controller AGC-100

Watchdog control

Behaviour of the system control (watchdog control) in the event of an error.

	Setting
WATCHDOG	
WATCHDOG AUTO	⇒ The system automatically acknowledges error messages of the watchdog after 2 s (factory setting)
WATCHDOG OFF	⇒ Error messages of the watchdog have to be acknowledged by the operator

Torr lock

The measurement units Torr and Micron can be suppressed in the corresponding parameter setting **UNIT TORR** (→ 42).

	Setting
TORR-LOCK	
TORR-LOCK OFF	⇒ Measurement units Torr and Micron available (factory setting)
TORR-LOCK ON	⇒ Measurement units Torr and Micron not available



Keylock

The keylock function prevents inadvertent entries in the parameter mode and thus malfunctions.

	Setting
KEY-LOCK	
KEY-LOCK OFF	⇒ Keylock function disabled (factory setting)
KEY-LOCK ON	⇒ Keylock function enabled

FLASH test

Test of the program memory.

	Test sequence
FLASH ▼+▲	Press   keys at the same time to start test
FLASH RUN	⇒ Test in progress (very briefly)
FLASH PASS	⇒ Test completed, no error found. After the test, an 8-digit checksum (e.g. FLASH 0x12345678) is displayed.
FLASH ERROR	⇒ Test completed, error found. After the test, an 8-digit checksum (e.g. FLASH 0x12345678) is displayed. If the error persists after repeating the test, please contact your nearest Agilent Vacuum service center.

Vacuum Gauge Controller AGC-100

EEPROM test

Test of the parameter memory.

	Test sequence
EEPROM ▼+▲	Press (▼) (▲) keys at the same time to start test
EEPROM RUN	⇒ Test in progress.
EEPROM PASS	⇒ Test completed, no error found.
EEPROM ERROR	⇒ Test completed, error found. If the error persists after repeating the test, please contact your nearest Agilent Vacuum service center.


Display test


Test of the display.

	Test sequence
DISPLAY ▼+▲	Press (▼) (▲) keys at the same time to start test
	⇒ After starting the test, all display elements are lit at the same time for 10 s.

I/O test

Test of the unit relays. The test program tests their switching function.


Caution



Caution: The relays switch irrespective of the pressure. Starting a test program may cause unwanted effects in connected control systems.

Disconnect all sensor and control system lines to ensure that no control commands or messages are triggered by mistake.

The relays switch on and off cyclically. The switching operations are indicated optically and are also clearly audible.

The switching function contacts are connected to the *CONTROL* connector on the rear of the unit (→ 20). Check their function with an ohmmeter.

	Test sequence
I/O ▼+▲	Press (▼) (▲) keys at the same time to start test
I/O OFF	⇒ All relays deactivated
I/O REL1 ON	⇒ Switching function relay 1
I/O REL1 OFF	⇒ Switching function relay 1
I/O REL2 ON	⇒ Switching function relay 2
⋮	

4.5.6 Data Logger Mode

DATA LOGGER

The data logger group is used for

- recording measurement data on a USB memory stick (interface type A on the front of the Center unit)
- deleting recorded measurement data from the USB memory stick



This group is only available when a USB memory stick formatted for the FAT file system (FAT32) is plugged in. Use a max. 32 GB memory stick.



Not all USB memory sticks are automatically recognized by the AGC-100, as they (in particular cheaper brands) do not always conform to USB standard requirements. Try a different memory stick before contacting your nearest Agilent Technologies service center.

Parameters in this group

DATE	Current date
TIME	Current time
INTERVAL	Display interval
DEC-SEPARATOR	Decimal separator
FILENAME	File name
START / STOP	Start / stop record
CLEAR	Deletion of files with displayed measurement data

Date

	Value
DATE	Current date in the format YYYY-MM-DD
e.g. DATE 2015-04-15	⇒ e.g. 2016-01-15

Time

	Value
TIME	Current time in the format hh:mm [24 h]
e.g. TIME 15:45	⇒ e.g. 15:45

Interval

Data logging interval.

	Value
INTERVAL	
INTERVAL 1s	⇒ Recording interval 1/s
INTERVAL 10s	⇒ Recording interval 1/10 s
INTERVAL 30s	⇒ Recording interval 1/30 s
INTERVAL 1min	⇒ Recording interval 1/60 s
INTERVAL 1%	⇒ Recording interval: in the event of measurement value changes ≥1%
INTERVAL 5%	⇒ Recording interval: in the event of measurement value changes ≥5%

Vacuum Gauge Controller AGC-100

Decimal separator

Decimal separator for measurement values in the measurement data file.

	Wert
DEC-SEPARATOR	
DEC-SEPARATOR ,	⇒ Decimal comma
DEC-SEPARATOR .	⇒ Decimal point

File name

	Value
FILENAME	Name of the measurement data file, max. 7 digits
e.g. FILENAME DATALOG	⇒ File ending: CSV

After entering the 7th digit the display stops flashing. The name of the data file is saved and the unit is in the read mode again.






Is the file name shorter than 7 digits, a blank space must be set to each remaining digit.

Start / Stop


Starting / stopping measurement value display.



 flashes during measurement data display (in the measurement mode only).



	Value
START	
START ▲	⇒ Press  key to start record: Data record is running, display has changed to STOP ▼ and the down arrow ▼ is blinking.
STOP ▼	⇒ Press  key to stop saving: Data record is stopped, display has changed to START ▲ and the up arrow ▲ is blinking.



The unit does not return automatically to the measurement mode, as long as the arrows ▼ or ▲ in the display are blinking. Press the  key to leave the editing mode. Then, after approx. 10 s, the unit returns automatically to the measurement mode.

Deletion

Deletion of all measurement data files (ending CSV) from USB memory stick.

	Value
CLEAR ▼+▲	Press   keys at the same time to delete files
CLEAR RUNNING	⇒ CSV files are being deleted
CLEAR DONE	⇒ CSV files have been deleted

4.5.7 Parameter Transfer Mode

SETUP

- This group is used for
- saving all parameters on a USB memory stick (interface type A on the front of the AGC-100)
 - loading all parameters from a USB memory stick onto the AGC-100
 - formatting a USB memory stick
 - deleting files with saved parameters from the USB memory stick



This group is only available when a USB memory stick formatted for the FAT file system (FAT32) is plugged in. Use a max. 32 GB memory stick.

Parameters in this group

SAVE

Saving all parameters

RESTORE

Loading all parameters onto the AGC-100

FORMAT

Formatting USB memory stick (FAT32)

CLEAR

Deletion of files with saved parameters

<

One level back

Saving a parameter

Saving all parameters of the AGC-100 to a USB memory stick (file ending: CSV).

	Value
SAVE	
SAVE SETUP	⇒ File name on the USB memory stick: SETUP01.CSV
:	
SAVE SETUP99	⇒ File name on the USB memory stick: SETUP99.CSV
SAVE RUNNING	⇒ CSV file is being saved
SAVE DONE	⇒ Saving completed

Loading parameters



Loading all parameters from a USB memory stick onto the AGC-100.

	Value
RESTORE	
RESTORE SETUP01	⇒ File name on the USB memory stick: SETUP.CSV
:	
RESTORE SETUP99	⇒ File name on the USB memory stick: SETUP99.CSV
RESTORE RUNNING	⇒ CSV file is being loaded
RESTORE DONE	⇒ Loading completed
RESTORE ERROR	⇒ Error occurred

Vacuum Gauge Controller AGC-100



Formatting

Formatting USB memory stick.

	Value
FORMAT ▼+▲	Press   keys at the same time to start formatting
FORMAT RUNNING	⇒ Formatting in progress
FORMAT DONE	⇒ Formatting completed

Deleting

Deleting all parameter files (ending CSV) from the USB memory stick.

	Value
CLEAR ▼+▲	Press   keys at the same time to delete files
CLEAR RUNNING	⇒ CSV files are being deleted
CLEAR DONE	⇒ CSV files have been deleted

5 Communication Protocol (Serial Interface)

The AGC-100 communicates with a computer via virtual serial interfaces (COM ports). Thus the user software can access the AGC-100 via USB Type B or via Ethernet interface.

Communication via USB Type B interface

The corresponding driver for the virtual COM port is installed automatically, when the AGC-100 is connected to a computer via the USB Type B interface. If the driver is not installed automatically, it can be downloaded from the FTDI website (www.ftdichip.com/Drivers/VCP.htm).

The installed virtual COM port appears as additional serial interface in the device manager of the computer.

Communication via Ethernet interface

With the Ethernet Configuration Tool a virtual serial interface (COM) can be assigned to an IP address. In addition, it allows configuration of the Ethernet interface via a computer (→ 86).

