



Agilent Technologies

Vacuum Products Division



Communications Protocol

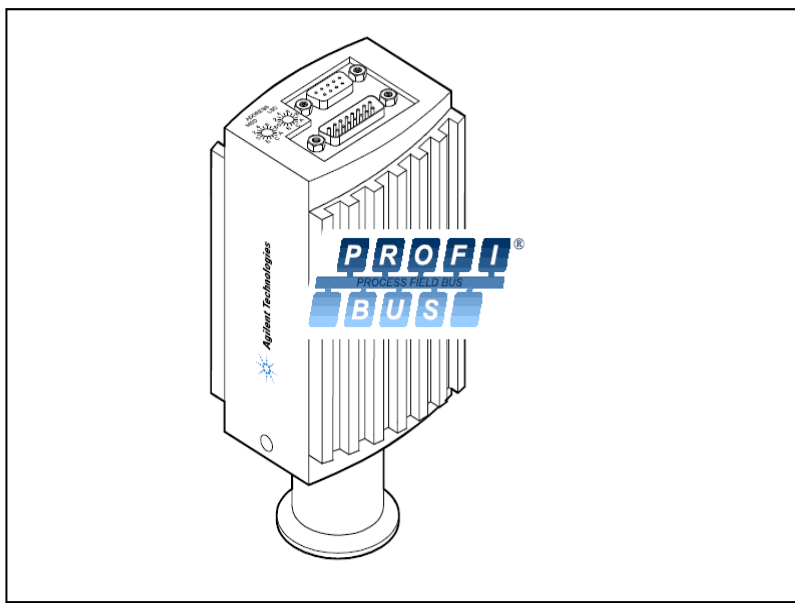
Profibus

***DP/V1 Interface for Bayard-Alpert Pirani
Gauge
FRG-730***

Manual No. TQRa73e1
Revision 3
March 2012

Profibus

DP/V1 Interface for Bayard-Alpert Pirani Gauge
FRG-730



Contents

About this Document	5
Product Identification	5
Validity	5
Intended Use	5
Functional Principle	5
Trademarks	5
1 General Data	6
1.1 Data Rate	6
1.2 Device Address	6
1.3 Ident Number	6
1.4 Configuration Data	7
1.5 User Parameter Data	7
1.6 Types of Communication	7
2 Data Exchange Mode	8
2.1 Acyclic Data Transmission with Profibus DPV1 Functionality	8
2.2 Structure of the Cyclic Data Telegrams in Data Exchange Mode	9
2.2.1 Parameter Channel	10
2.2.1.1 PKE Parameter Signature Value	10
2.2.1.2 PWE Parameter (Process Value)	11
2.2.1.3 Error Code (Error Message)	11
2.3 Cyclic Message Telegrams	12
3 Block Model	12
3.1 Device Block	14
3.1.1 Information on the Individual Indices	15
3.1.1.1 Block Type ID 16	15
3.1.1.2 Device Type ID 17	15
3.1.1.3 Standard Revision Level ID 18	15
3.1.1.4 Device Manufacturer Identifier ID 19	15
3.1.1.5 Manufacturer Model Number ID 20	15
3.1.1.6 Software or Firmware Revision Level ID 21	15
3.1.1.7 Hardware Revision Level ID 22	15
3.1.1.8 Device State ID 25	15
3.1.1.9 Exception Status ID 26	16
3.1.1.10 Exception Detail Alarm ID 27	17
3.1.1.11 Exception Detail Warning ID 28	19
3.1.1.12 Copy Common Exception Detail Alarm 0 ID 204	21
3.1.1.13 Copy Device Exception Detail Alarm 0 ... 3 ID 205	21
3.1.1.14 Copy Manufacturer Exception Detail Alarm 0 ID 207	21
3.1.1.15 Copy Common Exception Detail Warning 0 ID 208	21
3.1.1.16 Copy Device Exception Detail Warning 0 ... 3 ID 209	21
3.1.1.17 Copy Device Exception Detail Warning 4 ID 210	21
3.1.1.18 Copy Manufacturer Exception Detail Warning 0 ID 211	21
3.1.2 Device Block, Device Behavior	22
3.1.2.1 Device Block State Command	23
3.2 Analog Input Block	24
3.2.1 One Of N Analog Input Function Block / SLOT 1	24
3.2.1.1 AI Block Adjust Command ID 15	24
3.2.1.2 Block Type ID 16	24
3.2.1.3 Channel Instance Selector ID 46	24
3.2.1.3 PV Selector ID 47	24
3.2.2 Analog Sensor Input Function Block SLOT 1 / Instance 1	25
3.2.2.1 Process Value ID 19	25
3.2.2.2 Status ID 20	25
3.2.2.3 Data Type ID 21	25
3.2.2.4 Data Unit ID 22	26
3.2.2.5 Reading Valid ID 23	26
3.2.2.6 Full Scale ID 24	26
3.2.2.7 Safe State ID 39	27
3.2.2.8 Safe Value ID 40	27
3.2.2.9 Overrange ID 44	27
3.2.2.10 Underrange ID 45	27
3.2.2.11 AI Block Adjust Command (Pirani) ID 15	27

3.2.3	Analog Sensor Input Function Block SLOT 1 / Instance 2	28
3.2.3.1	Process Value ID 19	28
3.2.3.2	Status ID 20	28
3.2.3.3	Data Type ID 21	28
3.2.3.4	Data Unit ID 22	29
3.2.3.5	Reading Valid ID 23	29
3.2.3.6	Full Scale ID 24	29
3.2.3.7	Safe State ID 39	30
3.2.3.8	Safe Value ID 40	30
3.2.3.9	Overrange ID 44	30
3.2.3.10	Underrange ID 45	30
3.2.4	Analog Sensor Input Function Block SLOT 1 / Instance 3, 4	31
3.2.4.1	Process Value ID 19	31
3.2.4.2	Status ID 20	31
3.2.4.3	Data Type ID 21	31
3.2.4.4	Data Unit ID 22	32
3.2.4.5	Reading Valid ID 23	32
3.3	Transducer Block	33
3.3.1	One Of N Vacuum Gauge Transducer Block / SLOT 1	33
3.3.1.1	One Of N Status Extension	33
3.3.2	Heat Transfer Vacuum Gauge Transducer Block / SLOT 1 / Instance 1	33
3.3.2.1	Block Type ID 101	33
3.3.2.2	Status Extension ID 102	33
3.3.2.3	Sensor Alarm ID 103	33
3.3.2.4	Sensor Warning ID 104	34
3.3.3	Hot Cathode Ion Gauge Transducer Block / SLOT 1 / Instance 2	34
3.3.3.1	Block Type ID 101	34
3.3.3.2	Status Extension ID 102	34
3.3.3.3	Sensor Alarm ID 103	35
3.3.3.4	Sensor Warning ID 104	35
3.3.3.5	Emission Status ID 105	35
3.3.3.6	Emission Current ID 106	36
3.3.3.7	Active Filament ID 108	36
3.3.3.8	Degas Status ID 109	36
3.3.3.9	Active Degas Filament ID 116	36
3.3.3.10	Mode Filament Selection ID 119	36
3.3.3.11	Emission User Mode State ID 201	36
3.3.3.12	Hot Cathode Block State Command ID 14	37
Appendix A: Definitions		38
Appendix B: Block Type		41
Appendix C: Electrical Connections		42
Appendix D: Literature		44

For cross-references to other documents, the symbol (→  [XY]) is used.

About this Document

This document describes the functionality and programming of the Profibus interface of the FRG-730 gauge.



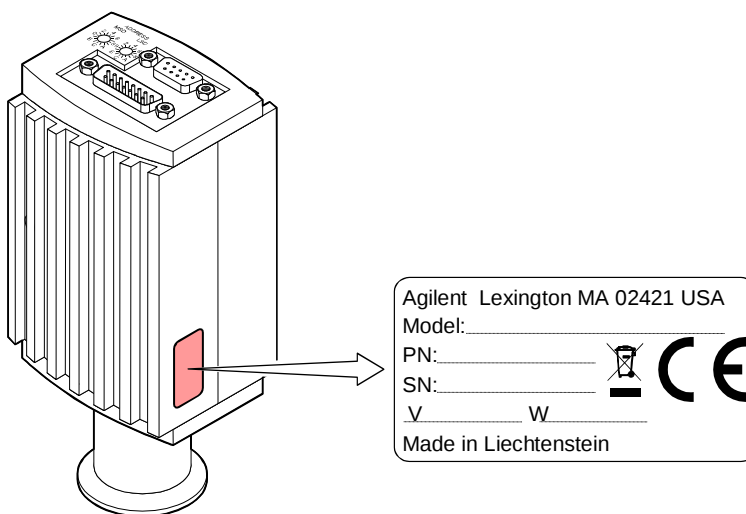
For safety information on and technical data of the gauges, please refer to the respective operating manuals (→ [1], [2]).

In information referring to the ionization vacuum measuring part of the gauge, the short designation "BA" (Bayard-Alpert measuring principle) is used.

The designation "Pirani" is used in information referring to the Pirani vacuum measuring part of the gauge.

Product Identification

In all communications with Agilent, please specify the information on the product nameplate. For convenient reference copy that information into the space provided below.



Validity

This document applies to products with part numbers

FRG-730 (with Profibus interface and switching functions)

FRG730KF25SP (DN 25 ISO-KF)

FRG730CF35SP (DN 40 CF-R)

The part number (PN) can be taken from the product nameplate.

Intended Use

The FRG-730 gauge has been designed for vacuum measurement of gases in the pressure range of 5×10^{-10} ... 1000 mbar.

It must not be used for measuring flammable or combustible gases in mixtures containing oxidants (e.g. atmospheric oxygen) within the explosion range.

The gauge can be operated in connection with an Agilent AGC-100 Vacuum Gauge Controller, an Agilent Turbo AG Rack Controller, or with another controller.

Functional Principle

The function of the gauge is described in the operating manual (→ [1], [2]).

The integrated Profibus interface allows operating the gauge in connection with other suitable devices in a Profibus network according to the standard described in [3], [4].

Trademarks

SEMI® Semiconductor Equipment and Materials International, California

Profibus This document contains Profibus-specific information described and defined in the Profibus Profiles (→ [3], [5], [6]).

1 General Data

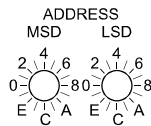
1.1 Data Rate

The gauge supports all data rates defined in the EN 50170 standard (→ [4]) up to 12 Mbaud. Automatic data rate setting is supported. Alternatively, a fixed data rate can be selected.

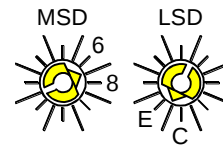
1.2 Device Address

The device address (= node address) must be set via two rotary switches when the gauge is installed.

For unambiguous identification of the gauge in a Profibus environment, a node address is required. The node address setting is made on the gauge.



The node address (0 ... 125_{dec}) is set in hexadecimal form (00 ... 7D_{hex}) (default value 5C_{hex}) via the "ADDRESS" switches. The "MSD" switch is used for setting the high-order address nibble and the "LSD" switch for defining the low-order address nibble.



