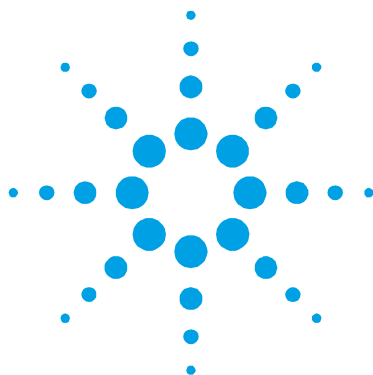




Ion Pumps Control Unit (No Battery)

**Model
829-5000**

**Manuale di Istruzioni
Bedienungshandbuch
Notice de Mode D'Emploi
User Manual**

**87-900-116-01 (E)
04/2014**



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: 87-900-116-01 (E)

Edition

Edition 04/2014

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

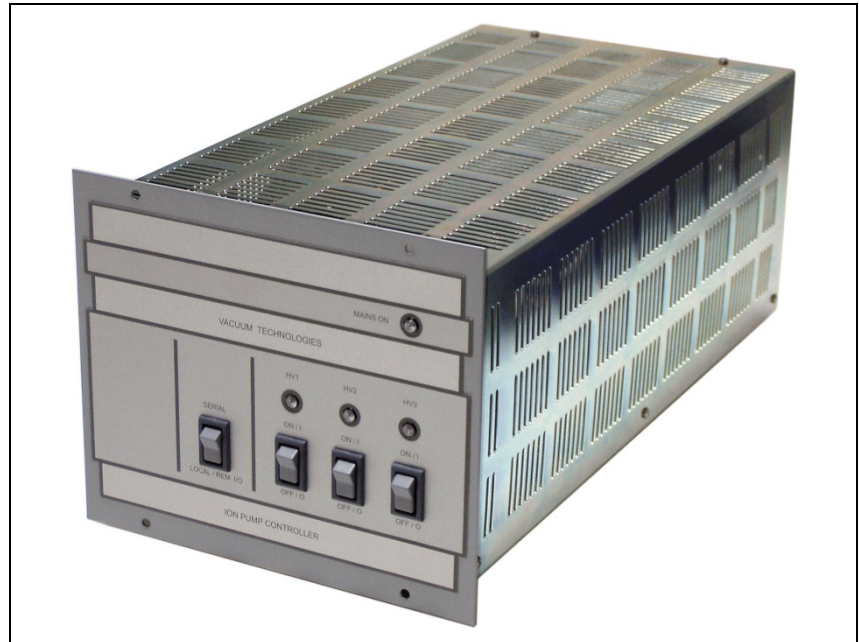
CAUTION

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 (No Battery)



Contents

1 **Procedura per l'installazione** 9

Informazioni Generali	10
Simboli utilizzati	11
Descrizione del controller	12
Immagazzinamento	12
Preparazione per l'installazione	13
Installazione	14
Uso	15
Procedure di Uso	20
Manutenzione	21
Modalità Operative	22
Smaltimento	24
Messaggi di Errore	25

2 **Anleitung zur Installation** 27

Allgemeine Hinweise	28
Verwendete Symbole	29
Beschreibung des Kontrollers	29
Lagerung	30
Installationsvorbereitung	31

Contents

Installation	32	
Benutzung	33	
Gebrauch	37	
Wartung	38	
Betriebsarten	38	
Entsorgung	41	
Fehlermeldungen	42	
3	Procédure pour l'installation	45
Indications Generales	46	
Symboles utilisés	47	
Description du Controleur	48	
Emmagasinage	48	
Préparation pour l'installation	49	
Installation	50	
Utilisation	51	
Procedures d'utilisation	55	
Maintenance	56	
Modalité Operationnelle	57	
Mise au rebut	59	
Messages d'erreur	60	

4 Installation Procedure 63

General Information 64

Symbols used 65

Controller Description 65

Storage 66

Preparation for Installation 67

Installation 68

Use 69

Usage 74

Maintenance 75

Operating Modes 76

Disposal 78

Error Messages 79

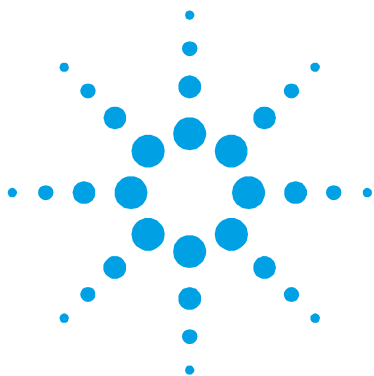
5 Technical Information 81

ION Pump Controller Description 82

Pin Functions 89

Procedure to Connect the Serial and I/O Ports
to an External Cable 90

Contents



1

Procedura per l'installazione

Informazioni Generali	10
Simboli utilizzati	11
Descrizione del controller	112
Immagazzinamento	12
Preparazione per l'installazione	13
Installazione	14
Uso	15
Comandi e indicatori del pannello anteriore dell'unità controller per pompe ioniche	16
Pannello posteriore dell'unità controller per pompe ioniche	18
Procedure di Uso	20
Manutenzione	21
Modalità Operative	22
Collegamento alla Linea di Alimentazione	22
Modalità operativa I/O LOCALE-REMOTA	22
Funzionamento SERIALE	23
Modalità operativa START/PROTECT	23 23
Smaltimento	24
Messaggi di Errore	25
Sul pannello anteriore	25
Sul connettore "Remote" nel pannello posteriore	26

Traduzione delle istruzioni originali



Informazioni Generali

Questa apparecchiatura è destinata ad uso professionale. L'utilizzatore deve leggere attentamente il presente Manuale di Istruzioni e ogni altra informazione aggiuntiva fornita dalla Agilent prima di utilizzare l'apparecchiatura. La Agilent si ritiene sollevata da eventuali responsabilità dovute all'inosservanza totale, o parziale, delle istruzioni, all'uso improprio da parte di personale non addestrato, ad interventi non autorizzati o ad un utilizzo non conforme alle normative nazionali specifiche.

Nei paragrafi seguenti sono riportate tutte le informazioni necessarie per garantire la sicurezza dell'operatore durante l'utilizzo dell'apparecchiatura. Nell'Appendice "Technical Information" sono forniti maggiori dettagli.

Il presente manuale utilizza il seguente protocollo standard:

AVVERTENZA!



I messaggi di avvertenza attirano l'attenzione dell'operatore su una procedura o una pratica specifica che, se non eseguita in modo corretto, potrebbe provocare gravi lesioni personali.

ATTENZIONE!

I messaggi di attenzione sono visualizzati prima di procedure che, se non osservate, potrebbero causare danni all'apparecchiatura.

NOTA

Le note contengono informazioni importanti estrapolate dal testo.

Simboli utilizzati

I simboli qui di seguito riportati vengono utilizzati coerentemente in tutte le immagini:

Simbolo	Descrizione
	On (Alimentazione)
	Off (Alimentazione)
	Attenzione, rischio di scossa elettrica
	Avvertenza, “consultare le istruzioni per l’uso/l’installazione”

Descrizione del controller

Controller della Agilent per pompe ioniche, in grado di azionare fino a 3 pompe simultaneamente e in modo indipendente. Questo controller è progettato per fornire l'alta tensione alle pompe ioniche mediante collegamento all'alimentazione principale (tensione in uscita max. = 5000 Vcc). (Per maggiori dettagli, vedere paragrafo "Technical Information").

Questo controller può essere azionato in modalità operativa I/O remota/locale oppure in modalità seriale attraverso la porta RS232. La selezione della modalità operativa è eseguita per mezzo dell'interruttore

remote-local/serial I/O presente sul pannello anteriore del controller.

Nella modalità I/O remota-locale, i comandi sono eseguiti agendo sul pannello anteriore oppure mediante collegamento al connettore "REMOTE" DB25 "REMOTE" presente nel pannello posteriore.

I dati inerenti le condizioni operative della pompa sono disponibili sulla linea seriale. Nella modalità seriale, i comandi sono eseguiti attraverso il collegamento seriale e tutti i dati relativi alle condizioni operative della pompa sono disponibili sul collegamento seriale.

Immagazzinamento

Durante il trasporto e l'immagazzinamento del controller, non devono essere superate le seguenti condizioni ambientali:

- temperatura: da -20° a +70 °C.
- umidità relativa: 0 – 95 % (non condensante)

Preparazione per l'installazione

Il controller è fornito all'interno di un apposito imballo protettivo. In caso di segni evidenti di danneggiamento accorsi durante il trasporto, contattare l'ufficio vendite locale. In fase di disimballaggio del controller, fare attenzione a non lasciarlo cadere ed evitare di sottoporlo ad urti o vibrazioni.

Trattare il materiale di imballaggio secondo le apposite procedure autorizzate. Tali materiali sono 100 % riciclabili e rispondono alla Direttiva EEC 85/399.

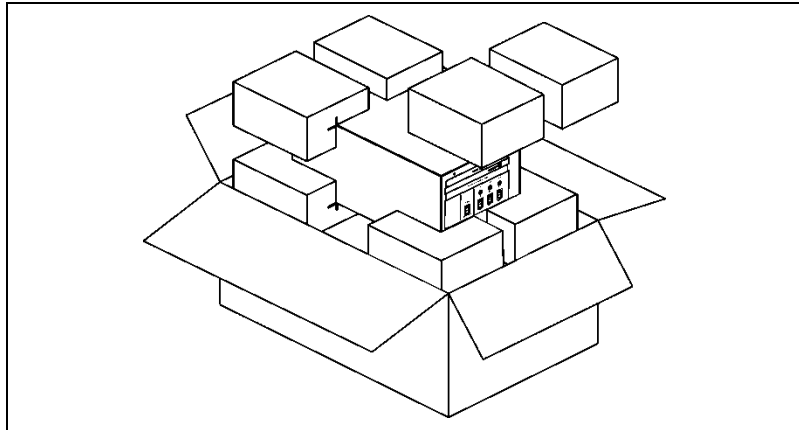


Figura 1 Disimballaggio

Installazione

AVVERTENZA!



Il controller è dotato di un cavo di alimentazione a 3 fili senza spina. Applicare una spina appropriata a seconda del paese di utilizzo del controller.

Si noti che il filo giallo/verde del cavo è il collegamento a massa del controller. Utilizzare sempre questo cavo di alimentazione ed inserire la spina in una presa con adeguato collegamento a massa onde evitare scariche elettriche. Verificare la correttezza dei collegamenti a massa. L'alta tensione generata nel controller può causare gravi danni o addirittura il decesso. Prima di eseguire qualsiasi intervento di manutenzione sull'unità, scollegare il cavo di alimentazione.

ATTENZIONE!

Il controller può essere utilizzato come unità da banco o modulo rack; in ogni caso deve essere posizionato in modo tale da consentire la circolazione d'aria attraverso le apposite aperture presenti sulla copertura superiore e laterale.

Se il controller è utilizzato come modulo rack, DEVE essere inserito in un adattatore per rack a 4 scomparti per evitare che il controller cada all'interno del rack. Il pannello anteriore del controller non è progettato per supportare il peso dell'unità. Non installare o utilizzare il controller in ambienti esposti ad agenti atmosferici (pioggia, gelo, neve), polveri, gas aggressivi, in ambienti esplosivi o ad elevato rischio di incendio.

NOTA

Se il controller è installato su un rack, rimuovere i quattro piedini e posizionarlo lasciando uno spazio minimo di 30 mm (1.2 pollici) in corrispondenza della superficie superiore e inferiore.

AVVERTENZA!



Il controller deve essere installato in modo tale da poter interrompere senza difficoltà la tensione (scollegamento della spina o interruzione mediante interruttore della linea generale del rack).

Durante il funzionamento dell'apparecchiatura è necessario che siano rispettate le seguenti condizioni ambientali:

- temperatura: da +5 °C a +40 °C
- umidità relativa: 0 – 95 % (non condensante)

Per collegare il controller alla pompa, utilizzare gli appositi cavi forniti.

Per le informazioni dettagliate circa i collegamenti sopra indicati o di altra natura, fare riferimento all'Appendice "Technical Information".

Uso

In questo paragrafo sono riportate le principali procedure operative. Per ulteriori dettagli e per le informazioni inerenti i collegamenti opzionali, fare riferimento al paragrafo "USO" dell'Appendice "Technical Information".

Prima di azionare il controller, eseguire tutti i collegamenti elettrici e verso la pompa e leggere il Manuale di Istruzioni della pompa collegata.

AVVERTENZA!



Per evitare danni personali o all'apparecchiatura, se la pompa è collocata su un tavolo, accertarsi che quest'ultimo sia stabile. Non azionare la pompa qualora la relativa presa sia scollegata dal sistema oppure sia ostruita.

Comandi e indicatori del pannello anteriore dell'unità controller per pompe ioniche

La figura e la tabella seguenti riportano i comandi e gli indicatori del pannello anteriore del controller, con le relative descrizioni e funzionalità.

NOTA

La funzionalità e il significato di alcuni comandi e indicatori cambia a seconda della posizione dell'interruttore HV ON OFF. La seguente tabella riporta tutte le funzionalità e tutti i significati dei segnali.