The installed virtual COM port appears as additional serial interface in the device manager of the computer.

When the AGC-100 is put into operation, it starts transmitting measured values in intervals of 1 s. As soon as the first character is transferred to the AGC-100, the automatic transmission of measured values stops. After the necessary inquiries or parameter modifications have been made, the transmission of measured values can be started again with the **COM** command (→ 58).

5.1 Data Transmission

The data transmission is bi-directional, i.e. data and control commands can be transmitted in either direction.

Data format

1 start bit, 8 data bits, no parity bit, 1 stop bit, no hardware handshake

Definitions

The following abbreviations and symbols are used:

Symbol	Meaning		
HOST	Computer or terminal		
[...]	Optional elements		
ASCII	American Standard Code for Information Interchange		
		Dez	Hex
<ETX>	END OF TEXT (CTRL C) Reset the interface	3	03
<CR>	CARRIAGE RETURN Go to beginning of line	13	0D
<LF>	LINE FEED Advance by one line	10	0A
<ENQ>	ENQUIRY (CTRL E) Request for data transmission	5	05
<ACK>	ACKNOWLEDGE Positive report signal	6	06
<NAK>	NEGATIVE ACKNOWLEDGE Negative report signal	21	15

"Transmit": Data transfer from HOST to AGC-100

"Receive": Data transfer from AGC-100 to HOST

Flow Control

After each ASCII string, the HOST must wait for a report signal (<ACK><CR><LF> or <NAK> <CR><LF>).

The input buffer of the HOST must have a capacity of at least 32 bytes.

5.2 Communication Protocol

Transmission format

Messages are transmitted to the AGC-100 as ASCII strings in the form of mnemonic operating codes and parameters. All mnemonics comprise three ASCII characters.

Spaces are ignored. <ETX> (CTRL C) clears the input buffer in the AGC-100.

Transmission protocol

HOST	AGC-100	Explanation
Mnemonics [and parameters]	—————>	Receives message with "end of message"
<CR>[<LF>]	—————>	
<—————	<ACK><CR><LF>	Positive acknowledgment of a received message

Reception format

When requested with a mnemonic instruction, the AGC-100 transmits the measurement data or parameters as ASCII strings to the HOST.

<ENQ> (CTRL E) must be transmitted to request the transmission of an ASCII string. Additional strings, according to the last selected mnemonic, are read out by repetitive transmission of <ENQ>.

If <ENQ> is received without a valid request, the ERROR word is transmitted.

Reception protocol

HOST	AGC-100	Explanation
Mnemonics [and parameters]	—————>	Receives message with "end of message"
<CR>[<LF>]	—————>	
<—————	<ACK><CR><LF>	Positive acknowledgment of a received message
<ENQ>	—————>	Requests to transmit data
<—————	Measurement values or parameters	Transmits data with "end of message"
<—————	<CR><LF>	
	:	:
<ENQ>	—————>	Requests to transmit data
<—————	Measurement values or parameters	Transmits data with "end of message"
<—————	<CR><LF>	


Error processing

The strings received are verified in the AGC-100. If an error is detected, a negative acknowledgment <NAK> is output.

Error recognition protocol

HOST	AGC-100	Explanation
Mnemonics [and parameters]	—————>	Receives message with "end of message"
<CR>[<LF>]	—————>	
***** Transmission or programming error *****		
<————	<NAK><CR><LF>	Negative acknowledgment of a received message
Mnemonics [and parameters]	—————>	Receives message with "end of message"
<CR>[<LF>]	—————>	
<—————	<ACK><CR><LF>	Positive acknowledgment of a received message

5.3 Mnemonics

		→ 
AYT	Are you there?	77
BAL	Backlight	67
BAU	Transmission rate (USB)	67
COM	Continuous mode of measurement values	58
COR	Calibration factor	61
DAT	Date	72
DCB	Display control bar graph	68
DCC	Display control contrast	69
DCD	Display resolution	62
DCS	Display control screensaver	69
DGS	Degas	62
ERA	Error relay allocation	69
ERR	Error status	58
ETH	Ethernet configuration	78
EUM	Emission user mode	62
EVA	Measurement range end value	70
FIL	Measurement value filter	63
FMT	Number format (measurement value)	70
FSR	Measurement range (linear gauges)	64
FUM	Filament user mode	63
GAS	Gas type correction	65
HDW	Hardware version	73
HVC	HV control, EMI on/off	73
ITR	Data output	65
LCM	Start / stop data logger	72
LNG	Language (display)	70
LOC	Keylock	73
MAC	Ethernet MAC address	74
OFS	Offset correction	66
PNR	Firmware version	74
PR1	Measurement data	59
PRE	Pirani range extension	71
RES	Reset	59
RHR	Operating hours	74
SAV	Save parameters (EEPROM)	71
SC1	Gauge control	66
SCM	Save / load parameters (USB)	73
SP1	Switching function 1	61
SP2	Switching function 2	61
SPS	Switching function status	60
TAD	A/D converter test	74
TAI	ID resistance test	74
TDI	Display test	75
TEE	EEPROM test	75
TEP	FLASH test	75
TID	Gauge identification	60
TIM	Time	72
TIO	I/O test	76

Vacuum Gauge Controller AGC-100

TKB	Operator key test	76
TLC	Torr lock	76
TMP	Inner temperature of the unit	77
TRS	Serial Interface test	77
UNI	Pressure unit	71
WDT	Watchdog control	77

5.4 Measurement Mode

COM - Continuous output of measurement values

Transmit: **COM** [,a] <CR>[<LF>]

	Description
a	Time interval, a = 0 → 100 ms 1 → 1 s (default) 2 → 1 minute

Receive: <ACK><CR><LF>
<ACK> is immediately followed by the continuous output of the measurement value in the desired interval.

Receive: b,sx.xxxxEsxx <CR><LF>

	Description
b	Status, b = 0 → Measurement data okay 1 → Underrange 2 → Overrange 3 → Measurement value error (sensor error) 4 → Sensor off 5 → No sensor 6 → Identification error 7 → Error FRG
sx.xxxxEsxx	Measurement value ¹⁾ [in current pressure unit] (s = sign)



¹⁾ Values always in exponential format.
For logarithmic gauges, the 3rd and 4th decimal are always 0.

ERR - Error status

Transmit: **ERR** <CR>[<LF>] Error status

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: aaaa <CR><LF>

	Description
aaaa	Error status, aaaa = 0000 → No error 1000 → ERROR (controller error (see display on front panel) 0100 → NO HWR (no hardware) 0010 → PAR (inadmissible parameter) 0001 → SYN (Syntax error)



The ERROR word is cancelled when read out. If the error persists, it is immediately set again.

Vacuum Gauge Controller AGC-100

PR1 - Measurement data gauge

Transmit: **PR1** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a,sx.xxxxEsxx <CR><LF>

	Description
a	Status, a = 0 → Measurement data okay 1 → Underrange 2 → Overrange 3 → Sensor error 4 → Sensor off 5 → No sensor 6 → Identification error 7 → Error FRG
sx.xxxxEsxx	Measurement value ¹⁾ [in current pressure unit] (s = sign)



¹⁾ Values always in exponential format.
 For logarithmic gauges, the 3rd and 4th decimal are always 0.

RES - Reset

Transmit: **RES** [,a] <CR>[<LF>]

	Description
a	a = 1 → Cancels currently active error and returns to measurement mode

Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: b[,b][,b][...] <CR><LF>

	Description
b	List of all present error messages, b = 0 → No error 1 → Watchdog has responded 2 → Task fail error 3 → FLASH error 4 → RAM error 5 → EEPROM error 6 → DISPLAY error 7 → A/D converter error 8 → UART error 9 → General error 10 → ID error

Vacuum Gauge Controller AGC-100

TID - Gauge identification

Transmit: TID <CR>[<LF>] Gauge identification
Receive: <ACK><CR><LF>
Transmit: <ENQ>
Receive: a <CR><LF>

	Description
a	Identification gauge 1, a = PVG5xx (Pirani Gauge) PCG75x (Pirani / Capacitance Gauge) IMG500 (Cold Cathode Gauge) FRG70x (Cold Cathode / Pirani Gauge) CDG500 (Capacitance Gauge) FRG720 (Hot Ionization / Pirani Gauge) FRG730 (Hot Ionization / Pirani Gauge) noSEn (No sensor) noid (No identifier)

5.5 Switching Function Parameters

SPS - Switching function status

Transmit: SPS <CR>[<LF>]
Receive: <ACK><CR><LF>
Transmit: <ENQ>
Receive: a,b <CR><LF>

	Description
a	Status switching function 1, a = 0 -> Off 1 -> On
b	Status switching function 2

Vacuum Gauge Controller AGC-100

SP1, SP2 - Switching function 1 and 2

Transmit: **SPx** [,a,x.xxxxEsxx,y.yyyyEsyy] <CR>[<LF>]

	Description
x	Switching function, x = 1 → Switching function 1 2 → Switching function 2
a	Switching function assignment, a = 0 → Turned off 1 → Turned on 2 → Measurement channel 1
x.xxxxEsxx	Lower threshold ¹⁾ [in current pressure unit] (default = depending on gauge) (s = sign)
y.yyyyEsyy	Upper threshold ¹⁾ [in current pressure unit] (default = depending on gauge) (s = sign)



¹⁾ Values can be entered in any format.
They are internally converted into the floating point format.