Example: Node address = 7D_{hex}:

The node address is polled by the firmware when the gauge is switched on. If the setting deviates from the stored value, the new value is taken over into the NVRAM. If a value >7D_{hex} (>125_{dec}) is entered, the node address setting currently stored in the device remains valid. However, the address can be set via the Profibus master with the "Set Slave Address" service. This address setting will be stored in the EEPROM of the gauge.

1.3 Ident Number

The ident number assigned to the gauge by the PNO (→ [3]) is:

Gauge	Ident number (hexadecimal)
FRG-730	0x09AA

1.4 Configuration Data

Depending on the standard telegrams used (→ section "Cyclic Message Telegrams"), the following configuration data have to be transmitted to the gauge during the configuration phase:

Standard telegram Master ⇒ Slave	Standard telegram Slave ⇒ Master	Configuration data
-	4	0x44, 0x84, 0x05, 0x05, 0x05, 0x03
-	5	0x44, 0x86, 0x05, 0x05, 0x05, 0x08
1	4	0xC6, 0x81, 0x84, 0x05, 0x05, 0x05, 0x05, 0x05, 0x05, 0x03
1	5	0xC6, 0x81, 0x86, 0x05, 0x05, 0x05, 0x05, 0x05, 0x05, 0x08
3	6	0xC6, 0x87, 0x8C, 0x0A, 0x0A, 0x05, 0x05, 0x05, 0x03
3	7	0xC6, 0x87, 0x8E, 0x0A, 0x0A, 0x05, 0x05, 0x05, 0x08
2	6	0xC8, 0x89, 0x8C, 0x0A, 0x05, 0x05, 0x0A, 0x05, 0x05, 0x05, 0x03
2	7	0xC8, 0x89, 0x8E, 0x0A, 0x05, 0x05, 0x0A, 0x05, 0x05, 0x05, 0x08

1.5 User Parameter Data

Depending on the pressure unit setting (= data unit), the following configuration string has to be transmitted to the gauge (parameter data in hexadecimal format):

Pressure unit	User parameter data string
COUNTS ¹⁾	00 00 00 03 E9
Torr	00 00 00 05 15
Micron	00 00 00 05 16
mbar	00 00 00 05 1C
Pascal	00 00 00 05 1D

¹⁾ If COUNTS is selected as pressure unit, a value is output, which can be converted into a corresponding pressure value by means of a formula (→ section "Analog Sensor Input Function Block" for more information).

1.6 Types of Communication

FRG-730 works according to the Profibus DPV1 specification and can be addressed in cyclic or acyclic data traffic (→ [3]).

Acyclic data traffic should be used to make device or process specific settings such as definition of the Safe Values, Safe States etc. or for reading or writing of rarely used attributes.

Cyclic data traffic is used for continuous exchange of the required process parameter values, i.e. pressure value and status indications. A number of standard telegrams are available for cyclic data traffic. They can be selected according to requirements (→ section "Cyclic Message Telegrams").

2 Data Exchange Mode

2.1 Acyclic Data Transmission with Profibus DPV1 Functionality

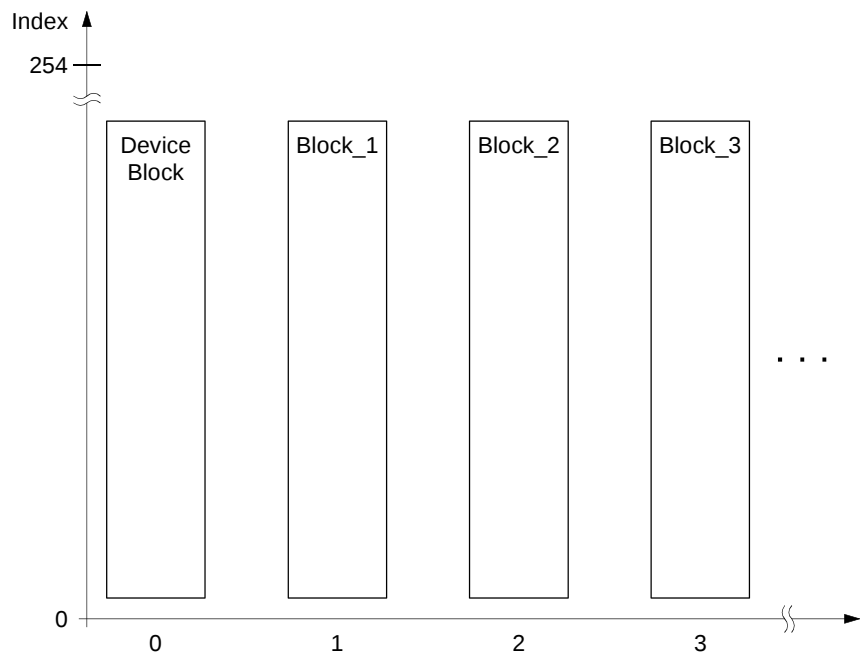
The reading and writing operations defined in Profibus are based on a slot index address scheme. In FRG-730, all device functions are organized in the following blocks:

- A device block describing all organizational parameters of the gauge (serial number, manufacturer, software version, ...)
- An Analog Sensor Function Block describing the function of the pressure presentation
- An Analog Sensor Transducer Block describing the physical interface between the gauge and the process (emission current, ion current, ...).

The block model is described in detail in section "Block Model".

Each block is assigned to a separate slot. The exact assignment Block $\Rightarrow$ Slot $\Rightarrow$ Index is described in section "Block Model". The Device Block is assigned to Slot 0, the transducer and functional blocks to Slot 1.

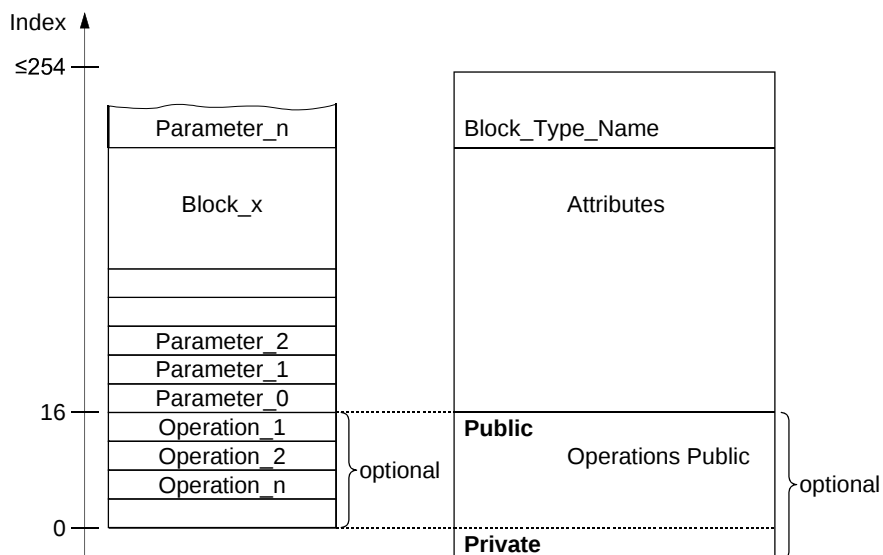
Block, slot and index assignment



There are 254 indices per slot. The indices can have a width of 255 bytes. All values that can be accessed via Profibus have to be mirrored to one of these slots/indices.

The parameters are generally numbered in ascending order, starting with index 16. Services such as "Degas On" or "Full scale" are numbered in descending order, starting with index 15.

Assignment of the block elements to the slot indices



The parameters of a block are numbered in an ascending order starting at index 16. Public operations are numbered in a descending order starting at index 15. Private block operations are hidden and therefore they will not be assigned to the indices of the corresponding block.

2.2 Structure of the Cyclic Data Telegrams in Data Exchange Mode

In Data Exchange mode, the DP master class 1 cyclically transmits data from and to all slaves that are connected to the bus.

In this document, data transmitted from the slave to the master are called "input data" and data transmitted from the master to the slave are called "output data".

The input and output data of the FRG-730 has two logic parts:

- 1) the parameter channel
- 2) the process data channel

There is a number of standard telegrams, consisting of:

- a) the parameter channel only
- b) the process data channel only
- c) both, the parameter and process data channel

The parameter channels allows masters without Profibus DPV1 to access device specific parameters that are not part of the normal cyclic data telegram. For masters with Profibus DPV1, no parameter channel is required.

Input data

The input data (transmitted by the FRG-730) consists of the 8 bytes of the parameter channel (if there is a parameter channel in the standard telegram) and of up to 15 bytes of process data depending on the selected standard telegram.

Byte								Byte				
1	2	3	4	5	6	7	8	9	...	23		
Parameter Channel								Process Data				
PKE		IND	res.	PWE								

Where: PKE = Parameter Signature Value = Reading or writing command and definition of the slot
IND = Sub Index = Index No. of the index to be read (→ "Block Model")
res. = reserved
PWE = Process Value = Value to be read or written

Output data

The output data (transmitted by the master) consist of 8 bytes of the parameter channel and up to 10 bytes of process data (control bytes).

Byte								Byte		
1	2	3	4	5	6	7	8	9	...	18
Parameter Channel								Process Data		
PKE	IND	res.	PWE							

2.2.1 Parameter Channel

The structure of the parameter channel is described in the table below.

The parameter channel (called PKW Interface hereinafter) consists of 8 bytes.

Byte							
1	2	3	4	5	6	7	8
PKE	IND	res.	PWE				

The PKW Interface allows reading and writing of slave parameters with a maximum data length of 4 bytes. Strings cannot be read.

The slave generates exactly one response per instruction transmitted by the master. The instruction and response cannot be blocked. This means that exactly one instruction per output telegram can be transmitted to the slave and that exactly one response per input telegram can be transmitted to the master. 4 bytes of actual data can thus be transmitted at a time.

2.2.1.1 PKE Parameter Signature Value

The instruction and response are represented in the first two bytes (PKE) of the parameter channel:

Bit position															
15	14	13	12	11	10	9	8	7	6	5	4	3	2	1	0
AK				res.				Slot							

Where:	Bits	Meaning
	15 ... 12	AK = Instruction/response signature
	11 ... 8	Reserved
	7 ... 0	Define the slot from which data are read or onto which a value is to be written

Instruction signature

In Master ⇒ Slave communication, the AK field contains the instruction signature of the master.

In Slave ⇒ Master communication, the AK field contains the response signature of the slave.

AK	Function Master ⇒ Slave (Instruction signature)	AK normal	Function Slave ⇒ Master (Response signature)	AK error
0	No instruction	0	No response	
1	Read parameter value	1 2 11	Transmit parameter value (word) Transmit parameter value (double word) Transmit parameter value (byte)	7 ¹⁾
2	Write parameter value (data type: word)	1	Transmit parameter value (word)	7 ¹⁾
3	Write parameter value (data type: double word)	2	Transmit parameter value (double word)	7 ¹⁾
10	Write parameter value (data type: byte)	11	Transmit parameter value (byte)	7 ¹⁾

¹⁾ Instruction cannot be executed (error code)

On the left of the table, the instruction signatures of the master are listed according to their function. On the right of the table, the corresponding normal responses (AK Normal) and error codes (AK Error) transmitted by the slave are listed.

Instruction – response sequence

- 1) The master transmits an instruction to the slave and repeats that instruction until it receives a response from the slave.
- 2) The slave keeps transmitting the response to the instruction until the master transmits a new instruction.
- 3) The master marks the end of the first instruction cycle by setting AK to zero. Only after that, a new instruction/response cycle may be started.