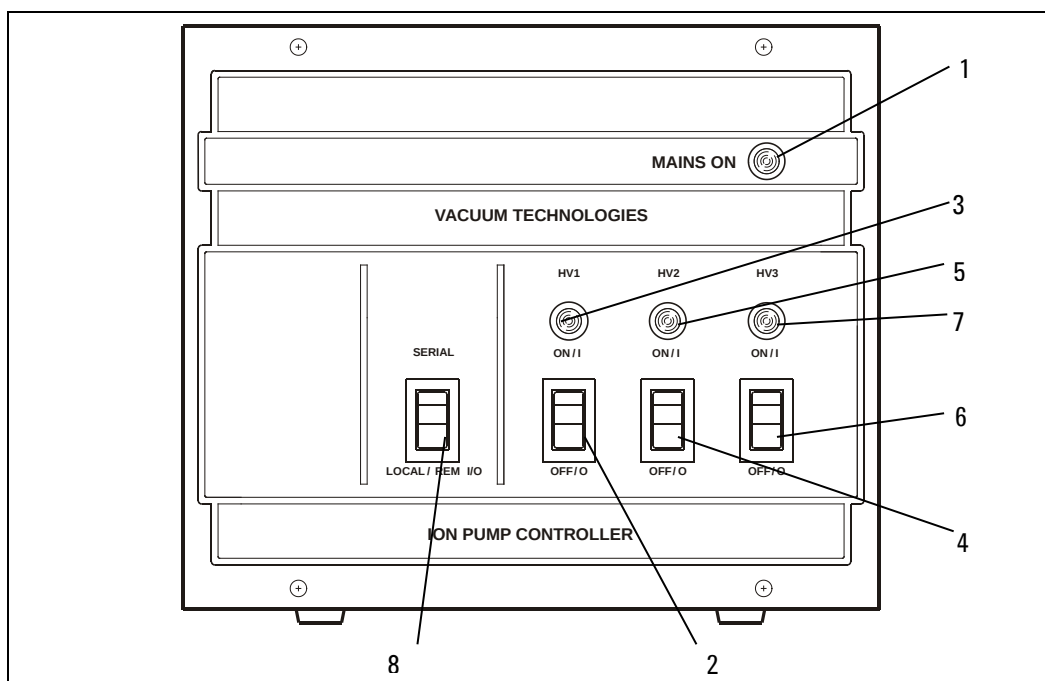


Figura 2 Comandi e indicatori del pannello anteriore dell'unità controller per pompe ioniche

Tab. 1

RIF.	SEGNALE	DESCRIZIONE
1	MAINS ON LED	Quando è On, significa che l'unità è collegata alla linea di alimentazione e che l'interruttore Line sul pannello posteriore è in posizione ON (I).
2	HV1 ON/OFF SWITCH	Quando è in posizione ON (I), abilita l'alimentazione ad alta tensione per la pompa 1 se il controller è impostato in modalità operativa I/O remota/locale e gli interlock del cavo esterno e del cavo alta tensione sono chiusi.
3	HV1 LED	È On quando il controller fornisce l'alta tensione per la pompa 1. Lampeggia in caso di anomalia (vedere paragrafo "Informazioni tecniche" per maggiori dettagli).
4	HV2 ON/OFF SWITCH	Quando è in posizione ON (I), abilita l'alimentazione ad alta tensione per la pompa 2 se il controller è impostato in modalità operativa I/O remota/locale e gli interlock del cavo esterno e del cavo alta tensione sono chiusi.
5	HV2 LED	È On quando il controller fornisce l'alta tensione per la pompa 2. Lampeggia in caso di anomalia (vedere paragrafo "Informazioni tecniche" per maggiori dettagli).
6	HV3 ON/OFF SWITCH	Quando è in posizione ON (I), abilita l'alimentazione ad alta tensione per la pompa 3 se il controller è impostato in modalità operativa I/O remota/locale e gli interlock del cavo esterno e del cavo alta tensione sono chiusi.
7	HV3 LED	È On quando il controller fornisce l'alta tensione per la pompa 3. Lampeggia in caso di anomalia (vedere paragrafo "Informazioni tecniche" per maggiori dettagli).
8	SERIAL/LOCAL-REM I/O SWITCH	Consente di selezionare la modalità operativa: SERIALE: I comandi sono forniti tutti attraverso la linea seriale. I/O remota/locale: I comandi sono forniti tutti agendo sul pannello anteriore dell'unità oppure attraverso il connettore REMOTE a 25 pin presente sul pannello posteriore.

Pannello posteriore dell'unità controller per pompe ioniche

La seguente figura illustra i comandi e gli indicatori del pannello posteriore del controller. Per le informazioni dettagliate circa i collegamenti, fare riferimento all'Appendice "Technical Information".

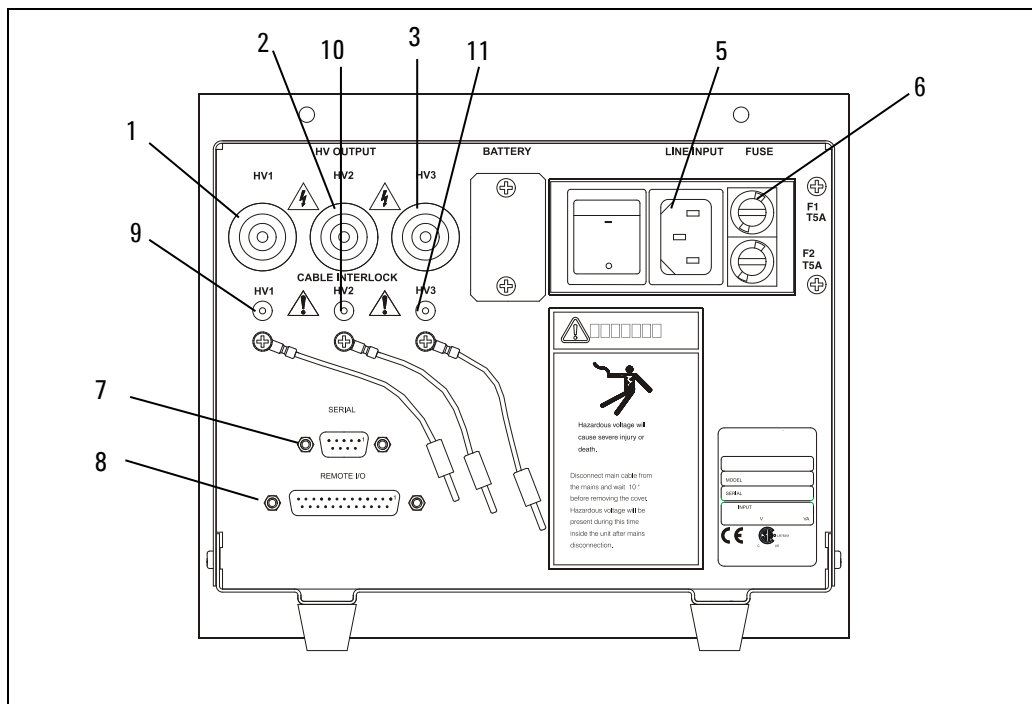


Figura 3 3 Pannello posteriore dell'unità controller per pompe ioniche

Tab. 2

RIF.	SEGNALE	DESCRIZIONE
1	HV1 Connector	Connettore HV per cavo di collegamento alla pompa ionica 1
2	HV2 Connector	Connettore HV per cavo di collegamento alla pompa ionica 2
3	HV3 Connector	Connettore HV per cavo di collegamento alla pompa ionica 3
5	LINE INPUT	Modulo contenente la spina del cavo di ingresso della linea e l'interruttore LINE
6	FUSE	Supporto fusibili (2 fusibili)
7	SERIAL Connector	Connettore DB9 per cavo seriale
8	REMOTE Connector	Connettore DB25 per collegamenti I/O remota
9	HV1 Cable Interlock	
10	HV2 Cable Interlock	
11	HV3 Cable Interlock	

Procedure di Uso

Questo paragrafo fornisce informazioni generiche inerenti l'uso del controller. Per i dettagli, fare riferimento al paragrafo "Technical Information". Collegando il cavo di alimentazione del controller alla linea di alimentazione e posizionando l'interruttore LINE presente nel pannello posteriore su ON (I), l'apparecchiatura è in grado di fornire le seguenti prestazioni:

- 3 canali alta tensione indipendenti
- tensione in uscita max. di -5000 Vcc
- potenza in uscita max. di 40 W per ogni canale
- corrente in uscita max. >30 mA (condizione di cortocircuito) per ogni canale
- modalità operativa seriale/I/O remota-locale
- funzionamento START/PROTECT

Manutenzione

Il Controller non richiede alcuna manutenzione. Qualsiasi intervento sul controller deve essere eseguito da personale autorizzato.

In caso di guasto, rivolgersi al servizio assistenza Agilent. I controller sostitutivi sono disponibili previo accordo mediante Agilent.

AVVERTENZA!

Prima di intervenire sul controller, scollegare l'alimentazione.



Modalità Operative

Collegamento alla Linea di Alimentazione

Modalità operativa I/O LOCALE-REMOTA

NOTA

I 3 interruttori alta tensione presenti sul pannello anteriore e i 3 ingressi On/Off HV sul connettore DB25 "Remote" sul pannello posteriore sono in serie.

L'alimentazione ad alta tensione alle 3 pompe può essere abilitata mediante i 3 interruttori HV (HV1, HV2, HV3) presenti sul pannello anteriore (in questo caso la spina DB25 con i ponticelli interni forniti con l'unità deve essere collegata e i 3 interlock del cavo alta tensione devono essere chiusi) oppure mediante la chiusura dei 3 ingressi On/Off HV sul connettore DB25 "Remote" presente sul pannello posteriore (in questo caso i 3 interruttori HV sul pannello anteriore devono essere in posizione ON).

Quando l'alta tensione è ON, il relativo led sul pannello anteriore si accende.

NOTA

Se il collegamento tra l'IPCU e la pompa è realizzato per mezzo di un cavo ad alta tensione con interlock di sicurezza, il cavo deve essere collegato all'F/T della pompa e il connettore unipolare di dimensioni ridotte deve essere inserito nella relativa spina "HV Cable Interlock" presente sul pannello posteriore dell'unità al fine di poter attivare l'alta tensione.

Se il collegamento è realizzato con un cavo alta tensione unipolare senza interlock di sicurezza, accertarsi che il ponticello fornito sul pannello posteriore dell'unità sia inserito nella spina "HV Cable Interlock".

Funzionamento SERIALE

NOTA

La modalità seriale è impostata al fine di utilizzare l'apparecchiatura collegando il controller ad un PC esterno che utilizza il programma "HYPER TERMINAL" di Windows. Per operare in modalità seriale, collegare la porta seriale al connettore "Serial" del controller sul pannello posteriore, quindi eseguire il programma "HYPER TERMINAL" con le seguenti impostazioni:

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

Questa modalità operativa consentirà di verificare la funzionalità dell'unità, lanciare i comandi e accedere a tutti i dati operativi.

Per maggiori dettagli, fare riferimento al paragrafo "Technical Information".

Modalità operativa START/PROTECT

Il controller può funzionare nelle due seguenti modalità: Start e Protect.

In modalità Start, i 3 canali in uscita possono fornire tutta la potenza in uscita a prescindere dalle condizioni della pompa, fino allo stato di cortocircuito.

Questa modalità viene utilizzata per azionare la pompa a pressione elevata.

In modalità Protect, il controller limita la corrente in uscita e disattiva l'alta tensione quando la corrente supera il valore di soglia (8 mA) per oltre 2 secondi.

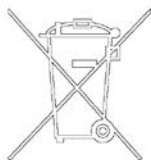
Smaltimento

Significato del logo "WEEE" presente sulle etichette.

Il simbolo qui sotto riportato applicato in ottemperanza alla direttiva CE denominata "WEEE".

Questo simbolo (**valido solo per i paesi della Comunità Europea**) indica che il prodotto sul quale è applicato, NON deve essere smaltito insieme ai comuni rifiuti domestici o industriali, ma deve essere avviato ad un sistema di raccolta differenziata.

Si invita pertanto l'utente finale a contattare il fornitore del dispositivo, sia esso la casa madre o un rivenditore, per avviare il processo di raccolta e smaltimento, dopo opportuna verifica dei termini e condizioni contrattuali di vendita.



Messaggi di Errore

Sul pannello anteriore

Durante il funzionamento del controller, se è rilevato un errore, i LED alta tensione sul pannello anteriore forniranno le seguenti indicazioni:

HV 1,2,3

LED ON: Funzionamento normale = HV ON

Led HV 1,2,3 lampeggianti: condizione di errore:

Lampeggio continuo = sovracorrente in Protect ($I > 8$ mA)

1 lampeggio ogni 2 sec = Interlock cavo HV

2 lampeggi ogni 2 sec = remote Interlock

3 lampeggi ogni 2 sec = HV Overoltage (Warning: non spegne l'HV)

4 lampeggi ogni 2 sec = errore interno ADC

5 lampeggi ogni 2 sec = sovracorrente sull'alta tensione (corto circuito) ($I > 30$ mA per oltre 30 sec)

6 lampeggi ogni 2 sec = Lo switch Local/Remote è stato spostato con HV ON

7 lampeggi ogni 2 sec = sovratemperatura interna all'unità

In modalità I/O remota-locale, la condizione di anomalia è resettata posizionando su Off (0) l'interruttore del canale corrispondente presente sul pannello anteriore.

In modalità seriale, la condizione di anomalia è resettata inviando il comando "FOn" attraverso il computer.

NOTA

Nel caso in cui si verifichi una delle seguenti condizioni mentre l'alta tensione del canale è attiva, l'uscita del canale è impostata immediatamente su Off, il LED sul pannello anteriore indica la condizione di anomalia corrispondente e l'informazione inerente l'anomalia è inviata all'uscita seriale.

