

Ion Pumps Control Unit

Model
529-5001R001

User Manual

87-900-128-01 (E)
03/2014



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2014

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

Manual Part Number

Publication Number: 87-900-128-01 (E)

Edition

Edition 03/2014

Printed in ITALY

Agilent Technologies Italia S.p.A.

Vacuum Products Division

Via F.lli Varian, 54

10040 Leini (TO)

ITALY

Warranty

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

Technology Licenses

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

Restricted Rights Legend

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or

contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

Trademarks

Windows and MS Windows are U.S. registered trademarks of Microsoft Corporation.

Safety Notices

CAUTION

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

WARNING

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

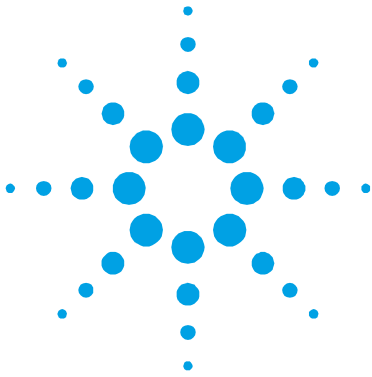
Ion Pumps Control Unit



Contents

1	Installation procedure	7
	General Information	8
	Symbols used	9
	Controller Description	10
	Storage	10
	Preparation for Installation	111
	Installation	13
	Use	14
	Usage	18
	Operating modes	19
	Error Messages	22
	Maintenance	24
	Disposal	25
2	Technical Information	26
	Ion Pump Controller Description	28
	PIN Functions	38

Contents



1 Installation Procedure

General Information	8
Symbols used	9
Controller Description	10
Storage	10
Preparation for installation	11
Installation	13
Use	14
Ion Pumps C.U. Front Panel Controls and Indicators	15
Ion Pumps C.U. Rear Panel	17
Usage	18
Operating modes	19
Power Line supply	19
LOCAL-REMOTE I/O Operation	19
SERIAL Operation	20
START/PROTECT Operation	21
Error Messages	22
On Front Panel	22
Maintenance	24
Disposal	25

Original Instructions



General Information

This equipment is destined for use by professionals. The user should read this instruction manual and any other additional information supplied by Agilent before operating the equipment. Agilent will not be held responsible for any events occurring due to non-compliance, even partial, with these instructions, improper use by untrained persons, non-authorized interference with the equipment or any action contrary to that provided for by specific national standards.

The following paragraphs contain all the information necessary to guarantee the safety of the operator when using the equipment. Detailed information is supplied in the appendix "Technical Information".

This manual uses the following standard protocol:

CAUTION!

The caution messages are displayed before procedures which, if not followed, could cause damage to the equipment.

WARNING!







The warning messages are for attracting the attention of the operator to a particular procedure or practice which, if not followed correctly, could lead to serious injury.

NOTE

The notes contain important information taken from the text.

Symbols used

The following symbols are used consistently throughout in all illustration:

Symbol	Description
	On (Power)
	Off (Power)
	Caution, risk of electric shock.
	Warning “see installation/operating mode instruction”

Controller description

Agilent's controller is an ion pumps controller.

It can drive up to 2 Ion Pumps simultaneously and independently. The controller is designed to give the HV to the Ion Pumps when it is connected to the Mains supply (Max. Output Voltage = 5000 Vdc). (See the paragraph "Technical Informations" for details).

Agilent's Ion Pump controller can be driven in Local/Remote I/O operating mode or in the Serial mode via the RS232 port. The selection of the operating mode is done by means of the SERIAL/LOCAL-Remote I/O switch on the controller front panel.

In the LOCAL-REMOTE I/O mode, all the commands are given by the front panel or by the connections to the DB25 "REMOTE" connector on the rear panel.

The pump operating conditions data are anyway available on the serial connection. In the SERIAL mode, all the commands are given through the serial connection and all the pump operating conditions data are available on the serial connection.

Storage

When transporting and storing the controller, the following environmental requirements should not be exceeded:

- temperature: from -20° to +70 °C
- relative humidity: 0 – 95 % (non-condensing)

Preparation for installation

The controller is supplied in a special protective packing. If this shows signs of damage which may have occurred during transport, contact your local sales office. When unpacking the controller, be sure not to drop it and avoid any kind of sudden impact or shock vibration to it.

Do not dispose of the packing materials in an unauthorized manner. The material is 100 % recyclable and complies with EEC Directive 85/399.

1 Installation Procedure
Preparation for installation

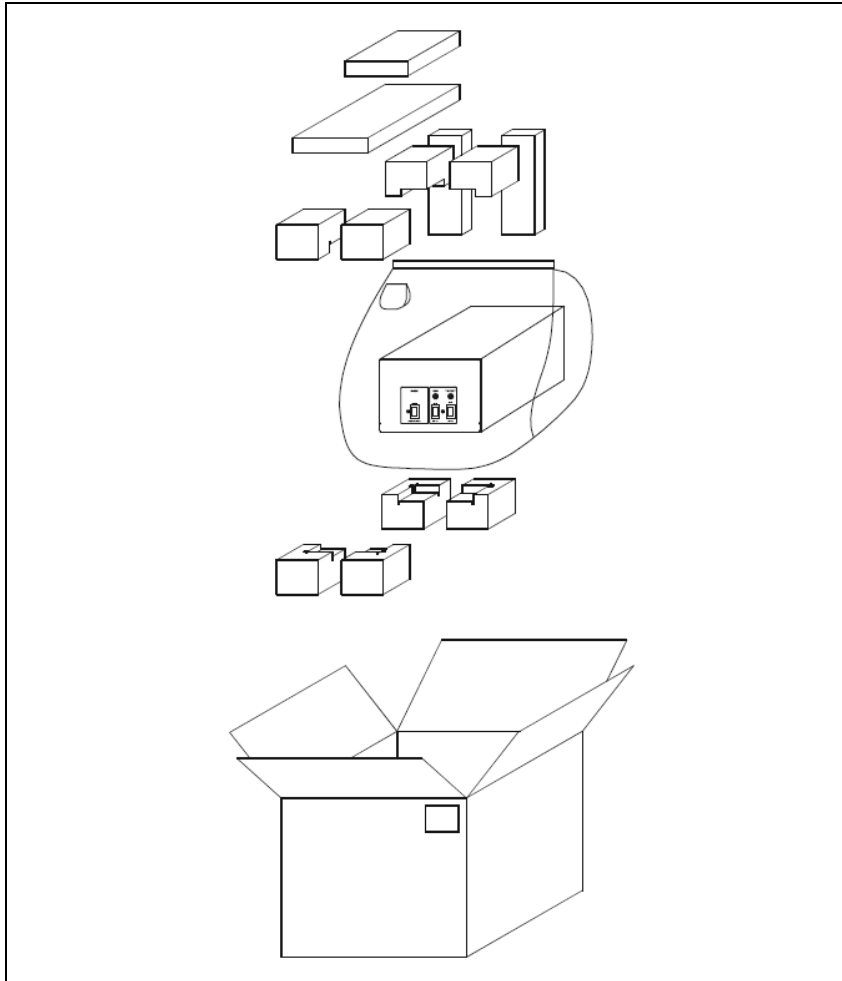


Figure 1 Packing

Installation

WARNING!



The controller is equipped with a 3-wire power cord without plug. Install the proper plug foreseen by the country where the controller will be used.

Note that the Yellow/Green wire of the cable is the Ground connection of the controller. Use this power cord and the installed plug in conjunction with a properly grounded power socket to avoid electrical shock. Verify that the ground connections are properly connected. High voltage developed in the controller can cause severe injury or death. Before servicing the unit, disconnect the power cable.

CAUTION!

The controller can be used as a bench unit or as a rack module, but it must be positioned so that free air can flow through the holes of the top and the side cover. If the controller is used as a rack module, it MUST be inserted in a 4 unit height rack adapter chassis to avoid the controller to fall inside the rack. The Controller front panel is not designed to support the unit weight. Do not install or use the controller in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

NOTE

If the controller is installed on a rack, remove the four feet and position it with at least 30 mm (1.2 inches) of clearance on top and bottom.

1 Installation Procedure

Use

WARNING!



The controller must be installed in a way that allows an easy interruption of the line voltage (disconnection of the line plug or interruption on the rack general line switch).

During operation, the following environmental conditions must be respected:

- temperature: from +5 °C to +40 °C
- relative humidity: 0 – 95 % (non-condensing)

To connect the controller to the pump use the specific cables supplied.

See the appendix “Technical Information” for detailed information about the above mentioned and the other connections.