2.2.1.2 PWE Parameter (Process Value)

The PWE represents the data element to be transmitted.

If a byte is to be transmitted, that byte has to be in position 8 of the parameter channel.

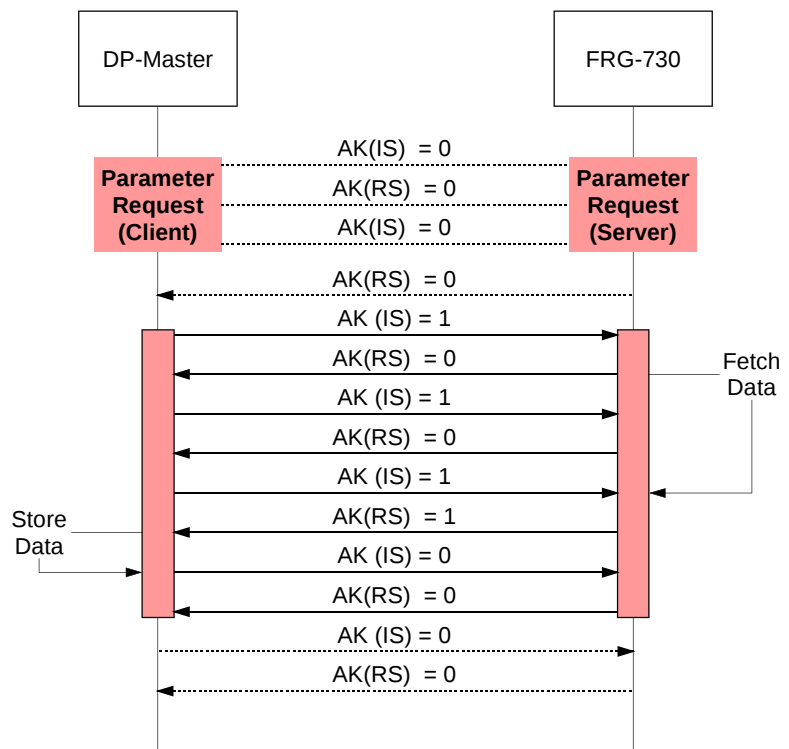
Integers are transmitted with bytes 7 and 8. Double integer and float values are transmitted with bytes 5 ... 8.

2.2.1.3 Error Code (Error Message)

In the event of a transmission error (AK response signature = 7), the slave transmits an error code in byte positions 7 and 8 (data type: INT16).

Error code	Meaning
0	Undefined slot
1	Parameter not changeable
2	Lower or upper value range limit overflow
3	Index error
5	Data type error
17	Instruction not allowed in this state
18	Other errors
201	Already in requested state
202	Object state conflict

The following diagram shows an example of a data request from a master to a FRG-730 via parameter channel.



2.3 Cyclic Message Telegrams

The message telegrams listed below are implemented in the gauge. They can be selected according to requirements. When selecting a message telegram, ascertain what output format of the measured value (integer/float) is required and whether a parameter channel is needed or not. The gauge can also be operated in such a way that the master does not transmit any output data to the slave.

Standard telegram	Master ↔ Slave	Byte	Meaning
1	M ⇒ S	0 1	Transition Command Transition Command Value for Hot Ion Gauges
2	M ⇒ S	0 ... 7 8 9	Parameter Channel Transition Command Transition Command Value for Hot Ion Gauges
3	M ⇒ S	0 ... 7	Parameter Channel
4	S ⇒ M	0 1 2 3 ... 4	Exception status One Of N status extension One Of N PV selector Process value UINT16
5	S ⇒ M	0 1 2 3 ... 6	Exception status One Of N status extension One Of N PV selector Process value float
6	S ⇒ M	0 ... 7 8 9 10 11 ... 12	Parameter channel Exception status One Of N status extension One Of N PV selector Process value UINT16
7	S ⇒ M	0 ... 7 8 9 10 11 ... 14	Parameter channel Exception status One Of N status extension One Of N PV selector Process value float

Configuration data

Depending on the standard telegrams used, the respective configuration data have to be transmitted to the gauge during the configuration phase (→ table on 7).

3 Block Model

Data to the FRG-730 can be transmitted by means of a number of communication protocols and corresponding masters. Profibus defines a master class 1 as normal control unit of the slave (typically a PLC) and a master class 2 as configuration and service unit. The following communication protocols are defined according to the Profibus DPV1 standard.

- MS0 Cyclic data traffic between master class 1 and slave
- MS1 Acyclic data traffic between master class 1 and slave
- MS2 Acyclic data traffic between master class 2 and slave

In the FRG-730, all functions that are made available by the gauge via Profibus are organized in blocks. Access to the individual parameters of the blocks is possible via acyclic services or, for byte, integer and float values, also in cyclic data traffic via the parameter channel.

Block types

The following block types are defined in the gauge.

Device Block	The Device Block contains all data that are required for describing the device and handling its state (status of Device State Machine).
Transducer Block	The physical, process specific functions or interfaces between the FRG-730 and the process such as current and voltage values are represented in transducer blocks. The following transducer blocks are implemented: • One of N Vacuum Gauge Transducer Block • Heat Transfer Vacuum Gauge Transducer Block (Pirani) • Hot Cathode Ion Gauge Transducer Block (BA)
Function Block	Application specific values such as pressure values that result from or can be calculated from the values of the transducer block are represented in the function blocks. • One of N Analog Input Function Block • Analog Input Function Block, Instance 1, Instance 2, Instance 3, Instance 4.

3.1 Device Block

The following table lists the services and parameters integrated in the Device Block (→ Appendix A for abbreviations).

ID	Name	Structure	Data type	Bytes	Access	Value	Store
15	Device Block State	Simple	UINT8	1	1/2_R/W		
16	Block Type	Simple	Octet string	4	1/2_R	1	N
17	Device Type	Simple	Visible string	8	1/2_R	CG	N
18	Standard Revision Level	Simple	Visible string	9	1/2_R	E54-0997	N
19	Device Manufacturer Identifier	Simple	Visible string	20	1/2_R	Agilent	N
20	Manufacturer Model Number	Simple	Visible string	20	1/2_R	e.g. FRG730KF25SP	N
21	Software or Firmware Revision Level	Simple	Visible string	8	1/2_R	e.g. 020_1.01	N
22	Hardware Revision Level	Simple	Visible string	8	1/2_R	e.g. 1.0	N
23	Serial Number	Simple	Visible string	30	1/2_R	e.g. 100	N
25	Device State	Simple	UINT8	1	1/2_R		V
26	Exception Status	Simple	UINT8	1	0_XI 1/2_R		V
27	Exception Detail Alarm	Record	→ below	-	1/2_R		V
28	Exception Detail Warning	Record	→ below	-	1/2_R		V
204	Common Exception Detail Alarm 0	Simple	UINT8	1	1/2_R		V
205	Device Exception Detail Alarm 0 ... 3	Struct	Array of 4 bytes	4	1/2_R		V
207	Manufacturer Exception Detail Alarm 0	Simple	UINT8	1	1/2_R		V
208	Common Exception Detail Warning 0	Simple	UINT8	1	1/2_R		V
209	Device Exception Detail Warning 0 ... 3	Struct	Array of 4 bytes	4	1/2_R		V
210	Device Exception Detail Warning 4	Simple	UINT8	1	1/2_R		V
211	Manufacturer Exception Detail Warning 0	Simple	UINT8	1	1/2_R		V

3.1.1 Information on the Individual Indices

3.1.1.1 Block Type ID 16

The Block Type Parameter contains an ID which describes the block type. The block type ID of the Device Block 1. The other defined block types are listed in Appendix B.

3.1.1.2 Device Type ID 17

The Device Type identifies the device type which is connected to the field bus via Profibus.
The Device Type of the BFRG-730 gauge is "CG", the abbreviation of Combination Gauge.

3.1.1.3 Standard Revision Level ID 18

This parameter describes the version of the "Sensor/Actuator Network Specific Device Model" published by the SEMI® (Semiconductor Equipment and Materials International, California), according to which the profile of this device has been developed.
The fixed setting of this parameter is "E54-0997".

3.1.1.4 Device Manufacturer Identifier ID 19

This parameter describes the manufacturer of the device, "Agilent".

3.1.1.5 Manufacturer Model Number ID 20

This parameter represents the part number of the gauge (→ section "Validity").
FRG-730 is available with different vacuum connection types:

Gauge	Vacuum connection	Part number
FRG-730	25 KF	FRG730KF25SP
	40 CF	FRG730CF35SP

3.1.1.6 Software or Firmware Revision Level ID 21

This parameter indicates the software version of the Profibus option in the following format: xxx.y.yy (where xxx is the version of the FRG-730 firmware and y.yy is the version of the Profibus board).

3.1.1.7 Hardware Revision Level ID 22

This parameter indicates the hardware version of the gauge in the format "1.0".

3.1.1.8 Device State ID 25

This parameter indicates the status of the gauge. Due to the structure of the Device State Machine, the following statuses are possible:

Parameter value	Status
0	Undefined
1	Self testing
2	Idle
3	Self test exception
4	Executing
5	Abort
6	Critical fault

The device statuses are described in detail in section "Device Block, Device Behavior".

3.1.1.9 Exception Status ID 26

The Exception Status describes the alarm and warning states of the gauge in an "Expanded error output format".

A difference is made between warnings and errors.

Alarms and errors are divided into three groups (→ sections "Exception Detail Alarm" and "Exception Detail Warning" for details):

- ALARM / Warning Device Common For errors that occur independently of the type of device used, e.g. supply error, RAM, ROM, or EEPROM error.
- ALARM / Warning Device Specific For device specific errors (e.g. filament rupture (Pirani) or cathode rupture (BA)).
- ALARM / Warning Manufacturer Specific For errors defined by the manufacturer that are not mentioned in the standard.

In each of the above groups, there are several error or warning conditions. The individual fields are presented in the "Exception Detail Alarm" and "Exception Detail Warning". If an error message occurs in "Exception Detail Alarm" or "Exception Detail Warning", the corresponding bit is set in the Exception Status. Therefore, if bits 0 ... 6 of the Exception Status are on "0" there is no warning message pending.

If a bit is set, the actual error can be read in the corresponding group.

The Exception Status is output in cyclic data and informs on the current error status using only one byte. If an error occurs, the current error status can be read via acyclic services or in cyclic data exchange via the parameter channel. This ensures that while the current error status is always available in the cyclic data, no unnecessary data overhead is transmitted.

Bit	Function	Meaning
0	ALARM, device common	The bit is set if an error of the Alarm Device Common group is detected.
1	ALARM, device specific	The bit is set if an error of the Alarm Device Specific group is detected.
2	ALARM, manufacturer specific	The bit is set if an error of the Alarm Manufacturer Specific group is detected.
3	-	-
4	WARNING, device common	The bit is set if an error of the Warning Device Common group is detected.
5	WARNING, device specific	The bit is set if an error of the Warning Device Specific group is detected.
6	WARNING, manufacturer specific	The bit is set if an error of the Warning Manufacturer Common group is detected.
7	Expanded Format	Is constantly on "1" and marks the use of the expanded error output format.

