

**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, unauthorized interference with the equipment or any action contrary to that provided for by specific national standards. The SQ405 is a high voltage feeder used to feed the ionic pumps.

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 section "Technical Information".

This manual uses the following standard protocol:

**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.

**CAUTION**

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

NOTE

The notes contain important information taken from the text.

STORAGE

When transporting and storing the Controller, the following environmental requirements should be satisfied:

- temperature: from -20 °C to + 70 °C
- relative humidity: 0 - 95%
(without condensation)

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, ensure that the module is not dropped or subjected to any form of impact. Do not dispose of the packing materials in an unauthorised manner. The material is 100% recyclable and complies with EEC Directive 85/399.

INSTALLATION**WARNING!**

The controller is equipped with a 3-wire power cord. Use this power cord in conjunction with a properly grounded power socket to avoid electrical shock.

High voltage developed in the controller can cause severe injury or death. Before servicing the unit, disconnect the input power cable.

NOTE

The controller must be installed inside a rack module, but it must be positioned so that free air can flow through the holes. 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.

During operation, the following environmental conditions must be respected:

- temperature: from 0 °C to +45 °C
- relative humidity: 0 - 95%
(without condensation)

To connect the controller to the pump use the specific cable supplied with the controller.

The following paragraph describes the fundamental operating procedures.

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

Power Failure

In the event of a power failure (momentary or long period) the controller is switched off. When power is restored, the controller will automatically restart.



WARNING!

High voltage in the controller can cause severe injury or death. Before servicing, turn power off and remove the power cable.

Technical Specifications	
Mains Supply:	
Voltage	200Vac $\pm$ 10% Single phase
Frequency	50 Hz
Power	800 VA max
Outputs:	
Voltage	7000 V max
Current	100 mA nominal – 200mA short circuit
Power	200 W max
Operating Temperature	From 0 to +40 °C
Storage Temperature	From -20 to+70 °C
Mains fuses	2 x T 6.3A

Connectors Description

The C.U. has connectors for the following functions:

- **Mains connection:** Plug IEC type with integrated fuses for 200V 50Hz supply
- **Connections to the Pump: J1,** Fisher D105A049 type connector

Connector I/O: J2, Delta type 25 pin female:

<i>N.Pin</i>	<i>Description</i>	<i>Type</i>
14	Output +24Vcc	Supply voltage
1	Input HV ON/OFF positive	Optically Insulated Input
2	Input HV ON/OFF negative	
15	GND	Supply voltage
5	FAULT n.o. output	RELAY output
18	FAULT n.o. output	
6	HV ON n.o. output	
19	HV ON n.o. output	
7	PROTECT/START n.o. output	
20	PROTECT/START n.o. output	
8	LOCAL/REMOTE n.o. output	
21	LOCAL/REMOTE n.o. output	
3	CURRENT analog output positive	Logarithmic analog output
16	CURRENT analog output GND	
4	PRESSURE analog output positive	
17	PRESSURE analog output GND	
9	RESERVED	Analog Output

- **Serial Connector: J3,** Delta type 9 pin female for RS422

<i>Pin</i>	<i>Description</i>	<i>Type</i>
6	TX-A (+)	RS 422 output
8	TX-B (-)	
4	RX-A (+)	RS 422 input
7	RX-B (-)	
5	GND	Supply
9	RESERVED	Input

The Front Panel has a 16 x 2 alphanumeric display and a 4 keys keyboard:

<i>Tasto</i>	<i>Description</i>
H.V. ON/OFF	Pump supply ON/OFF
MODE	Working mode configuration
MEASURE	Selection of the displayed measurement
CHANGE	Current parameter value change

Display messages related to the pump working condition

When the C.U. is switched on, the display shows:

				A	U	T	O	T	E	S	T			
							O	K						

Main page: if in Local mode

P	U	M	P		R	E	A	D	Y		F	O	R	
L	O	C	A	L		S	T	A	R	T				

If in Remote mode

P	U	M	P		R	E	A	D	Y		F	O	R	
R	E	M	O	T	E		S	T	A	R	T			

From this page it is possible to Start the pump or to go to the C.U. configuration.

Configuration: In this condition it is possible to change the operating mode.

The configuration is just possible with the HV off. To enter in the Configuration, press the Mode push button

L	O	C	A	L	/	R	E	M	O	T	E			
S	E	L	E	C	T	I	O	N	:		X	X	X	

Where XXX can be LOC, REM or SERIAL.

The corresponding output on the rear panel connector will be: open contact for LOC or SERIAL and closed contact for REM.