Use

This paragraph describes the fundamental operating procedures. Detailed information and operating procedures that involve optional connections or options are supplied in the paragraph “USE” of the appendix “Technical Information”.

Make all vacuum manifold and electrical connections and refer to the connected pump instruction manual prior to operating the controller..

WARNING!



To avoid injury to personnel and damage to the equipment, if the pump is laying on a table make sure it is steady. Never operate the pump if the pump inlet is not connected to the system or blanked off.

Ion Pumps C.U. Front Panel Controls and Indicators

The figure and the table of the following page illustrate the Controller front panel controls and indicators with relevant description and function.

NOTE

Some controls and indicators change their function and meaning according to the position of HV ON OFF switch. The table describes all function and meaning.

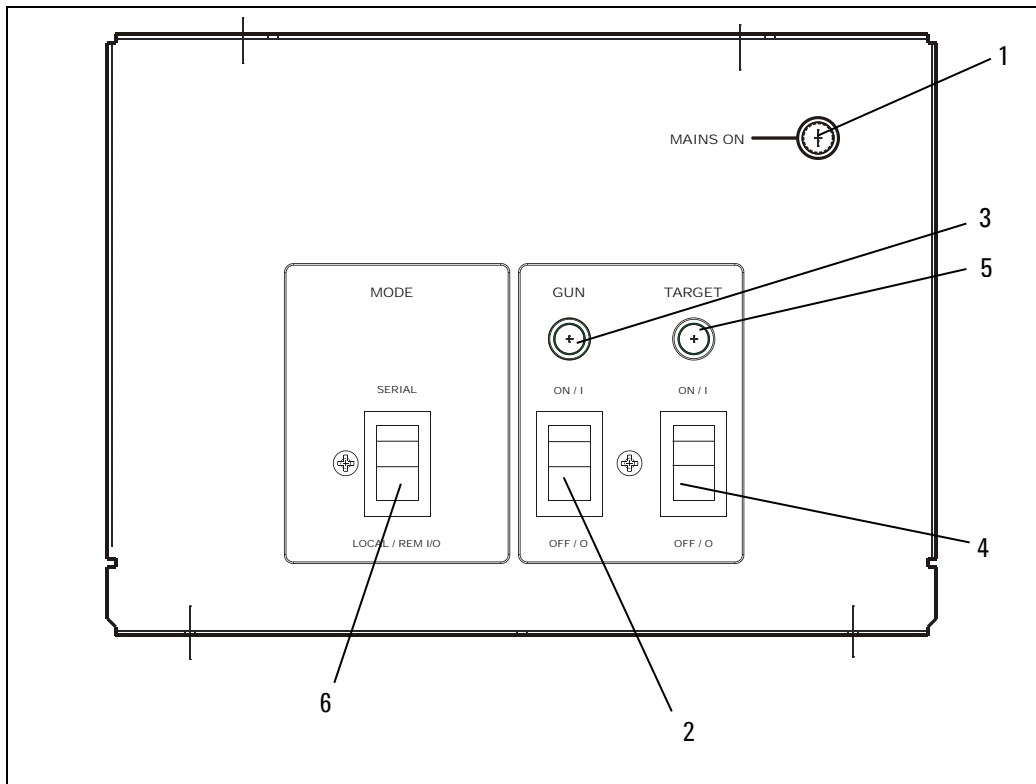


Figure 2 Ion Pumps C.U. Front Panel Controls and Indicators

1 Installation Procedure

Use

REF.	SIGNAL	DESCRIPTION
1	MAINS ON LED	When On it means that the unit is connected to the Power Line and the Line Switch on the rear panel is in ON (I) position.
2	GUN ON/OFF SWITCH	When in ON (I) position, it switches On the HV to the GUN PUMP if the controller is set for LOCAL/REMOTE I/O operation and the external and HV Cable interlocks are closed.
3	GUN LED	It is On bold when the HV GUN PUMP is on and the voltage is over 1250 V (+/- 250 V hysteresys) volt. It is slow flashing (1 flash at second) when the voltage on HV GUN PUMP is lower than 1250 V (+/- 250 V hysteresys) volt. It is fast flashing in case of Fault (see paragraph "Technical Information" for details)
4	TARGET ON/OFF SWITCH	When in ON (I) position, it switches On the HV to the TARGET PUMP if the controller is set for LOCAL/REMOTE I/O operation and the external and HV Cable interlocks are closed.
5	TARGET LED	It is On bold when the HV TARGET PUMP is on and the voltage is over 1250 V (+/- 250 V hysteresys). It is slow flashing (1 flash at second) when the voltage on HV TARGET PUMP is lower than 1250 V (+/- 250 V hysteresys). It is fast flashing in case of Fault (see paragraph "Technical Information" for details).
6	SERIAL/LOCAL-REM I/O SWITCH	It allows to select the operating mode: SERIAL: all the commands are given through the serial line. LOCAL-REM I/O: all the commands are given through the unit front panel or through the 25 pin REMOTE connector on the rear panel.

Ion Pumps C.U. Rear Panel

The controller rear panel controls and connections are shown in the following figure. For detailed information about the connections see the appendix “Technical Information”.

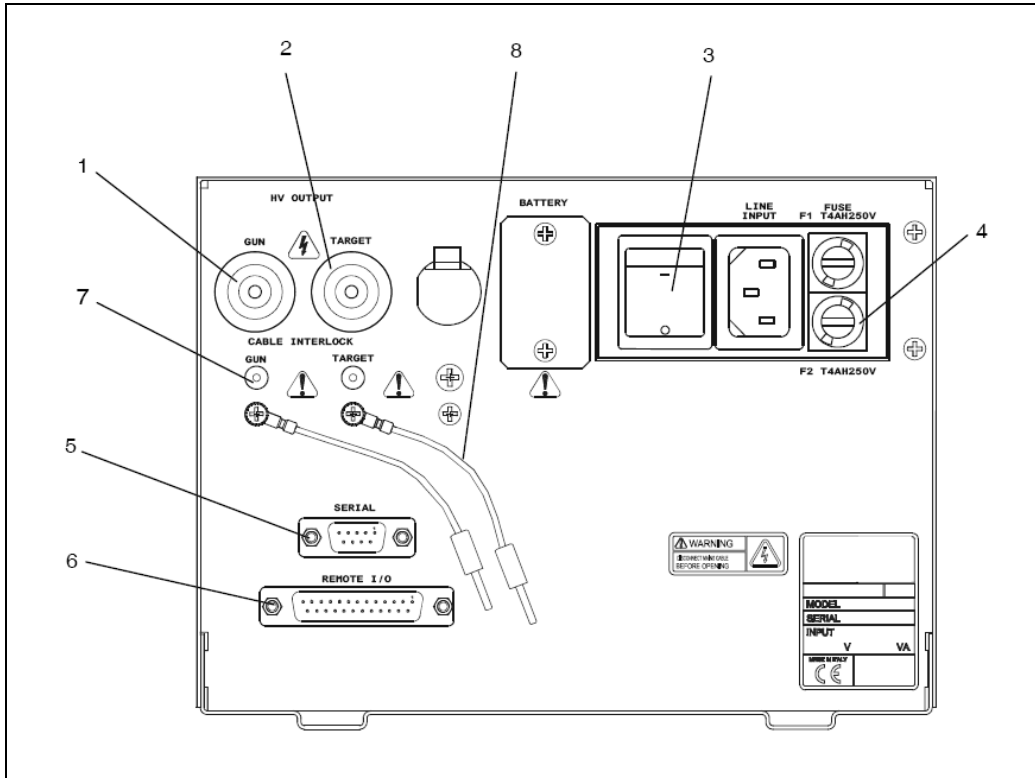


Figure 3 Ion Pumps C.U. Rear Panel

1 Installation Procedure

Usage

REF.	SIGNAL	DESCRIPTION
1	GUN Connector	HV Connector for Cable to GUN Pump
2	TARGET Connector	HV Connector for Cable to TARGET Pump
3	LINE INPUT	Module containing the Line Input cable plug and the Line Switch
4	FUSE	Protection Fuse holder (2 fuses)
5	SERIAL Connector	DB9 connector for the Serial cable
6	REMOTE Connector	DB25 connector for the REMOTE I/O connections
7	GUN Cable Interlock	
8	TARGET Cable Interlock	

Usage

This paragraph gives a general information on the use of the controller. For details refer to the “Technical Information” paragraph. When the controller is plugged to the Power Line and the Line Switch on the rear panel is ON (I), it is able to operate with its full set of features:

- 2 independent HV channels
- +5000 Vdc maximum output Voltage
- 40 W maximum output Power for channel 1
- 80 W maximum output Power for channel 2
- 30 mA maximum output current (short circuit condition) for gun
- 50 mA maximum output current (short circuit condition) for target
- LOCAL-REMOTE I/O/SERIAL operating mode
- START/PROTECT operation

Operating modes

Power Line supply

LOCAL-REMOTE I/O Operation

NOTE

The 2 HV switches on the front panel and the 2 HV On/Off inputs on the DB25 “Remote” connector on the rear panel are functionally in series.