3.1.1.10 Exception Detail Alarm ID 27

If, in the Exception Status, one of the bits 0 ... 2 is set, the current error can be read in the "Exception Detail Alarm" parameter. The "Exception Detail Alarm" parameter consists of a total of 10 bytes that inform on the error status of the gauge. Due to the use of the expanded error output format, these bytes have the following structure:

Byte No	Name	Description	Value
0	Common Exception Detail Size	Indicates the number of subsequent bytes that contain the Common Exception Detail Alarm.	2
1	Common Exception Detail 0	Contains current error messages from the Common Exception Detail Alarm group.	Depending on error status
2	Common Exception Detail 1	not supported	0
3	Device Exception Detail Size	Indicates the number of subsequent bytes that contain the Device Exception Detail Alarm.	4
4	Device Exception Detail 0 (Pirani error data)	This error information from the Device Exception Detail Alarm group refers to Pirani.	Depending on error status
5	Device Exception Detail 1 (Pirani error data)	This error information from the Device Exception Detail Alarm group refers to Pirani.	Depending on error status
6	Device Exception Detail 2 (BA error data)	This error information from the Device Exception Detail Alarm group refers to BA.	Depending on error status
7	Device Exception Detail 3 (BA error data)	This error information from the Device Exception Detail Alarm group refers to BA.	Depending on error status
8	Manufacturer Exception Detail Size	Indicates the number of subsequent bytes that contain the Manufacturer Exception Detail Alarm.	1
9	Manufacturer Exception Detail 0	Contains current error messages from the Manufacturer Exception Detail Alarm group.	Depending on error status

Common Exception Detail Alarm

Bit	Common Exception Detail 0
0	0
1	0
2	EPROM exception
3	EEPROM exception
4	RAM exception
5	0
6	0
7	0

Bit	Common Exception Detail 1
0	0
1	0
2	0
3	0
4	0
5	0
6	0
7	0

Device Exception
Detail Alarm

Bit	Device Exception Detail 0 Referring to Pirani
0	0
1	0
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Alarm byte 0 of Pirani Transducer Block.	

Bit	Device Exception Detail 1 Referring to Pirani
0	0
1	Electronics / sensor error
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Alarm byte 1 of Pirani Transducer Block.	

Bit	Device Exception Detail 2 Referring to BA
0	Sensor filament 1 Alarm
1	Sensor filament 2 Alarm
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Alarm byte 0 of BA Transducer Block.	

Bit	Device Exception Detail 3 Referring to BA
0	0
1	Electronics / sensor error
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Alarm byte 1 of BA Transducer Block.	

Manufacturer Exception
Detail Alarm 0

Manufacturer Exception Detail 0 is set to "1", if there is an internal communication error in the gauge.

3.1.1.11 Exception Detail Warning ID 28

If, in the Exception Status, one of bits 4 ... 6 is set, the current warning can be read in the parameter "Exception Detail Warning". The Exception Detail Warning parameter consists of a total of 11 bytes that inform on the error status of the gauge.

Due to the use of the expanded error output format, these bytes have the following structure:

Byte No	Name	Description	Value
0	Common Exception Detail Size	Indicates the number of subsequent bytes that contain the Common Exception Detail Warning.	2
1	Common Exception Detail 0	Contains current error messages from the Common Exception Detail Warning group.	Depending on warning status
2	Common Exception Detail 1	Contains current error messages from the Common Exception Detail Warning group.	Depending on warning status
3	Device Exception Detail Size	Indicates the number of subsequent bytes that contain the Device Exception Detail Warning.	5
4	Device Exception Detail 0	One of N Status Extension.	Depending on warning status
5	Device Exception Detail 1 (Pirani)	This error information from the Device Exception Detail Warning group refers to Pirani.	Depending on warning status
6	Device Exception Detail 2 (Pirani)	This error information from the Device Exception Detail Warning group refers to Pirani.	Depending on warning status
7	Device Exception Detail 3 (BA)	This error information from the Device Exception Detail Warning group refers to BA.	Depending on warning status
8	Device Exception Detail 4 (BA)	This error information from the Device Exception Detail Warning group refers to BA.	Depending on warning status
9	Manufacturer Exception Detail Size	Indicates the number of subsequent bytes that contain the Manufacturer Exception Detail Warning.	1
10	Manufacturer Exception Detail	Contains current error messages from the Manufacturer Exception Detail Warning group.	Depending on warning status

Common Exception
Detail Warning

Bit	Common Exception Detail 0
0	0
1	0
2	EPROM exception
3	EEPROM exception
4	RAM exception
5	0
6	0
7	0

Bit	Common Exception Detail 1
0	0
1	0
2	0
3	0
4	0
5	0
6	0
7	0

The warning bits are set in the same way as the error bits because here, warnings have the same meaning as errors.

Device Exception
Detail Warning

Bit	Device Exception Detail 0
0	Bit set if Reading Invalid
1	Bit set if Device Overrange
2	Bit set if Device Underrange
3	0
4	0
5	0
6	0
7	0
This byte is a copy of the One Of N Vacuum Gauge Transducer Block.	

Bit	Device Exception Detail 1 Referring to Pirani
0	0
1	0
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Warning byte 0 of the Pirani Transducer Block.	

Bit	Device Exception Detail 3 Referring to BA
0	Sensor filament 1 Warning
1	Sensor filament 2 Warning
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Warning byte 0 of the BA Transducer Block.	

Bit	Device Exception Detail 2 Referring to Pirani
0	0
1	Electronics / sensor warning
2	0
3	0
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Warning byte 1 of the Pirani Transducer Block.	

Bit	Device Exception Detail 4 Referring to BA
0	0
1	Electronics / sensor warning
2	0
3	Pressure too high for degas
4	0
5	0
6	0
7	0
This byte is a copy of Sensor Warning byte 0 of the BA Transducer Block.	

Manufacturer Exception
Detail Warning 0

Manufacturer Exception Detail 0 is set to "1", if there is an internal communication error in the gauge.



The following parameters are copies of the ID 27 and ID 28. They are used only if you want to access these parameters by the parameter channel.

**3.1.1.12 Copy Common Exception
Detail Alarm 0
ID 204**

This parameter corresponds to the Common Exception Detail Alarm of the Exception Detail Alarm (ID 27).

**3.1.1.13 Copy Device Exception
Detail Alarm 0 ... 3
ID 205**

This parameter corresponds to the Device Exception Detail Alarm Byte 0 ... 3 of the Exception Detail Alarm (ID 27).

**3.1.1.14 Copy Manufacturer
Exception Detail Alarm 0
ID 207**

This parameter corresponds to the Manufacturer Exception Detail Alarm Byte 0 of the Exception Detail Alarm (ID 27).

**3.1.1.15 Copy Common Exception
Detail Warning 0
ID 208**

This parameter corresponds to the Common Exception Detail Warning Byte 0 of the Exception Detail Warning (ID 28).

**3.1.1.16 Copy Device Exception
Detail Warning 0 ... 3
ID 209**

This parameter corresponds to the Device Exception Detail Warning Byte 0 ... 3 of the Exception Detail Warning (ID 28).

**3.1.1.17 Copy Device Exception
Detail Warning 4
ID 210**

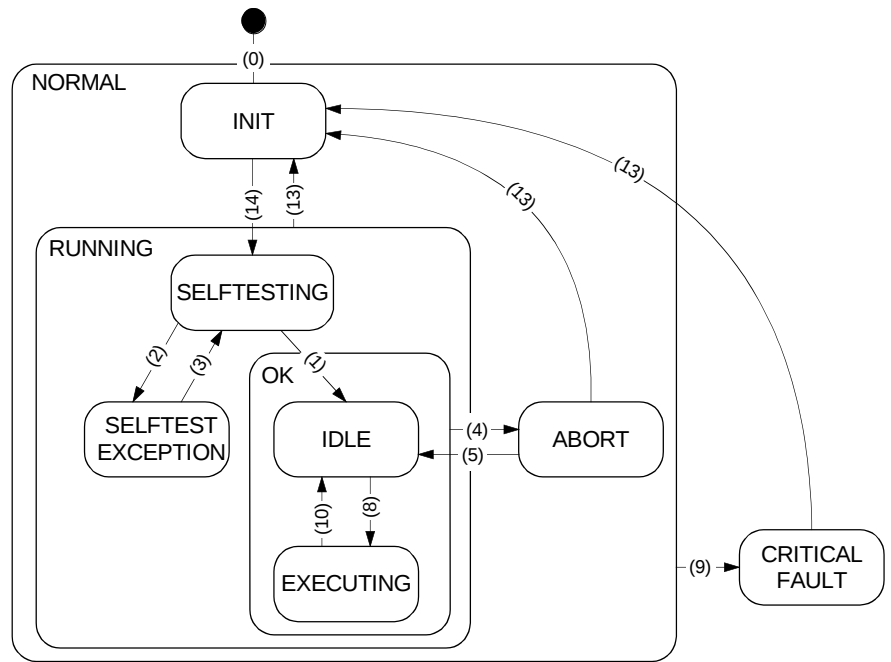
This parameter corresponds to the Device Exception Detail Warning Byte 4 of the Exception Detail Warning (ID 28).

**3.1.1.18 Copy Manufacturer
Exception Detail
Warning 0
ID 211**

This parameter corresponds to the Manufacturer Exception Detail Warning Byte 0 of the Exception Detail Warning (ID 28).

3.1.2 Device Block, Device Behavior

The FRG-730 behaves as described in the status diagram below.



After the start, the gauge independently runs through the INIT and SELFTESTING status and eventually changes to the IDLE status (if there is no error) or to the SELFTEST_EXCEPTION status (if there is a gauge error).

When data traffic is taken up, a difference has to be made between cyclic and acyclic data traffic.

Cyclic data traffic

As soon as cyclic data interchange is taken up, the gauge automatically changes to the EXECUTING status.

Acyclic data traffic

In acyclic data traffic, a START service has to be transmitted to bring the gauge to the EXECUTING status.

Device statuses

Status name	Description
NORMAL	The communication interface can respond to commands. All defined blocks of the gauge exist.
RUNNING	All block instances are initialized and the individual parameters have their initial or default values.
SELFTESTING	In this status, the gauge performs a selftest.
IDLE	All blocks defined in the gauge are initialized and the complete hardware has been tested and found free of errors. The value defined in the Safe State is output as measured value.
SELFTEST EXCEPTION	During the self test, an error has been detected. The gauge has changed to the SELFTEST EXCEPTION status. The cause of the error can be found in the attributes Device Exception Detail Alarm/Warning of the Device Block. The value defined in the Safe State is output as measured value.
EXECUTING	The gauge functions normally and outputs the measured value.
ABORT	The Device Block is in the ABORT status. The gauge specific measurement functions are not carried out.
CRITICAL FAULT	A device error has occurred. The gauge specific measurement functions are not carried out. The value defined in the Safe State is output as measured value.

3.1.2.1 Device Block State Command

There are a number of special commands for bringing the gauge into a status it does not automatically go to.