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a,x.xxxxEsxx,y.yyyyEsyy <CR><LF>

	Description
a	Switching function assignment
x.xxxxEsxx	Lower threshold [in current pressure unit] (s = sign)
y.yyyyEsyy	Upper threshold [in current pressure unit] (s = sign)

5.6 Gauge Parameters

COR - Calibration factor

Transmit: **COR** [,a.aaa] <CR>[<LF>]

	Description
a.aaa	Calibration factor, 0.100 ... 10.000 (default = 1.000)

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a.aaa,b.bbb,c.ccc <CR><LF>

	Description
a.aaa	Calibration factor

Vacuum Gauge Controller AGC-100

DCD - Display resolution

Transmit: **DCD** [,a] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Resolution a = 0 → AUTO (default) 1 → One digit 2 → Two digits 3 → Three digits 4 → Four digits

When the PrE (→ 71) is ON and the pressure is in the range $p < 1.0E-4$ mbar the display resolution of the PCG75x gauges is reduced by one decimal digit.

DGS - Degas

Transmit: **DGS** [,a] <CR>[<LF>]

	Description
a	Degas, a = 0 → Degas off (default) 1 → Degas on (3 minutes)

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Degas status

EUM - Emission user mode

Transmit: **EUM** [,a] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Emission, a = 0 → Manually 1 → Automatic (default)

Vacuum Gauge Controller AGC-100

FIL - Measurement value filter

Transmit: **FIL** [,a] <CR>[<LF>]

	Description
a	Filter, a = 0 → Filter off 1 → Fast 2 → Normal 3 → Slow

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Filter time constant

FUM - Filament user mode
FRG730

Transmit: **FUM** [,a] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Filament, a = 0 → Automatic (default) 1 → Filament 1 2 → Filament 2

Vacuum Gauge Controller AGC-100

FSR - Measurement range
(linear gauges)



The full scale value of the measurement range (Full Scale) of linear analog gauges has to be defined by the user. The full scale value of linear digital gauges and logarithmic gauges is automatically recognized.

Transmit: FSR [,a] <CR><LF>

	Description
a	Full scale value, a = 0 → 0.01 mbar 1 → 0.01 Torr 2 → 0.02 mbar 3 → 0.02 Torr 4 → 0.05 mbar 5 → 0.05 Torr 6 → 0.10 mbar 7 → 0.10 Torr 8 → 0.25 mbar 9 → 0.25 Torr 10 → 0.50 mbar 11 → 0.50 Torr 12 → 1 mbar 13 → 1 Torr 14 → 2 mbar 15 → 2 Torr 16 → 5 mbar 17 → 5 Torr 18 → 10 mbar 19 → 10 Torr 20 → 20 mbar 21 → 20 Torr 22 → 50 mbar 23 → 50 Torr 24 → 100 Torr 25 → 100 mbar 26 → 200 mbar 27 → 200 Torr 28 → 500 mbar 29 → 500 Torr 30 → 1000 mbar 31 → 1100 mbar 32 → 1000 Torr 33 → 2 bar 34 → 5 bar 35 → 10 bar 36 → 50 bar

Receive: <ACK><CR><LF>
Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Full scale value

Vacuum Gauge Controller AGC-100

GAS - Gas type correction

Transmit: **GAS** [,a] <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a <CR><LF>

	Description
a	Gas type correction, a = 0 → nitrogen / air (default) 1 → Argon 2 → Hydrogen 3 → Helium 4 → Neon 5 → Krypton 6 → Xenon 7 → Other gases

HVC - HV control, EMI on / off

Transmit: **HVC** [,a] <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a <CR><LF>

	Description
a	Mode, a = 0 → Off 1 → On

ITR - Data output FRG-720, FRG-730, CDG

Transmit: **ITR** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: aa,aa,aa,aa,aa,aa,aa,aa <CR><LF>

	Description
aa,aa,aa,aa,aa,aa,aa,aa	Data string (byte 0 ... 7 in hex format)

Vacuum Gauge Controller AGC-100

OFS - Offset correction
(linear gauges)

Transmit: **OFS** [,a,sx.xxxxEsxx] <CR>[<LF>]

	Description
a	Mode, a = 0 → Off (default) No offset value needs to be entered 1 → On If no offset value has been entered, the previously defined offset value is taken over 2 → Auto (offset measurement) No offset value needs to be entered 3 → Zero adjustment CDG No offset value needs to be entered
sx.xxxxEsxx	Offset ¹⁾ , [in current pressure unit] (default = 0.0000E+00) s = sign



¹⁾ Values can be entered in any format. They are internally converted into the floating point format.

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a,sx.xxxxEsxx <CR><LF>

	Description
a	Mode
sx.xxxxEsxx	Offset ¹⁾ , [in current pressure unit] s = sign

5.7 Gauge Control

SC1 - Gauge control

Transmit: **SC1** [,a,b,c.ccEscc,d.ddEsdd] <CR>[<LF>]

	Description
a	Gauge activation, a = 0 → Manual (default) 1 → Hot start 3 → Via measurement channel 1
b	Gauge deactivation, b = 0 → Manual (default) 1 → Self control 3 → Via measurement channel 1
c.ccEscc	ON threshold (s = sign)
d.ddEsdd	OFF threshold (s = sign)

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a,b,c.ccEscc,d.ddEsdd <CR><LF>

	Description
a	Gauge activation
b	Gauge deactivation
c.ccEscc	ON threshold (s = sign)
d.ddEsdd	OFF threshold (s = sign)

5.8 General Parameters

BAL - Backlight

Transmit: **BAL** [,a] <CR>[<LF>]

	Description
a	Backlight in percent, a = 0 ... 100 100% is full brightness

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Backlight

BAU - Transmission rate (USB)

Transmit: **BAU** [,a] <CR>[<LF>]

	Description
a	Transmission rate, a = 0 → 9600 Baud 1 → 19200 Baud 2 → 38400 Baud 3 → 57600 Baud 4 → 115200 Baud (default)



As soon as the new baud rate has been entered, the report signal is transmitted at the new transmission rate.

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: x <CR><LF>

	Description
a	Transmission rate

Vacuum Gauge Controller AGC-100

DCB - Display control bar graph Transmit: DCB [,a] <CR><LF>

	Description
a	Bar graph display, a = 0 -> Off (default) 1 -> Bar graph covering full scale range 2 -> Bar graph covering full scale range, high-level presentation 3 -> Bar graph covering full scale range and setpoint threshold 4 -> Bar graph covering a decade according to current measurement value 5 -> Bar graph covering a decade according to current measurement value, high-level presentation 6 -> Bar graph covering a decade according to current measurement value and setpoint threshold 7 -> $p = f(t)$, autoscaled, 0.2 seconds / pixel For each measurement every 200 ms a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 20 seconds. 8 -> $p = f(t)$, autoscaled, 1 second / pixel For each measurement every second a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 100 seconds. 9 -> $p = f(t)$, autoscaled, 6 seconds / pixel For each measurement every 6 seconds a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 10 minutes. 10 -> $p = f(t)$, autoscaled, 1 minute / pixel For each measurement every minute a measurement value is saved in tabular form and the last 100 measurement values (=100 pixel) are shown autoscaled. The represented data string corresponds to a logging duration of 100 minutes.