HV On/Off gun by front panel:

Connect together pin 8 and pin 21 of the remote I/O connector and move the gun switch on the front panel.

HV On/Off target by front panel:

Connect together pin 9 and pin 22 of the Remote I/O connector and move the gun switch on the front panel.

The correspondent HV cable Interlocks must be closed.

When the HV is ON, the corresponding HV LED on the front panel, will be ON.

WARNING!



Auto – start function:ON.

In case of power loss with GUN and/or TARGET ON, at next power-on the unit will switch on the HV (previously on) automatically.

Auto power-on function.

With front panel switches in ON position and remote I/O jumpers, the unit will switch on both channels as soon as the mains is on.

1 Installation Procedure

Operating modes

NOTE

If the connection between the IPCU and the pump is done by means of an HV Cable with Safety Interlock, the cable must be connected to the pump F/T and the small “banana” connector on the cable controller side must be inserted in the corresponding “HV Cable Interlock” plug on the unit rear panel in order to be able to switch on the HV.

If the connection is done with a single pole HV Cable without the Safety Interlock, ensure that the jumper provided on the unit rear panel is inserted in the “HV Cable Interlock” plug.

SERIAL Operation

NOTE

The SERIAL is set in order to operate by connecting the controller to an external PC that uses the “HYPER TERMINAL” or equivalent program (for example a proprietary user program) running under Windows Operating System or other. To operate in serial mode, connect the serial port to the controller “Serial” connector on the rear panel, run the Windows “HYPER TERMINAL” program on the PC with the following default settings:

- COM1
- 9600 baud
- No parity
- data bits
- 1 stop bit

This way of operation will allow to verify the functionality of the unit in the “SERIAL” operating mode, give the commands and get all the operating data.

NOTE

In the serial operation, depending on the user environment, the O.S. used must have real time capability.

See the “Technical Information” paragraph for details.

START/PROTECT Operation

The controller can operate in two modes: Start and Protect.

In Start, the 2 output channels can provide all the output power regardless of the pump condition up to the short circuit condition.

This operating mode must be used to start the pump at high pressure.

In Protect, the controller limits the output current and switches off the HV when the current exceeds the threshold value (8 mA for 40 W channel and 16 mA for 80 W channel) for more than 2 seconds.

The default mode is “Start” mode.

Error Messages

On Front Panel

During the controller operation, if an error condition is detected, the HV LEDs on the front panel will give the indication:

GUN, TARGET

LEDS ON: Normal Operation = HV ON

Led GUN, TARGET flashing: error condition:

- 1 flash every 4 seconds = Interlock HV cable
- 2 flashes every 4 seconds = remote Interlock
- 3 flashes every 4 seconds = HV Over voltage (This Fault doesn' t switch-off the HV)
- 4 flashes every 4 seconds = error inside ADC
- 5 flashes every 4 seconds = overcurrent on HV (short circuit)
- 6 flashes every 4 seconds = the Local/Remote switch was moved with HV ON
- 7 flashes every 4 seconds = over temperature inside unit
- 8 flashes every 4 seconds = overcurrent in Protect (8 mA for 40W channel and 16 mA for 80W channel)
- 9 flashes every 4 seconds = HV Under voltage

In Local-Remote I/O mode, the Reset of the Fault condition is accomplished by switching to Off (0) the front panel switch of the corresponding channel.

In Serial mode, the Reset of the Fault condition is accomplished by sending the "F0n" command by the computer.

NOTE

If one of the following conditions happens while the HV of the channel is On, the output of the channel is immediately set to Off, the LED on the front panel indicates the corresponding Fault condition and the Fault information is sent to the Serial output.

Auto retries function. If an under-voltage fail occur, for instance on GUN while both channels were on the unit do:

- Distinguish between a real low voltage and thermal switch (TS) intervention.
- Keep the Target ON, GUN switched off, no fail condition activated (fault relay remains open), current output recorder goes to 5 V.
- Start time out cont (30 min.)
- If the TS is restarted before the time out the gun goes on again and time out count resetted.
- If the TS is not restarted after the time out the IPCU goes into fail and both GUN and TARGET are off.
- On “Remote” connector on Rear Panel.
- “Fault” output: N.C. contact (idle state) =
- The contact is Open in “Fault” condition (relay Off)
- The contact is Close in normal condition (relay On)
- “HV 1,2” output: N.O. contact (idle state) =
- The contact is Open with HV in Off state (relay Off)
- The contact is Closed with HV in On state (relay On)
- “HV 1,2” output current set point: N.O. contact (idle state) =
- The contact is Open with output current value is below the current set point (relay Off)
- The contact is Closed with output current value above the current set point (relay On)

Maintenance

The Controller does not require any maintenance. Any work performed on the controller must be carried out by authorized personnel.

When a fault has occurred it is possible to use the Agilent repair service. Replacement controllers are available on an advance exchange basis through Agilent.

WARNING!

Before carrying out any work on the controller, disconnect it from the supply.



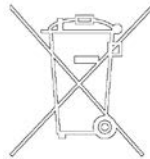
Disposal

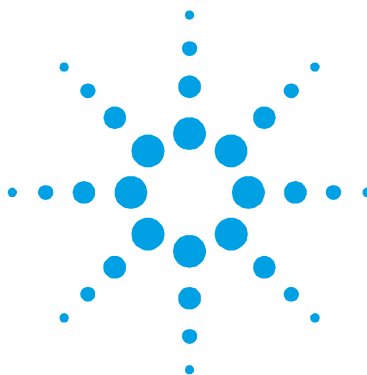
Meaning of the "WEEE" logo found in labels

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

This symbol (**valid only in countries of the European Community**) indicates that the product it applies to must NOT be disposed of together with ordinary domestic or industrial waste but must be sent to a differentiated waste collection system.

The end user is therefore invited to contact the supplier of the device, whether the Parent Company or a retailer, to initiate the collection and disposal process after checking the contractual terms and conditions of sale.





2 Technical Information

Ion Pump Controller Description	28
Controller Specifications	28
Remote I/O Connector Pinout	31
Input Commands Connection	33
Output Signals Connection	33
Analog Recorder Outputs	34
Serial Operation	36
Controller RS232 Standard Serial Pin outs (DB9 Female Connector)	37
PIN Functions	38
Procedure to Connect the Serial and I/O Ports to an External Cable	39



Ion Pump Control Unit - Operator / Host Serial
Communication 41

- Controller Report Mode 42
- Controller Commands Mode 42
- About Report 46
- Enable – Disable Remark Lines 47
- Continuous and NON Continuous Report Mode 47
- Enable - Disable Input ECO Characters 48
- About Commands 48
- Report TYPE 0...6 49
- User Operator Report Type 0 and User
Technician Report Type 1 49
- User Operator Report Type 2 and User
Technician Report Type 3 50
- User Operator Report Type 4 and User
Technician Report Type 5 51
- Old IPCU Releases - Compatibility Report Type 6 52
- Floating Point Notation 53
- Cleaning 54

Original Instructions

Ion Pump Controller Description

The Ion Pump controller is composed of a number of printed circuit boards:

- 1 board: (Power Supply), it converts the Line Voltage to the 24 Vdc that is used to generate the HV.
- 1 board: (Auxiliary Power Supply), it generates the + and – 12 Vdc for the analog circuitry, the 5 Vdc for the digital circuitry.
- 2 boards: (HV generation), they generate the HV for the 2 output channels.
- 1 board: (CPU), it contains all the digital circuitry and the microprocessor.
- 1 internal fan (OFF if for each channel the output current is <2 mA).