ID	Name	Structure	Data type	Bytes	Access	Store	Range
15	Device Block State Command	Simple	UINT8	1	1/2_R/W	N	→ below

ID value	Name	Description
0	Inactive	No action.
1	Reset	Used for reinitializing the device.
2	Abort	Brings the device to the ABORT status.
3	Recover	Used for bringing the device from the ABORT status into the Recovered State = IDLE.
4	Execute	Brings the unit to the EXECUTING status, in which the gauge functions normally. As soon as cyclic data traffic is initialized, this status command is executed automatically.
5	Stop	Brings the gauge to the IDLE status.
6	Perform Diagnostic	Stops the running activity and starts SELFTEST.

3.2 Analog Input Block

All gauge functions of the FRG-730 are described in the Analog Input Block.

Since the gauge includes two measuring systems, there are also two Analog Input Block Instances representing the Pirani and the ionization measuring part (BA) respectively.

3.2.1 One Of N Analog Input Function Block / SLOT 1

ID	Name ¹⁾	Structure	Data type	Bytes	Access	Store
15	AI Block Adjust Command	Simple	UINT8	1	1/2_W	-
16	Block Type	Simple	Octet string	4	2_R	N
46	Channel Instance Selector	Simple	UINT8	1	1/2_R/W	N
47	PV Selector	Simple	UINT8	1	0_XI 1/2_R	V

¹⁾ The above parameters are described below.

3.2.1.1 AI Block Adjust Command ID 15

To start "Full Scale Adjust" of the gauge, the value "1" has to be entered.

3.2.1.2 Block Type ID 16

The value of the Block Type ID is "3" (→ table in Appendix A).

3.2.1.3 Channel Instance Selector ID 46

The gauge has two instances of the Analog Input Function Block and two instances of the Vacuum Gauge Transducer Block, or one instance of each block per measuring system (Pirani and BA). There are two additional instances for describing the thresholds of the switching functions (→ section "Analog Sensor Input Function Block", Instances 3 and 4).

However, there is only one address range for both, querying and setting the corresponding parameters. The Parameter Channel Instance Selector is used for defining the parameters that are written into the address range.

If the Parameter Channel Instance Selector is set to "1", the attributes of Instance 1 are mapped into that address range and can be written or read by addressing Slot ⇒ Instance ⇒ Parameter-ID.

3.2.1.3 PV Selector ID 47

The PV Selector is determined by the gauge and defines the Analog Input Function Block Instance from which the measured value is copied into the cyclic output data telegrams. Therefore, the value output in the cyclic data is always the measured value of the active instance, i.e. the one that is currently measuring. While the gauge is measuring in the BA range, the measured value of the ionization vacuum meter is output in the cyclic data and while the gauge is measuring in the Pirani range, the measured value of the Pirani is output.



The measured values of Instances 3 and 4 (thresholds of the switching functions) are not output in the cyclic data.

The pressure ranges, in which measurement is performed either by the Pirani or by the BA measuring part, and are thus called "active", are indicated below:

Measuring range

Pressure [mbar]	PV Selector
$5.5 \times 10^{-3} < p \leq 1000$	Pirani
$5 \times 10^{-10} < p < 2 \times 10^{-2}$	BA

In the pressure range $5.5 \times 10^{-3} \dots 2 \times 10^{-2}$ mbar an averaging of the two corresponding measurement principles is done. The active instance in this pressure range is Pirani.

3.2.2 Analog Sensor Input Function Block SLOT 1 / Instance 1

Instance 1 of the Analog Sensor Input Function Blocks describes the functionality of the Pirani measuring part of the gauge.

The following attributes are supported:

ID	Name	Structure	Data type	Bytes	Access	Store
15	AI Block Adjust Command	Record		-	2_W	
19	Process Value (PV)	Simple	According to Data Type value (Parameter 21)	-	0_XI 1/2_R	V
20	Status	Simple	UINT8	1	0_XI 1/2_R	V
21	Data Type	Simple	UINT8	1	2_R/W	N
22	Data Units	Simple	UINT16	2	2_R/W	N
23	Reading Valid	Simple	Boolean	1	1/2_R	V
24	Full Scale	Simple	According to Data Type value (Parameter 21)	-	1/2_R	N
39	Safe State	Simple	Unsigned8	1	1/2_R/W	N
40	Safe Value	Simple	According to Data Type value (Parameter 21)	-	1/2_R/W	N
44	Overrange	Simple	According to Data Type value (Parameter 21)	-	1/2_R	N
45	Underrange	Simple	According to Data Type value (Parameter 21)	-	1/2_R	N

3.2.2.1 Process Value ID 19

The Process Value contains the measured value of the Pirani Device Instance in the currently selected data unit (ID 22) and in the selected data type (ID 21).

If the device is not in the EXECUTING State (ID 25, Device Block), the value defined in the Safe State is output.

Values output in the data unit COUNTS can be converted into a pressure value by means of the following formulas:

Calculation of the pressure (PV is the abbreviation used for Process Value):

$$PV_{\text{mbar}} = 10^{(\text{COUNTS} / 2000) - 12.5}$$

$$PV_{\text{Torr}} = 0.75006168 \times PV_{\text{mbar}}$$

$$PV_{\text{Micron}} = 10^{-3} \times PV_{\text{Torr}}$$

$$PV_{\text{Pa}} = 100 \times PV_{\text{mbar}}$$

3.2.2.2 Status ID 20

This parameter remains on "0".

3.2.2.3 Data Type ID 21

Two data types are supported: Float and Integer16.

In cyclic data exchange, the data type cannot be modified. After the gauge is switched on, the value stored in the EEPROM is loaded. The data type setting can only be modified when the gauge is in the IDLE status. By defining the configuration data for cyclic data exchange (selection of standard telegrams), the data type used in the selected standard telegram is taken over. All settings previously made in acyclic data traffic are thus overwritten.

If the data type is set in one instance, that data type setting applies to all instances. Likewise, when a standard telegram is selected, the data type used by that standard telegram will be valid for all instances:

Coding	Data type
3	Integer16
8	Float (default)

3.2.2.4 Data Unit ID 22

The gauge supports the following pressure units:

Coding (dec)	Data type
1001	COUNTS
1301	Torr
1302	mTorr (Micron)
1308	mbar
1309	Pascal



For safety reasons, it is not possible to change the pressure unit while the gauge is cyclically interchanging data with a DP/V0 master.

The data unit setting can only be modified when the gauge is in the IDLE status.

In cyclic data traffic, the data unit must be set in the User Parameter Data. All settings previously made in acyclic data traffic are overwritten (→ section "User Parameter Data").

If the data unit is set in one instance, that data unit setting applies to all instances. Likewise, the data unit setting made in the User Parameter Data is valid for all instances.

3.2.2.5 Reading Valid ID 23

This parameter indicates that the pressure reading is within a valid range. It is set to "1" if:

- The gauge is in the EXECUTING status.
- There is no error
- The measured value is lower than the overrange value and higher than the underrange value.



If this value is set to zero, the pressure reading is not valid. In such a case, either check Exception Status (ID 26, Device Block) to find out whether there is an error or check One Of N Status Extension (ID 120, One Of N Vacuum Gauge Transducer Block) to find out whether the measured value is out of the specified measuring range (overrange or underrange mode).

3.2.2.6 Full Scale ID 24

This parameter contains the valid maximum value of the pressure reading in the currently selected data unit and data type.

Pressure unit	Full scale
COUNTS	31000
Torr	750.06168
mTorr (Micron)	750061.68
mbar	1000
Pascal	100000

3.2.2.7 Safe State ID 39

When the gauge is not in the EXECUTING status (ID 25, Device Block) or if there is a device error, a value defined by Safe State is output as pressure value:

Safe State	Coding	PV behavior
Zero	0	The Process Value (measured value ID 19) is set to 0.
Full Scale	1	The Process Value (measured value ID 19) is set to the full scale value (ID 24).
Hold Last Value	2	The Process Value is set to the last valid value obtained in the EXECUTING status.
Use Safe Value	3	The Process Value (measured value ID 19) is set to the Safe Value (ID 40).

3.2.2.8 Safe Value ID 40

The Safe Value is the value output with the Process Value parameter (ID 19) when an error occurs or the gauge goes to the NOT EXECUTING status. If this value is set to zero, it will remain on zero when the data unit is changed.

3.2.2.9 Overrange ID 44

Overrange is the highest valid measured value at which Reading Valid is still on "1".

Pressure unit	Overrange
COUNTS	31000
Torr	750.06168
mTorr (Micron)	750061.68
mbar	1000
Pascal	100000

3.2.2.10 Underrange ID 45

Underrange is the lowest valid measured value at which Reading Valid is still on "1".

Pressure unit	Underrange
COUNTS	20480
Torr	4.12534×10^{-3}
mTorr (Micron)	4.12534
mbar	5.5×10^{-3}
Pascal	0.55

3.2.2.11 AI Block Adjust Command (Pirani) ID 15

The parameter controls zero point and fullscale adjustments of the Pirani.

Byte	Name	Structure	Data type	Bytes	Access
0	State Command	Simple	UINT8	1	1/2_W

State Command	Name	Meaning
0	Zero Adjust	Pirani zero point adjustment
1	Gain Adjust	Pirani fullscale adjustment

3.2.3 Analog Sensor Input Function Block SLOT 1 / Instance 2

Instance 2 of the Analog Sensor Input Function Block describes the functionality of the BA measuring part of the gauge.

In Instance 2, the same attributes as in Instance 1 are supported:

ID	Name	Structure	Data type	Bytes	Access	Store
19	Process Value (PV)	Simple	According to Data Type value (parameter 21)	-	0_XI 1/2_R	V
20	Status	Simple	UINT8	1	0_XI 1/2_R	V
21	Data Type	Simple	UINT8	1	2_R/W	N
22	Data Units	Simple	UINT16	2	2_R/W	N
23	Reading Valid	Simple	Boolean	1	1/2_R	V
24	Full Scale	Simple	According to Data Type value (parameter 21)	-	1/2_R	N
39	Safe State	Simple	UINT8	1	1/2_R/W	N
40	Safe Value	Simple	According to Data Type value (parameter 21)	-	1/2_R/W	N
44	Overrange	Simple	According to Data Type value (parameter 21)	-	1/2_R	N
45	Underrange	Simple	According to Data Type value (parameter 21)	-	1/2_R	N

3.2.3.1 Process Value ID 19

The Process Value contains the measured value of the BA Device Instance in the currently selected data unit (ID 22) and in the selected data type (ID 21).

If the device is not in the EXECUTING status (ID 25, Device Block), the value defined in the Safe State is output.

Values output in the data unit COUNTS can be converted into a pressure value by means of the following formulas:

Calculation of the pressure (PV is the abbreviation used for Process Value):

$$PV_{\text{mbar}} = 10^{(\text{COUNTS} / 2000) - 12.5}$$

$$PV_{\text{Torr}} = 0.75006168 \times PV_{\text{mbar}}$$

$$PV_{\text{Micron}} = 10^{-3} \times PV_{\text{Torr}}$$

$$PV_{\text{Pa}} = 100 \times PV_{\text{mbar}}$$

3.2.3.2 Status ID 20

This parameter is remains on "0".

3.2.3.3 Data Type ID 21

Two data types are supported: Float and Integer16.

In cyclic data traffic, the data type cannot be modified. After the gauge is switched on, the value stored in the EEPROM is loaded. The data type setting can only be modified when the gauge is in the IDLE status. By defining the configuration data for cyclic data traffic (selection of standard telegram), the data type used in the selected standard telegram is taken over. All settings previously made in acyclic data traffic are thus overwritten.