Receive: <ACK><CR><LF>
Transmit: <ENQ>
Receive: a <CR><LF>

	Description
a	Bar graph display

Vacuum Gauge Controller AGC-100

DCC - Display control contrast

Transmit: **DCC** [,a] <CR><LF>

	Description
a	Contrast in percent, a = 0 ... 100 100% = full contrast

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Contrast

DCS - Display control screensaver

Transmit: **DCS** [,a] <CR><LF>

	Description
a	Screensaver, a = 0 → Off (default) 1 → After 10 minutes 2 → After 30 minutes 3 → After 1 hour 4 → After 2 hours 5 → After 8 hours 6 → The backlight is switched off completely after 1 minute

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Screensaver

ERA - Error relay allocation

Transmit: **ERA** [,a] <CR><LF>

	Description
a	Switching behaviour error relay, a = 0 → Switches for all errors (default) 1 → Only unit errors 2 → Error sensor and unit error

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Switching behaviour error relay

Vacuum Gauge Controller AGC-100

EVA - Measurement range end value

Transmit: **EVA** [,a] <CR>[<LF>]

	Description
a	Measurement range end value, a = 0 → UR or OR is displayed (default) when an underrange or overrange occurs 1 → The measurement range end value is displayed when an underrange or overrange occurs

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Measurement range end value

FMT - Number format (measurement value)

Transmit: **FMT** [,a] <CR>[<LF>]

	Description
a	Number format (measurement value), a = 0 → Floating point format, if possible (default) 1 → Exponential format

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Number format

LNG - Language (display)

Transmit: **LNG** [,a] <CR>[<LF>]

	Description
a	Language, a = 0 → English (default) 1 → German 2 → French

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Language

Vacuum Gauge Controller AGC-100

PRE - Pirani range extension

Transmit: **PRE** [,a] <CR>[<LF>]

	Description
a	Pirani range extension, a = 0 → Disabled (default) 1 → Enabled

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Pirani range extension



PCG-75x gauges only, measurement range up to 5×10^{-5} mbar.

SAV - Save parameters (EEPROM)

Transmit: **SAV** [,a] <CR>[<LF>]

	Description
a	Save parameters to EEPROM, a = 0 → Save default parameters (default) 1 → Save user parameters

Receive: <ACK><CR><LF>

UNI - Pressure unit

Transmit: **UNI** [,a] <CR>[<LF>]

	Description
a	Pressure unit, a = 0 → mbar/bar 1 → Torr 2 → Pascal 3 → Micron 4 → hPascal (default) 5 → Volt

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Pressure unit

5.9 Data Logger Parameters



The group is only available when a USB memory stick formatted for the the FAT file system (FAT32) is plugged in. Use a max. 32 GB memory stick.

DAT - Date

Transmit: **DAT** [,yyyy-mm-dd] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: yyyy-mm-dd <CR><LF>

	Description
yyyy-mm-dd	Current date in the format yyyy-mm-dd

LCM - Start / stop data logger

Transmit: **LCM** [,a,b,c,ddddddd] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a,b,c,ddddddd <CR><LF>

	Description
a	Data logger command, a = 0 → Stop / data logging stopped 1 → Start / data logging started 2 → Clear / deletion of measurement data file (ending CSV) from USB memory stick
b	Data logging interval, b = 0 → Logging interval 1/s 1 → Logging interval 1/10 s 2 → Logging interval 1/30 s 3 → Logging interval 1/60 s 4 → Logging interval in the event of measurement value changes ≥1% 5 → Logging interval in the event of measurement value changes ≥5%
c	Decimal separator, c = 0 → , (decimal comma) 1 → . (decimal point)
ddddddd	File name (max. 7 digits)

TIM - Time

Transmit: **TIM** [,hh:mm] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: hh:mm <CR><LF>

	Description
hh:mm	Current time in the format hh:mm [24 h]

5.10 Parameter Transfer



The group is only available when a USB memory stick formatted for the the FAT file system (FAT32) is plugged in. Use a max. 32 GB memory stick.

SCM - Save / load parameters (USB)

Transmit: **SCM** [,a,bb] <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Setup parameters, a = 0 → Saving completed (read only) 1 → CSV file is being saved (read only) 2 → Loading all parameters from the USB memory stick onto the AGC-100 3 → Formatting USB memory stick (FAT32) 4 → Deleting parameter files (ending CSV) from the USB memory stick
bb	Number in the file name (0 ... 99)

5.11 Test Parameters

(For service personnel)

HDW - Hardware version

Transmit: **HDW** <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a.a <CR><LF>

	Description
a.a	Hardware version, e.g. 1.0

LOC - Keylock

Transmit: **LOC** [,a] <CR>[<LF>]

	Description
a	Keylock, a = 0 → Off (default) 1 → On

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Keylock status

Vacuum Gauge Controller AGC-100

MAC - Ethernet MAC address

Transmit: **MAC** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: aa-aa-aa-aa-aa-aa <CR><LF>

	Description
aa-aa-aa-aa-aa-aa	Ethernet MAC address of the unit: 00-A0-41-0A-00-00 ... 00-A0-41-0B-FF-FF

PNR - Firmware version

Transmit: **PNR** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a.aa <CR><LF>

	Description
a.aa	Firmware version, e.g. 1.00

RHR - Operating hours

Transmit: **RHR** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a <CR><LF>

	Description
a	Run (operating) hours, e.g. 24 [hours]

TAD - A/D converter test

Transmit: **TAD** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: aa.aaaa <CR><LF>

	Description
aa.aaaa	A/D converter Measurement signal [0.0000 ... 11.0000 V]

TAI – ID resistance test

Transmit: **TAI** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ> starts the test (very brief)
 Receive: a.aa <CR><LF>

	Description
a.aa	Identification [kOhm]

Vacuum Gauge Controller AGC-100

TDI - Display test

Transmit: **TDI** [,a] <CR>[<LF>]

	Description
a	Display test, a = 0 → Stops the test - display according to current operating mode (default) 1 → Starts the test - all segments on

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: x <CR><LF>

	Description
x	Display test status

TEE - EEPROM test

Test of the parameter memory.

Transmit: **TEE** <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ> Starts the test (duration <1 s)



Do not keep repeating the test (EEPROM life).

Receive: aaaa <CR><LF>

	Description
aaaa	Error word

TEP - FLASH test

Test of the program memory.

Transmit: **TEP** <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ> Starts the test (very brief)

Receive: aaaa,bbbbbbbb <CR><LF>

	Description
aaaa	Error word
bbbbbbbbbb	Check sum (hex)

TIO - I/O test



Caution



Caution: The relays switch irrespective of the pressure.

Starting a test program may cause unwanted effects in connected control systems.

Disconnect the sensor cable and control system lines to ensure that no control commands or messages are triggered by mistake.

Vacuum Gauge Controller AGC-100

Transmit: **TIO** [,a,b] <CR>[<LF>]

	Description
a	Test status, a = 0 → Off 1 → On
b	Relay status (in hex format), bb = 00 → All relays deactivated 01 → Switching function relay 1 activated 02 → Switching function relay 2 activated 40 → Error relay activated 4F → All relays activated

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a,b <CR><LF>

	Description
a	I/O test status
b	Relay status

TKB - Operator key test

Transmit: **TKB** <CR>[<LF>]

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: abcd <CR><LF>

	Description
a	Key 1, a = 0 → Not pushed 1 → Pushed
b	Key 2, b = 0 → Not pushed 1 → Pushed
c	Key 3, c = 0 → Not pushed 1 → Pushed
d	Key 4, d = 0 → Not pushed 1 → Pushed

TLC - Torr lock

Transmit: **TLC** [,a] <CR>[<LF>]

	Description
a	Torr lock, a = 0 → Off (default) 1 → On

Receive: <ACK><CR><LF>

Transmit: <ENQ>

Receive: a <CR><LF>

	Description
a	Torr lock status

Vacuum Gauge Controller AGC-100

TMP - Inner Temperature of the Unit

Transmit: **TMP** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: aa <CR><LF>

	Description
aa	Temperature (± 2 °C) [°C]

TRS - Serial Interface test

Transmit: **TRS** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ> Starts the test (repeats each character, test is interrupted with <CTRL> C).

WDT - Watchdog control

Transmit: **WDT** [,a] <CR>[<LF>]

	Description
a	Watchdog control, a = 0 → Manual error acknowledgement 1 → Automatic error acknowledgement ¹⁾ (default)



¹⁾ If the watchdog has responded, the error is automatically acknowledged and cancelled after 2 s.

Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a <CR><LF>

	Description
a	Watchdog control

5.12 Further

AYT - Are you there?