Controller Specifications

Tab. 1

Line:	
Voltage	100 - 240 Vac \pm 10 %
Frequency	50 / 60 Hz \pm 5 %
Max Power	200 VA
Operating conditions:	
Internal use only	
Max Altitude:	3000 m
Max Line Voltage variation:	\pm 10 %
Installation category:	II
Pollution degree:	2

Output for the pumps:	Default value =+5000 V +/-5 %, settable from +3000 to
Output Voltage (each channel)	+5000 Vdc at steps of +/- 50 V by Serial Line command
Output Current (each channel)	30 mA ± 10 % for 40 W output, 50 mA +/- 10 % for 80 W
Maximum Output Power GUN	output
Maximum Output Power TARGET	40 W
	80 W
Current Recorder Output signal	0 to 5 Vdc logarithmic, proportional to 1.0e-8 to 1.0e-1 A 1 recorder output for each pump. Minimum recorder input impedance: 1MΩ Precision: < than 1% Current Measurement Resolution: 10nA under 100μA; 1μA over 100μA
Operating Temperature	10 °C to 40 °C
Relative Humidity	0 – 95 % (non-condensing)
Storage Temperature	-20 °C to 70 °C
Compliance to Norms	Radio Interference: EN55011 Class A Group 1 Harmonics: EN61000/3/2 Flickers: EN61000/3/3 ESD: EN61000/4/2 4 kV Contact 8kV AIR Burst: EN61000/4/4 Vac Signal Radiated RF Immunity: EN61000/4/3 Surge: EN61000/4/5 Conducted noise imm.: EN61000/4/6 Line supply interruptions: EN61000/4/11 Compliance Safety: EN61010-1
Line fuse	Two T4AH250V
High Voltage Connection	2 High Voltage Fischer connectors

2 Technical Information

Ion Pump Controller Description

Installation according to international rules, only of the following HV cables Part No. may be used:

HV Cables P/N

929-0705	HV bakeable cable, radiation resistant, 4 meter
929-0707	HV bakeable cable, radiation resistant, 7 meter
929-0708	HV bakeable cable, radiation resistant, 10 meter
929-0709	HV bakeable cable, radiation resistant, 20 meter

NOTE

All connecting cables for remote I/O and serial connections, must be of the shielded type.

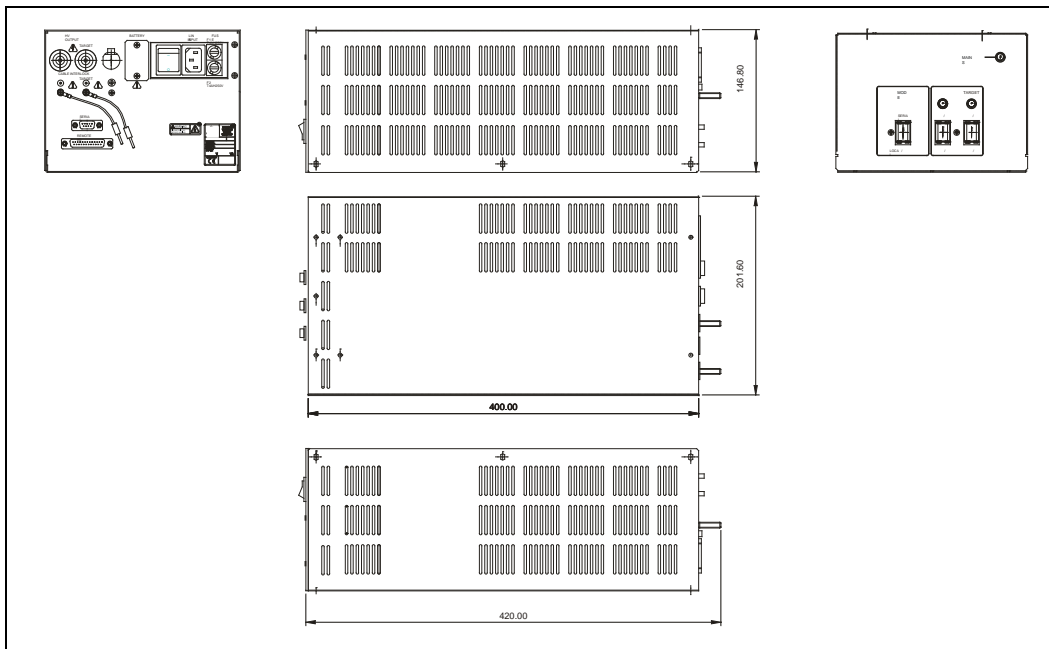


Figure 4 Controller Outline

Remote I/O Connector Pinout

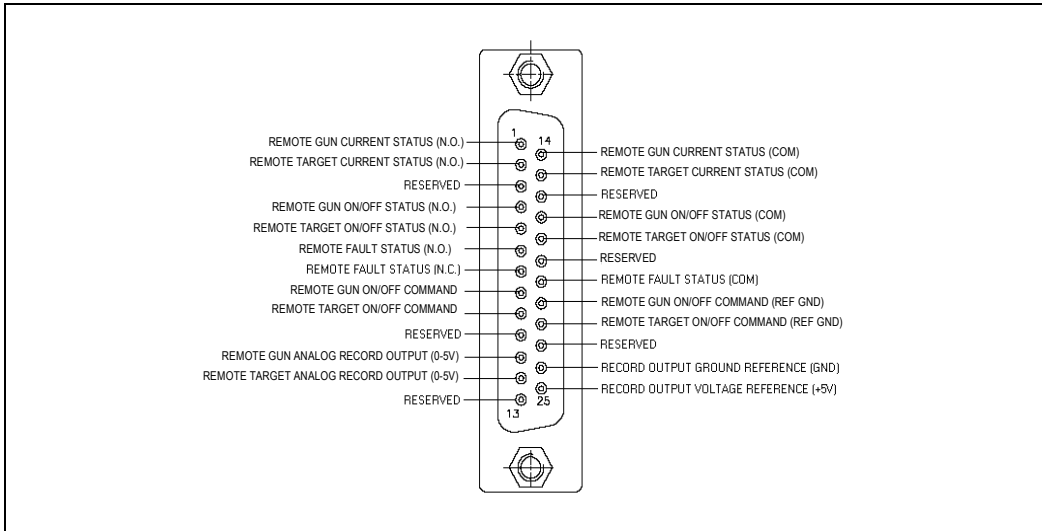


Figure 5 Remote I/O Pinout Rear Panel D-Shell Female 25P

Installation according to international rules.

All the input/output remote signals to/from the controller must be connected at mating connector. With the provided mating connector make the connection with AWG 24 (0.25 mm²) or smaller wire to the pins indicated in the figure to obtain the desired capability. It is a 15-pins D type connector; the available signals are detailed in the table, the following paragraphs describe the signal characteristics and use.

In case there is the need to use the Remote Interlock function or to switch On and Off the HV on the 2 pumps by Remote, remove the two jumpers and follow the indications given in the above input commands connection pin out drawing.

2 Technical Information

Ion Pump Controller Description

Tab. 2 Remote I/ O connector signals

25-Pin No.	IN - OUT	Name	Description
1-14	OUT	Remote GUN current status N.O.	Closed when Iout < Iset-point (2mA) Open when Iout > Iset-point (2mA)
2-15	OUT	Remote TARGET current status N.O.	Closed when Iout < Iset-point (2mA) Open when Iout > Iset-point (2mA)
4-17	OUT	Remote GUN status N.O.	Open: HV off Closed: HV on
5-18	OUT	Remote TARGET status N.O.	Open: HV off Closed: HV on
6-20	OUT	Remote FAULT status N.O.	Open: FAIL Closed: OK
7-20	OUT	Remote FAULT status N.C.	Open: OK Closed: FAIL
8-21	IN	Remote GUN ON/OFF command	Open: OFF Closed: ON
9-22	IN	Remote TARGET ON/OFF command	Open: OFF Closed: ON
11-24 (GND)	OUT	Remote GUN current analog record	0 to 5 volt (see below)
12-24 (GND)	OUT	Remote TARGET current analog record	0 to 5 volt (see below)

Input Commands Connection

Input signals: Optically Insulated

Insulated 24 Vdc (+10 %, -50 %) supply for input signal: Internally generated.

Output Signals Connection

Relay outputs.

Max voltage: 24 V

Max current: 200 mA

Min current: 10 μ A

Bouncing Time: NO=1msec; NC=3msec

Analog Recorder Outputs

The analog recorder outputs are two independent outputs for each channel.

The voltage is related to the log10 of the current absorbed by the pump.