If the data type is set in one instance, that data type setting applies to all instances. Likewise, when a standard telegram is selected, the data type used by that standard telegram will be valid for all instances.

Coding	Data type
3	Integer16
8	Float (default)

3.2.3.4 Data Unit ID 22

The gauge supports the following pressure units:

Coding (dec)	Pressure unit
1001	COUNTS
1301	Torr
1302	mTorr (Micron)
1308	mbar
1309	Pascal



For safety reasons, it is not possible to change the pressure unit while the gauge is cyclically interchanging data with a DP/V0 master.

The data unit setting can only be modified when the gauge is in the IDLE status.

In cyclic data traffic, the data unit must be set in the User Parameter Data. All settings previously made in acyclic data traffic are overwritten (→ section "User Parameter Data").

If the data unit is set in one instance, that data unit setting applies to all instances. Likewise, the data unit setting made in the User Parameter Data is valid for all instances.

3.2.3.5 Reading Valid ID 23

This parameter indicates that the pressure reading is within a valid range. It is set to "1" if:

- The gauge is in the EXECUTING status.
- There is no error
- The measured value is lower than the overrange value and higher than the underrange value.



If this value is set to zero, the pressure reading is not valid. In such a case, either check Exception Status (ID 26, Device Block) to find out whether there is an error or check One Of N Status Extension (ID 120, One Of N Vacuum Gauge Transducer Block) to find out whether the measured value is out of the specified measuring range (overrange or underrange mode).

3.2.3.6 Full Scale ID 24

This parameter contains the valid maximum value of the pressure reading in the currently selected data unit and data type.

Pressure unit	Full scale
COUNTS	21602
Torr	1.5×10^{-2}
mTorr (Micron)	15
mbar	2×10^{-2}
Pascal	2

3.2.3.7 Safe State ID 39

When the gauge is not in the EXECUTING state (ID 25, Device Block) or if there is a device error, a value defined by Safe State is output as pressure value:

Safe State	Coding	PV behavior
Zero	0	The Process Value (measured value ID 19) is set to 0.
Full Scale	1	The Process Value (measured value ID 19) is set to the full scale value (ID 24).
Hold Last Value	2	The Process Value is set to the last valid value obtained in the EXECUTING status.
Use Safe Value	3	The Process Value (measured value ID 19) is set to the Safe Value (ID 40).

3.2.3.8 Safe Value ID 40

The Safe Value is the value output with the Process Value Parameter (ID 19) when an error occurs or the gauge goes to the NOT EXECUTING status. If this value is set to zero, it will remain on zero when the data unit is changed.

3.2.3.9 Overrange ID 44

Overrange is the highest valid measured value at which Reading Valid is still on "1".

Pressure unit	Overrange
COUNTS	21602
Torr	1.5×10^{-2}
mTorr (Micron)	15
mbar	2×10^{-2}
Pascal	2

3.2.3.10 Underrange ID 45

Underrange is the lowest valid measured value at which Reading Valid is still on "1".

Pressure unit	Underrange
COUNTS	6398
Torr	3.7503×10^{-10}
mTorr (Micron)	3.7503×10^{-7}
mbar	5.0×10^{-10}
Pascal	5.0×10^{-8}

3.2.4 Analog Sensor Input Function Block SLOT 1 / Instance 3, 4

Instance 3 and 4 of the Analog Sensor Input Function Block describe the functionality of the threshold potentiometers used to set the gauges switching functions ("Setpoint A and B, → [2]).

Instance	Switching Function ("Setpoint")
3	A
4	B

ID	Name	Structure	Data type	Bytes	Access	Store
19	Process Value (PV)	Simple	According to Parameter Data Type	-	1/2_R	V
20	Status	Simple	UINT8	1	1/2_R	V
21	Data Type	Simple	UINT8	1	2_R/W	N
22	Data Units	Simple	UINT16	2	2_R/W	N
23	Reading Valid	Simple	Boolean	1	1/2_R	V

3.2.4.1 Process Value ID 19

The Process Value contains the current setting of the threshold potentiometers for "Setpoint A" and "Setpoint B" in the currently selected data unit (ID 22) and data type (ID 21).

If the pressure drops below the set threshold, the relay is closed. If after that, the pressure rises above that threshold with a hysteresis of 10%, the relay is opened again (→ [2]).

Values output in the pressure unit COUNTS can be converted into a pressure value by means of the following formulas:

$$p_{\text{mbar}} = 10^{(\text{COUNTS} / 2000) - 12.5}$$

The threshold voltages of the Setpoint potentiometer can be converted into a pressure value by means of the following formulas:

$$p_{\text{mbar}} = 10^{(U - 7.75) / 0.75}$$

Where: U is the threshold voltage (0 ... +10 V) of the corresponding Setpoint (A, B), measured at the D-Sub connector (sensor cable connector) of the gauge (→ [2]).

3.2.4.3 Status ID 20

This parameter is remains on "0".

3.2.4.4 Data Type ID 21

Two data types are supported: Float and Integer16.

In cyclic data traffic, the data type cannot be modified. After the gauge is switched on, the value stored in the EEPROM is loaded. The data type setting can only be modified when the gauge is in the IDLE status. By defining the configuration data for cyclic data traffic (selection of standard telegram), the data type used in the selected standard telegram is taken over. All settings previously made in acyclic data traffic are thus overwritten.

If the data type is set in one instance, that data type setting applies to all instances. Likewise, when a standard telegram is selected, the data type used by that standard telegram will be valid for all instances.

Coding	Data type
3	Integer16
8	Float (default)

3.2.4.5 Data Unit ID 22

The gauge supports the following pressure units:

Coding (dec)	Pressure unit
1001	COUNTS
1301	Torr
1302	mTorr (Micron)
1308	mbar
1309	Pascal



For safety reasons, it is not possible to change the pressure unit while the gauge is cyclically interchanging data with a DP/V0 master.

The data unit setting can only be modified when the gauge is in the IDLE status.

In cyclic data traffic, the data unit must be set in the User Parameter Data. All settings previously made in acyclic data traffic are overwritten (→ section "User Parameter Data").

If the data unit is set in one instance, that data unit setting applies to all instances. Likewise, the data unit setting made in the User Parameter Data is valid for all instances.

3.2.4.6 Reading Valid ID 23

This parameter is set to "1".

3.3 Transducer Block

3.3.1 One Of N Vacuum Gauge Transducer Block / SLOT 1

ID	Name	Structure	Data type	Bytes	Access	Store
120	One Of N Status Extension	Simple	UINT8	1	1/2_R	V

3.3.1.1 One Of N Status Extension

This parameter indicates whether the overrange or underrange of the gauge is exceeded.

Bit 7	Bit 6	Bit 5	Bit 4	Bit 3	Bit 2	Bit 1	Bit 0
0	0	0	0	0	Underrange Exceeded	Overrange Exceeded	Reading Invalid

If the gauge is operated in its overrange or underrange ($p < 5 \times 10^{-10}$ mbar or $p > 1000$ mbar), the corresponding bit and additionally the bit "Reading Invalid" is set.

If an error occurs, the bit "Reading Invalid" as well as the corresponding error bits in Device Block (ID 26, 27, 28 Device Block) are set.

3.3.2 Heat Transfer Vacuum Gauge Transducer Block / SLOT 1 / Instance 1

ID	Name	Structure	Data type	Bytes	Access	Store
101	Block Type	Simple	Octet string	4	2_R	N
102	Status Extension	Simple	UINT8	1	1/2_R	V
103	Sensor Alarm	Struct	Array of 2 bytes	2	1/2_R	V
104	Sensor Warning	Struct	Array of 2 bytes	2	1/2_R	V

3.3.2.1 Block Type ID 101

According to the table in Appendix A, the Block Type ID has the value "13".

3.3.2.2 Status Extension ID 102

This parameter indicates whether the overrange or underrange of the Heat Transfer Vacuum Gauge device instance (Pirani) is exceeded.

Bit 7	Bit 6	Bit 5	Bit 4	Bit 3	Bit 2	Bit 1	Bit 0
Reserved					Underrange Exceeded ¹⁾	Overrange Exceeded ²⁾	Reading Invalid ³⁾

¹⁾ Underrange: The parameter is set to "1", if the pressure value is below the value defined in ID 45 (Analog Input Function Block Instance 1).

²⁾ Overrange: The parameter is set to "1", if the pressure value is above the value defined in ID 44 (Analog Input Function Block Instance 1).

³⁾ Reading Invalid: The parameter is set to "1", if "Reading Valid" ID 23 (Analog Input Function Block Instance 1) is set to "0".

If the instance is operated in its overrange or underrange, the corresponding bit and additionally the bit "Reading Invalid" is set.

If an error occurs, the bit "Reading Invalid" as well as the corresponding error bits in Device Block (ID 26, 27, 28 Device Block) are set.

3.3.2.3 Sensor Alarm ID 103

This parameter indicates the detectable errors occurring in connection with the Pirani measuring part.

	Bit 7	Bit 6	Bit 5	Bit 4	Bit 3	Bit 2	Bit 1	Bit 0
Byte 0	0	0	0	0	0	0	0	0
Byte 1	0	0	0	0	0	0	Electronics Failure ¹⁾	0

¹⁾ Electronics Failure includes a sensor error.

The Sensor Alarm bits defined here are copied into the Device Block ID 27 in the "Device Exception Detail Alarm" range of the Pirani measuring part.

3.3.2.4 Sensor Warning ID 104

This parameter indicates the detectable warnings occurring in connection with the Pirani measuring part.

	Bit 7	Bit 6	Bit 5	Bit 4	Bit 3	Bit 2	Bit 1	Bit 0
Byte 0	0	0	0	0	0	0	0	0
Byte 1	0	0	0	0	0	0	Electronics Warning	0

The Sensor Warning bits defined here are copied into the Device Block ID 28 in the "Device Exception Detail Warning" range of the Pirani measuring part.

3.3.3 Hot Cathode Ion Gauge Transducer Block / SLOT 1 / Instance 2

ID	Name	Structure	Data type	Bytes	Access	Store
14	Hot Cathode Block State Command	Rec		1	1/2_W	
101	Block Type	Simple	Octet string	4	2_R	N
102	Status Extension	Simple	UINT8	1	1/2_R	V
103	Sensor Alarm	Struct	Array of 2 bytes	2	1/2_R	V
104	Sensor Warning	Struct	Array of 2 bytes	2	1/2_R	V
105	Emission Status	Simple	Boolean	1	1/2_R/W	V
106	Emission Current	Simple	FLOAT	4	1/2_R/W	N
108	Active Filament	Simple	Boolean	1	1/2_R/W	N
109	Degas Status	Simple	Boolean	1	1/2_R	V
116	Active Degas Filament	Simple	UINT8	1	1/2_R	V
119	Mode Filament Selection (Filament User Mode)	Simple	Boolean	1	2_R/W	N
201	Emission User Mode State	Simple	Boolean	1	1/2_R	V

3.3.3.1 Block Type ID 101

According to the table in Appendix A, the Block Type ID has the value "16".

3.3.3.2 Status Extension ID 102

This parameter indicates that the overrange or underrange of the Hot Cathode Ion Gauge instance is exceeded.

