

# Turbo-V Remote Panel

Model

969-8850

**Manuale di istruzioni**  
**Bedienungshandbuch**  
**Notice de mode d'emploi**  
**Manual de instrucciones**  
**Manual de instruções**  
**Bedrijfshandleiding**  
**Istruktionsbog**  
**Bruksanvisning**  
**Instruksjon manual**

**Ohjekäsikirja**  
**Felhasználói kézikönyv**  
**Podrecznik instrukcji**  
**Návod k použití**  
**Návod na obsluhu**  
**Priročnik za navodila**  
**User Manual**

87-900-987-01 (E)

11/2011



**Agilent Technologies**

## Notices

© Agilent Technologies, Inc. 2011

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: 87-900-987-01 (E)

### Edition

Edition 11/2011

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.

# Turbo-V Remote Panel



## **Turbo-V Remote Panel**

# Contents

## 1 Istruzioni per l'uso 13

Informazioni Generali	14
Immagazzinamento	16
Preparazione per l'installazione	16
Installazione	17
Uso	19
Manutenzione	23
Smaltimento	23
Messaggi di errore	24

## 2 Gebrauchsanleitung 27

Allgemeine Informationen	28
Lagerung	30
Installationsvorbereitung	30
Installation	31
Anwendung	33
Wartung	37
Entsorgung	37
Fehlermeldungen	38

## Contents

<b>3</b>	<b>Mode d'emploi</b>	<b>41</b>
	Indications generales	42
	Emmagasinage	44
	Preparation pour l'installation	44
	Installation	45
	Utilisation	47
	Entretien	51
	Mise au rebut	51
	Messages d'erreur	52
<b>4</b>	<b>Manual de instrucciones</b>	<b>55</b>
	Información general	56
	Almacenamiento	58
	Preparación para la instalación	58
	Instalación	59
	Uso	61
	Mantenimiento	65
	Eliminación	65
	Mensajes de error	66
<b>5</b>	<b>Manual de Instruções</b>	<b>69</b>
	Informações gerais	70
	Armazenagem	72
	Preparação para a instalação	72

Instalação	73
Utilização	75
Manutenção	79
Eliminação	79
Mensagens de erro	80

## **6 Bedrijfshandleiding 83**

Algemene Informatie	84
Opslag	86
Vorbereiding voor installatie	86
Installatie	87
Gebruik	89
Onderhoud	93
Afvalverwerking	93
Foutmeldingen	94

## **7 Istruktionsbog 97**

Generel information	98
Opbevaring	100
Forberedelser før installation	100
Installation	101
Anvendelse	103
Vedligeholdelse	107
Fejlmeddelelser	108

## Contents

### **8 Bruksanvisning 111**

- Allmän information 112
- Förvaring 114
- Förberedelser för installation 114
- Installation 115
- Användning 117
- Underhåll 121
- Bortskaffning 121
- Felmeddelanden 122

### **9 Instruksjon Manual 125**

- Generell informasjon 126
- Lagring 128
- Forberede installasjonen 128
- Installasjon 129
- Bruk 131
- Vedlikehold 135
- Eliminering 135
- Feilmeldinger 136

### **10 Ohjekäsikirja 139**

- Yleistä 140
- Säilytys 142
- Asennuksen valmistelu 142

Asennus	143
Käyttö	145
Huolto	149
Hävittäminen	149
Vianetsintä	150

## **11 Felhasználói Kézikönyv 153**

Általános információk	154
Tárolás	156
A telepítésre való előkészítés	156
Telepítés	157
Használat	159
Karbantartás	163
Megsemmisítés	163
Hibaüzenetek	164

## **12 Podrecznik Instrukcji 167**

Informacje ogólne	168
Magazynowanie	170
Przygotowanie do instalacji	170
Instalacja	171
Użytkowanie	173
Konserwacja	177
Przetworstwo odpadów	177

## Contents

	Bledne informacije	178
<b>13</b>	<b>Návod k Použití</b>	<b>181</b>
	Všeobecné informace	182
	Uskladnění	184
	Příprava k instalaci	184
	Instalace	185
	Použití	187
	Údržba	191
	Likvidace	191
	Chybné zprávy	192
<b>14</b>	<b>Návod na Obsluhu</b>	<b>195</b>
	Všeobecné informácie	196
	Uskladňovanie	198
	Príprava pre inštaláciu	198
	Inštalácia	199
	Použitie	201
	Údržba	205
	Likvidácia	205
	Oznamy vád	206
<b>15</b>	<b>Priročnik za Navodila</b>	<b>209</b>
	Splošne informacije	210
	Skladiščenje	212