Modifica dello stato dell'interruttore Local-Remote I/O
Interruttore HV del pannello anteriore posizionato su Off (0)
Interlock del cavo HV e/o esterno aperto

Sul connettore “Remote” nel pannello posteriore

Uscita “Fault”: contatto N.C. =

Il contatto è aperto durante il normale funzionamento (relè ON)

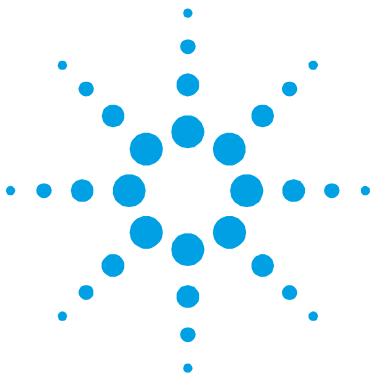
Il contatto è chiuso in condizione di anomalia (relè OFF)

Uscita “HV 1,2,3”: contatto N.O. =

Il contatto è aperto con HV Off (relè Off)

Il contatto è chiuso con HV On (relè On)

Rate dei relè delle uscite: 1 A 48 Vca/cc



2

Anleitung zur Installation

Allgemeine Hinweise	28
Verwendete Symbole	29
Beschreibung des Kontrollers	29
Lagerung	30
Installationsvorbereitung	31
Installation	32
Benutzung	33
Steuerungen und Anzeigen auf dem vorderen Bedienfeld des Ionenpumpen-Kontrollers	34
Rückseite des Ionenpumpen-Kontrollers	36
Gebrauch	37
Wartung	38
Betriebsarten	39
Netzversorgung	39
Betrieb LOCAL-REMOTE I/O	39
SERIELLER Betrieb	40
START/PROTECT Betrieb	40
Entsorgung	41
Fehlermeldungen	42
Auf dem vorderen Bedienfeld	42
Auf dem "Remote" Stecker auf der Rückseite	43

Übersetzung der Originalanleitungen



Allgemeine Hinweise

Dieser Apparat ist für Fachbetriebe bestimmt. Vor Gebrauch sollte der Benutzer dieses Handbuch sowie alle weiteren von Agilent mitgelieferten Zusatzdokumente genau lesen. Bei - auch nur teilweise - Nicht-beachtung der enthaltenen Hinweise, unsachgemäßem Gebrauch durch ungeschultes Personal, nicht autorisierten Eingriffen und Missachtung der einheimischen, hier zur Geltung kommenden Bestimmungen übernimmt die Fa Agilent keinerlei Haftung. In den folgenden Abschnitten sind alle erforderlichen Informationen für die Sicherheit des Bedieners bei Anwendung des Geräts angeführt.

Ins einzelne gehende Informationen über das Gerät finden sich im Anhang " Technical Information ".

In dieser Gebrauchsanleitung werden Sicherheitshinweise folgendermaßen hervorgehoben:

WARNUNG!



Die Warnhinweise lenken die Aufmerksamkeit des Bedieners auf bestimmte Vorgänge oder Praktiken, die bei falscher Ausführung zu schweren Verletzungen führen können.

VORSICHT!

Die Vorsichtshinweise vor bestimmten Prozeduren machen den Bediener darauf aufmerksam, daß bei Nichteinhaltung Schäden am Gerät entstehen können.

HINWEIS

Die Hinweise enthalten wichtige Informationen, die aus dem Text hervorgehoben werden.

Verwendete Symbole

Folgende Symbol werden einheitlich auf allen Darstellungen verwendet:

Symbol	Beschreibung
	Ein (Netzversorgung)
	Aus (Netzversorgung)
	Vorsicht, Stromschlaggefahr.
	Warnung "siehe Installations-/Betriebsanleitung"

Beschreibung des Kontrollers

Der Agilent Controller dient zur Steuerung von Ionen-pumpen. Er kann gleichzeitig bis zu 3 Ionenpumpen unabhängig voneinander bedienen und ist so entworfen, dass er die Ionenpumpen mit HV versorgt, wenn er an das Netz angeschlossen ist (max. Ausgangsspannung = 5000 V=). Weitere Einzelheiten finden sich im Anhang "Technical Information.

Der Agilent Ionenpumpen-Controller kann in den Betriebsarten LOCAL-REMOTE I/O oder im seriellen Modus über eine serielle RS232 Schnittstelle benutzt werden. Die Betriebsart wird mit dem Schalter SERIAL/LOCAL-Remote I/O auf dem vorderen Bedienfeld des Controllers gewählt.

In der Betriebsart LOCAL-REMOTE I/O werden alle Befehle auf dem vorderen Bedienfeld oder durch Anschlüsse an den Stecker DB25 "REMOTE" auf der Rückseite gegeben.

Daten über die Betriebsbedingungen der Pumpen sind in jedem Fall über den seriellen Anschluss verfügbar. Im SERIELLEN Modus kommen die Befehle über die serielle Verbindung, die auch die Daten über die Betriebsbedingungen der Pumpen liefert.

Lagerung

Bei Transport und Lagerung des Controllers müssen die folgenden Umgebungsbedingungen eingehalten werden:

- Temperatur: -20° bis +70 °C
- rel. Luftfeuchtigkeit: 0 – 95 % (nicht kondensierend)

Installationsvorbereitung

Der Controller wird in einer speziellen Schutz-verpackung geliefert. Wenn diese Spuren von Beschädigung zeigt, die bei der Spedition erfolgt sein kann, kontaktieren Sie Ihr lokales Verkaufsbüro. Beim Auspacken des Controllers darauf achten, dass er nicht herunterfällt oder Stößen oder Vibrationen ausgesetzt wird.

Das Verpackungsmaterial umweltfreundlich entsorgen. Das Material kann zu 100 % recycelt werden und entspricht der EU Richtlinie 85/399.

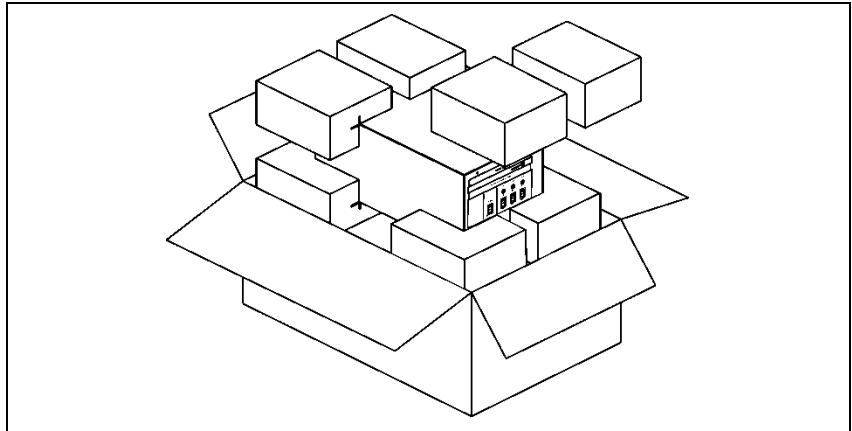


Abbildung 1 Verpackung

Installation

WARNUNG!



Der Controller hat ein dreiadriges Netzkabel ohne Stecker. Bringen Sie den richtigen Stecker für das Land an, in dem der Controller eingesetzt wird.

Der gelbgrüne Leiter des Kabels ist der Erdungsanschluss des Controllers. Benutzen Sie diesen Leiter und den angeschlossenen Stecker zusammen mit einer richtig geerdeten Steckdose, um elektrische Schläge zu verhindern. Prüfen Sie, dass die Erdungsverbindung richtig funktioniert. Die im Controller erzeugte Hochspannung kann zu schweren Verletzungen und zum Tode führen. Vor Wartungsarbeiten am Controller, das Netzkabel abziehen.

VORSICHT!

Der Controller kann als Tisch- oder als Gestellmodul benutzt werden. Er muss aber so positioniert sein, dass Luft frei durch die Löcher in der Seitenwand und der oberen Abdeckung fließen kann. Wenn der Controller in ein Gestell eingebaut ist, MUSS er in ein vier Einheiten hohes Adapterchassis eingebaut werden, damit er nicht in das Gestell kippen kann. Das vordere Bedienfeld ist nicht dafür ausgelegt, das Gewicht der Einheit zu halten. Den Controller nicht in einer Umgebung installieren, die Regen, Schnee, Eis, Staub, aggressiven Gasen usw. ausgesetzt, explosionsgefährdet oder stark feuergefährdet ist.

HINWEIS

Wenn der Controller in einem Gestell installiert wird, die vier Füße entfernen und oben und unten mindestens 30 mm Raum lassen.

WARNUNG!



Der Controller muss so installiert werden, dass er leicht von der Netzspannung getrennt werden kann (Abziehen des Netzsteckers oder Hauptschalter des Gestells).

Beim Betrieb müssen die folgenden Umgebungsbedingungen eingehalten werden:

- Temperatur: +5 °C bis +40 °C
- rel. Luftfeuchtigkeit: 0 – 95 % (nicht kondensierend)

Zum Anschluss des Controllers an die Pumpen die mitgelieferten Kabel benutzen.

Für Einzelheiten über die erwähnten und andere Anschlüsse den Anhang "Technical Information" zu Rate ziehen.

Benutzung

Dieser Abschnitt beschreibt die fundamentalen Betriebsprozeduren. Detaillierte Informationen und Betriebsprozeduren mit optionalen Anschlüssen oder Optionen werden im Abschnitt "GEBRAUCH" im Anhang "Technical Information" behandelt.

Vor Inbetriebnahme des Controllers alle elektrischen Anschlüsse und diejenigen an das Vakuumsammelrohr vornehmen und das Anleitungshandbuch für die angeschlossenen Pumpen zu Rate ziehen.

WARNUNG!



Wenn die Pumpe auf einem Tisch liegt, sicherstellen, dass sie ruhig liegt, um Verletzungen von Personen und Beschädigung des Geräts zu verhindern. Die Pumpe nie betreiben, wenn deren Einlass nicht am System angeschlossen oder verschlossen ist.

Steuerungen und Anzeigen auf dem vorderen Bedienfeld des Ionenpumpen-Kontrollers

Die Abbildung und die Tabelle auf der folgenden Seite zeigen die Steuerungen und Anzeigen auf dem vorderen Bedienfeld des Ionenpumpen-Kontrollers mit Beschreibung und Funktion.

NOTA

Bei einige Steuerungen und Anzeigen hängt die Funktion von der Stellung des Schalters HV ON OFF ab. Die Tabelle enthält all diese Funktionen und ihre Bedeutung.

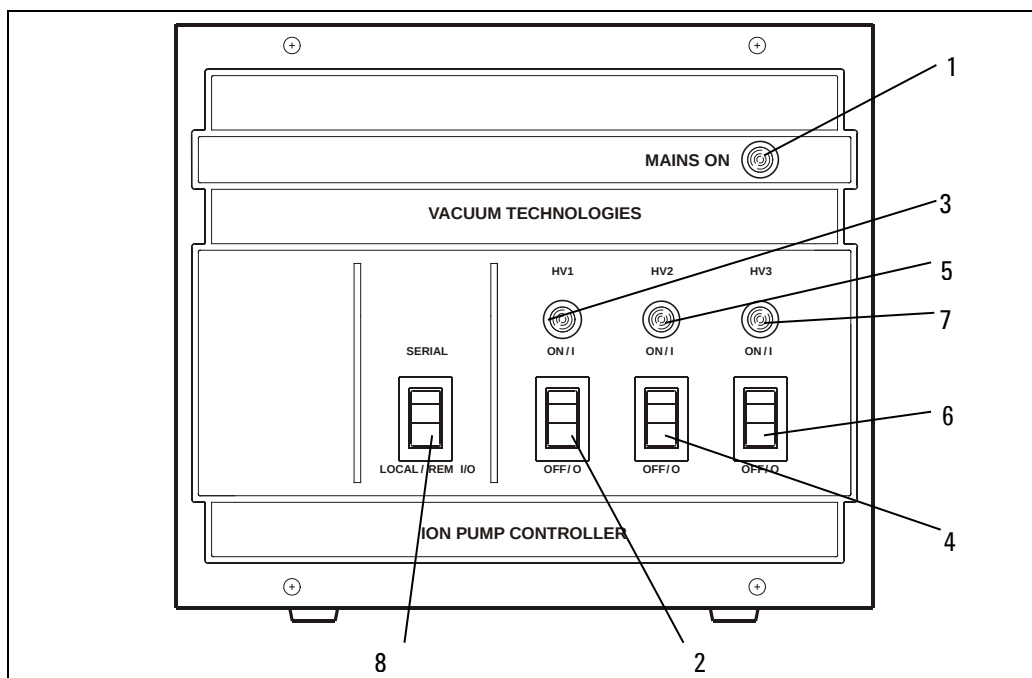


Abbildung 2 Steuerungen und Anzeigen auf dem vorderen Bedienfeld des Ionenpumpen-Kontrollers

Tab. 1

RIF.	SIGNAL	BESCHREIBUNG
1	MAINS ON LED	Wenn an, ist die Einheit ans Netz angeschlossen und der Netzschalter auf der Rückseite steht auf ON (I).
2	HV1 ON/OFF SWITCH	Wenn auf ON (I), wird die HV für die Pumpe 1 geschaltet, sofern der Controller auf LOCAL/REMOTE I/O Betrieb gesetzt ist und die externen wie die HV Kabel-Verriegelungen geschlossen sind.
3	HV1 LED	Wenn an, versorgt der Controller die Pumpe 1 mit HV. Wenn blinkend, liegt ein Fehler vor (Einzelheiten finden sich im Anhang "Technische Informationen").
4	HV2 ON/OFF SWITCH	Wenn auf ON (I), wird die HV für die Pumpe 2 geschaltet, sofern der Controller auf LOCAL/REMOTE I/O Betrieb gesetzt ist und die externen wie die HV Kabel-Verriegelungen geschlossen sind.
5	HV2 LED	Wenn an, versorgt der Controller die Pumpe 2 mit HV. Wenn blinkend, liegt ein Fehler vor (Einzelheiten finden sich im Anhang "Technische Informationen").
6	HV3 ON/OFF SWITCH	Wenn auf ON (I), wird die HV für die Pumpe 3 geschaltet, sofern der Controller auf LOCAL/REMOTE I/O Betrieb gesetzt ist und die externen wie die HV Kabel-Verriegelungen geschlossen sind.
7	HV3 LED	Wenn an, versorgt der Controller die Pumpe 3 mit HV. Wenn blinkend, liegt ein Fehler vor (Einzelheiten finden sich im Anhang "Technische Informationen").
8	SERIAL/LOCAL-REM I/O SWITCH	Wählt die Betriebsart: SERIAL: alle Befehle kommen über die serielle Leitung. LOCAL-REM I/O: alle Befehle werden über das vordere Bedienfeld der Einheit oder über den 25-stiftigen Stecker REMOTE auf der Rückseite gegeben.