Bit 7	Bit 6	Bit 5	Bit 4	Bit 3	Bit 2	Bit 1	Bit 0
Reserved					Underrange Exceeded ¹⁾	Overrange Exceeded ²⁾	Reading Invalid ³⁾

- ¹⁾ Underrange: The parameter is set to "1", if the pressure value is below the value defined in ID 45 (Analog Input Function Block Instance 2).
- ²⁾ Overrange: The parameter is set to "1", if the pressure value is above the value defined in ID 44 (Analog Input Function Block Instance 2).
- ³⁾ Reading Invalid: The parameter is set to "1", if "Reading Valid" ID 23 (Analog Input Function Block Instance 2) is set to "0".

If the instance is operated in its overrange or underrange, the corresponding bit and additionally the bit "Reading Invalid" is set.

If an error occurs, the bit "Reading Invalid" as well as the corresponding error bits in Device Block (ID 26, 27, 28 Device Block) are set.

3.3.3.3 Sensor Alarm ID 103

This parameter indicates the detectable errors occurring in connection with the BA measuring part.

	Bit7	Bit6	Bit5	Bit4	Bit3	Bit2	Bit 1	Bit 0
Byte 0	0	0	0	0	0	0	Sensor Filament 2 Alarm ¹⁾	Sensor Filament 1 Alarm ¹⁾
Byte 1	0	0	0	0	0	0	Electronics Failure ²⁾	0

¹⁾ Meaning → below, table Filament status.

²⁾ Electronics Failure includes a sensor error.

The Sensor Alarm bits defined here are copied into the Device Block ID 27 in the "Device Exception Detail Alarm" range of the BA measuring part.

3.3.3.4 Sensor Warning ID 104

This parameter indicates the detectable warnings occurring in connection with the BA measuring part.

	Bit 7 ... 4	Bit 3	Bit 2	Bit 1	Bit 0
Byte 0	0	0	0	Sensor Filament 2 Warning ¹⁾	Sensor Filament 1 Warning ¹⁾
Byte 1	0	Pressure too high for degas ²⁾	0	Electronics Warning	0

¹⁾ Meaning → below, table Filament status.

²⁾ Pressure too high for degas: the bit is set if the pressure value is $>7.2 \times 10^{-6}$ mbar during a degas command. The bit is reset if the pressure value drops below 7.2×10^{-6} mbar.

The Sensor Warning bits defined here are copied into the Device Block ID 28 in the "Device Exception Detail Warning" range of the BA measuring part.

Filament status

Byte 0 of Sensor Alarm and Sensor Warning represents the filament status of the gauge.

Bit 1	Bit 0	Filament status	Sensor Filament 1 Warning	Sensor Filament 2 Warning	Sensor Filament 1 Alarm	Sensor Filament 2 Alarm
0	0	both filaments OK	0	0	0	0
0	1	filament 1 interrupted, filament 2 OK	1	0	0	0
1	0	filament 1 OK, Filament 2 interrupted	0	1	0	0
1	1	Filament 1 interrupted, Filament 2 interrupted	1	1	1	1

3.3.3.5 Emission Status ID 105

This parameter indicates the emission status of the gauge.

Emission status	Meaning
FALSE	Emission Off
TRUE	Emission On

3.3.3.6 Emission Current ID 106

This parameter indicates the value of the emission current in mA.

Emission currents [mA]
0
0.025
5
20

3.3.3.7 Active Filament ID 108

The Parameter defines / indicates the active filament (→ Mode Filament Selection).

Active Filament	Byte value	Meaning
1	1	filament 1 active
2	2	filament 2 active

3.3.3.8 Degas Status ID 109

This parameter indicates whether the gauge is in Degas mode.

Degas Status	Meaning
FALSE	Degas Off
TRUE	Degas On

3.3.3.9 Active Degas Filament ID 116

For the FRG-730, the value of the parameter is the same as the parameter value of Active Filament (ID 108).

3.3.3.10 Mode Filament Selection ID 119

The parameter defines the operation mode of the filaments:

Mode Filament Selection (Filament User Mode)	Byte value	Meaning
Auto	0	Automatic Mode (default): The gauge alternates between filaments every time the emission is switched on.
Manual	1	Manual Mode: The gauge uses the filament selected by the attribute Active Filament (ID 108).

3.3.3.11 Emission User Mode State ID 201

This parameter indicates whether the gauge is in Automatic or Manual Emission Mode. The mode is set by Transition command 18 (ID 14). Automatic Mode is selected by default.

Emission User Mode State	Byte value	Meaning
FALSE	0	Manual Mode
TRUE	1	Automatic Mode

3.3.3.12 Hot Cathode Block State Command ID 14

This service is used for activating degas and emission modes via Profibus.

ID	Name	Structure	Data type	Bytes	Access	Range
14	Hot Cathode Block State Command	Rec	→ below	2	1/2_W	

Byte 0	Transition Command	Simple	UINT8	1		→ below
Byte 1	Value	Simple	UINT8	1		→ below

Transition Command	Name	Description
0	Inactive	No action
1	Set Degas State	Activates/deactivates the Degas mode This service is used for activating/deactivating the Degas mode ($p < 7.2 \times 10^{-6}$ mbar). If the Degas mode has not been deactivated with Degas Off before, it is automatically turned off after 3 minutes. The Degas State can only be changed if $p < 7.2 \times 10^{-6}$ mbar, otherwise the command returns the error "Object State Conflict". The Transition Command Data Field Value can have the following values: 0 ⇒ Degas Off 1 ⇒ Degas On
3	Set Emission State	Turns the emission On or Off according to the target value 0 ⇒ Switch emission Off 1 ⇒ Switch emission On
128	Emission User Mode	The device supports two modes: Automatic Mode: Emission is switched on or off by the integrated Pirani automatically. If the emission is switched on and is then switched off manually, the emission will be switched on again only if the pressure rises above the 3.2×10^{-2} mbar and then drops below 2.4×10^{-2} mbar. Manual Mode: Emission is switched on or off by the command "Set Emission State" - If during run time the pressure exceeds 3.2×10^{-2} mbar, the emission is switched off automatically. The measurement value equals the Pirani value. - If the emission is switched off by the user while the pressure is still in BA range, the measurement value equals the Pirani value. There is no automatic change to the BA range by the FRG-730. 0 ⇒ sets Manual Mode 1 ⇒ sets Automatic Mode

Appendix A: Definitions

Data types

Abbreviation	Range	Data type
INT8	$-2^7 \dots (2^7 - 1)$	Integer 1 byte
INT16	$-2^{15} \dots (2^{15} - 1)$	Integer 2 byte
INT32	$-2^{31} \dots (2^{31} - 1)$	Integer 4 byte
UINT8	$0 \dots (2^8 - 1)$	Unsigned integer 1 byte
UINT16	$0 \dots (2^{16} - 1)$	Unsigned integer 2 byte
UINT32	$0 \dots (2^{32} - 1)$	Unsigned integer 4 byte
FLOAT	$\pm 3.402 \times 10^{38}$	Floating Point, IEEE 754 Short Real Number, 4 byte
VSTRING(n)		ISO 646 and ISO 2375
OSTRING(n)		Octet string

Definitions

Term	Meaning
Store	This parameter defines whether the values are stored in non-volatile memory (→ store characteristics)
Default	Manufacturer-defined value

Store characteristics

Abbreviation	Meaning
V	"Volatile": Value is not saved to the RAM or EEPROM and is lost in the event of a power failure
N	"Nonvolatile": Value is saved to the RAM or EEPROM and is not lost in the event of a power failure

Data access

Abbreviation	Meaning
1_R/W	Acyclically readable and writeable by a Master Class 1
2_R/W	Acyclically readable and writeable by a Master Class 2
1/2_R/W	Acyclically readable and writeable by a Master Class 1 and 2
1_R	Acyclically readable by a master Class 1
2_R	Acyclically readable by a master Class 2
1/2_R	Acyclically readable by a master Class 1 and 2
1_W	Acyclically writeable by a master Class 1
2_W	Acyclically writeable by a master Class 2
1/2_W	Acyclically writeable by a master Class 1 and 2
0_XI	Cyclic output data with master Class 1

Excerpts from:
"PROFIBUS Profile for SEMI"
(→ [3], [5], [6])

The following table explains terms used in connection with the Profibus.

Term	Meaning
Alert Elements	Alert Elements are used to communicate notification messages from slave to master when warnings, alarms or events are detected.
Application	A <i>software functional unit</i> consisting of an interconnected aggregation of <i>function blocks</i> , <i>events</i> and <i>objects</i> , which may be distributed and which may have <i>interfaces</i> with other <i>applications</i> .
Characteristic	An characteristic is a property or characteristic of an <i>entity</i> . (Au) In block applications a block interface is defined by input/output parameters. These parameters have characteristics called parameter characteristics. Examples are access rights and identification names. (IT) The UML defines characteristics as a feature within a classifier that describes a range of values that instances of the classifier may hold. It is a property of a class instance (object).
Block (Block Instance)	A logical processing unit of software comprising an individual, named copy of the block and associated parameters specified by a block type, which persists from one invocation of the block to the next. Concept similar to the class/object approach, but well suited to the automation requirements.
Class	(IT) A class represents a template for several objects and describes how these objects are structured internally. Objects of the same class have the same definition both for their operations and for their information structures.
Configuration (of a system/device)	A step in system design: selecting functional units, assigning their locations and identifiers and defining their interconnections.
Data Structure	An <i>aggregate</i> whose elements need not be of the same <i>data type</i> , and each of them is uniquely referenced by an <i>offset identifier</i> .
Data Type	A data item with certain characteristics and permissible operations on that data, e.g. INT8.
Device	A physical entity capable of performing one or more specified functions in a particular context and delimited by its interfaces.
Direction of Data	<i>Input data</i> are transmitted from the device to the bus. <i>Output data</i> are transmitted from the bus to the device.
Direction of Flow	A positive set point causes a flow from P to A.
Entity	A particular thing, such as a person, place, <i>process</i> , object, concept, association or <i>event</i> .
Function	(1) A specific purpose of an entity. (2) One of a group of actions performed by an entity.
Function Block	A named <i>block</i> consisting of one or more input, output and contained parameters. Function blocks represent the basic automation functions performed by an application which is as independent as possible from the specifics of I/O devices and the network. Each function block processes input parameters according to a specified algorithm and an internal set of contained parameters. They produce output parameters that are available for use within the same function block application or by other function block applications.

Excerpts from:
"PROFIBUS Profile for SEMI"
(cont.)