Priprava za montažo 212

Montaža 213

Uporaba 215

Vzdrževanje 219

Odlaganje odpadkov 219

Obvestilo o napaki 220

## **16 Instructions for Use 223**

General Information 224

Storage 226

Preparation for Installation 226

Installation 227

Use 229

Maintenance 233

Disposal 233

Error Messages 234

## **17 Technical Information 237**

Specifications 239

Controller Outline 240

Controller Connection 242

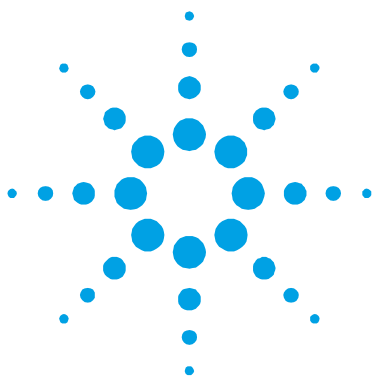
J5 – Gauge Connector 251

J2 – Serial Connector 252

Window Protocol Description 255

## Contents

How to use by Front Panel	266
Use	272
Programming	278
How to use by Remote I/O	292
How to use in Serial Mode	292
Error Messages	294
Orderable Parts	297



# 1

## Istruzioni per l'uso

Informazioni Generali	14
Immagazzinamento	16
Preparazione per l'installazione	16
Installazione	17
Uso	19
Comandi, indicatori e connettori	19
Descrizione pannello frontale	20
Accensione del Controller	22
Avvio della pompa	22
Arresto della Pompa	22
Manutenzione	23
Smaltimento	23
Messaggi di errore	24

Traduzione delle istruzioni originali



## Informazioni Generali

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

Il Turbo-V Remote Panel è controllato da un microprocessore e consente all'utente di comandare una pompa di tipo "controller on-board" mediante un pannello frontale:

TV551 con Navigator controller

TV701 con Navigator controller

TV1001 con Navigator controller

TV2k-G con controller on-board

TV3k-T con controller on-board

TV3k-G con controller on-board

TV70 con Navigator controller

TV81 con Navigator controller

TV301 con Navigator controller

TV303 con Navigator controller

Nei paragrafi seguenti sono riportate tutte le informazioni necessarie a garantire la sicurezza dell'operatore durante l'utilizzo dell'apparecchiatura. Informazioni dettagliate sono fornite nell'appendice "Technical Information".

**Questo manuale utilizza le seguenti convenzioni:**

**ATTENZIONE!**

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

---

**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.

---

**NOTA**

Le note contengono informazioni importanti estrapolate dal testo.

---

## Immagazzinamento

Durante il trasporto e l'immagazzinamento dei controller devono essere soddisfatte le seguenti condizioni ambientali:

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

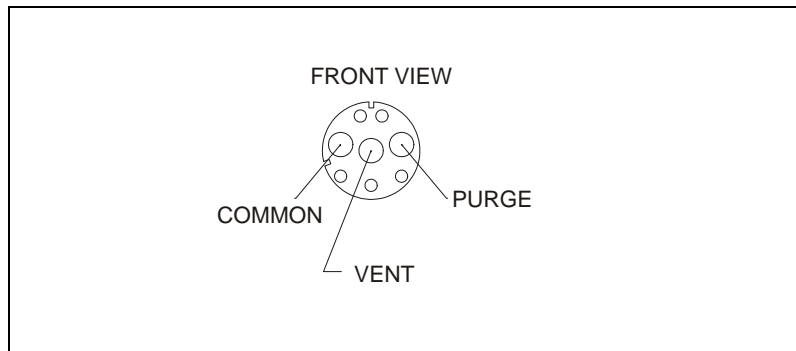
## Preparazione per l'installazione

Il Turbo-V Remote Panel viene fornito in un imballo protettivo speciale; se si presentano segni di danni, che potrebbero essersi verificati durante il trasporto, contattare l'ufficio vendite locale. Durante l'operazione di disimballaggio, prestare particolare attenzione a non lasciar cadere il controller e a non sottoporlo ad urti. Non disperdere l'imballo nell'ambiente. Il materiale è completamente riciclabile e risponde alla direttiva CEE 85/399 per la tutela dell'ambiente.