The CHANGE push button allows to change the present selection.

To confirm the selection and proceed to the following page, press the MODE push button.

P	R	O	T	E	C	T	/	S	T	A	R	T		
S	E	L	E	C	T	I	O	N	:		X	X	X	X

Where XXXXX can be PROT or START. The selection is done by means of the CHANGE button.

The corresponding output on the I/O rear panel connector will be: open contact for PROT and closed contact for START.

Press the MODE push button again to go to the Serial communication Baude Rate selection.

S	E	R	I	A	L		B	A	U	D		R	A	T	E
			X	X	X	X		B	A	U	D				

XXXX can be 9600, 4800, 1200 or 600 selectable by means of the CHANGE button.

Press the MODE button again to go to the SERIAL address selection page.

S	E	R	I	A	L		A	D	D	R	E	S	S	:	
							X	X							

XX is the address from 1 to 32 that can be modified by means of the CHANGE button.

To pass from the tens to the units press the MODE button.

Press the MODE button again to go back to the main menu.

Pump in Start: In the following pages the operating mode and the analog values are displayed.

The pump is switched on by pressing the HV ON/OFF push button in Local mode or through the HV ON/OFF input in REMOTE mode.

If the HV cable is disconnected from the pump, the Error (Fault) condition "Cable Interlock" will be generated. When the HV cable is connected, the display shows:

H	V		O	N				P	R	O	T	E	C	T
1	.	3	E	-	7		P	a				X	X	X

Where the Protect indication is shown only if the Protect mode has been selected.

The HV ON output on the Remote I/O connector will be: open contact with HV OFF and closed contact with HV ON.

The MEASURE push button allows to change the page on the display.

H	V		O	N				P	R	O	T	E	C	T
1	.	3	E	-	6		A					X	X	X

Error Condition:

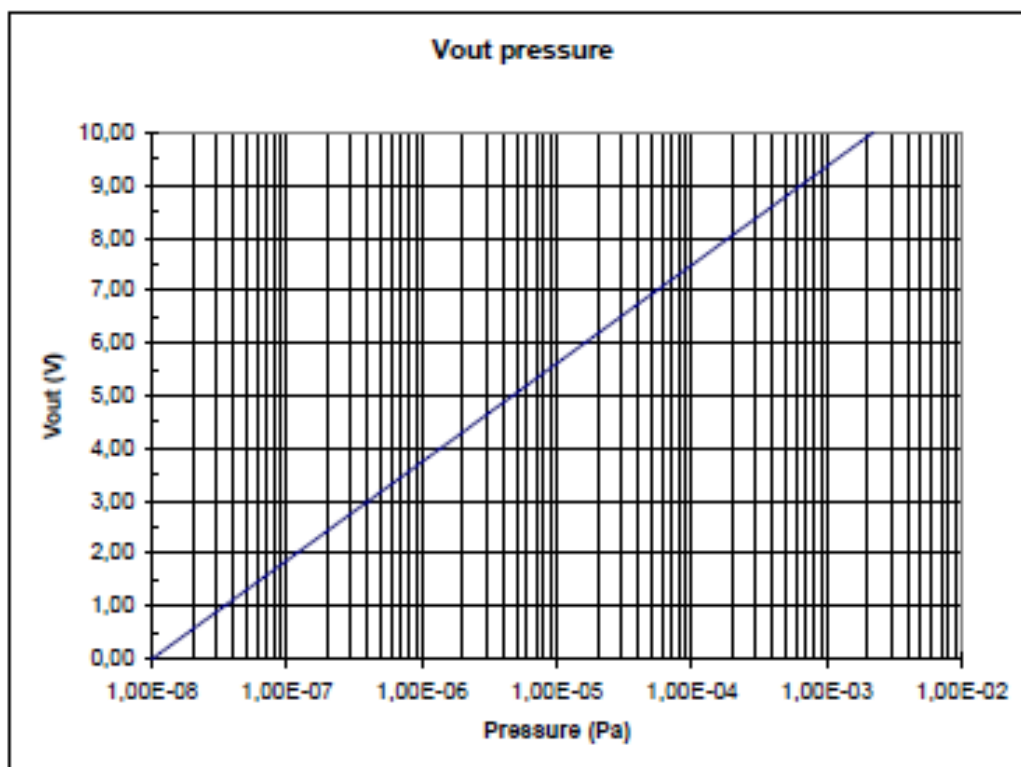
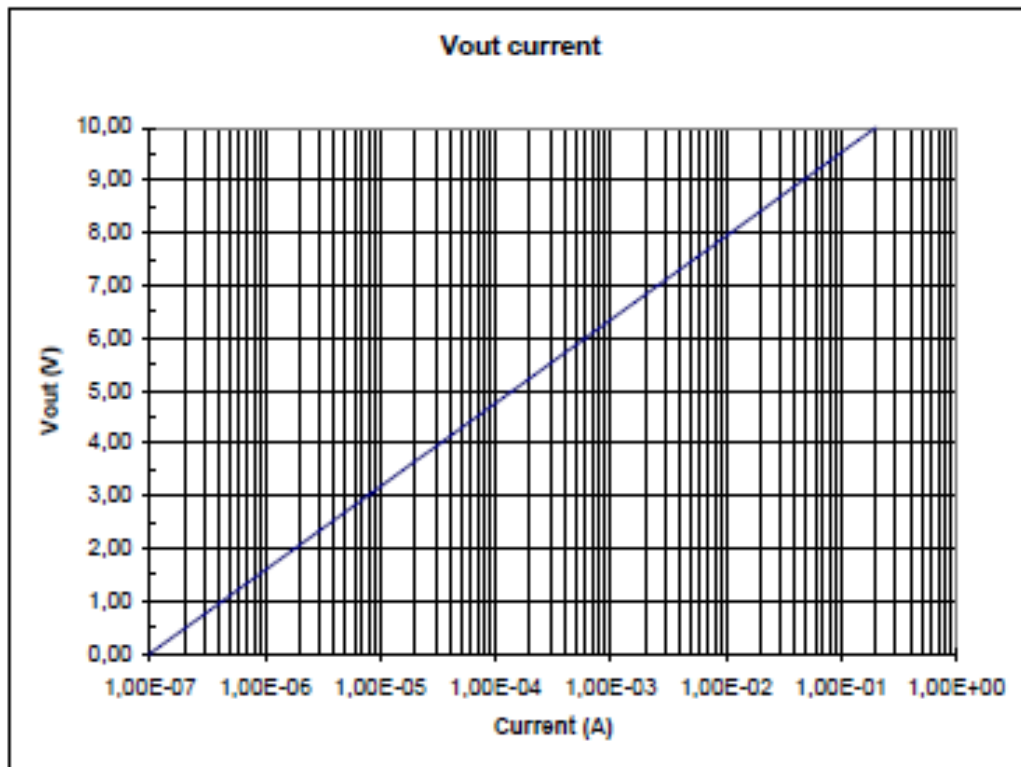
If an Error condition is detected, the FAULT contact on the I/O connector will close .