Rückseite des Ionenpumpen-Kontrollers

Die folgende Abbildung zeigt die Steuerungen und Stecker auf der Rückseite des Ionenpumpen-Kontrollers. Einzelheiten über die Anschlüsse finden sich im Anhang "Technical Information".

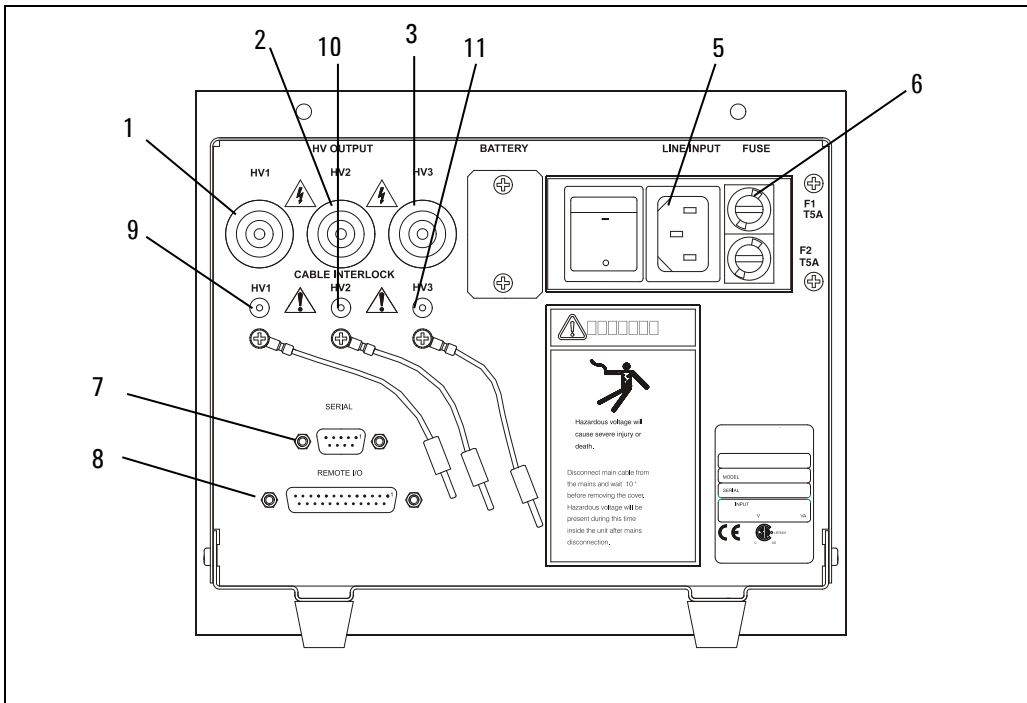


Abbildung 3 Rückseite des Ionenpumpen-Kontrollers

Tab. 2

RIF.	SIGNAL	BESCHREIBUNG
1	HV1 Connector	HV-Stecker für das Kabel zur Ionenpumpe 1
2	HV2 Connector	HV-Stecker für das Kabel zur Ionenpumpe 2
3	HV3 Connector	HV-Stecker für das Kabel zur Ionenpumpe 3
5	LINE INPUT	Modul mit dem Stecker für das Netzkabel und dem Netzschalter
6	FUSE	Halter für die Schmelzsicherungen (2 Sicherungen)
7	SERIAL Connector	DB9 Stecker für das serielle Kabel
8	REMOTE Connector	DB25 Stecker für die REMOTE I/O Anschlüsse
9	HV1 Cable Interlock	
10	HV2 Cable Interlock	
11	HV3 Cable Interlock	

Gebrauch

Dieser Abschnitt enthält allgemeine Informationen über den Gebrauch des Controllers. Einzelheiten finden sich im Anhang “Technische Informationen”. Wenn der Controller ans Netz angeschlossen ist und der Netzschalter auf der Rückseite auf ON (I) steht, können alle seine Funktionen genutzt werden:

- 3 unabhängige HV Kanäle
- max. -5000 V= Ausgangsspannung
- max. 40 W Ausgangsleistung für jeden Kanal
- max. >30 mA Ausgangsstrom (Kurzschlussbedingung) für jeden Kanal
- Betriebsarten LOCAL-REMOTE I/O/SERIAL
- START/PROTECT Betrieb

Wartung

Der Controller braucht nicht gewartet zu werden. Jede Maßnahme am Controller muss durch genehmigtes Fachpersonal erfolgen.

Beim Auftreten eines Fehlers kann der Agilent Reparaturdienst in Anspruch genommen werden. Ersatz-Controller stehen über ein Agilent Austauschsystem im Vorfeld zur Verfügung.

WARNUNG!

Vor jeglichen Arbeiten am Controller die Netzversorgung ausschalten.



Betriebsarten

Netzversorgung

Betrieb LOCAL-REMOTE I/O

HINWEIS

Die 3 HV Schalter auf dem vorderen Bedienfeld und die 3 HV On/Off Eingänge auf dem Stecker DB25 "Remote" auf der Rückseite sind in Serie.

Die Hochspannung für die 3 Pumpen können mit den 3 HV Schaltern (HV1, HV2, HV3) auf dem vorderen Bedienfeld eingeschaltet werden (in diesem Fall müssen der DB25 Stecker mit den internen Brücken, die mit der Einheit geliefert werden, überbrückt und die 3 HV-Kabelverriegelungen geschlossen sein) oder bei Schließen der 3 HV Ein/Auseingänge auf dem Stecker DB25 "Remote" auf der Rückseite (in diesem Fall müssen die 3 HV-Schalter auf dem vorderen Bedienfeld in der AN Position sein).

Wenn die Hochspannung AN ist, ist die entsprechende HV-LED auf dem vorderen Bedienfeld erleuchtet.

HINWEIS

Wenn die Verbindung zwischen dem Controller und der Pumpe mit einem HV-Kabel mit Sicherungs-verriegelung realisiert ist, muss das Kabel mit dem F/T der Pumpe verbunden und der kleine Bananenstecker auf der Controllerseite des Kabels in die entsprechenden HV-Kabel-Verriegelungsbuchse auf der Rückseite der Einheit eingesteckt sein, um die Hochspannung einschalten zu können.

Erfolgt der Anschluss mit einem einadrigen HV-Kabel ohne Sicherheitsverriegelung, sicherstellen, dass auf der Rückseite der Einheit die vorgesehene Brücke in die Buchse für HV-Kabelverriegelung eingesteckt ist.

SERIELLER Betrieb

HINWEIS

SERIAL ist eingestellt, um den Controller von einem externen PC aus zu steuern, auf dem das Windowsprogramm "HYPER TERMINAL" läuft. Um im seriellen Modus zu arbeiten, die serielle Schnittstelle mit dem seriellen Stecker auf der Rückseite verbinden und das Windowsprogramm "HYPER TERMINAL" auf dem PC mit den folgenden Einstellungen starten:

- COM1
- 9600 Baud
- keine Parität
- 8 Datenbits
- 1 Stopbit

Bei dieser Betriebsart können die Funktion der Einheit im seriellen Betrieb getestet, die Befehle gegeben und alle Betriebsdaten empfangen werden.

Weitere Einzelheiten finden sich im Anhang "Technical Information".

START/PROTECT Betrieb

Der Controller kann in zwei Betriebsarten arbeiten: Start und Protect.

In Start können die 3 Ausgangskanäle die ganze Ausgangsleistung unabhängig von dem Pumpen-zustand bis zur Kurzschlussleistung liefern. Diese Betriebsart muss zum Starten der Pumpen bei hohem Druck benutzt werden.

In Protect begrenzt der Controller den Ausgangsstrom und schaltet die Hochspannung ab, wenn der Strom den Schwellwert (8 mA) für mehr als 2 Sek. übersteigt.

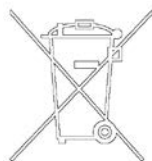
Entsorgung

Bedeutung des "WEEE" Logos auf den Etiketten.

Das folgende Symbol ist in Übereinstimmung mit der EU-Richtlinie WEEE (Waste Electrical and Electronic Equipment) angebracht.

Dieses Symbol (**nur in den EU-Ländern gültig**) zeigt an, dass das betreffende Produkt nicht zusammen mit Haushaltsmüll entsorgt werden darf sondern einem speziellen Sammelsystem zugeführt werden muss.

Der Endabnehmer sollte daher den Lieferanten des Geräts - d.h. die Muttergesellschaft oder den Wiederverkäufer - kontaktieren, um den Entsorgungsprozess zu starten, nachdem er die Verkaufsbedingungen geprüft hat.



Fehlermeldungen

Auf dem vorderen Bedienfeld

Wenn während des Controllerbetriebs ein Fehler entdeckt wird, wird dieser von den HV-LEDs auf dem vorderen Bedienfeld angezeigt:

HV 1,2,3

LEDS AN: Normalbetrieb = HV AN

LED HV 1,2,3 blinkend: Fehlerbedingung:

Dauernd blinkend = Überstrom in Protect ($I > 8 \text{ mA}$)

1 Blinken alle 2 Sek = Interlock HV-Kabel

2 Blinken alle 2 Sek = remote Interlock

3 Blinken alle 2 Sek = HV Überspannung

(Warnung: HV nicht abschalten)

4 Blinken alle 2 Sek = interner ADC Fehler

5 Blinken alle 2 Sek = Überstrom der Hochspannung

(Kurzschluss)

($I > 30 \text{ mA}$ für länger als 30 Sek)

6 Blinken alle 2 Sek = der Local/Remote Schalter wurde mit HV
ON umgeschaltet

7 Blinken alle 2 Sek = interne in der Einheit

In der Betriebsart Local-Remote I/O werden Fehlerbedingungen durch Schalten des betreffenden Kanals auf dem vorderen Bedienfeld in OFF (0) zurückgesetzt.

In der seriellen Betriebsart werden Fehlerbedingungen durch Übersenden des Befehls "F0n" vom PC zurückgesetzt.

HINWEIS

Wenn eine der folgenden Bedingungen auftritt, während die Hochspannung des Kanals AN ist, wird der Ausgang dieses Kanals sofort auf AUS gesetzt. Die LED auf dem vorderen Bedienfeld zeigt die Fehlerbedingung an und Fehlerinformationen werden zum seriellen Ausgang gesendet.

Statusänderung des Schalters Local-Remote I/O –Serial.
HV Schalter auf dem vorderen Bedienfeld in Off (0) gesetzt.
Externe und/oder HV-Kabelverriegelung offen

Auf dem “Remote” Stecker auf der Rückseite

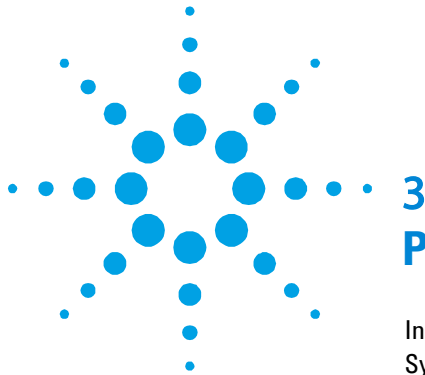
Fehlerausgang: Öffnerkontakt =
der Kontakt ist offen bei Normalbetrieb (Relais AN)
der Kontakt ist geschlossen bei Fehlerbedingung (Relais aus)

“HV 1,2,3” Ausgangs-Schließerkontakt =
der Kontakt ist offen bei HV AUS (Relais aus)
der Kontakt ist geschlossen bei HV AN (Relais an)

Ausgangsrelais: 1 A 48 V~/V=

2 Anleitung zur Installation

Fehlermeldungen



3

Procédure pour l'installation

Indications Generales	46
Symboles utilisés	47
Description du Controleur	48
Emmagasinage	48
Préparation pour l'installation	49
Installation	50
Utilisation	51
Commandes et Indicateurs du Panneau Antérieur de l'unité	
Contrôleur pour Pompes Ioniques	52
Panneau postérieur de l'unité contrôleur pour Pompes	
Ioniques	54
Procedures d'utilisation	55
Maintenance	56
Modalite Operationnelle	57
Connexion à la ligne d'alimentation	57
Modalité opérationnelle I/O LOCALE-A DISTANCE	57
Fonctionnement SERIEL	58
Modalité opérationnelle START/PROTECT	58
Mise au rebut	59
Messages d'erreur	60
Sur le panneau antérieur	60
Sur le connecteur "Remote" dans le panneau postérieur	61

Traduction de la mode d'emploi originale



Indications Generales

Cet appareillage a été conçu en vue d'une utilisation professionnelle. Il est conseillé à l'utilisateur de lire attentivement cette notice d'instructions ainsi que toute autre indication supplémentaire fournie par Agilent, avant l'utilisation de l'appareillage. Agilent décline toute responsabilité en cas d'inobservation totale ou partielle des instructions données, d'utilisation incorrecte de la part d'un personnel non formé, d'opérations non autorisées ou d'un emploi contraire aux réglementations nationales spécifiques.