## Installazione

Il Turbo-V Remote Panel deve essere collegato al controller mediante l'apposito cavo (vedere tabella delle parti ordinabili), come illustrato nella seguente figura. Poiché la potenza/i segnali richiesti sono forniti da questo cavo, non è necessaria nessun'altra alimentazione esterna.

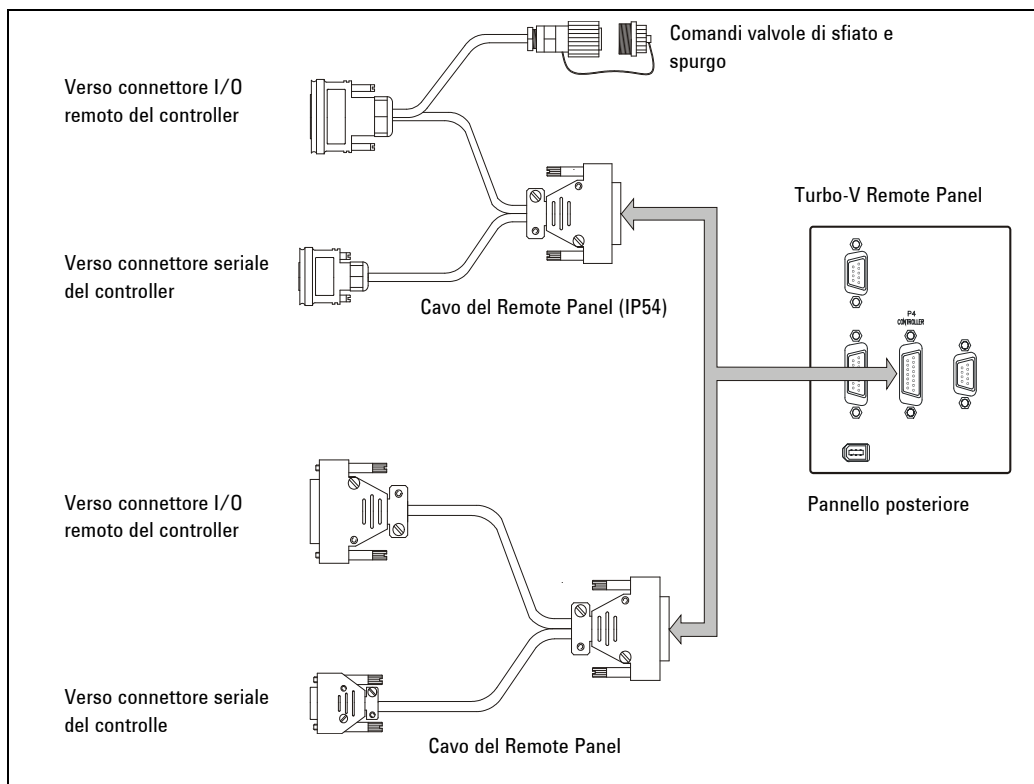
Se l'operatore desidera comandare le valvole di sfiato e spurgo utilizzando i comandi esterni, utilizzare la versione con cavo IP54, impostare il controller su "valve setting" "on command" e collegare i cavi come illustrato nella figura. Le valvole sono normalmente chiuse; fornendo 24 V tra la valvola di sfiato e spurgo e il terminale comune, le valvole si apriranno.



**Figura 1**

# 1 Istruzioni per l'uso

## Installazione



**Figura 2** Collegamenti

### NOTA

Il Turbo-V Remote Panel può essere installato su di un tavolo o all'interno di un apposito rack. In ogni caso occorre che l'aria di raffreddamento possa circolare liberamente intorno all'apparato. Non installare e/o utilizzare il controller in ambienti esposti ad agenti atmosferici (pioggia, gelo, neve), polveri, gas aggressivi, in ambienti esplosivi o con elevato rischio di incendio.

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

- temperatura: da 5 °C a +45 °C
- umidità relativa: 0 – 95 % (non condensante).

Per gli altri collegamenti e l'installazione degli accessori opzionali, vedere la sezione "Technical Information".