Transmit: **AYT** <CR>[<LF>]
 Receive: <ACK><CR><LF>
 Transmit: <ENQ>
 Receive: a,b,c,d,e <CR><LF>

	Description
a	Type of the unit, e.g. AGC-100
b	Model No. of the unit, e.g. AGC100
c	Serial No. of the unit, e.g. LI1605F001
d	Firmware version of the unit, e.g. 1.00
e	Hardware version of the unit, e.g. 1.0

Vacuum Gauge Controller AGC-100

ETH - Ethernet configuration

Transmit: **ETH** [,a,bbb.bbb.bbb.bbb,ccc.ccc.ccc.ccc,ddd.ddd.ddd.ddd]
<CR>[<LF>]
Receive: <ACK><CR><LF>
Transmit: <ENQ>
Receive: a,bbb.bbb.bbb.bbb,ccc.ccc.ccc.ccc,ddd.ddd.ddd.ddd <CR><LF>

	Description
a	DHCP (dynamic host configuration protocol), a = 0 -> Statically 1 -> Dynamically
bbb.bbb.bbb.bbb	IP address
ccc.ccc.ccc.ccc	Subnet address
ddd.ddd.ddd.ddd	Gateway address

5.13 Example



"Transmit (T)" and "Receive (R)" are related to host.

T: TID <CR> [<LF>]	Request for gauge identification
R: <ACK> <CR> <LF>	Positive acknowledgement
T: <ENQ>	Request for data transmission
R: PVG <CR> <LF>	Gauge identification
T: SP1 <CR> [<LF>]	Request for parameters of switching function 1 (setpoint 1)
R: <ACK> <CR> <LF>	Positive acknowledgement
T: <ENQ>	Request for data transmission
R: 1,1.0000E-09,9.0000E-07 <CR> <LF>	Thresholds
T: SP1 ,1,6.80E-3,9.80E-3 <CR> [<LF>]	Modification of parameters of switching function 1 (setpoint 1)
R: <ACK> <CR> <LF>	Positive acknowledgement
T: FOL ,2 <CR> [<LF>]	Modification of filter time constant (syntax error)
R: <NAK> <CR> <LF>	Negative acknowledgement
T: <ENQ>	Request for data transmission
R: 0001 <CR> <LF>	ERROR word
T: FIL ,2 <CR> [<LF>]	Modification of filter time constant
R: <ACK> <CR> <LF>	Positive acknowledgement
T: <ENQ>	Request for data transmission
R: 2 <CR> <LF>	Filter time constants
T: PR1 <CR> [<LF>]	Request for measurement data
R: <ACK> <CR> <LF>	Positive acknowledgement
T: <ENQ>	Request for data transmission
R: 0,8.3400E-03 <CR> <LF>	Status and pressure
T: <ENQ>	Request for data transmission
R: 1,8.0000E-04 <CR> <LF>	Status and pressure

6 Maintenance

Cleaning the AGC-100

For cleaning the outside of the unit a slightly moist cloth will usually do. Do not use any aggressive or scouring cleaning agents.

DANGER

DANGER: mains voltage
 Contact with live parts is extremely hazardous when liquids penetrate into the unit.

Make sure no liquids penetrate into the equipment.

Battery replacement





The product contains a battery (type CR2032, service life >10 years) in order to maintain the data integrity of the real-time clock. Battery replacement is necessary if the real-time clock repeatedly shows an incorrect date. Please contact your local Agilent Technologies service center.

7 Troubleshooting

Signalization of errors

The error is shown in the dot matrix and the error relay opens (→ 20).

Error messages

	Possible cause and remedy/acknowledgement
SENSOR ERROR	<p>Interruption or instability in sensor line or connector (Sensor error).</p> <p>⇒ Acknowledge with the  key.</p>
WATCHDOG ERROR	<p>The AGC-100 has been turned on too fast after power off.</p> <p>⇒ Acknowledge with the  key.</p> <p>If the watchdog is set to Auto, the AGC-100 acknowledges the message automatically after 2 s (→ 48).</p> <p>The watchdog has tripped because of a severe electric disturbance or an operating system error.</p> <p>⇒ Acknowledge with the  key.</p> <p>If the watchdog is set to WATCHDOG AUTO, the AGC-100 acknowledges the message automatically after 2 s (→ 48).</p>
DATA CORRUPTED	<p>Parameter memory error (EEPROM).</p> <p>⇒ Acknowledge with the  key.</p>

Technical support



If the problem persists after the message has been acknowledged several times and/or the gauge has been exchanged, please contact your nearest Agilent Technologies service center.

8 Repair


Return defective products to your nearest Agilent Technologies service center for repair.


Agilent Technologies assumes no liability and the warranty is rendered null and void if repair work is carried out by the end-user or by third parties.

9 Accessories

	Ordering number
Adapter panel for installation into a 19" rack chassis adapter, height 3 U	AGC100ADPT

10 Storage


Caution





Caution: electronic components.

Inappropriate storage (static electricity, humidity etc.) may damage electronic components.

Store the product in an antistatic bag or container. Observe the relevant specifications under Technical Data (→ 10).

11 Disposal


WARNING



WARNING: substances detrimental to the environment.

Products or parts thereof (mechanical and electric components, operating fluids etc.) may be detrimental to the environment.

Please dispose of such materials in accordance with the relevant local regulations.

Separating the components

Electronic and non-electronic components

After disassembling the product, separate its components in accordance with the following criteria:

Such components must be separated according to their materials and recycled.

Appendix

A: Conversion Tables

Weights

	kg	lb	slug	oz
kg	1	2.205	68.522×10^{-3}	35.274
lb	0.454	1	31.081×10^{-3}	16
slug	14.594	32.174	1	514.785
oz	28.349×10^{-3}	62.5×10^{-3}	1.943×10^{-3}	1

Pressures

	N/m ² , Pa	Bar	mBar, hPa	Torr	at
N/m ² , Pa	1	10×10^{-6}	10×10^{-3}	7.5×10^{-3}	9.869×10^{-6}
Bar	100×10^3	1	10^3	750.062	0.987
mBar, hPa	100	10^{-3}	1	750.062×10^{-3}	0.987×10^{-3}
Torr	133.322	1.333×10^{-3}	1.333	1	1.316×10^{-3}
at	101.325×10^3	1.013	1.013×10^3	760	1

Pressure units used in the vacuum technology

	mBar	Bar	Pa	hPa	kPa	Torr mm HG
mBar	1	1×10^{-3}	100	1	0.1	0.75
Bar	1×10^3	1	1×10^5	1×10^3	100	750
Pa	0.01	1×10^{-8}	1	0.01	1×10^{-3}	7.5×10^{-3}
hPa	1	1×10^{-3}	100	1	0.1	0.75
kPa	10	0.01	1×10^3	10	1	7.5
Torr mm HG	1.332	1.332×10^{-3}	133.32	1.3332	0.1332	1

$$1 \text{ Pa} = 1 \text{ N/m}^2$$

Linear measurements

	mm	m	inch	ft
mm	1	10^{-3}	39.37×10^{-3}	3.281×10^{-3}
m	10^3	1	39.37	3.281
inch	25.4	25.4×10^{-3}	1	8.333×10^{-2}
ft	304.8	0.305	12	1

Temperature

	Kelvin	Celsius	Fahrenheit
Kelvin	1	$^{\circ}\text{C} + 273.15$	$(^{\circ}\text{F} + 459.67) \times 5/9$
Celsius	$\text{K} - 273.15$	1	$5/9 \times ^{\circ}\text{F} - 17.778$
Fahrenheit	$9/5 \times \text{K} - 459.67$	$9/5 \times (^{\circ}\text{C} + 17.778)$	1

B: Firmware Update



If your AGC-100 firmware needs updating, e.g. for implementing a new gauge type, please contact your nearest Agilent Technologies service center.

A firmware update is possible

- via a USB memory stick (type A connector on the front of the unit), or
- with the USB Update Tool via the USB type B connector on the rear of the unit.

User Parameters

Most of the settings you may have made in the Parameter mode will not be affected by a firmware update. However, we recommend that you save the parameters before an update (→ 52).

Firmware update with a USB memory stick (type A)



Not all USB memory sticks are automatically recognized by the AGC-100, as they (particularly cheaper brands) do not always conform to USB standard specifications. Try a different memory stick before contacting your nearest Agilent Technologies service center.

1 Download two files with the ending ".S19" and ".CNF" from our website "www.agilent.com" to a USB memory stick.

2 Switch off the unit.

3 Plug in the memory stick and then turn on the unit.

4 The update occurs automatically in the following steps:

BOOTING	Very brief.
BOOTLOADER V1.x	Very brief.
ERASING FW...	Old firmware is being deleted from the unit.
UPDATING FW...	New firmware is being loaded onto the unit.
UPDATE COMPLETE	Update completed.