Term	Meaning
Function Block Application	Application of an automation system performed by a Device Block, Function Block, Transducer Block and accompanied elements.
Instance	A set of data related to an invocation of a function block or a class.
Internal Resolution (ir)	The internal resolution is 16383 (3FFF _{hex}) for 100% and -16384 (C000 _{hex}) for -100% of the range.
Mode	Determines the block operating mode and available modes for a block instance.
Object	(IT) A software entity having identity, attributes and behavior.
Parameter	A <i>variable</i> that is given a constant value for a specified <i>application</i> and that may denote the <i>application</i> .
Device Block	A Device Block is a named block. Hardware specific parameters of a field device, which are associated with a resource, are made visible through the Device Block. Similar to transducer blocks, they insulate function blocks from the physical hardware by a set of implementation independent hardware parameters.
Record	A set of <i>data items</i> of different data types treated as a unit.
Resource	A resource is considered to be a logical subdivision within the software (and possibly hardware) structure of a device. Resources have independent control of their operation. The definition of a resource may be modified without affecting other resources within a device. A resource accepts and processes data and/or events from the process and/or communication interfaces and returns data and/or events to the process and/or communication interfaces, as specified by the applications utilizing the resource. An interoperable network view of applications is provided through device resources. Each resource specifies the network visible aspects of one or more local applications (or parts of distributed applications).
Simple Variable	A single variable which is characterized by a defined Data Type.
Substitute Value	In case an optional parameter has not been implemented, the device behaves according to the substitute value for this parameter.
Transducer Block	Transducer Block is a named block. Transducer blocks insulate function blocks from the specifics of I/O devices, such as sensors, actuators, and switches. Transducer blocks control access to I/O devices through a device independent interface defined for use by function blocks. Transducer blocks also perform functions, such as calibration and linearization, on I/O data to convert it to a device independent representation. Their interface to function blocks is defined as one or more implementation independent I/O channels.
Variable	A <i>software</i> entity that may assume any one of a set of values. The values of a variable are usually restricted to a certain data type.

Appendix B: Block Type

Currently defined Block Type IDs

Block Name	Block Type ID
Device Block	1
Sensor Analog Input Function Block	2
One of N Channel Sensor Analog Input Function Block	3
Multi Channel Sensor Analog Input Function Block	4
Discrete Input Function Block	5
Actuation Analog Output Function Block	6
Discrete Output Function Block	7
Analog Output Function Block	8
Single Stage Controller Function Block	9
Gas Calibration Transducer Block	10
Flow Transducer Block	11
Sensor Analog Input Ambient Temperature Transducer Block	12
Heat Transfer Vacuum Gauge... ¹⁾	13
Diaphragm Gauge	14
Cold Cathode Ion Gauge	15
Hot Cathode Ion Gauge ²⁾	16
Trip Point Function Block	17
Reserved	18 ... $2^8 - 1$
Manufacturer-specific	$2^8 \dots 2^{16} - 1$

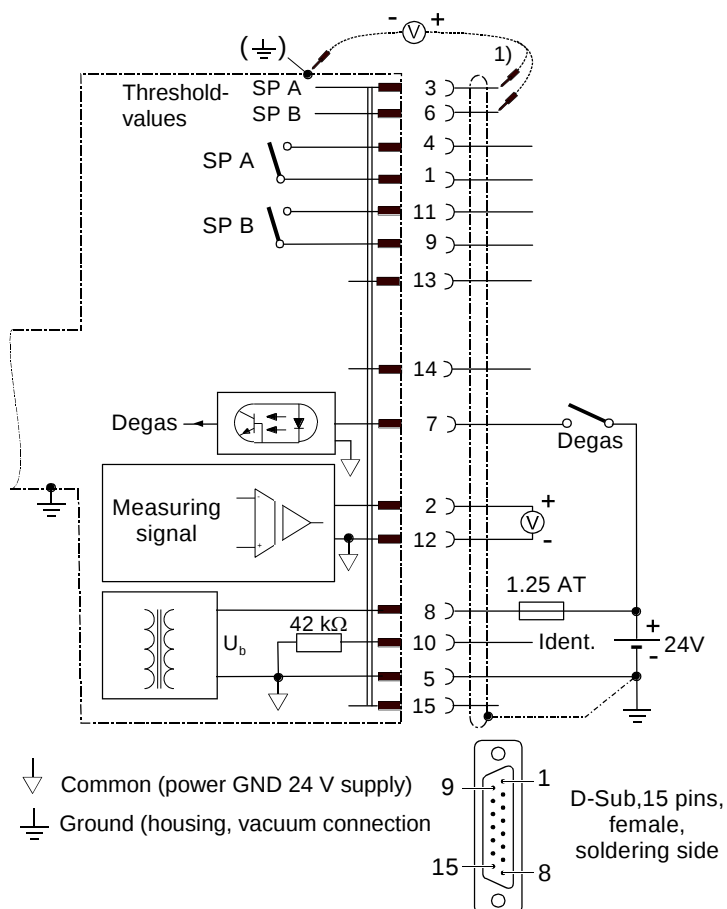
¹⁾ Pirani measuring system

²⁾ BA measuring system

Appendix C: Electrical Connections

Technical data of FRG-730 gauges → [1], [2].

Sensor cable connection

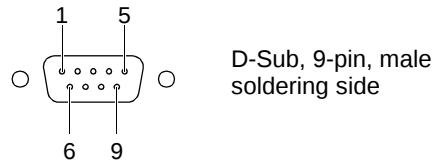


Electrical connection

Pin 1	Relay switching function A, com contact	
Pin 2	Measuring signal output	0 ... +10 V
Pin 3	Threshold (setpoint) A	1) 0 ... +10 V
Pin 4	Relay switching function A, n.o. contact	
Pin 5	Supply common	0 V
Pin 6	Threshold (setpoint) B	1) 0 ... +10 V
Pin 7	Degas on, active high	0 V/+24 V
Pin 8	Supply	+24 V
Pin 9	Relay switching function B, com contact	
Pin 10	Gauge identification	
Pin 11	Relay switching function B, n.o. contact	
Pin 12	Measuring signal common	
Pin 13	Do not connect	
Pin 14	Do not connect	
Pin 15	Do not connect	

¹⁾ Do not connect pin 3 and pin 6 for normal operation of the gauge. These pins are reserved for adjustment of the setpoint potentiometers.

Profibus connection



Pin 1 not connected

Pin 2 not connected

Pin 3 RxD/TxD-P

Pin 4 CNTR-P ¹⁾

Pin 5 DGND ²⁾

Pin 6 VP ²⁾

Pin 7 not connected

Pin 8 RxD/TxD-N

Pin 9 not connected

¹⁾ Only to be connected if an *optical link* module is used.

²⁾ Only required as line termination for devices at both ends of bus cable
(→ [3]).

Appendix D: Literature

-  [1] www.agilent.com
Short Operating Instructions
FRG-730
tqma72e1
Agilent Technologies, Lexington, MA 02421, USA
-  [2] www.agilent.com
Operating Manual
FRG-730
tqna72e1
Agilent Technologies, Lexington, MA 02421, USA
-  [3] www.profibus.com
(Profibus user organization)
-  [4] European Standard for Profibus EN 50170
-  [5] Profile for Semi-Conductor Devices; Part 1
Common Definitions v1.0; 8/2002
-  [6] Profile for Semi-Conductor Devices; Part 3
Vacuum Pressure Gauges v1.0; 8/2002

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>



Agilent Technologies

**Vacuum Products Division
Request for Return Form
(Health and Safety Certification)**

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: ☐ Taxable ☐ 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
- Does not start - Does not spin freely - Does not reach full speed - Mechanical Contact - Cooling defective	- Noise - Vibrations - Leak - Overtemperature - Clogging	- Vertical - Horizontal - Upside-down - Other: _____
		Power: _____ Rotational Speed: _____ Current: _____ Inlet Pressure: _____ Temp 1: _____ Foreline Pressure: _____ Temp 2: _____ Purge flow: _____ OPERATING TIME: _____

ION PUMPS/CONTROLLERS

- Bad feedthrough - Vacuum leak - Error code on display	- Poor vacuum - High voltage problem - Other
---	--

VALVES/COMPONENTS

- Main seal leak - Solenoid failure - Damaged sealing area	- Bellows leak - Damaged flange - Other
--	---

LEAK DETECTORS

- Cannot calibrate - Vacuum system unstable - Failed to start	- No zero/high background - Cannot reach test mode - Other
---	--

INSTRUMENTS

- Gauge tube not working - Communication failure - Error code on display	- Display problem - Degas not working - Other
--	---

SCROLL AND ROTARY VANE PUMPS

- Pump doesn't start - Doesn't reach vacuum - Pump seized	- Noisy pump (describe) - Over temperature - Other
---	--

DIFFUSION PUMPS

- Heater failure - Doesn't reach vacuum - Vacuum leak	- 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
121 Hartwell Avenue
Lexington, MA 02421 USA
Tel.: +1 781 861 7200
Toll-Free: +1 800 882 7426
Fax: +1 781 860 5437
vpl-customerservice@agilent.com

Benelux

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

China

Agilent Technologies (China) Co. Ltd
No.3, Wang Jing Bei Lu,
Chao Yang District, Beijing,
100102 China Tel.: +86
(10) 6439 7888
Fax: +86 (10) 6439 1318
Toll-Free: 800 820 8266
vpc-customerservice@agilent.com

France

Agilent Technologies France
7 avenue des Tropiques
Z.A. de Courtaboeuf - B.P. 12
91941 Les Ulis cedex France
Tel.: +33 (0) 1 69 86 38 84
Fax: +33 (0) 1 69 86 29 88
Toll free: 00 800 234 234 00
vpf.sales@agilent.com

Germany & Austria

Agilent Technologies
Lyoner Str. 20
60 528 Frankfurt am Main Germany
Tel.: +49 69 6773 43 2230
Fax: +49 69 6773 43 2250
Toll free: 00 800 234 234 00

India

Agilent Technologies India Pvt. Ltd.
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
cag_india@agilent.com

Italy

Agilent Technologies Italia S.p.A.
via F.lli Varian 54
10040 Leini, (Torino) Italy
Tel.: +39 011 997 9111
Fax: +39 011 997 9350
Toll-Free: 00 800 234 234 00
vpt.sales@agilent.com
vpt-customerservice@agilent.com

Japan

Agilent Technologies Japan, Ltd.
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
vpj-customerservice@agilent.com

Korea

Agilent Technologies Korea Ltd.
Shinsa 2nd Bldg. 2F
966-5 Daechi-dong
Kangnam-gu, Seoul Korea 135-280
Tel.: +82 2 3452 2455
Toll-Free: 080 222 2452
Fax: +82 2 3452 2451
vpk-customerservice@agilent.com

Singapore

Agilent Technologies Singapore Pte Ltd
No.1 Yishun Avenue 7 Singapore 768923
Tel: +65 6215 8045
Fax: +65 6754 0574
Toll-Free: 1 800 2762622
vps-customerservice@agilent.com

South East Asia

Agilent Technologies Sales Sdn Bhd
Unit 201, Level 2 uptown 2,
2 Jalan SS21/37, Damansara Uptown
47400 Petaling Jaya, Selangor, Malaysia
Tel: +603 7712 6106
Fax: +603 6733 8121
Toll-Free: 1 800 880 805
vps-customerservice@agilent.com

Taiwan

Agilent Technologies Taiwan Limited
20 Kao-Shuang Rd.,
Pin-Chen City, 324
Taoyuan Hsien, Taiwan, R.O.C.
Tel. +886 34959281
Toll Free: 0800 051 342
vpw-customerservice@agilent.com

UK & Ireland

Agilent Technologies UK Ltd
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
vpt-customerservice@agilent.com

Learn more:

www.agilent.com/chem/vacuum

This information is subject to change
without notice.

© Agilent Technologies, Inc., 2011
Published in USA, October, 2011



Agilent Technologies