F	A	U	L	T	:										
I	N	T	E	R	L	O	C	K							

This message is displayed when the HV Cable is disconnected

Analog Outputs

The following graphs give the relationship between the Current and the Pressure versus the Recorder Outputs values.



Serial Communication Description

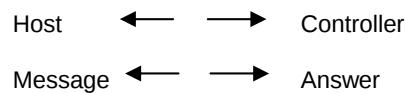
Communication Format

- 8 Data bits
- No Parity
- 1 Stop bit
- Baud Rate: 600/1200/2400/4800/9600 baud, programmable

Communication Protocol

- Master/Slave type (Computer = Host, Controller = Slave)
- Max. number of peripheral units = 32

The communication takes place according to the following scheme:



Message Format

<ADR> + <LDAT> + <COM> + <CHAN> + <DATA> + <CRC>

Where:

- <ADR>: 0x80 + Peripheral unit address (from 0x1 to 0x20)
- <LDAT>: COM, CHAN and DATA fields length (coded in decimal on 2 characters) (00 to 99)
- <COM>: Capital or non capital letter identifying the command (see the "Command Table") followed by the SubCommand (always 0x30 equivalent to the 0n character)
- <CHAN>: Channel number (it is always 0x30, equivalent to 0, for this type of controller)
- <DATA>: String of length and content that depend from the command.

If the command is a request of reading of a data, this field will contain the "?" (0x3f hexadecimal) character.

If the command is a data setting command, or if the message is the answer from the controller to a reading command, this field will contain a data string like shown in the following table.