Les paragraphes suivants donnent toutes les indications nécessaires à garantir la sécurité de l'opérateur pendant l'utilisation de l'appareillage. Dans l'Annexe "Technical Information" sont fournis de plus amples détails.

Le présent manuel utilise les signes conventionnels suivants:

AVERTISSEMENT!



Les messages d'avertissement attirent l'attention de l'opérateur sur une procédure ou une manoeuvre spéciale qui, si elle n'est pas effectuée correctement, risque de provoquer de graves lésions personnelles.

ATTENTION!

Les messages d'attention sont affichés avant les procédures dont le non respect pourrait endommager sérieusement l'appareillage.

NOTE

Les notes contiennent des renseignements importants, extrapolés du texte.

Symboles utilisés

Les symboles qui suivent sont très souvent utilisés dans les illustrations :

Symbole	Description
	Allumé (Tension)
	Éteint (Tension)
	Attention, risque d'électrocution.
	Avertissement « voir notice d'installation/mode de fonctionnement »

Description du Contrôleur

Contrôleur de Agilent pour pompes ioniques en mesure d'actionner jusqu'à 3 pompes simultanément et de manière autonome. Ce contrôleur est conçu pour fournir la haute tension aux pompes ioniques par l'intermédiaire d'une connexion à l'alimentation principale (tension en sortie max. = 5000 Vcc). Pour de plus amples détails, voir le paragraphe "Technical Information".

Ce contrôleur peut être actionné en modalité opérationnelle I/O à distance/locale ou bien en modalité sérieuse à travers le port série RS232. La sélection de la modalité opérationnelle est effectuée au moyen de l'interrupteur remote-local/serial I/O présent sur le panneau antérieur du contrôleur.

En modalité I/O à distance/locale, les commandes sont effectuées en agissant sur le panneau antérieur ou bien au moyen d'une connexion au connecteur "REMOTE" DB25 "REMOTE" présent sur le panneau postérieur.

Les données inhérentes aux conditions opérationnelles de la pompe sont disponibles sur la connexion sérieuse. En modalité sérieuse, les commandes sont effectuées à travers la connexion sérieuse et toutes les données relatives aux conditions opérationnelles de la pompe sont disponibles sur la connexion sérieuse.

Emmagasinage

Durant le transport et l'emmagasinage du contrôleur, il faudra veiller à respecter les conditions environnementales suivantes:

- température: de -20 °C à +70 °C.
- humidité relative: 0 – 95 % (non condensante)

Préparation pour l'installation

Le contrôleur est fourni dans un emballage de protection. En cas de signes évidents d'endommagements qui ont eu lieu durant le transport, contacter le bureau de vente local. Pendant la phase d'ouverture de l'emballage du contrôleur, veiller à ne pas le laisser tomber et éviter de le soumettre à des chocs ou des vibrations.

Traiter le matériel d'emballage suivant les procédures spéciales autorisées. Ces matériaux sont 100 % recyclables et sont conformes à la Réglementation EEC 85/399.

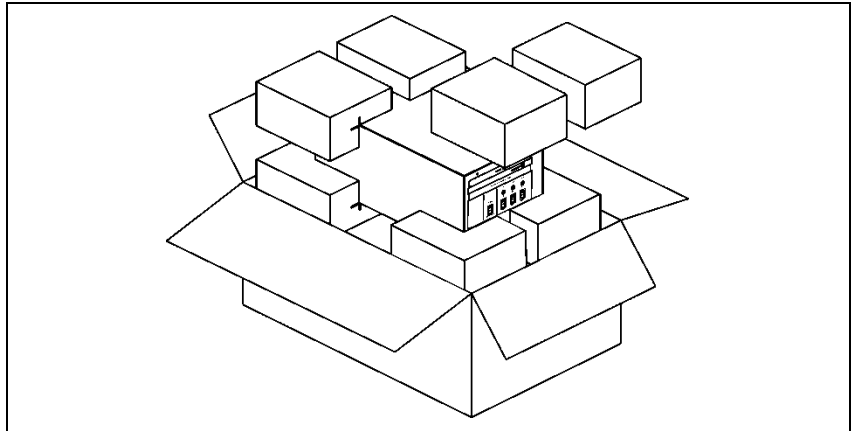


Figure 1 Ouverture de l'emballage

Installation

AVERTISSEMENT!



Le contrôleur est pourvu d'un câble d'alimentation à 3 fils sans fiche. Appliquer une fiche appropriée selon le pays d'utilisation du contrôleur. A remarquer que le fil jaune/vert du câble est la connexion à masse du contrôleur. Utiliser toujours ce câble d'alimentation et introduire la fiche dans une prise avec la connexion appropriée à masse afin d'éviter les décharges électriques. Vérifier le bon usage des connexions à masse. La haute tension générée dans le contrôleur peut provoquer de graves lésions voire le décès. Avant de procéder à toute opération de maintenance sur l'unité, débrancher le câble de l'alimentation.

ATTENTION!

Le contrôleur peut être utilisé comme unité de banc ou module rack; en tout cas, il doit être positionné de manière à permettre la circulation de l'air à travers les ouvertures appropriées présentes sur le cache supérieur et latéral. Si le contrôleur est utilisé comme module rack, DIL DOIT ETRE introduit dans un adaptateur pour rack à 4 compartiments afin d'éviter que le contrôleur ne tombe à l'intérieur du rack. Le panneau antérieur du contrôleur n'est pas conçu pour supporter le poids de l'unité. Ne pas installer ou utiliser le contrôleur dans des milieux exposés aux agents atmosphériques (pluie, gelée, neige), aux poussières, aux gaz agressifs, dans des milieux à risque d'explosion ou d'incendie.

NOTE

Si le contrôleur est installé sur un rack, enlever les quatre pieds et le positionner en laissant un espace minimum de 30 mm (1.2 pouces) en correspondance de la superficie supérieure ou inférieure.

AVERTISSEMENT!



Le contrôleur doit être installé de manière à pouvoir interrompre sans problèmes la tension (débranchement de la fiche ou interruption au moyen de l'interrupteur de la ligne générale du rack).

Le fonctionnement de la pompe n'est optimisé que si celle-ci est utilisée avec l'une des unités de contrôle Agilent spécifiques (Dual, MidiVac ou MiniVac).

ATTENTION!

Lors de l'utilisation des pompes, le respect des normes de sécurité est impérativement subordonné à l'emploi des unités de contrôle Agilent.

Durant le fonctionnement de l'appareillage, il est nécessaire que soient respectées les conditions environnementales suivantes:

- température: de +5 °C à +40 °C
- humidité relative: 0 – 95 % (non condensante)

Pour brancher le contrôleur à la pompe, utiliser les câbles appropriés fournis. Pour les informations détaillées à propos des connexions ci-dessus indiquées ou d'une autre nature, se référer à l'annexe "Technical Information".

Utilisation

Dans ce paragraphe sont reportées les principales procédures opérationnelles. Pour de plus amples détails et pour les informations inhérentes aux connexions facultatives, se référer au paragraphe "UTILISATION" de l'Annexe "Technical Information". Avant d'actionner le contrôleur, effectuer toutes les connexions électriques ainsi que vers la pompe et lire le Manuel d'Instructions de la pompe branchée.

AVERTISSEMENT!



Afin d'éviter des lésions personnelles ou des endommagements à l'appareillage, si la pompe est placée sur une table, vérifier que cette dernière soit stable. Ne pas actionner la pompe au cas où la relative fiche est débranchée du système ou bien obstruée.

Commandes et Indicateurs du Panneau Antérieur de l'unité Contrôleur pour Pompes Ioniques

La figure et le tableau suivants indiquent les commandes et les indicateurs du panneau antérieur du contrôleur avec les relatives descriptions et fonctions.

NOTA

La fonctionnalité et la signification de certaines commandes et indicateurs changent selon la position de l'interrupteur HV ON OFF. Le tableau suivant montre toutes les fonctions et toutes les significations des signaux.

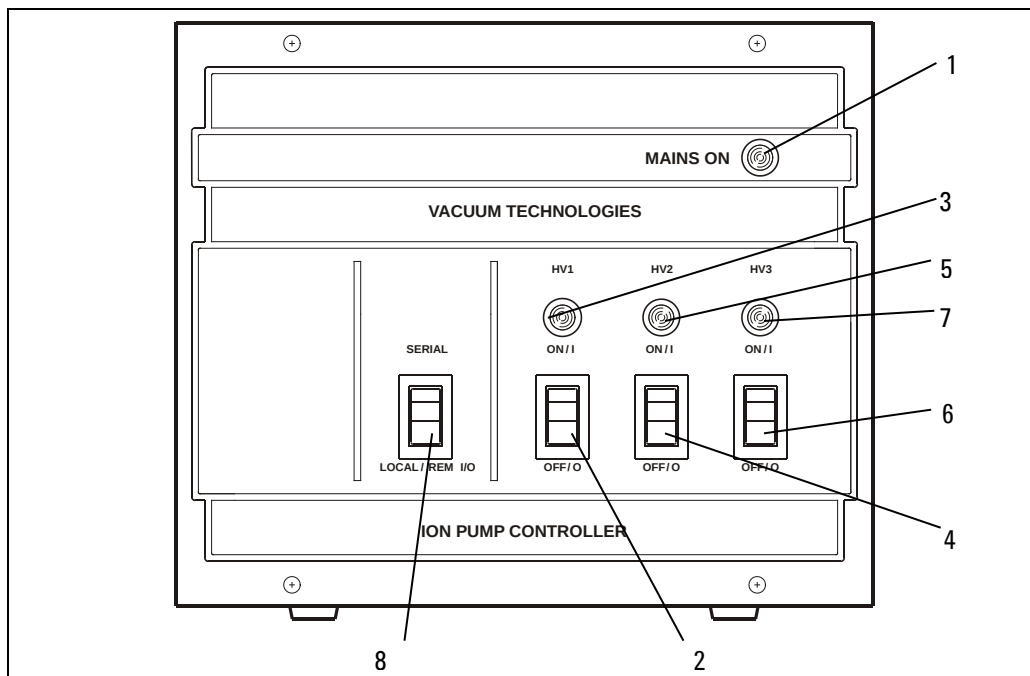


Figure 2 Commandes et indicateurs du panneau antérieur de l'unité contrôleur pour pompes ioniques

Tab. 1

RIF.	SIGNAL	DESCRIPTION
1	MAINS ON LED	Lorsqu'il y a une condition de On, cela veut dire que l'unité est branchée à la ligne de l'alimentation et que l'interrupteur Line sur le panneau postérieur est en position ON (I).
2	HV1 ON/OFF SWITCH	Lorsqu'il est en position ON (I), il habilite l'alimentation à haute tension pour la pompe 1 si le contrôleur est sur modalité opérationnelle I/O à distance/locale et que les interlock du câble extérieur et du câble haute tension sont fermés.
3	HV1 LED	Il est sur On lorsque le contrôleur fournit la haute tension pour la pompe 1. Il clignote en cas d'anomalie (voir paragraphe "Informations techniques" pour de plus amples détails).
4	HV2 ON/OFF SWITCH	Lorsqu'il est sur ON (I), il habilite l'alimentation à haute tension pour la pompe 2 si le contrôleur est sur modalité opérationnelle I/O à distance/locale et que les interlock du câble extérieur et du câble à haute tension sont fermés.
5	HV2 LED	Il est sur On lorsque le contrôleur fournit la haute tension pour la pompe 2. Il clignote en cas d'anomalie (voir paragraphe "Informations techniques" pour de plus amples détails).
6	HV3 ON/OFF SWITCH	Lorsqu'il est en position ON (I), il habilite l'alimentation à haute tension pour la pompe 3 si le contrôleur est sur la modalité opérationnelle I/O à distance /locale et que les interlock du câble extérieur et du câble haute tension sont fermés.
7	HV3 LED	Il est sur On lorsque le contrôleur fournit la haute tension pour la pompe 3. Il clignote en cas d'anomalie (voir paragraphe "Informations techniques" pour de plus amples détails).
8	SERIAL/LOCAL-REM I/O SWITCH	Il permet de sélectionner la modalité opérationnelle: SÉRIELLE: Les commandes sont fournies à travers la ligne série. I/O à distance/locale: Les commandes sont fournies en agissant sur le panneau antérieur de l'unité ou bien à travers le connecteur REMOTE à 25 pin présent sur le panneau postérieur.

Panneau postérieur de l'unité contrôleur pour Pompes Ioniques

La figure suivante montre les commandes et les indicateurs du panneau postérieur du contrôleur. Pour de plus amples renseignements détaillés à propos des connexions, se référer à l'annexe "Technical Information".