The exact formula is:

$$V_{rec-out} = 1.35 \cdot \log_{10} I_{out} + 6,77$$

$$[V_{rec-out}] = \text{Volt}; [I_{out}] = \text{A}$$

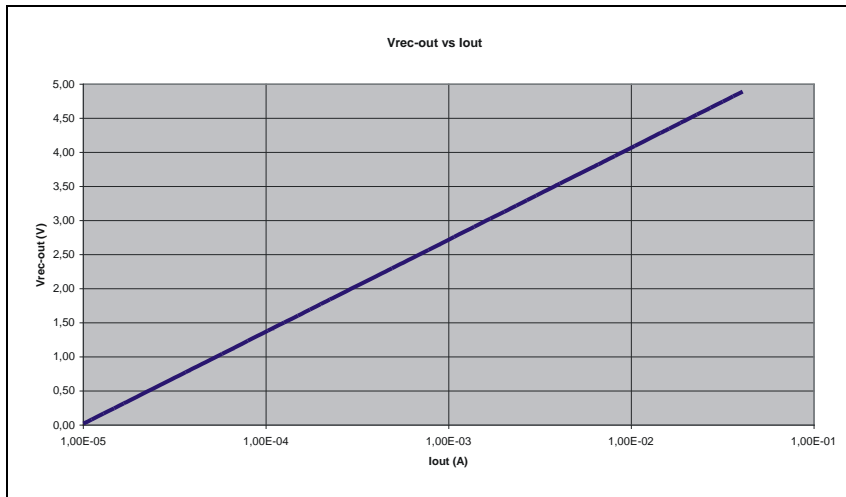


Figure 6

If the pump connected to the output is a 35l/s SEM the relation between Vrec-out and pressure is:

$$P[\text{mbar}] = 0,003 \cdot 10^{\frac{V_{\text{recout}} \cdot 6,77}{1,35}} + 5 \cdot 10^{-8}$$

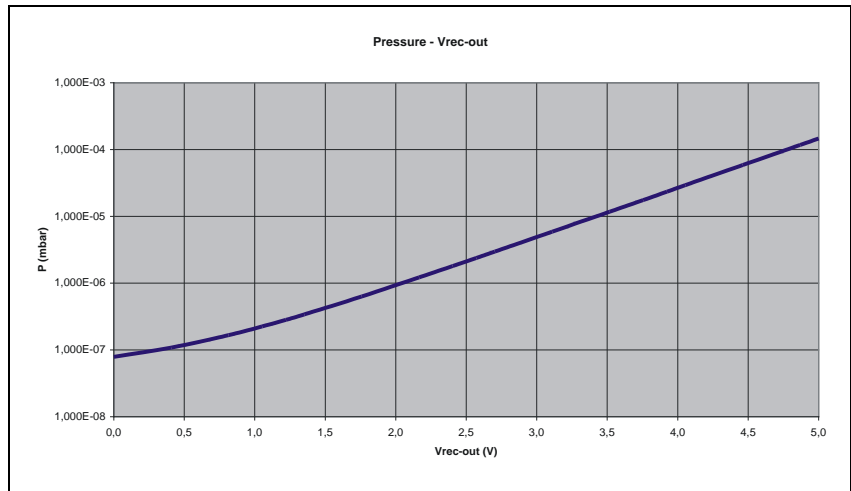


Figure 7

NOTE The current reading (using the analog recorder output or serial line) is fixed to the max current (30 mA for GUN, 50 mA for TARGET) until the HV voltage cross the 1.5 KV.

2 Technical Information

Ion Pump Controller Description

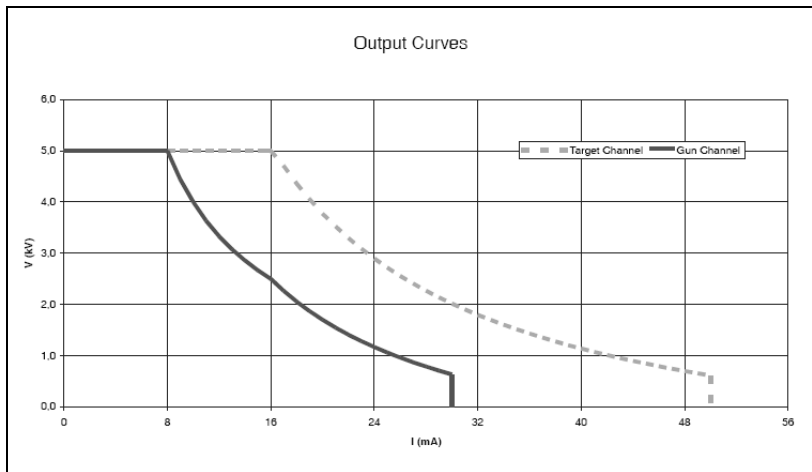


Figure 8

NOTE

At Pin 25 a reference voltage corresponding to the 2 recorder outputs full-scale value (5 V +/- 5 %) is given. This reference allows getting a more precise reading from the 2 recorder outputs. It allows the scaling of the actual reading to the real value of the full scale.

Serial Operation

The Controller works on RS232 DCE Mode (like a Modem with DB9 FEMALE connector).

The Host/Terminal/Personal Computer works always on RS232 DTE Mode (like a PC with DB9 MALE connector).

The Cable Interconnections are 1 to 1 signals (DB9 Male/Female Standard CABLE).

Controller RS232 Standard Serial Pin outs (DB9 Female Connector)

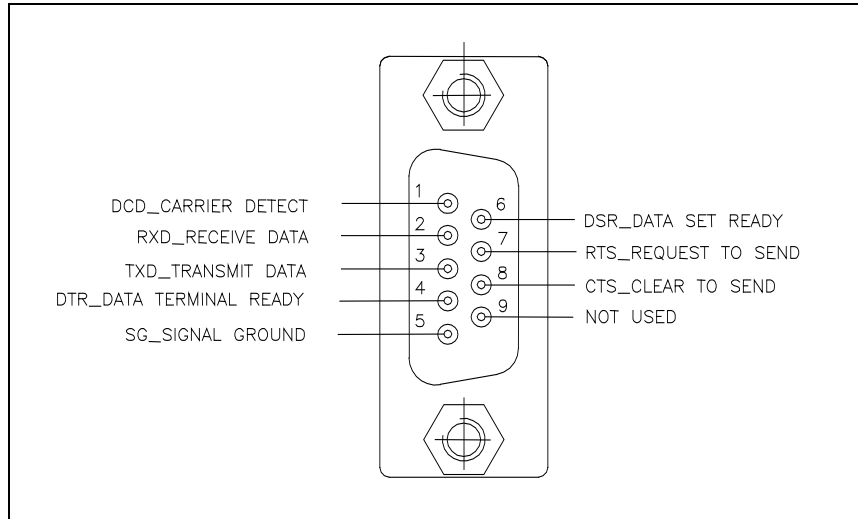


Figure 9 RS232 D-Shell Female 9P

Installation according to international rules.

Tab. 3 D Type 9 Pin FEMALE Connector

D-Type-9 Pin No.	Abbreviation	Full Name
Pin 3	TXD	Transmit Data
Pin 2	RXD	Receive Data
Pin 7	RTS	Request To Send
Pin 8	CTS	Clear To Send
Pin 6	DSR	Data Set Ready
Pin 5	SG	Signal Ground
Pin 1	DCD	Carrier Detect
Pin 4	DTR	Data Terminal Ready
Pin 9	RI	Ring Indicator

PIN Functions

Tab. 4

Abbreviation	Full Name	Function
TXD	Transmit Data	DTE Serial Data Output (TXD), IPCU Serial Data Input
RXD	Receive Data	DTE Serial Data Input (RXD), IPCU Serial Data Output
CTS	Clear to Send	This line indicates that the Modem (Controller) is ready to exchange data.
DCD	Data Carrier Detect	When the Modem (IPCU) detects a "Carrier" from the other end of the line, this line becomes active.
DSR	Data Set Ready	This tells the DTE that the Modem (IPCU) is ready to establish a link.
DTR	Data Terminal Ready	This is the opposite of DSR. This tells the Modem (Controller) that the DTE is ready to link.
RTS	Request To Send	This line informs the Modem (IPCU) is ready to exchange data.
RI	Ring Indicator	Goes active when Modem detects a ringing signal from the PSTN. Not used on the Controller.

RTS / CTS are jumpered together on the Controller connector and DTR is jumpered to DSR / DCD on IPCU connector. The DTE sees the IPCU like a Standard MODEM.

Procedure to Connect the Serial and I/O Ports to an External Cable

The following picture shows the right procedure to connect a cable to the I/O or to the serial port connector.

A shielded cable of 30 m maximum length has to be utilized for both serial and I/O port connections.

NOTE

Take care to have a good contact (soldered) between the metallic connector case and the external shield of the cable. Moreover, this connection has to be assured at least on the controller side.

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

In this way, you will be sure to reduce the influence of the external noise and to accomplish the EMC requests. In picture d is showed the cable assembled.