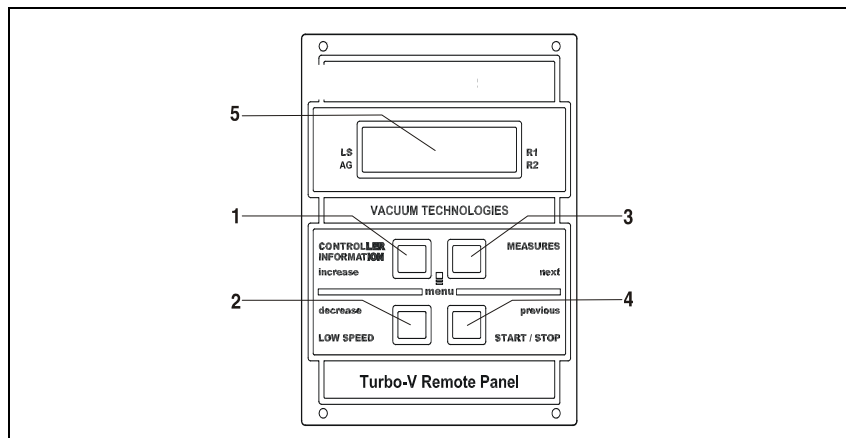
## Uso

In questo paragrafo sono riportate le principali procedure operative. Per ulteriori dettagli e per procedure che coinvolgono collegamenti o particolari opzionali, fare riferimento al paragrafo "Use" dell'appendice "Technical Information".

## Comandi, indicatori e connettori

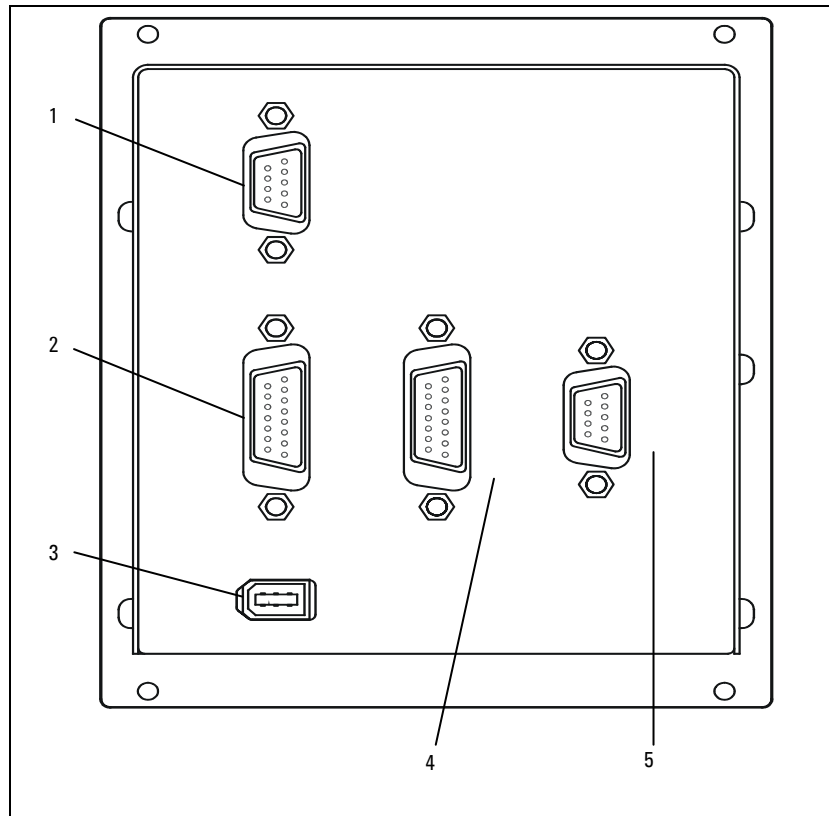
Di seguito sono illustrati il pannello frontale e posteriore del controller. Per maggiori dettagli fare riferimento alla sezione "Technical Information".