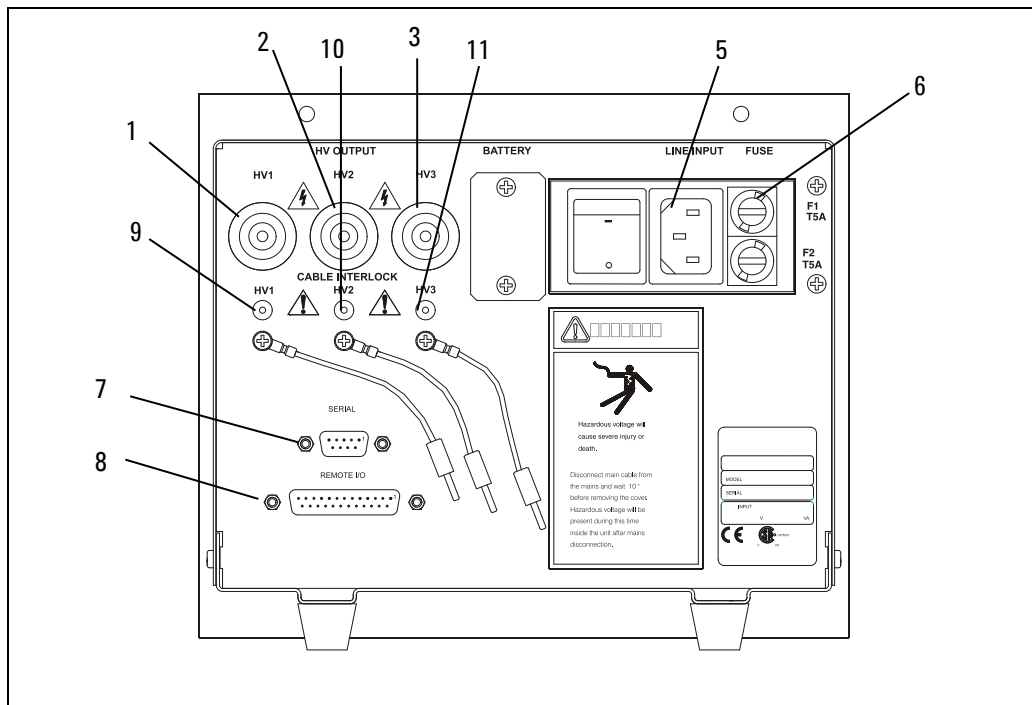


Figure 2 Commandes et indicateurs du panneau antérieur de l'unité contrôleur pour pompes ioniques

Tab. 2

RIF.	SIGNAL	DESCRIPTION
1	HV1 Connector	Connecteur HV pour câble de connexion à la pompe 1
2	HV2 Connector	Connecteur HV pour câble de connexion à la pompe inique 2
3	HV3 Connector	Connecteur HV pour câble de connexion à la pompe inique 3
5	LINE INPUT	Module contenant la fiche du câble d'entrée de la ligne et l'interrupteur LINE
6	FUSE	Support fusibles (2 fusibles)
7	SERIAL Connector	Connecteur DB9 pour câble sériel
8	REMOTE Connector	Connecteur DB25 pour connexion I/O à distance
9	HV1 Cable Interlock	
10	HV2 Cable Interlock	
11	HV3 Cable Interlock	

Procédures d'utilisation

Ce paragraphe fournit les informations générales relatives à l'utilisation du contrôleur. Pour les détails, se référer au paragraphe "Informations techniques" en branchant le câble d'alimentation du contrôleur à la ligne d'alimentation et en positionnant l'interrupteur LINE présent dans le panneau postérieur sur ON(I), l'appareillage est en mesure de fournir les prestations suivantes:

- 3 canaux haute tension indépendants
- tension en sortie max. di -5000 Vcc
- puissance en sortie max. di 40 W pour chaque canal
- courant en sortie max. >30 mA (condition de court circuit) pour chaque canal
- modalité opérationnelle sérielle I/O à distance-locale
- fonctionnement START/PROTECT

Maintenance

Le Contrôleur ne nécessite aucune maintenance. Tout travail réalisé sur le contrôleur doit être effectué par un personnel autorisé.

En cas de panne, il est possible de contacter le service de réparation Agilent. Il est également possible de commander des contrôleurs d'avance par le biais d'Agilent.

AVERTISSEMENT!



Avant de procéder à une quelconque intervention sur le contrôleur, ne pas oublier de débrancher l'appareil du courant.

Modalite Operationnelle

Connexion à la ligne d'alimentation

Modalité opérationnelle I/O LOCALE-A DISTANCE

NOTE

Les 3 interrupteurs haute tension présents sur le panneau antérieur et les 3 entrées On/Off HV sur le connecteur DB25 "Remote" sur le panneau postérieur sont en série.

L'alimentation à haute tension aux 3 pompes peut être habilitée au moyen de 3 interrupteurs HV (HV1, HV2, HV3) présents sur le panneau antérieur (dans ce cas la fiche DB25 avec les shunts intérieurs fournis avec l'unité doit être connectée et les 3 interlock du câble haute tension doivent être fermés) ou bien au moyen de la fermeture des 3 entrées On/Off HV sur le connecteur DB25 "Remote" présent sur le panneau postérieur (dans ce cas les 3 interrupteurs HV sur le panneau antérieur doivent être en position ON).

Lorsque la haute tension est sur ON, le relatif voyant LED sur le panneau antérieur s'allume.

NOTE

Si la connexion entre le IPCU et la pompe se fait par l'intermédiaire d'un câble haute tension avec interlock de sécurité, le câble doit être connecté au F/T de la pompe et le connecteur unipolaire aux dimensions réduites doit être introduit dans la relative fiche "HV Cable Interlock" présente sur le panneau postérieur de l'unité afin de pouvoir activer la haute tension.

Si la connexion est réalisée avec un câble haute tension unipolaire sans interlock de sécurité, vérifier que le shunt fourni sur le panneau postérieur de l'unité soit inséré dans la fiche "HV Cable Interlock".

Fonctionnement SERIEL

NOTE

La modalité sériele est programmée de manière à utiliser l'appareillage en tranchant le contrôleur à un PC externe qui utilise le programme "HYPER TERMINAL" de Windows. Pour travailler en modalité sériele, connecter le port série au connecteur "Serial" du contrôleur sur le panneau postérieur, ensuite effectuer le programme "HYPER TERMINAL" avec les suivantes attributions:

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

Cette modalité opérationnelle permetta de vérifier la fonctionnalité de l'unité, lancer les commandes et accéder à toutes les données opérationnelles.

Pour de plus amples détails, se référer au paragraphe "Technical Information".

Modalité opérationnelle START/PROTECT

Le contrôleur peut fonctionner dans les deux modalités suivantes: Start et Protect.

En modalité Start, les 3 canaux en sortie peuvent fournir toute la puissance nécessaire en sortie sans tenir compte des conditions de la pompe jusqu'à la condition de court circuit.

Cette modalité est utilisée pour actionner la pompe à une pression élevée.

En modalité Protect, le contrôleur limite le courant en sortie et désactive la haute tension lorsque le courant dépasse la valeur de seuil (8 mA) pour plus de 2 secondes.

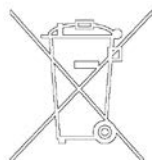
Mise au rebut

Signification du logo "WEEE" imprimé sur les étiquettes.

Le symbole ci-dessous est appliqué conformément à la directive CE nommée "WEEE".

Ce symbole (**uniquement valide pour les pays de la Communauté européenne**) indique que le produit sur lequel il est appliqué NE doit PAS être mis au rebut avec les ordures ménagères ou les déchets industriels ordinaires, mais passer par un système de collecte sélective.

Après avoir vérifié les termes et conditions du contrat de vente, l'utilisateur final est donc prié de contacter le fournisseur du dispositif, maison mère ou revendeur, pour mettre en œuvre le processus de collecte et mise au rebut.



Messages d'erreur

Sur le panneau antérieur

Durant le fonctionnement du contrôleur, on a trouvé une erreur, les voyants LED haute tension sur le panneau antérieur fourniront les indications nécessaires suivantes:

HV 1,2,3

LED ON: Fonctionnement normal = HV ON

Led HV 1,2,3 clignotants: condition d'erreur:

Clignotement continu = trop de jus en Protect ($I > 8$ mA)

1 clignotement toutes les 2 secondes = Interlock câble HV

2 clignotements toutes les 2 secondes = remote Interlock

3 clignotements toutes les 2 secondes =

HV survoltage (attention: ne pas éteindre l'HV)

4 clignotements toutes les 2 secondes = erreur interne ADC

5 clignotements toutes les 2 secondes =

rop de jus sur la haute tension (court circuit)

($I > 30$ mA pendant plus de 30 secondes)

6 clignotements toutes les 2 secondes =

Le commutateur local/Remote, à distance,
a été déplacé avec HV ON

7 clignotements toutes les 2 secondes = surchauffe interne à
unité

En modalité I/O à distance-locale, la condition d'anomalie est remise à zéro en positionnant sur Off (0) l'interrupteur du canal correspondant présent sur le panneau antérieur.

En modalité série, la condition d'anomalie est remise à zéro en envoyant la commande "FOn" à travers l'ordinateur.

NOTE

Au cas où une des conditions suivantes s'avèrent tandis que la haute tension du canal est active, la sortie du canal est programmée immédiatement sur OFF, le voyant LED sur le panneau antérieur inique la condition d'anomalie correspondante et l'information inhérente à cette dernière est envoyée à la sortie sérielle.

Modification de la condition de l'interrupteur Local-Remote I/O
Interrupteur HV du panneau antérieur positionné sur Off (0)
Interlock du câble HV et/ou externe ouvert

Sur le connecteur "Remote" dans le panneau postérieur

Sortie "Fault": contact N.C. =

Le contact est ouvert durant le fonctionnement normal (relais ON)

Le contact est fermé en condition d'anomalie (relais OFF)

Sortie "HV 1,2,3": contact N.O. =

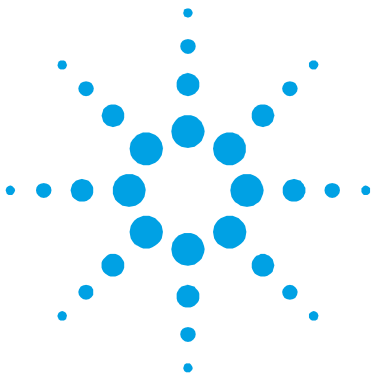
Le contact est ouvert avec HV Off (relais Off)

Le contact est fermé avec HV On (relais On)

Tranche des relais des sorties: 1 A 48 Vca/cc

3 Procédure pour l'installation

Messages d'erreur



4

Installation Procedure

General Information	64
Symbols used	65
Controller Description	65
Storage	66
Preparation for Installation	67
Installation	68
Use	69
Ion Pumps C.U. Front Panel Controls and Indicators	70
Ion Pumps C.U. Rear Panel	72
Usage	74
Maintenance	75
Operating Modes	76
Power Line Supply	76
LOCAL-REMOTE I/O Operation	76
SERIAL Operation	77
START/PROTECT Operation	77
Disposal	78
Error Messages	79
On Front Panel	79
On "Remote" Connector on Rear Panel	80

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-authorised 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:

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.

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

It can drive up to 3 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 Information” 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.

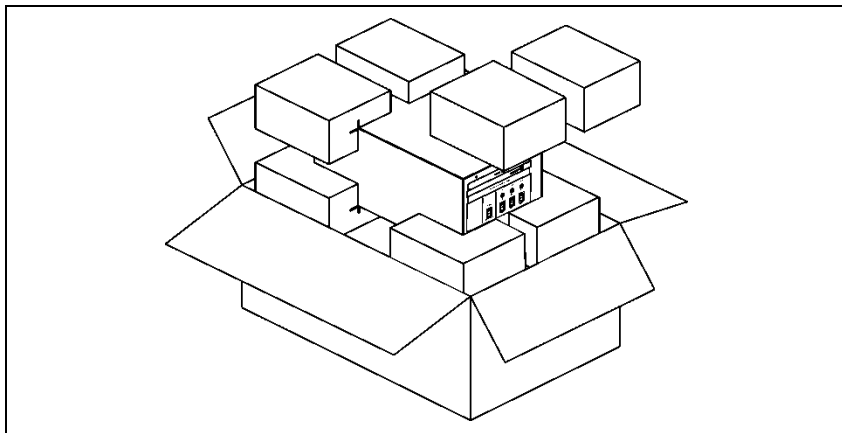


Figure 1 Packing

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.

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.

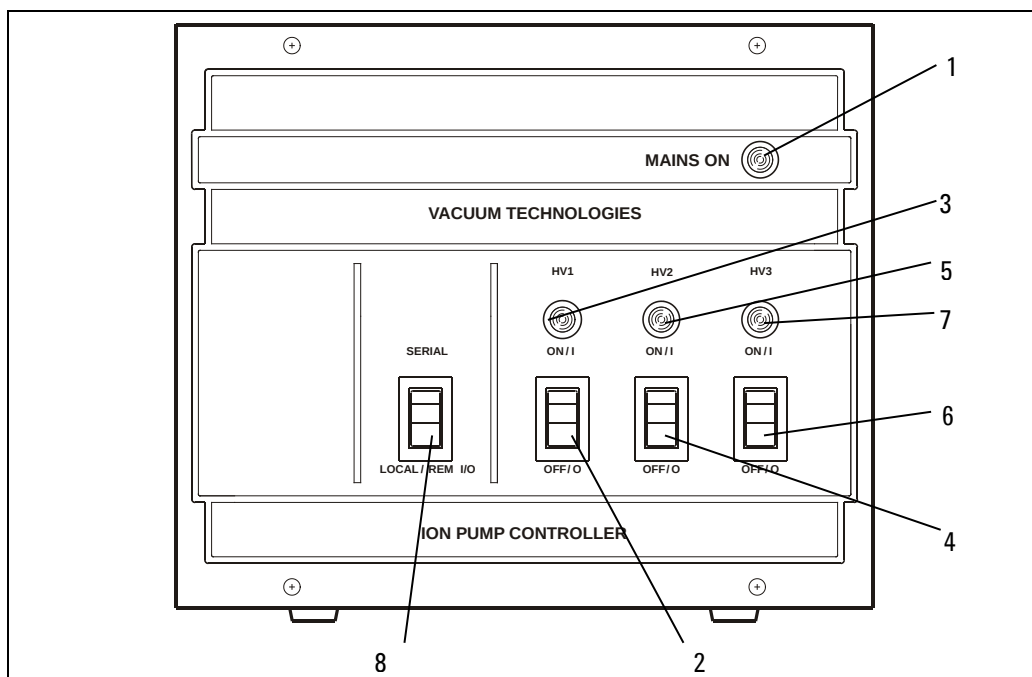


Figure 2 Ion Pumps C.U. Front Panel Controls and Indicators

Tab. 1

RIF.	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	HV1 ON/OFF SWITCH	When in ON (I) position, it switches On the HV to the pump 1 if the controller is set for LOCAL/REMOTE I/O operation and the external and HV Cable interlocks are closed
3	HV1 LED	It is On when the controller is supplying the HV to the pump 1. It is flashing in case of Fault (see paragraph "Technical Information" for details)
4	HV2 ON/OFF SWITCH	When in ON (I) position, it switches On the HV to the pump 2 if the controller is set for LOCAL/REMOTE I/O operation and the external and HV Cable interlocks are closed
5	HV2 LED	It is On when the controller is supplying the HV to the pump 2. It is flashing in case of Fault (see paragraph "Technical Information" for details)
6	HV3 ON/OFF SWITCH	When in ON (I) position, it switches On the HV to the pump 3 if the controller is set for LOCAL/REMOTE I/O operation and the external and HV Cable interlocks are closed
7	HV3 LED	It is On when the controller is supplying the HV to the pump 3. It is flashing in case of Fault (see paragraph "Technical Information" for details)
8	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”.