5 Remove the memory stick and the unit will restart automatically.

6 If necessary, customer-specific settings saved before the update may now be resaved to the unit (→ 52).

Firmware update with USB Update Tool (USB type B)

Precondition: Microsoft Windows XP, 7, 8 or 10 operating system



Your operating system should be updated first. Additionally administrator rights are required.



During firmware update, no USB memory stick should be connected on the front of the unit.

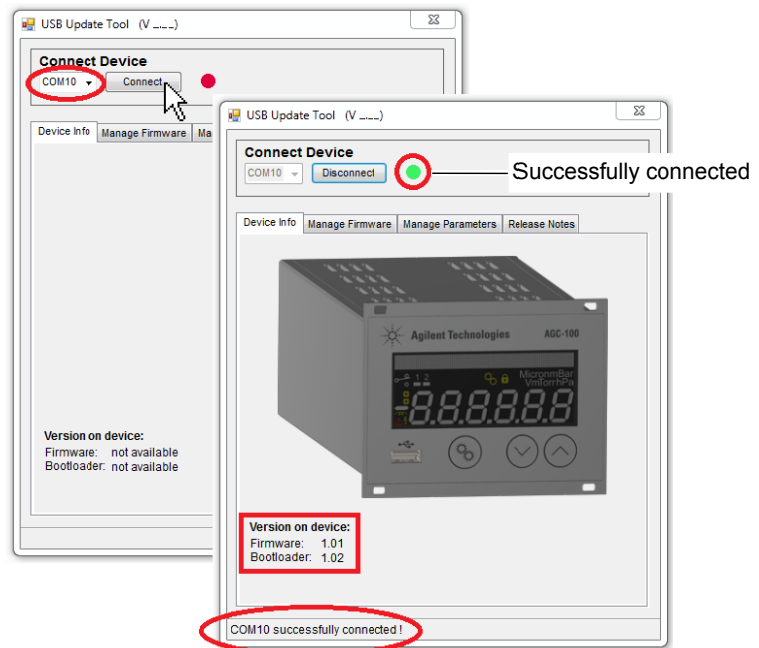


If a virtual serial interface (COM) is not automatically established, you may download and then install the driver from the website "www.ftdichip.com/drivers/vcp.htm".

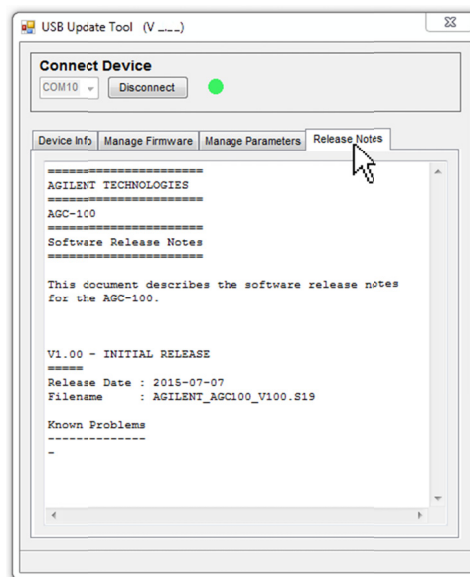
1 Download the USB UpdateTool from the CD ROM or from our website "www.agilent.com".

Vacuum Gauge Controller AGC-100

- 2 Using a USB cable type A/B connect the unit to the computer.
- 3 Start USB UpdateTool, select the COM interface from the menu and click on <Connect>.



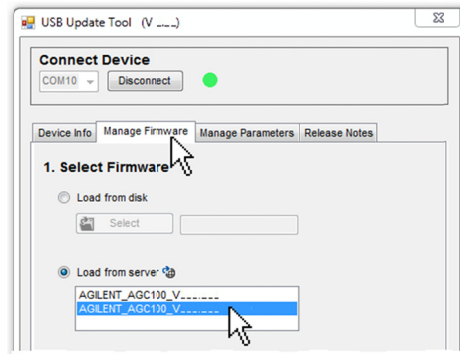
- 4 Click on <Release Notes> to view the software release notes.



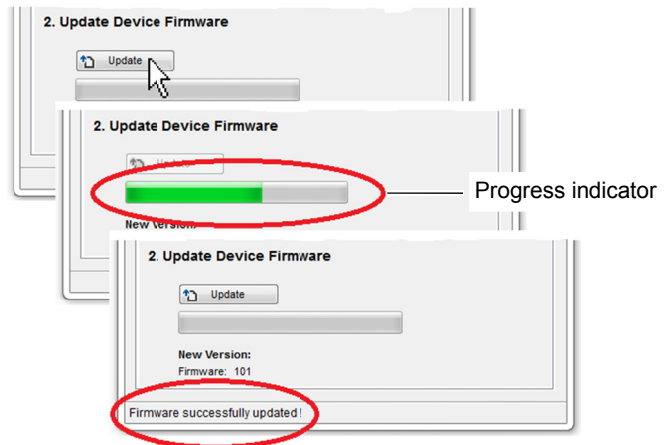
Vacuum Gauge Controller AGC-100

5 Click on <Manage Firmware>, select firmware ...

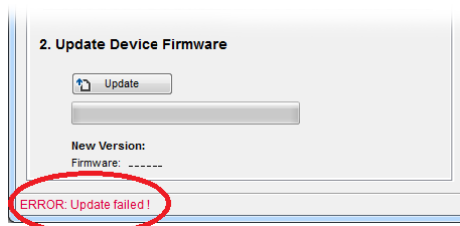
- Option <Load from disk>: Download a copy of the firmware from our website "www.agilent.com". Then, select the appropriate folder.
- Option <Load from server>: The update tool connects to the internet. Select the desired firmware version from the selection list.



... and click <Update>: The firmware is updated.



If the update was not successful, try again.



C: Ethernet Configuration

The user program (e.g. terminal program, LabView, etc.) must support serial interfaces. Under Microsoft Windows operating systems the AGC-100 is listed as a virtual COM interface.



Please contact your network administrator, before starting Ethernet configuration.



Your operating system should be updated first. Additionally administrator rights are required.

C 1: Connect the AGC-100 to a Network

With registration

1

Readout the MAC address of the AGC-100 (→ 47).

2

The AGC-100 should be registered in the network by the network administrator. After registration ask him for the Ethernet parameters (IP ADDRESS, GATEWAY, NETMASK and DHCP).

3

Configuring the AGC-100:

- Save all AGC-100 parameters on a USB memory stick ("SAVE SETUP", → 52).
- Set the Ethernet parameters (IP ADDRESS, GATEWAY, NETMASK and DHCP) in the saved CSV file on the memory stick.
- Load the modified parameters onto the AGC-100 ("RESTORE SETUP", → 52).
- Connect the AGC-100 with an Ethernet patch cable to the network.

4

Search for the AGC-100 in the network using the Ethernet Configuration Tool and assign it to a virtual COM interface (→ 87).

5

Start the program for communication with the AGC-100 and connect it to the assigned COM interface.

Without registration

1

If unknown, ask the network administrator for the Ethernet parameters (IP ADDRESS, GATEWAY, NETMASK and DHCP).

2

Configuring the AGC-100:

- Save all AGC-100 parameters on a USB memory stick ("SAVE SETUP", → 52).
- Set the Ethernet parameters (IP ADDRESS, GATEWAY, NETMASK and DHCP) in the saved CSV file on the memory stick.
- Load the modified parameters onto the AGC-100 ("RESTORE SETUP", → 52).
- Connect the AGC-100 with an Ethernet patch cable to the network.

3

Search for the AGC-100 in the network using the Ethernet Configuration Tool and assign it to a virtual COM interface (→ 87).

4

Start the program for communication with the AGC-100 and connect it to the assigned COM interface.

C 2: Connect the AGC-100 to a Computer

Computer with DHCP server

- ➊ Connect the AGC-100 to a computer ...
 - with a crossover Ethernet cable,
 - via a switch, or
 - with an Ethernet patch cable (precondition: the interface is auto MDI-X capable).
- ➋ The DHCP server assigns automatically an IP address.
Precondition: DHCP = ON (standard)
- ➌ Search for the AGC-100 in the network using the Ethernet Configuration Tool and assign it to a virtual COM interface (→ 87).
- ➍ Start the program for communication with the AGC-100 and connect it to the assigned COM interface.

Computer without DHCP server

- ➊ Save all AGC-100 parameters on a USB memory stick ("SAVE SETUP", → 52).
- ➋ Set the following Ethernet parameters in the saved CSV file on the memory stick:
 IP ADDRESS: 192.168.0.1 (192.168.0.2 for a second unit, and so on)
 NETMASK: 255.255.0.0
 DHCP: OFF
- ➌ Load the modified parameters onto the AGC-100 ("RESTORE SETUP", → 52).
- ➍ Connect the AGC-100 to a computer ...
 - with a crossover Ethernet cable,
 - via a switch, or
 - with an Ethernet patch cable (precondition: the interface is auto MDI-X capable).
- ➎ Search for the AGC-100 in the network using the Ethernet Configuration Tool and assign it to a virtual COM interface (→ 87).
- ➏ Start the program for communication with the AGC-100 and connect it to the assigned COM interface.