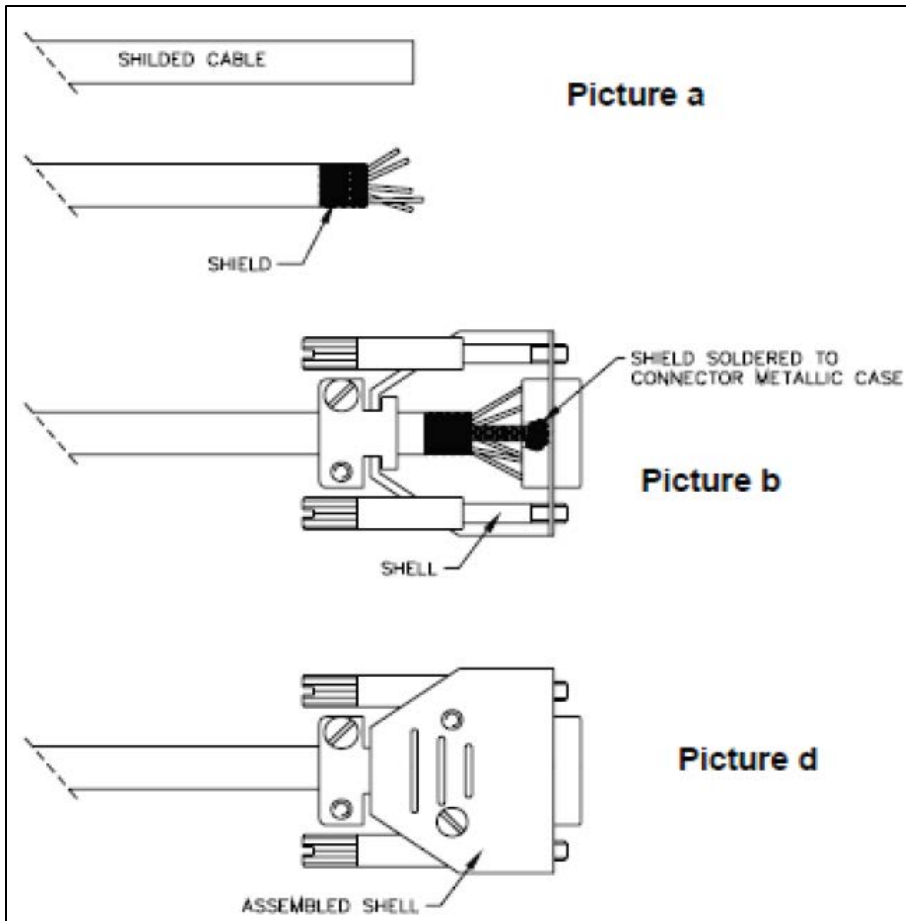


Figure 10

Ion Pump Control Unit - Operator / Host Serial Communication

The Controller Report and Operator/Host Command are ASCII strings CR (0D hexadecimal, 13 decimal) terminated.

Default Communication Parameters are 9600 bits/sec, no parity, 8 bits data, 1 stop bit.

Usually the human interface of the controller is a terminal emulator (Windows "HyperTerminal" communication program for example) running on Personal Computer.

Software program on PC or PLC can use the same ASCII protocol described below for human interface for doing automatic control and monitoring of the controller.

Every 100 (min) to 1000 (max) msec, if the controller is in Report mode (default), all the 2 Report strings made of 30 characters each, corresponding to the 2 HV channels are sent by the IPCU.

The string is made of different length depending on the report type selected for each HV channel (independently by the channel status).

A transmission from the external PC, turns the IPCU to the Command Mode.

To go back to the Report Mode complete the command string with CR (0x0D) or ESC (0x1B).

Every single received character is Echoed back in order to allow a control by the operator.

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

Controller Report Mode

Tab. 5 " HVn status fault start/ protect current voltage"

HV	[HV]	2 characters 'HV'
n	[1, 2, 3]	1 character
' '	[0x20]	1 character SPACE
status	[OFF_, FAULT, ON_]	5 characters
' '	[0x20]	1 character SPACE
fault	[00, ..., FF]	2 characters
' '	[0x20]	1 character SPACE
start/protect	[0, 1] 0=Start(default), 1=Protect	1 character
' '	[0x20]	1 character SPACE
current	[10 ... 99999 nA, 100 ... 60000 uA]	7 characters
' '	[0x20]	1 character SPACE
voltage	[0 ... 5999V]	6 characters 'CR'
' '	[0x0D]	1 character CR

Controller Commands Mode

Tab. 6 HV Channel On/ Off Setting

"A0ns"[Enter] where :	
A0	HV Channel Serial Command for Status Set
N	HV Channel Number 1 or 2
s	0 for OFF, 1 for ON

Tab. 7 HV Channel Start/ Protect Mode Setting

"C0ns"[Enter] where :	
C0	HV Channel Serial Command for Start/Protect Setting
n	HV Channel Number 1 or 2
s	0 for START mode, 1 for PROTECT mode

Tab. 8 Clear Fault (Reset)

"F0n"[Enter] where:	
F0	HV Channel Serial Command for Fault reset
n	HV Channel Number 1 or 2

Procedure to Connect the Serial and I/O Ports to an External Cable

Tab. 9 Bit positional Fault table when HV ON:

Err #	Description	Front Panel LED
0001]	Cable Interlock when HV On	1 blink of 200 ms every 4 s
[0002]	Remote Interlock when HV On	2 blink of 200 ms every 4 s
[0004]	HV Over Voltage when HV On	3 blink of 200 ms every 4 s
[0008]	Bad Current Offset starting HV On	4 blink of 200 ms every 4 s
[0010]	HV Over Power when HV On	5 blink of 200 ms every 4 s
[0020]	Panel Switch Moved when HV On	6 blink of 200 ms every 4 s
[0040]	HV Over Temperature when HV On	7 blink of 200 ms every 4 s
[0080]	HV Over Current – Protect=Enable	8 blink of 200 ms every 4 s
[0100]	HV Under Voltage when HV On	9 blink of 200 ms every 4 s

Tab. 10 Bit positional Event table when HV ON or OFF:

Err #	Description	Front Panel LED
[0001]	Cable Interlock when HV Off	No blink
[0002]	Remote Interlock when HV Off	No blink
[0004]	HV Over Voltage when HV On	No blink
[0008]	Bad Current Offset when HV Off	No blink
[0010]	HV Over Power when HV On	No blink
[0020]	Spare	No blink.
[0040]	Over Temperature	No blink
[0080]	HV Over Current – Protect=Disable	No blink
[0100]	Current Below Threshold when HV On	No blink
[0200]	HV Under Voltage when HV On	No blink

Tab. 11 Max Output Voltage setting

"H0nxxx"[Enter] where:	
H0	HV Channel Serial Command for Voltage Target
n	HV Channel number 1 or 2
xxxx	= Voltage Target (min 3000V, max 5000V in 50 V steps)

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

Tab. 12 Set-Point Current Threshold Setting (Minimum Access Level Required: "0")

"WR[space]Y[space]xxxxx"[Enter]	
WR	Write
Y	14 for channel 1 15 for channel 2
xxxxx = Current Set-Point value in nA (from 0 to 100000000)	

Tab. 13 Set-Point Current Threshold reading

"RD[space]Y"[Enter]	
RD	Read
Y	14 for channel 1 15 for channel 2

Tab. 14 Report Setting

"RT[space]X"	
or	
"RT[space]X[space]Y"[Enter]	
or	
"RT[space]X[space]Y[space]Z"[Enter]	
X	type [0, ..., 5] (see detailed description on "About Report" chapter)
Y	time rate[0, ..., 255] based on 100 msec tick, default is "3". For time rate = 0, to stop continuous reporting mode and start report on request (see Request Report Command)
Z = mode[0, ..., 4]	
where:	
'0' only Remarks, no channel reports	
'1' reports only channel 1	
'2' reports only channel 2	
'4' reports all channels	

Tab. 15 Report Type Read

"RT"[Enter]	
-------------	--

Procedure to Connect the Serial and I/O Ports to an External Cable

Tab. 16 Request Report (in NON Continuous Report Mode)

"RR"[Enter]

Tab. 17 Enable, Dis able Receiving Input ECO

(Access Level 1 Required)
"WR[space]Y[space]X"[Enter]

WR Write

Y 29 (parameter number)

X 0 for No input ECO 1 for input ECO (default value)

Tab. 18 Enable, Disable Remark Line Reporting

(Access Level 1 Required)
"WR[space]Y[space]X"[Enter]

WR Write

Y 30 (parameter number)

X 0 for No remark lines (except the first three)1 for Remark Line (default value)

Tab. 19 Enable, Disable Value in Floating Point Report Mode

(Access Level 1 Required)
"WR[space]Y[space]X"[Enter]

WR Write

Y 31 (parameter number)

X 0 for Standard Report Mode (default value) 1 for Floating Point Report Mode

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

About Report

The continuous report sent by the IPCU protocol communication is structured as follow:

At each main power on start up six remarking lines are sent.