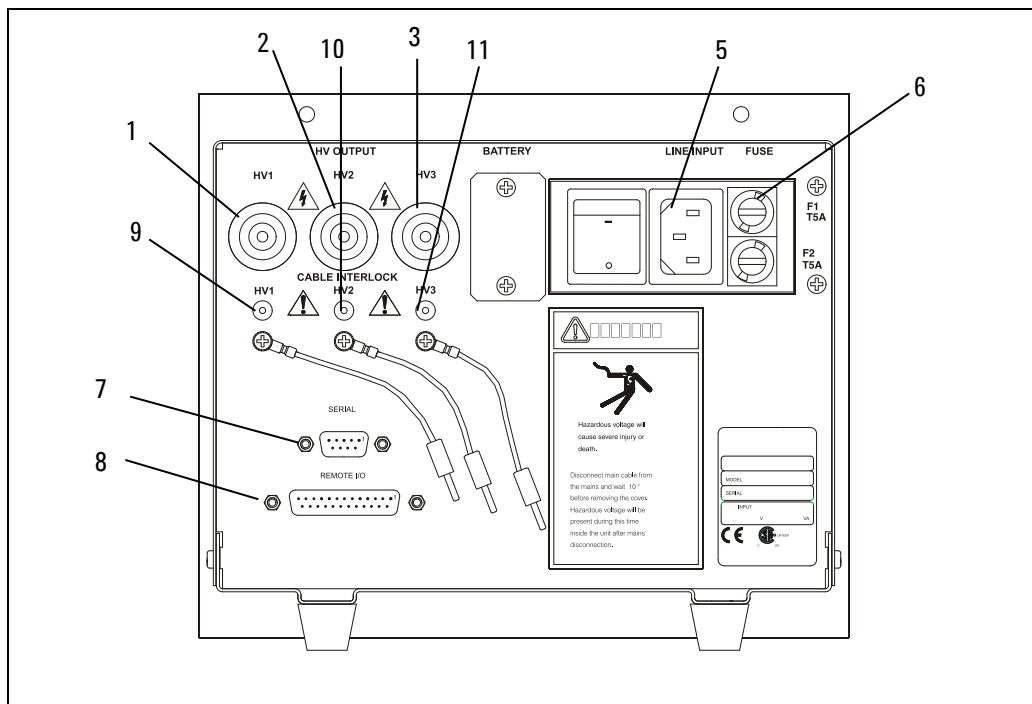


Figure 3 Ion Pumps C.U. Rear Panel

Tab. 2

RIF.	SIGNAL	DESCRIPTION
1	HV1 Connector	HV Connector for Cable to Ion Pump 1
2	HV2 Connector	HV Connector for Cable to Ion Pump 2
3	HV3 Connector	HV Connector for Cable to Ion Pump 3
5	LINE INPUT	Module containing the Line Input cable plug and the Line Switch
6	FUSE	Protection Fuse holder (2 fuses)
7	SERIAL Connector	DB9 connector for the Serial cable
8	REMOTE Connector	DB25 connector for the REMOTE I/O connections
9	HV1 Cable Interlock	
10	HV2 Cable Interlock	
11	HV3 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:

- 3 independent HV channels
- -5000 Vdc maximum output Voltage
- 40 W maximum output Power for each channel
- >30 mA maximum output current (short circuit condition) for each channel
- LOCAL-REMOTE I/O/SERIAL operating mode
- START/PROTECT operation

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.

Operating Modes

Power Line Supply

LOCAL-REMOTE I/O Operation

NOTE

The 3 HV switches on the front panel and the 3 HV On/Off inputs on the DB25 “Remote” connector on the rear panel are in series.

The HV to the 3 pumps can be switched ON by the 3 HV (HV1, HV2, HV3) switches on the front panel (in this case the DB25 plug with the internal jumpers provided with the unit must be connected and the 3 HV Cable Interlocks must be closed) or by the closure of the 3 HV On/Off inputs on the DB25 “Remote” connector on the rear panel (in this case the 3 HV switches on the front panel must be in the ON position).

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

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” program of Windows. 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 settings:

- COM1
- 9600 baud
- No parity
- 8 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.

See the “Technical Information” paragraph for details.

START/PROTECT Operation

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

In Start, the 3 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 more than 2 seconds.

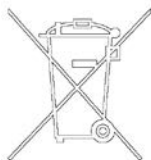
Disposal

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.



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:

HV 1,2,3

LEDS ON: Normal Operation = HV ON

Led HV 1,2,3 flashing: error condition:

continuous flashing = overcurrent in Protect ($I > 8 \text{ mA}$)

1 flash every 2 seconds = Interlock HV cable

2 flashes every 2 seconds = remote Interlock

3 flashes every 2 seconds = HV Overvoltage (Warning: do not switch off HV)

4 flashes every 2 seconds = error inside ADC

5 flashes every 2 seconds = overcurrent on HV (short circuit) ($I > 30 \text{ mA}$ for more than 30 seconds)

6 flashes every 2 seconds = the Local/Remote switch was moved with HV ON

7 flashes every 2 seconds = overtemperature inside unit

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.

Change of the Local-Remote I/O switch status

Front panel HV switch set to Off (0)

External and/or HV Cable Interlock open

On “Remote” Connector on Rear Panel

“Fault” output: N.C. contact =

 The contact is Open in normal operation (relay ON)

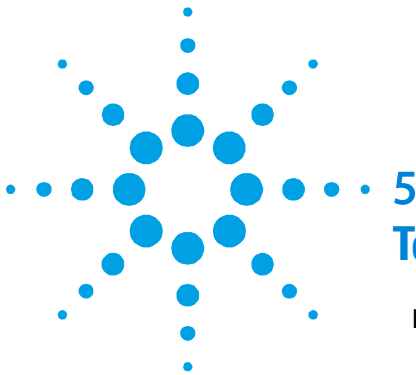
 The contact is Closed in “ Fault” condition (relay OFF)

“HV 1,2,3” output: N.O. contact =

 The contact is Open with HV Off (relay Off)

 The contact is Closed with HV On (relay On)

Outputs relays rating: 1 A 48 Vac/dc



5 Technical Information

ION Pump Controller Description	82
Controller Specifications	82
Remote I/O Connector Pinout	85
Input Commands Connection	86
Output Signals Connection	87
SERIAL Operation	87
Controller RS232 Standard Serial Pinouts (DB9 Female Connector)	88
Pin Functions	89
Procedure to Connect the Serial and I/O Ports to an External Cable	90
Ion Pump Control Unit - Operator / Host Serial Communication	92
Controller Report Mode	93
Controller Commands Mode	93
Clear Fault (Reset)	94
Fault Table	94
Cleaning	95

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.
- 3 boards: (HV generation), they generate the HV for the 3 output channels
- 1 board: (CPU), it contains all the digital circuitry and the microprocessor.

Controller Specifications

Tab. 3

Line:	
Voltage	100 - 240 Vac
Frequency	50 / 60 Hz
Power	200 VA
Operating conditions:	
Internal use only	
Max Altitude:	3000 m
Max Line Voltage variation:	± 10 %
Installation category:	II
Pollution degree:	2
External Battery specifications	12 Vdc 24 Ah Lead Acid Type
Output for the pumps:	
With Line Supply:	
Output Voltage (each channel)	-5000 Vdc +1, -5 %
Output Current (each channel)	>30 mA ± 10 %
Maximum Output Power (each channel)	40 W

With Battery supply:	
Output Voltage	-3000 Vdc +20, - 5 %
Output Current (total)	0.3 mA
Output Power	1 W
Operating Temperature	0 °C to 40 °C
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 Burst: EN61000/4/4 Radiated RF Immunity: EN61000/4/3 Surge: EN61000/4/5 Conducted noise imm.: EN61000/4/6 Line supply interruptions: EN61000/4/11 Safety: EN61010-1
Line Cable	3 m long cable (3 wires, 3 m long), no plug
Line fuse	Two T5A
High Voltage Connection	3 High Voltage Fischer connectors
Battery Connection	3 pin connector (for Lead Acid Battery 12 Vdc 24 Ah)

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

5 Technical Information

ION Pump Controller Description

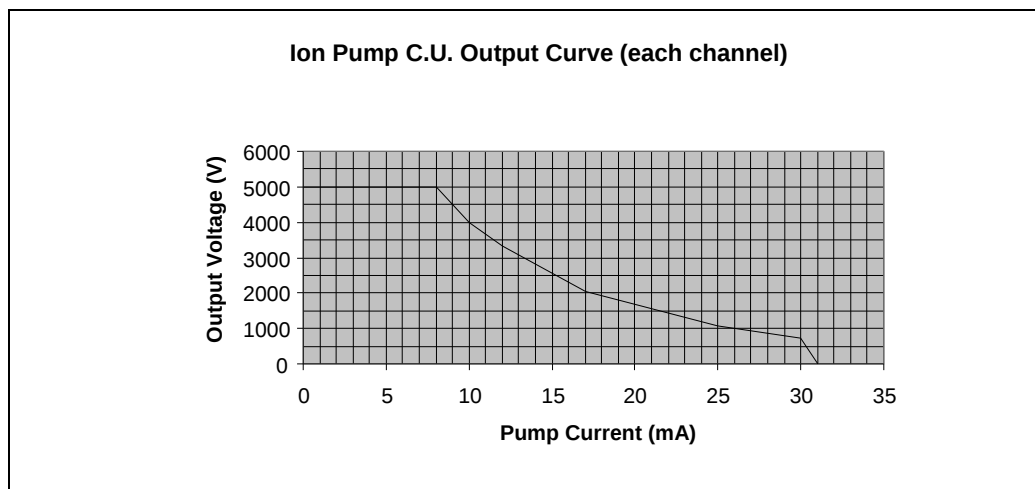


Figure 4 Output Voltage Curve for Line Supply

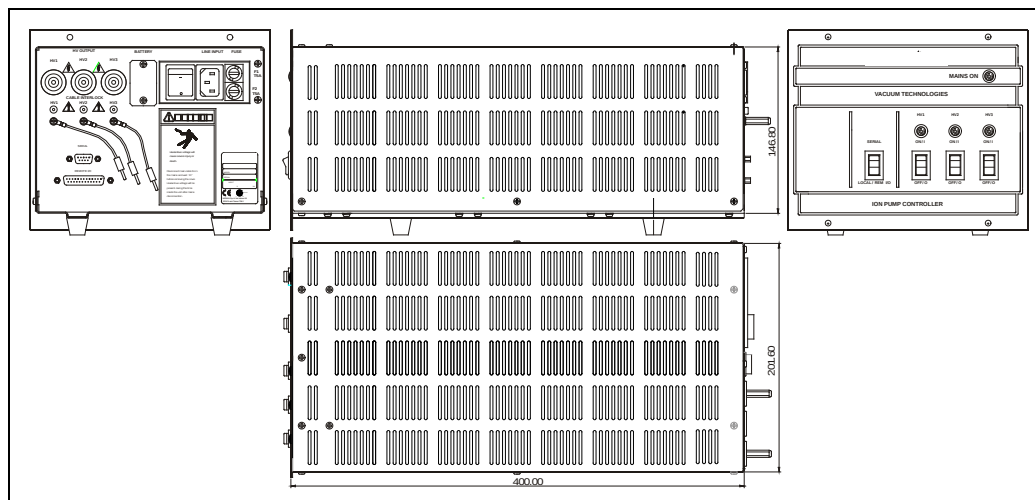


Figure 5 Controller Outline

Remote I/O Connector Pinout

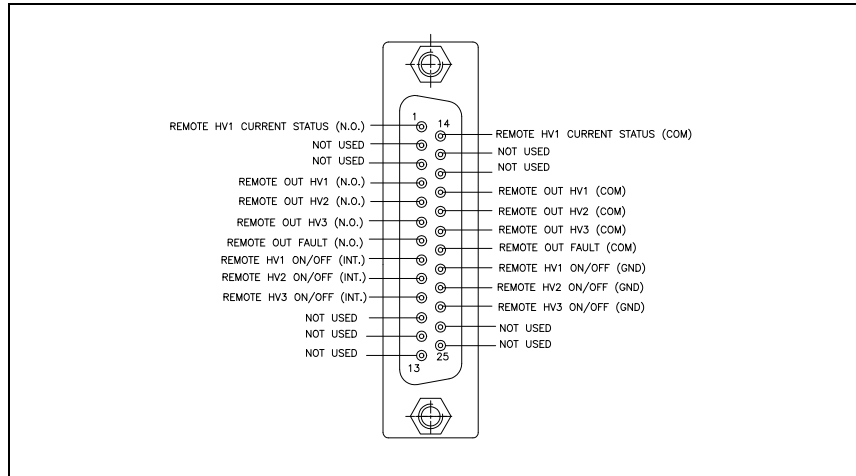


Figure 6 Remote I/O Pinout Rear Panel D-Shell Female 25P

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.