C 3: Ethernet Configuration Tool

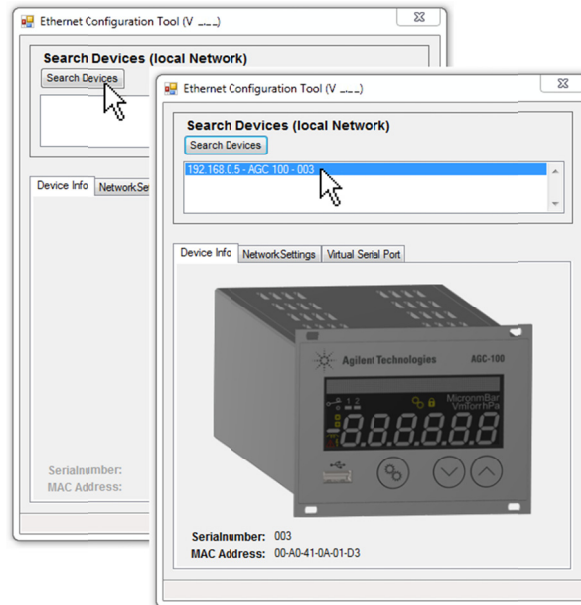
With the Ethernet Configuration Tool a virtual serial interface (COM) can be assigned to an IP address. In addition, it allows configuration of the Ethernet interface via a computer.

Precondition: Windows 7 or 8 operating system (does not work under Windows XP)

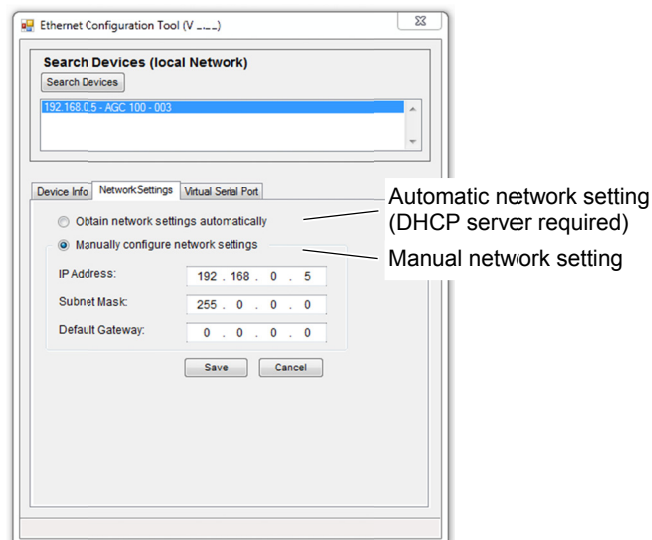
- ➊ Download the Ethernet Configuration Tool from the CD ROM or from our website "www.agilent.com".

Vacuum Gauge Controller AGC-100

- 2 Start the Ethernet Configuration Tool and click on <Search Devices>: the Tool searches the local network for connected devices and lists the devices thus found in the selection window. The <Device Info> register shows basic information about the selected device.

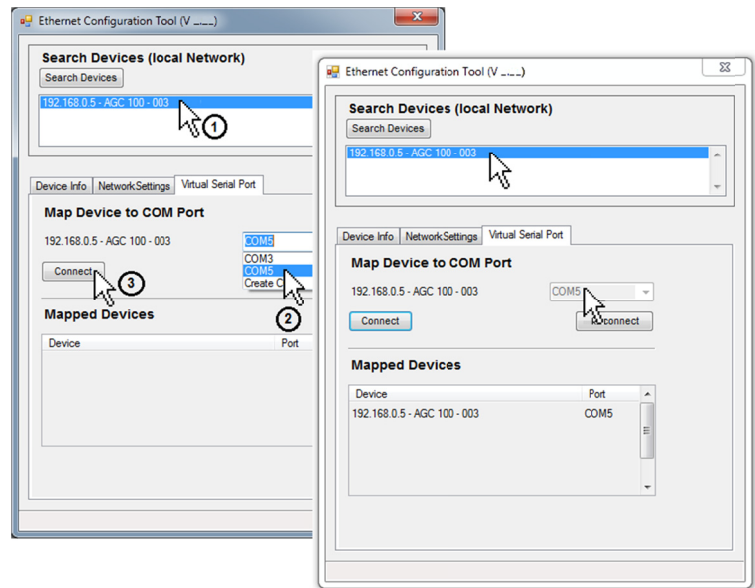


- 3 Automatic or manual network setting occurs in the <Network Settings> register.

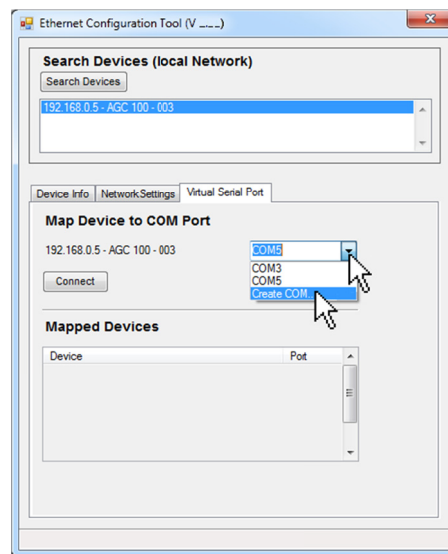


Vacuum Gauge Controller AGC-100

- 4 In the <Virtual Serial Port> register a specific COM Port can be assigned to each device, and/or ...











... a new COM Port can be created.



The new created virtual interface (COM) appears in the list box and in the Windows Device Manager.

D: Literature

-  [1] www.agilent.com
Operating Instructions
Inverted Magnetron Pirani Gauge FRG-700, FRG-702
tqna74e1
Agilent Technologies, Lexington, MA 02421, USA
-  [2] www.agilent.com
Operating Instructions
Bayard-Alpert Pirani Gauge FRG-720
tqna70e1
Agilent Technologies, Lexington, MA 02421, USA
-  [3] www.agilent.com
Operating Instructions
Bayard-Alpert Pirani Gauge FRG-730
tqna72e1
Agilent Technologies, Lexington, MA 02421, USA
-  [4] www.agilent.com
Operating Instructions
Pirani Standard Gauge PVG-500, PVG-502
tqna69e1
Agilent Technologies, Lexington, MA 02421, USA
-  [5] www.agilent.com
Operating Instructions
Pirani Capacitance Diaphragm Gauge PCG-750, PCG-752
tqna77e1
Agilent Technologies, Lexington, MA 02421, USA
-  [6] www.agilent.com
Operating Instructions
Pirani Standard Gauge PVG-550, PVG-552
tqna79e1
Agilent Technologies, Lexington, MA 02421, USA
-  [7] www.agilent.com
Operating Instructions
Capacitance Diaphragm Gauge CDG-500
tqna76e1
Agilent Technologies, Lexington, MA 02421, USA
-  [8] www.agilent.com
Operating Instructions
Compact Cold Catode Gauge IMG-500
tqna75e1
Agilent Technologies, Lexington, MA 02421, USA

ETL Certification



ETL LISTED

The product AGC-100

- conforms to the UL Standards UL 61010-1 and UL 61010-2-030
- is certified to the CSA Standards CSA C22.2 # 61010-1 and CSA C22.2 # 61010-2-030

Notes

Vacuum Products Division

Instructions for returning products

Dear Customer:

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

- 1) Complete the attached Request for Return form and send it to Agilent Technologies (see below), taking particular care to identify all products that have pumped or been exposed to any toxic or hazardous materials.
- 2) After evaluating the information, Agilent Technologies will provide you with a Return Authorization (RA) number via email or fax, as requested.
Note: Depending on the type of return, a Purchase Order may be required at the time the Request for Return is submitted. We will quote any necessary services (evaluation, repair, special cleaning, eg).
- 3) Important steps for the shipment of returning product:
 - Remove all accessories from the core product (e.g. inlet screens, vent valves).
 - Prior to shipment, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
 - If ordering an Advance Exchange product, 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.
 - Agilent Technologies is not responsible for returning customer provided packaging or containers.
 - Clearly label package with RA number. Using the shipping label provided will ensure the proper address and RA number are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will be returned.
- 4) Return only products for which the RA was issued.
- 5) Product being returned under a RA must be received within 15 business days.
- 6) Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information. Customer is responsible for freight charges on returning product.
- 7) Return shipments must comply with all applicable Shipping Regulations (IATA, DOT, etc.) and carrier requirements.