- The first line sent is the firmware release, for instance:
“* REL [20051117 ICPU]”.
- The second line is the notification that the EEPROM's parameter are wrote into RAM:
*** EVT [LOAD_EEPROM_PARAMETERS_INTO_RAM]**
- The third line is the notification that the calibration of the Analog to Digital Convert used for the measurement of output high voltage and current are done:
*** EVT [ADC_CALIBRATION]**
- The fourth line is reporting that the internal I/O BUS are initialized:
*** EVT [IO_BUS_INIT]**
- The fifth line is the notification that a main power on cold system startup reset is happened:
*** EVT [COLD_RESET_SYSTEM_STARTUP]**
- The sixth and last line reports the number “n” of Power On Cold Reset happened up to that moment:
*** POR [n]**

These six lines are sent also if the EEPROM's parameter n. 30 are set to “0” = Disable Remark Lines.

After these six lines, the normal report sequence starts, accordingly to the report type chosen notifying the user about the values of voltage, current and power, faults and events status for each channel.

Procedure to Connect the Serial and I/O Ports to an External Cable

When an ion pump discharge happens, the unit performs a warm system start reset automatically bringing to 0 V the HV to stop the discharge, then it sends again the first five remarking lines with the fifth changed to:

* EVT [WARM_RESET_SYSTEM_RESTART]

to inform that a warm reset system restart is happened and the HV output, that was in the ON state before the discharge, is automatically restarted.

After these five lines, the normal report sequence starts, accordingly to the report type chosen.

Enable – Disable Remark Lines

These five lines are sent also if the EEPROM's parameter n. 30 are set to "0" = Disable Remark Lines.

If the EEPROM's parameter n. 30 is set to "1", all the remark lines, the lines that begin with the symbol "*", are sent.

For EEPROM's parameter n. 30 = "0", all the remark lines are NOT sent, except the first six lines at the main power on reset (cold reset) and the first five lines after an ion pump discharge reset (warm reset) sent anyway.

Continuous and NON Continuous Report Mode

IPCU is able to send report only on request.

To do this, it is necessary to set the report mode with the time rate = 0.

After the "enter", the continuous report mode stops immediately.

According to the report type and mode, the unit sends one instantaneous report each time a "Request Report" (RR) command is received.

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

To return in the continuous report mode it is necessary to send a new “Report Setting” (RT) command with time rate $\neq 0$.

After the “enter” the continuous report mode starts immediately.

Enable - Disable Input ECO Characters

By default, the unit transmits the ECO of all characters received from the host.

If required, It is possible to disable this feature.

See the command “Enable, Disable Receiving Input ECO” to do this.

The ECO of input characters may be used to understand that the physical communication between the unit and host is good so, disabling the ECO this feature is lost.

About Commands

When the unit receives a command, it transmits the plus symbol: “+” to confirm the command is executable.

If the command is not correct, the unit transmits the minus symbol: “-”.

When the parameter of a command is out of range, the unit sends:

“- [PARAMETER_ERROR]”.

The unit receives continuously input characters from host but they are hold only if the panel switch “Local-Remote/Serial” is on Serial.

If this switch is on Local-Remote and a character is received, the unit sends:

Report TYPE 0...6

The User Operator Reports Type 0, 2, 4 and 6 display the ‘continuous average value’ for the HV Channel Current, Voltage and Power, both in normal or scientific notation and the ‘instantaneous measured or control values’ in the hexadecimal format.

Technical Reports Type 1, 3 and 5 report the ‘instantaneous measured values’ for the HV Channel Current, Voltage and Power, both in normal or scientific notation and the ‘continuous average measured values’ in the hexadecimal format.

User Operator Report Type 0 and User Technician Report Type 1

Report Format: “HVC sssss iiiixA vvvvV F=ffff E=eeee <CR>”

Report Fields:

- HVC HV Channel c= 1 or 2
- sssss HV Channel Status [‘OFF__’, ‘ON___’, ‘FAULT’]
- iiiix HV Channel Current in xA
- xA nA or uA (nano Ampere or micro Ampere)
- vvvv HV Channel Voltage in V (Volts)
- ffff Hexadecimal representation of HV Channel Faults
- eeee Hexadecimal representation of HV Channel Events

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

Example:

RT 0[Enter]

+

HV2 OFF 0uA 0V F=0000 E=0000

HV1 OFF 0uA 0V F=0000 E=0000

HV2 OFF 0uA 0V F=0000 E=0000

User Operator Report Type 2 and User Technician Report Type 3

Report Format:

“HVc sssss iiiixA vvvvV F=ffff E=eeee pppppmW i=yyyy v=zzz
c=www <CR>”

Report Fields:

HVc HV Channel c= 1 or 2

“- [LOCAL_MODE]”

to inform that a character is received but it is not possible to execute commands coming from the serial line.

When a command received is formally correct but can no be executed because the status of the unit does not allow this, (for instance a command: “A011” is sent but the cable interlock is not OK) the unit sends:

“- [COMMAND_UNEXECUTABLE]”

to inform that the request is not coherent with the status of the unit.

- sssss HV Channel Status [‘OFF__’, ‘ON___’, ‘FAULT’]
- iiiii HV Channel Current in nA or uA
- vvvv HV Channel Voltage in V

Procedure to Connect the Serial and I/O Ports to an External Cable

- ffff Hexadecimal representation of HV Channel Faults
- eeee Hexadecimal representation of HV Channel Events
- ppppp HV Channel Power in mW
- yyyy 16 Bits hexadecimal Current
- zzz 10 Bits hexadecimal Voltage
- www 12 Bits hexadecimal Control Output

Example:

RT 2[Enter]

+

```

HV2 OFF  0uA  0V F=0000 E=0000  0mW i=0001 v=000 c=D80
HV1 OFF  0uA  0V F=0000 E=0000  0mW i=0003 v=000 c=D80
HV2 OFF  0uA  0V F=0000 E=0000  0mW i=0001 v=000 c=D80
    
```

User Operator Report Type 4 and User Technician Report Type 5

Report Format:

“Hvc sssss iiiixA vvvvV F=ffff E=eeee pppppmW i=yyyy v=zzz
c=www in=hhhh dac=dd <CR>”

Report Fields:

- Hvc HV Channel c= 1 or 2
- sssss HV Channel Status [‘OFF__’, ‘ON___’, ‘FAULT’]
- iiiix HV Channel Current in nA or uA
- vvvv HV Channel Voltage in V
- ffff hexadecimal representation of HV Channel Faults
- eeee hexadecimal representation of HV Channel Events

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

- ppppp HV Channel Power in mW
- yyyy 16 Bits hexadecimal Current
- zzz 10 Bits hexadecimal Voltage
- www 12 Bits hexadecimal Control Output
- hhhh 16 Bits hexadecimal Digital Debounced Inputs
- dd 8 Bits hexadecimal DAC Output

Example:

RT 4[Enter]

+

HV2 OFF 0uA 0V F=0000 E=0000 0mW i=0001 v=000 c=D80
d=7407 r=00

HV1 OFF 0uA 0V F=0000 E=0000 0mW i=0003 v=000 c=D80
d=7407 r=00

HV2 OFF 0uA 0V F=0000 E=0000 0mW i=0001 v=000 c=D80
d=7407 r=00

HV1 OFF 0uA 0V F=0000 E=0000 0mW i=0003 v=000 c=D80
d=7407 r=00

HV2 OFF 0uA 0V F=0000 E=0000 0mW i=0001 v=000 c=D80
d=7407 r=00

Old IPCU Releases - Compatibility Report Type 6

Report Format: "HVc sssss ff p iiiixA vvvvV <CR>"

Report Fields:

HVc HV Channel c= 1 or 2

Floating Point Notation

The floating point notation follows the ANSI/IEEE Standard 754-1985, for Binary Floating Point Arithmetic.