The controller is shipped with the mating (male) connector with the 3 jumpers on the HV 1,2,3 On/Off (External Interlock) inputs in order to allow to switch the HV On and Off by the unit front panel.

In case there is the need to use the Remote Interlock function or to switch On and Off the HV on the 3 pumps by Remote, remove the 3 jumpers and connect the appropriate connections.

5 Technical Information

ION Pump Controller Description

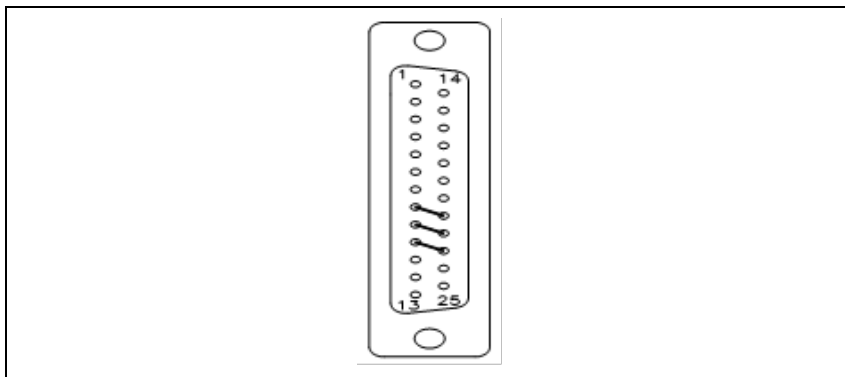


Figure 7 Remote I/O Jumpers Conn. D-Shell Male 25 Pin

Input Commands Connection

Input signals: Optically Insulated

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

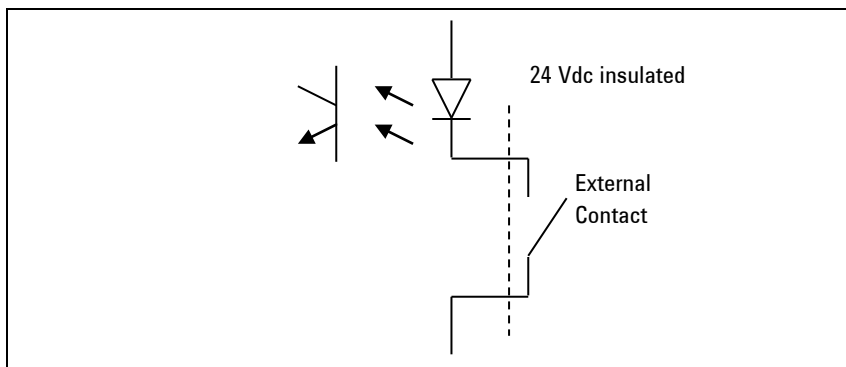


Figure 8

Output Signals Connection

Relay outputs: Single contact (Rating: 1A 48Vac/dc)

The HV1,2,3 output contacts will be Closed when the HV is ON (relay ON) and Open when the HV is OFF (relay OFF). The Fault output contact (N.C.) will be Open in Fault condition (relay OFF) and Closed in normal operation (relay ON)

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 Pinouts (DB9 Female Connector)

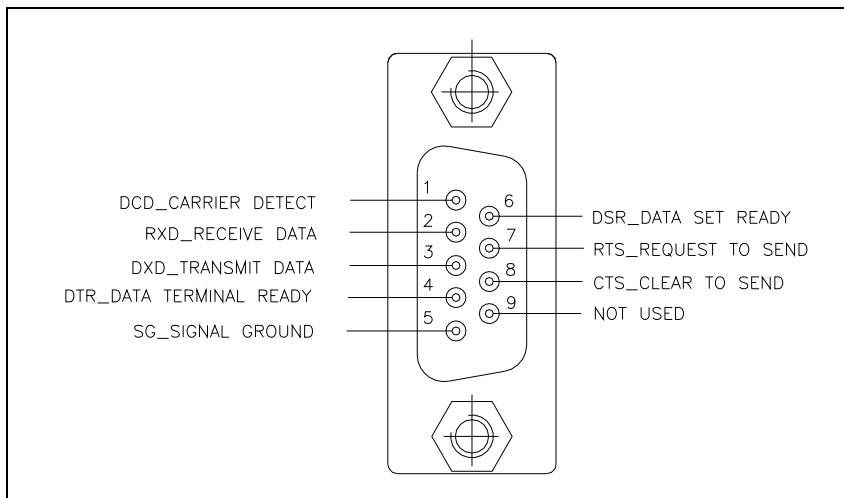


Figure 9 RS232 D-Shell Female 9P

Installation according to international rules

Tab. 4 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. 5

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 to 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 connected on the Controller connector and DTR is connected to DSR / DCD on IPCU connector then the DTE see 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.

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.

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

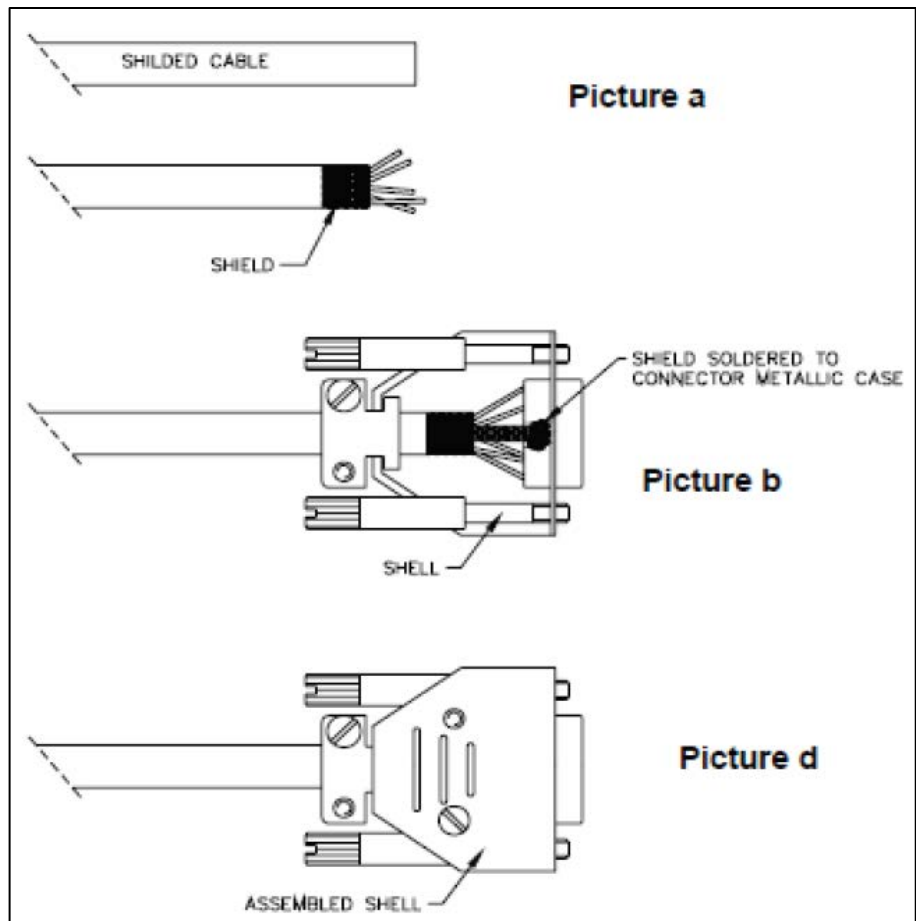


Figure 10

Ion Pump Control Unit - Operator / Host Serial Communication

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

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 500 msec, if the controller is in Report mode (default), all the 3 Report strings made of 30 characters each, corresponding to the 3 HV channels are sent by the IPCU.

The string is always made of 30 characters 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.

Controller Report Mode

"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 ... 99999nA, 100 ... 50000uA]	7 characters
' '	[0x20]	1 character SPACE
voltage	[0 ... 5999V]	6 characters 'CR '
	[0x0D]	1 character CR
Total 30 characters		

Controller Commands Mode

HV Channel Status Setting

"A0ns"

where :

A0 = HV Channel Serial Command for Status Set

n = HV Channel Number 1, 2 or 3

s = 0 for OFF, 1 for ON

5 Technical Information

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

HV Channel Start/Protect Mode Setting

"C0ns"

where :

C0 = HV Channel Serial Command for Start/Protect Setting

n = HV Channel Number 1, 2 or 3

s = 0 for START mode, 1 for PROTECT mode

Clear Fault (Reset)

"F0n"

where:

F0 = HV Channel Serial Command for Fault reset

N = HV Channel number 1,2,3

Fault Table

Err. nb	Description	Front Panel Led
[01]	Cable Interlock	1 blink every 2 sec.
[02]	Remote Interlock	2 blink every 2 sec.
[04]	HV Overvoltage	3 blink every 2 sec.
[08]	ADC Offset high	4 blink every 2 sec.
[10]	Output short circuit	5 blink every 2 sec.
[20]	Loc/Ser switch moved	6 blink every 2 sec.
[40]	Overtemperature	7 blink every 2 sec.
[80]	Overcurrent in Protect	8 blink every 2 sec. (Continuous)

Cleaning

For safety reasons, before cleaning the controller:
Turn the controller off;
disconnect the controller power plug from the electrical outlet
disconnect all cables.
If the exterior of the controller becomes dirty, use a dry soft cloth.
The internal cleaning must be done only by Agilent operator.

sssss 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", with a small flourish at the end.

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.

CUSTOMER REQUEST FOR CORRECTIVE / PREVENTIVE / IMPROVEMENT ACTION

TO: AGILENT VACUUM PRODUCTS DIVISION TORINO – QUALITY ASSURANCE FAX

N°: XXXX-011-9979350

ADDRESS: AGILENT TECHNOLOGIES ITALIA S.p.A. – Vacuum Products Division –
Via F.lli Varian, 54 – 10040 Leini (TO) – Italy

E-MAIL: vpd-qualityassurance_pdl-ext@agilent.com

NAME	COMPANY	FUNCTION
ADDRESS:		
TEL. N° : FAX N° :		
E-MAIL:		
PROBLEM / SUGGESTION :		
REFERENCE INFORMATION (model n°, serial n°, ordering information, time to failure after installation, etc.):		
DATE		
CORRECTIVE ACTION PLAN / ACTUATION (by AGILENT VPD)		LOG N°

XXX = Code for dialing Italy from your country (es. 01139 from USA; 00139 from Japan, etc.)

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

- o Remove all accessories from the core product (e.g. inlet screens, vent valves).
- o 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.
- o If ordering an Advance Exchange product, please use the packaging from the Advance Exchange to return the defective product.
- o 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.
- o Include a copy of the Health and Safety Declaration in the shipping documentation on the outside of the shipping box of your returning product.
- o 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.
- o Return only products for which the RA was issued.

Shipping

- o 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.
- o 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.

- o 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.
- o Agilent Technologies is not responsible for returning customer provided packaging or containers.
- o 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.
- o 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.
- o 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.
- o Products returned that have not been drained from oil will be disposed.
- o A Special Cleaning fee will apply to all exposed products.
- o If requesting a calibration service, units must be functionally capable of being calibrated.



Customer information		
Company :		Contact Name:
Address:		Tel: Fax:
		Email:

Equipment			
Product description	Agilent PartNo	Agilent Serial No	Original Purchasing Reference
Failure description		Type of process (for which the equipment was used)	

Type of return	
☐ Non Billable ☐ Billable  New PO # (hard copy must be submitted with this form): _____	
☐ Exchange ☐ Repair ☐ Upgrade ☐ Demo ☐ Calibration ☐ Evaluation ☐ Return for Credit	

Health and safety		Substances (please refer to MSDS forms)			
The product has been exposed to the following substances: (by selecting 'YES' you MUST complete the table to the right)		* Agilent will not accept delivery of any product that is exposed to radioactive, biological, explosive substances or dioxins, PCB's without written evidence of decontamination.			
		Trade name	Chemical name	Chemical Symbol	CAS Number
Toxic	☐ YES ☐ NO				
Harmful	☐ YES ☐ NO				
Corrosive	☐ YES ☐ NO				
Reactive	☐ YES ☐ NO				
Flammable	☐ YES ☐ NO				
Explosive (*)	☐ YES ☐ NO				
Radioactive (*)	☐ YES ☐ NO				
Biological (*)	☐ YES ☐ NO				
Oxidizing	☐ YES ☐ NO				
Sensitizer	☐ YES ☐ NO				
Other dangerous substances	☐ YES ☐ NO				

Goods preparation	
If you have replied YES to one of the above questions. Has the product been purged? ☐ YES ☐ NO	
If yes, which cleaning agent/method:	
Has the product been drained from oil? ☐ YES ☐ NOT APPLICABLE	
I confirm to place this declaration on the outside of the shipping box. ☐	

I declare that the above information is true and complete to the best of my knowledge and belief. I understand and agree to the terms and conditions on page 2 of this document.	
Name:	Authorized Signature:
Position:	
Date:	
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.	

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 Penteado de Ulhoa
Rodrigues, 939 - 6º 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-customercare@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

04/2014

Publication Number: 87-900-116-01 (E)



Agilent Technologies



Agilent Technologies