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

EUROPE:
Fax: 00 39 011 9979 330
Fax Free: 00 800 345 345 00
Toll Free: 00 800 234 234 00
vpt-customer-care@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>



Please read important policy information on Page 3 that applies to all returns.

1) CUSTOMER INFORMATION

Company Name:		Contact Name:	
Tel:	Email:	Fax:	
Customer Ship To:		Customer Bill To:	
Europe only: VAT reg. Number:		USA/Canada only: <input type="checkbox"/> Taxable <input type="checkbox"/> Non-taxable	

2) PRODUCT IDENTIFICATION

Product Description	Agilent P/N	Agilent S/N	Original Purchasing Reference

3) TYPE OF RETURN (Choose one from each row and supply Purchase Order if requesting a billable service)

- 3A. ☐ Non-Billable ☐ Billable ☒ New PO # (hard copy must be submitted with this form):
- 3B. ☐ Exchange ☐ Repair ☐ Upgrade ☐ Consignment/Demo ☐ Calibration ☐ Evaluation ☐ Return for Credit

4) HEALTH and SAFETY CERTIFICATION

AGILENT TECHNOLOGIES CANNOT ACCEPT ANY PRODUCTS CONTAMINATED WITH BIOLOGICAL OR EXPLOSIVE HAZARDS, RADIOACTIVE MATERIAL, OR MERCURY AT ITS FACILITY.

Call Agilent Technologies to discuss alternatives if this requirement presents a problem.

The equipment listed above (check one):

- ☐ HAS NOT pumped or been exposed to any toxic or hazardous materials. OR
- ☐ HAS pumped or been exposed to the following toxic or hazardous materials. If this box is checked, the following information must also be filled out. Check boxes for all materials to which product(s) pumped or was exposed:

☐ Toxic ☐ Corrosive ☐ Reactive ☐ Flammable ☐ Explosive ☐ Biological ☐ Radioactive

List all toxic/hazardous materials. Include product name, chemical name, and chemical symbol or formula:

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.

Print Name: _____ Authorized Signature: _____ Date: _____

5) FAILURE INFORMATION:

Failure Mode (REQUIRED FIELD. See next page for suggestions of failure terms):

Detailed Description of Malfunction: (Please provide the error message)

Application (system and model):

I understand and agree to the terms of Section 6, Page 3/3.

Print Name: _____ Authorized Signature: _____ Date: _____



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

TURBO PUMPS and TURBO CONTROLLERS

APPARENT DEFECT/MALFUNCTION	POSITION	PARAMETERS
<ul style="list-style-type: none"> - Does not start - Does not spin freely - Does not reach full speed - Mechanical Contact - Cooling defective 	<ul style="list-style-type: none"> - Noise - Vibrations - Leak - Overtemperature - Clogging 	<ul style="list-style-type: none"> - Vertical - Horizontal - Upside-down - Other: _____
		Power: _____ Rotational Speed: _____ Current: _____ Inlet Pressure: _____ Temp 1: _____ Foreline Pressure: _____ Temp 2: _____ Purge flow: _____ OPERATING TIME: _____

ION PUMPS/CONTROLLERS

<ul style="list-style-type: none"> - Bad feedthrough - Vacuum leak - Error code on display 	<ul style="list-style-type: none"> - Poor vacuum - High voltage problem - Other
---	--

VALVES/COMPONENTS

<ul style="list-style-type: none"> - Main seal leak - Solenoid failure - Damaged sealing area 	<ul style="list-style-type: none"> - Bellows leak - Damaged flange - Other
--	---

LEAK DETECTORS

<ul style="list-style-type: none"> - Cannot calibrate - Vacuum system unstable - Failed to start 	<ul style="list-style-type: none"> - No zero/high background - Cannot reach test mode - Other
---	--

INSTRUMENTS

<ul style="list-style-type: none"> - Gauge tube not working - Communication failure - Error code on display 	<ul style="list-style-type: none"> - Display problem - Degas not working - Other
--	---

SCROLL AND ROTARY VANE PUMPS

<ul style="list-style-type: none"> - Pump doesn't start - Doesn't reach vacuum - Pump seized 	<ul style="list-style-type: none"> - Noisy pump (describe) - Over temperature - Other
---	--

DIFFUSION PUMPS

<ul style="list-style-type: none"> - Heater failure - Doesn't reach vacuum - Vacuum leak 	<ul style="list-style-type: none"> - Electrical problem - Cooling coil damage - Other
---	--

Section 6) ADDITIONAL TERMS

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

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

Service & Support

North America

Agilent Technologies
Vacuum Products Division
121 Hartwell Avenue
Lexington, MA 02421 USA
Tel.: +1 781 881 7200
Toll-Free: +1 800 882 7426
Fax: +1 781 880 5437

Benelux

Agilent Technologies Netherlands B.V.
Vacuum Products Division
Herculesweg 8
4338 PL Middelburg
The Netherlands
Tel: +31 118 671570
Fax: +31 118 671569
Toll free: 00 800 234 234 00

Canada

Central coordination through
Agilent Technologies
Vacuum Products Division
121 Hartwell Avenue
Lexington, MA 02421 USA
Tel.: +1 781 881 7200
Toll-Free: +1 800 882 7426
Fax: +1 781 880 5437

China

Agilent Technologies (China) Co. Ltd
Vacuum Products Division
No.3, Wang Jing Bei Lu, Chao Yang
District
Beijing, 100102 China
Tel.: +86 (0)10 64397888
Fax: +86 (0)10 64391318
Toll-Free: 800 820 3278

France

Agilent Technologies France
Vacuum Products Division
3 avenue du Canada, Parc Technopolis
Z.A. Courtaboeuf - CS 90263
91978 Les Ulis cedex
France
Tel.: +33 (0)1 64 53 61 15
Fax: +33 (0)1 64 53 50 01
Toll free: 00 800 234 234 00
e-Mail: vpf.sales@agilent.com

Germany & Austria

Agilent Technologies
Sales & Services GmbH & Co. KG
Vacuum Products Division
Lyoner Str. 20
60 528 Frankfurt am Main
Germany
Tel.: +49 69 6773 43 2230
Fax: +49 69 6773 43 2250

This information is subject to change
without notice.

© Agilent Technologies, Inc., 2016

India

Agilent Technologies India Pvt. Ltd.
Vacuum Products Division
G01, Prime corporate Park,
230/231, Sahar Road,
Opp. Blue Dart Centre,
Andheri (East), Mumbai – 400 099, India
Tel: +91 22 30648287/8200
Fax: +91 22 30648250
Toll Free: 1800 113037

Italy

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

Japan

Agilent Technologies Japan, Ltd.
Vacuum Products Division
8th Floor, Sumitomo Shibaura Building
4-16-36 Shibaura Minato-ku
Tokyo 108-0023 JAPAN
Tel.: +81 3 5232 1253
Toll-Free: 0120 655 040
Fax: +81 3 5232 1710

Korea

Agilent Technologies
Vacuum Products Division
Shinsa 2nd Bldg. 2F
986-5 Daechi-dong
Kangnam-gu, Seoul KOREA 135-280
Tel.: +82 (0)2 3452 2455
Toll-Free: 080 222 2452
Fax: +82 (0)2 3452 3947

Singapore

Agilent Technologies Singapore Pte. Ltd
Vacuum Products Division
Agilent Technologies Building,
No.1 Yishun Avenue 7
Singapore 768923
Tel: +65 6215 8045
Fax: +65 6754 0574

Southeast Asia

Agilent Technologies Sales Sdn Bhd
Vacuum Products Division
Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya,
Selangor, Malaysia
Tel : +603 7712 6106
Fax: +603 6733 8121

Taiwan

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

UK & Ireland

Agilent Technologies
Vacuum Products Division
6 Mead Road
Oxford Industrial Park
Yarnton, Oxford OX5 1QU UK
Tel.: +44 (0) 1865 291570
Fax: +44 (0) 1865 291571
Toll free: 00 800 234 234 00

Other Countries

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

Customer Service and Support

North America
Toll Free: 800 882 7426, Option 3
vpl-ra@agilent.com

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

Pacific Rim:
Please visit our website for individual
office information <http://www.agilent.com>

Worldwide Web Site, Catalog and Order

On-line:
www.agilent.com
Representatives in most countries



Agilent Technologies



1998061