Data Type	Valid Character
Logical	0 1
Numerical	5 character string completed with "0" characters
Exponential	String with the following format x.xEsxx

- <CRC>: XOR of all the characters of the message, excluding the CRC itself and with the MSB set to 0

The answer structure from the addressed Slave unit, will depend on the received command message:

- No characters if the message has a wrong CRC or a wrong Slave address.
- ACK (0x6 hexadecimal) to confirm the setting of the parameter associated to the command sent by Master, if the command is a writing command
- A message with the same structure as the message described above, but with the <DATA> field that contains the Parameter requested by the Master and with the <ADR> with the MSB set to 0 when the command is a reading command.
- If the message is wrong, the answer will contain the error code identified by the "!" character followed by the "Error Code" coded in decimal on a single character:

- **2: Not existing command**
- **4: Not a reading command**
- **5: Data not valid**
- **6: Out of range value**

Command Table

Following is the command table with the description of the allowed operations and the associated data type:

Command	Channel	Description	Type	Read/Write	Data
L0	0	LOC/REM Configuration	Numerical	R/W	'0':Loc '1':Rem '2':Serial
R0	0	PROTECT/START Configuration	Logical	R/W	'0':Protect '1':Start
A0	0	Address	Numerical	R/W	01 – 32
O0	0	HV On/Off	Logical	R/W	'0':Off '1':On
B0	0	Baud Rate	Numerical	R/W	'0': 600baud, '1': 1200baud, '2': 2400baud, '3': 4800baud, '4': 9600baud
I0	0	Current Measurement	Exponential	R	'x.xE-x'
P0	0	Pressure Measurement	Exponential	R	'x.xE-x'
S0	0	Status	Numerical	R	'0':Stop '1':Start '2':Fault
E0	0	Error Code	Numerical	R	'0':No err '1':Overcurrent '2':Overtemp '3': interlock
f0	0	Crc memory	Numerical	R	Flash memory Crc16

Data type:

- Logical: 1 character data field (0/1)
- Exponential: Numerical data field (x.xEsxx)
- Numerical: Numerical data field (xxxxx)

Serial communication examples

HV ON command by Serial line with controller in Serial mode

0x81 0x30 0x34 0x4F 0x30 0x30 0x31 0x7B

Answer from the control unit:

0x06

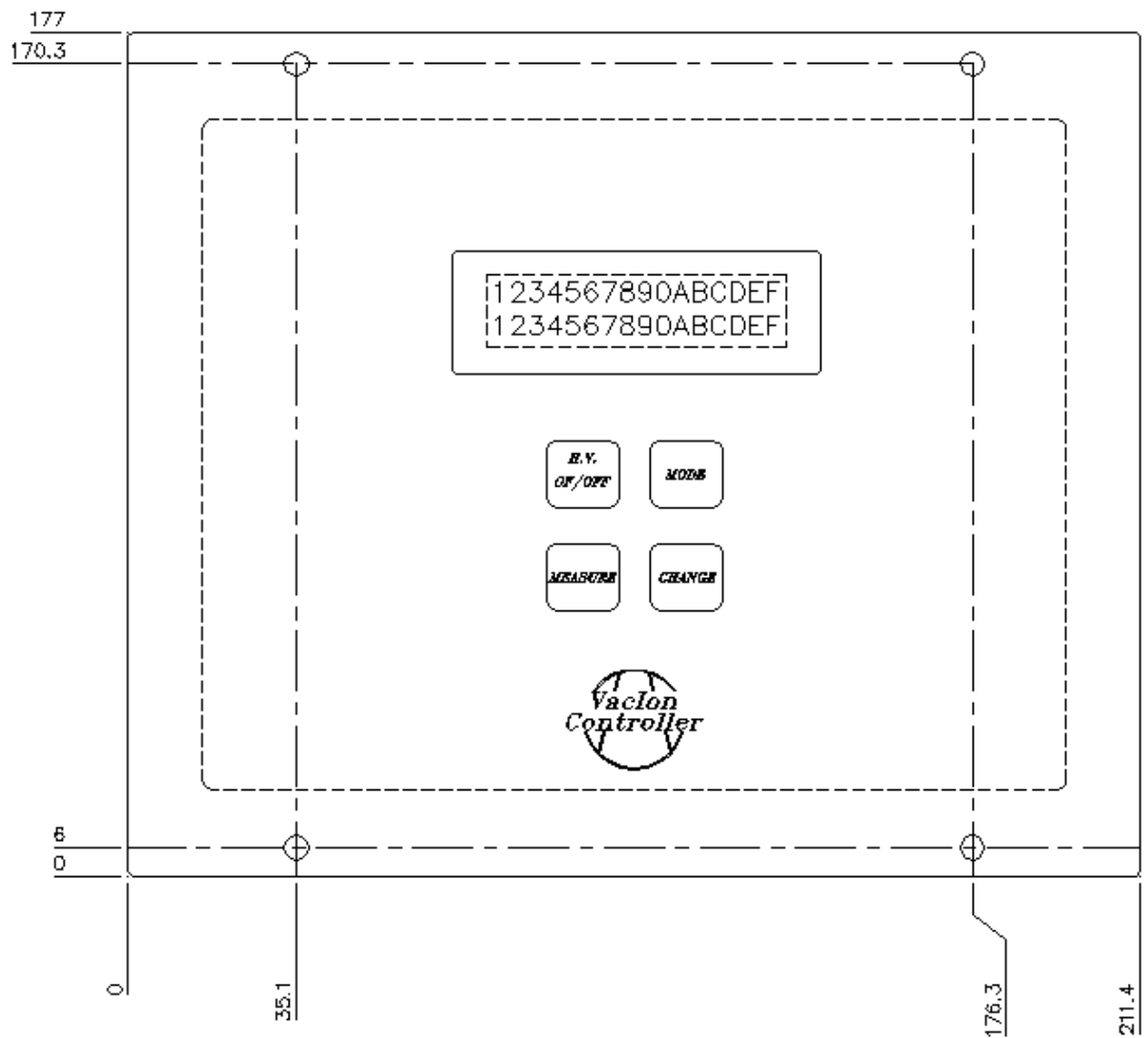
Pressure reading request:

0x81 0x30 0x34 0x50 0x30 0x30 0x3F 0x6A

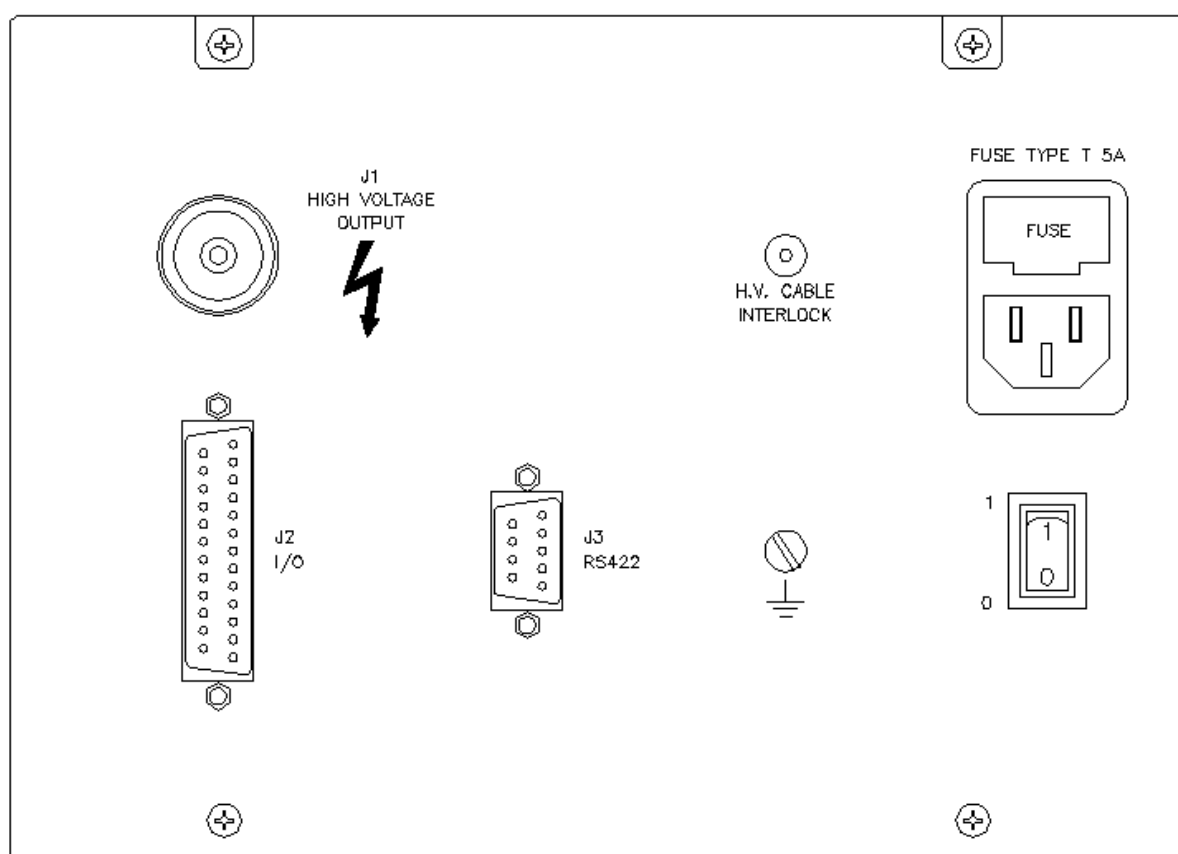
Answer from the control unit:

0x01 0x31 0x30 0x50 0x30 0x30 0x34 0x2E 0x31 0x45 0x2D 0x30 0x35 0x16

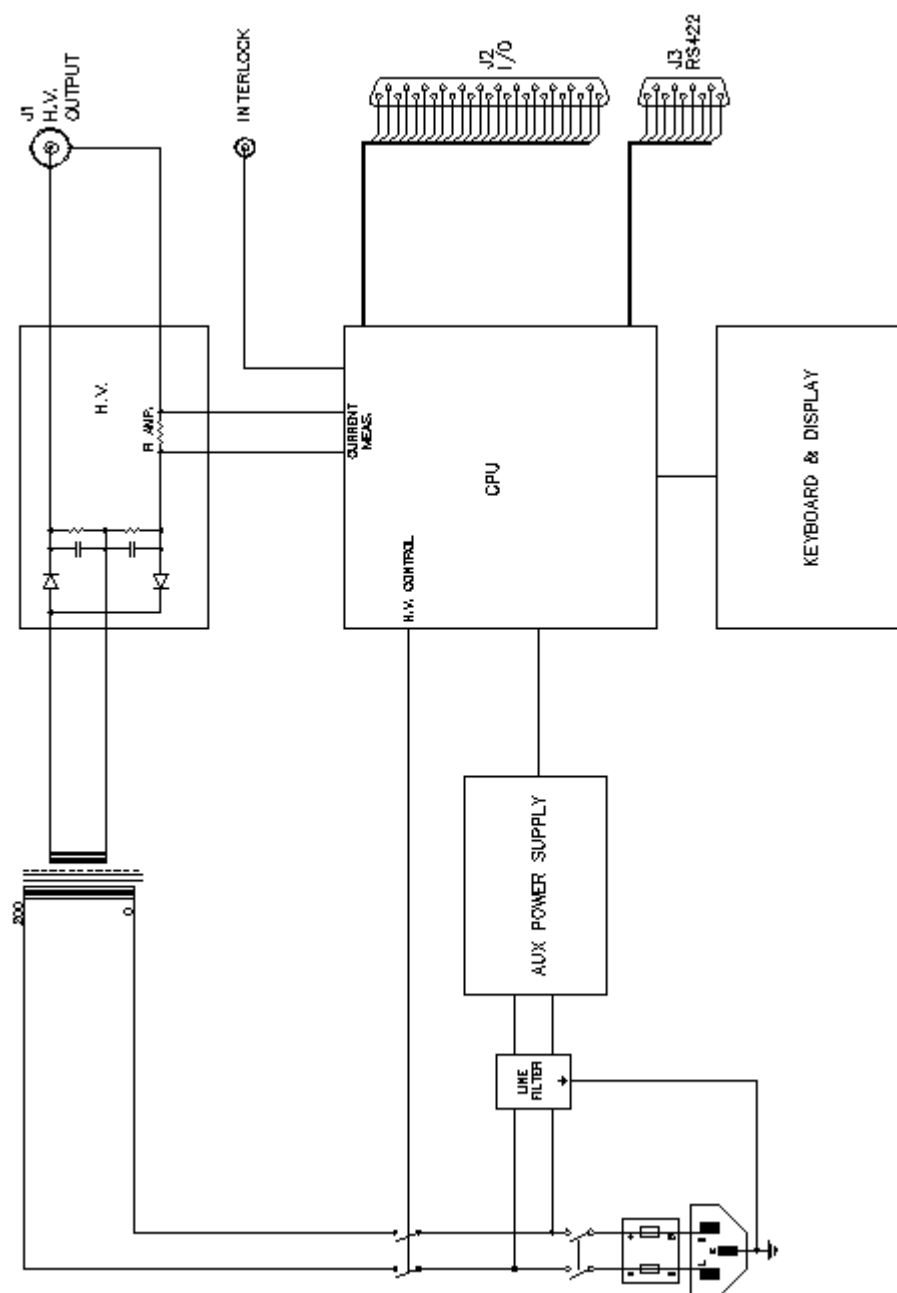
Front Panel



Rear panel



Block Diagram



Interconnections schematic