Example:

AL 1 11111111[Enter] (to grant access level 1)

1

wr 31 1[Enter] (set parameter n. 31 to 1)

+

RT 0 (select report type "0")

+

HV2 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV1 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV2 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV1 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV2 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV1 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV2 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV1 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

HV2 OFF 0.0E-9A 0.0E+0V F=0000 E=0000

2 Technical Information

Procedure to Connect the Serial and I/O Ports to an External Cable

Cleaning

For safety reasons, before cleaning the controller:

- 1 Turn the controller off;
- 2 disconnect the controller power plug from the electrical outlet
- 3 disconnect all cables.
- 4 If the exterior of the controller becomes dirty, use a dry soft cloth.

The internal cleaning must be done only by Agilent operator.

- ssss HV Channel Status ['OFF__', 'ON__', 'FAULT']
- ff 8 bits HEX representation of HV Channel Faults
- p '0'=Start, '1'=Protect
- iiii HV Channel Current in xA
- xA nA or uA (nano Ampere or micro Ampere)
- vvvv HV Channel Voltage in V (Volts)

Example:

RT 6

+

HV2 OFF 00 0 0uA 0V

HV1 OFF 00 0 0uA 0V

HV2 OFF 00 0 0uA 0V

HV1 OFF 00 0 0uA 0V

HV2 OFF 00 0 0uA 0V



Vacuum Products Division

Dear Customer,

Thank you for purchasing an Agilent vacuum product. At Agilent Vacuum Products Division we make every effort to ensure that you will be satisfied with the product and/or service you have purchased.

As part of our Continuous Improvement effort, we ask that you report to us any problem you may have had with the purchase or operation of our products. On the back side you find a Corrective Action request form that you may fill out in the first part and return to us.

This form is intended to supplement normal lines of communications and to resolve problems that existing systems are not addressing in an adequate or timely manner.

Upon receipt of your Corrective Action Request we will determine the Root Cause of the problem and take the necessary actions to eliminate it. You will be contacted by one of our employees who will review the problem with you and update you, with the second part of the same form, on our actions.

Your business is very important to us. Please, take the time and let us know how we can improve.

Sincerely,

A handwritten signature in black ink, appearing to read "Giampaolo LEVI".

Giampaolo LEVI

*Vice President and General Manager
Agilent Vacuum Products Division*

Note: Fax or mail the Customer Request for Action (see backside page) to Agilent Vacuum Products Division (Torino) – Quality Assurance or to your nearest Agilent representative for onward transmission to the same address.

Vacuum Products Division Instructions for returning products

Dear Customer,

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

Complete the attached **Request for Return form** and send it to Agilent Technologies (see below), taking particular care to include the completed **Health and Safety** declaration Section. No work can be started on your unit until we receive a completed copy of this form.

After evaluating the information, Agilent Technologies will provide you with a **Return Authorization (RA) number** via email or fax, as requested. Note: Depending on the type of return, a Purchase Order may be required at the time **the Request for Return is submitted**. We will quote any necessary services (evaluation, repair, special cleaning, eg).

Product preparation

- Remove all accessories from the core product (e.g. inlet screens, vent valves).
- Prior to shipment and if applicable for your product, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
- If ordering an Advance Exchange product, please use the packaging from the Advance Exchange to return the defective product.
- Seal the product in a plastic bag, and package product carefully to avoid damage in transit. You are responsible for loss or damage in transit.
- Include a copy of the Health and Safety Declaration in the shipping documentation on the outside of the shipping box of your returning product.
- Clearly label package with RA number. Using the shipping label provided will ensure the proper address and RA number are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will be returned.
- Return only products for which the RA was issued.

Shipping

- Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information. Customer is responsible for freight charges on returning product.
- Return shipments must comply with all applicable Shipping Regulations (IATA, DOT, ADR, etc.) and carrier requirements.

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

EUROPE:

Fax: 00 39 011 9979 330
Fax Free: 00 800 345 345 00
Toll Free: 00 800 234 234 00

NORTH AMERICA:

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

PACIFIC RIM:

Please visit our website for individual office information

vpt-customer@agilent.com

vpl-ra@agilent.com

<http://www.agilent.com>



TERMS AND CONDITIONS

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

- Unless otherwise pre-negotiated, customer is responsible for the freight charges for the returning product. Return shipments must comply with all applicable **Shipping Regulations** (IATA, DOT, etc.) and carrier requirements.
- Agilent Technologies is not responsible for returning customer provided packaging or containers.
- Customers receiving an Advance Exchange product agree to return the defective, rebuildable part to Agilent Technologies **within 15 business days**. Failure to do so, or returning a non-rebuildable part (crashed), will result in an invoice for the non-returned/non-rebuildable part.
- Returns for credit toward the purchase of new or refurbished Products are subject to prior Agilent approval and may incur a restocking fee. Please reference the original purchase order number.
- Units returned for evaluation will be evaluated, and a quote for repair will be issued. If you choose to have the unit repaired, the cost of the evaluation will be deducted from the final repair pricing. A Purchase Order for the final repair price should be issued within 3 weeks of quotation date. Units without a Purchase Order for repair will be returned to the customer, and the evaluation fee will be invoiced.
- Products returned that have not been drained from oil will be disposed.
- A Special Cleaning fee will apply to all exposed products.
- If requesting a calibration service, units must be functionally capable of being calibrated.

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies

121 Hartwell Avenue
Lexington, MA 02421 - USA
Tel.: +1 781 861 7200
Fax: +1 781 860 5437
Toll-Free: +1 800 882 7426
vpl-customerservice@agilent.com

Benelux

Agilent Technologies Netherlands B.V.

Groenelaan 5, 1186 AA Amstelveen
The Netherlands
Tel: +31 20 547 2000
Fax: +31 20 547 2093

Brazil

Agilent Technologies Brasil

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

China

Agilent Technologies (China) Co. Ltd

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
vacuum.cnmarketing@agilent.com
vpc-customerservice@agilent.com

France

Agilent Technologies

Parc Technopolis - Z.A. de Courtaboeuf
3, avenue du Canada - CS 90263
91978 Les Ulis cedex, France
Tel: +33 (0) 1 64 53 61 15
Fax: +33 (0) 1 64 53 50 01
vpf.sales@agilent.com

Germany and Austria

Agilent Technologies

Sales & Services GmbH & Co. KG

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

India

Agilent Technologies India Pvt. Ltd.

Unit Nos 105-116
First Floor, Splendor Forum,
Plot No.-3, District Centre, Jasola
New Delhi-110025
Ph: +91 11 4623 7100
Fax: +91 4623 7105
Toll Free: 18001801517

Italy

Agilent Technologies Italia S.p.A.

Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel: +39 011 9979 111
Fax: +39 011 9979 350
Toll free: 00 800 234 234 00
vpt-customerservice@agilent.com

Japan

Agilent Technologies Japan, Ltd.

8th Floor Sumitomo Shibaura Building
4-16-36 Shibaura Minato-ku
Tokyo 108-0023 - Japan
Tel.: +81 3 5232 1253
Fax: +81 3 5232 1710
Toll-Free: 0120 655 040
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 (0)2 2194 9449
Fax: +82 (0)2 3452 3947
Toll free: 080 222 2452
vpk-customerservice@agilent.com

Mexico

Agilent Technologies

Concepcion Beistegui No 109
Col Del Valle
C.P. 03100 – Mexico, D.F.
Tel.: +52 5 523 9465
Fax: +52 5 523 9472

Singapore

Agilent Technologies Singapore Pte. Ltd.

1 Yishun Avenue 7,
Singapore 768923
Tel : (65) 6215 8045
Fax : (65) 6754 0574
Toll free: 1 800 2762622
vps-customerservice@agilent.com

Southeast Asia

Agilent Technologies Sales Sdn Bhd

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

Taiwan

Agilent Technologies Taiwan Limited

20 Kao-Shuang Road Ping-Chen City
Tao-Yuan Hsien,
32450 Taiwan, R.O.C.
Tel: +886 3 4959004
Toll free: 0800 018 768
vpw-customerservice@agilent.com

UK and 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

Other Countries

Agilent Technologies Italia S.p.A.

Via F.lli Varian, 54
10040 Leini, (Torino) - Italy
Tel.: +39 011 997 9111
Fax: +39 011 997 9350
Toll-Free: 00 800 234 234 00

Customer Support & Service

NORTH AMERICA:

Toll Free: 800 882 7426
vpl-ra@agilent.com

EUROPE:

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

PACIFIC RIM:

please visit our website for individual office
information

<http://www.agilent.com/chem/vacuum>

Worldwide Web Site, Catalog and Order On-line:

www.agilent.com/chem/vacuum

Representatives in most countries

12/13

© Agilent Technologies, Inc. 2014
Printed in ITALY
03/2014
Publication Number: 87-900-128-01 (E)



Agilent Technologies



Agilent Technologies