## Descrizione pannello frontale



**Figura 3** Turbo-V Remote Panel

1	Pulsante per richiamare sul display i parametri cycle number, cycle time e pump life.
2	Pulsante per la selezione del modo LOW SPEED. È attivo solo quando è selezionato il modo di comando dal pannello frontale. Premendolo una volta, la pompa ruota a velocità "stand-by". Premendolo ancora si disattiva il modo LOW SPEED.
3	Pulsante per richiamare sul display i parametri pump current, pump temperature, pump power e rotational speed. È sempre attivo indipendentemente dal modo di funzionamento scelto. Premendo assieme i pulsanti 3 e 4 per almeno 2 secondi viene attivato un programma con il quale è possibile programmare alcuni parametri operativi.
4	Pulsante per inviare i comandi di START, STOP/RESET. È attivo solo quando è selezionato il modo di comando dal pannello frontale. Premendolo una volta si attiva la fase di avvio; premendolo nuovamente si arresta la pompa. Se la pompa si è fermata automaticamente a causa di un guasto, occorre premere questo pulsante una prima volta per eseguire il reset del controller ed una seconda volta per riavviare la pompa.
5	Display alfanumerico a cristalli liquidi: matrice di punti, 2 linee x 16 caratteri.



**Figura 4** Pannello posteriore del Turbo-V Remote Panel

- |   |                                                                                                                                 |
|---|---------------------------------------------------------------------------------------------------------------------------------|
| 1 | Connettori di ingresso dei segnali logici (il connettore di accoppiamento viene fornito con l'apposito ponticello di chiusura). |
| 2 | Connettore segnali logici in uscita e monitoraggio frequenza pompa più uscita analogica programmabile.                          |
| 3 | Connettore di uscita alimentazione (24 Vcc per calibro).                                                                        |
| 4 | Connettore per ingresso alimentazione dal controller della pompa e comunicazione seriale con il controller della pompa.         |
| 5 | Connettore seriale per controllo remoto.                                                                                        |

## Accensione del Controller

Per accendere il controller è sufficiente inserire il cavo di alimentazione nella presa di rete.

## Avvio della pompa

Per avviare la pompa occorre abilitare l'ingresso di interlock e premere il pulsante START del pannello frontale.

## Arresto della Pompa

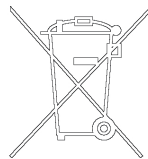
Per arrestare la pompa occorre premere il pulsante STOP del pannello frontale.

## Manutenzione

I controller della serie Turbo-V 301 Rack non richiedono alcuna manutenzione. Qualsiasi intervento deve essere eseguito da personale autorizzato. In caso di guasto è possibile usufruire del servizio di riparazione Agilent o del "Agilent advance exchange service", che permette di ottenere un controller rigenerato in sostituzione di quello guasto. Qualora un controller dovesse essere rottamato, procedere alla sua eliminazione nel rispetto delle normative nazionali specifiche.

## Smaltimento

**Significato del logo "WEEE" presente sulle etichette.** Il simbolo qui sotto riportato è 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

In alcuni casi di guasto la circuiteria di autodiagnosi del controller presenta alcuni messaggi di errore elencati nella tabella seguente.

<b>MESSAGGIO</b>	<b>DESCRIZIONE</b>	<b>AZIONE CORRETTIVA</b>
COMMUNICATION ERROR	Il Remote Panel non riesce ad avviare la comunicazione con il controller.	Verificare la connessione seriale tra il Remote Panel e il controller.
COMMUNICATION BROKEN	Il Remote Panel non riesce a continuare la comunicazione con il controller.	Verificare la connessione seriale tra il Remote Panel e il controller.
XXXXXXX MODEL	Controller non supportato. Il Remote Panel funziona soltanto con le seguenti pompe: TV551 con navigator controller, TV701 con navigator controller, TV1001 con navigator controller, TV2k-G e TV3k-T.	Collegare il Remote Panel al controller supportato.
CHECK CONNECTION TO PUMP	Malfunzionamento nel collegamento tra pompa e controller. Oppure La pompa ha una temperatura inferiore a 0 °C.	Verificare che il cavo di collegamento tra pompa e controller sia ben fissato da entrambe le estremità e non sia interrotto. Premere due volte il pulsante START per riavviare la pompa.
WAITING INTERLOCK	È attivo il segnale di interlock presente sul connettore P1 a causa dell'interruzione del corto circuito tra il pin 3 ed il pin 8 del connettore J1, o a causa dell'apertura del segnale di interlock esterno.	Ripristinare il corto circuito tra il pin 3 ed il pin8 del connettore J1, o chiudere il segnale di interlock esterno.
PUMP OVERTEMP.	La temperatura del cuscinetto superiore o della pompa ha superato i 60 °C.	Attendere che la temperatura ritorni al di sotto della soglia. Premere due volte il pulsante START per riavviare la pompa.
CONTROLLER OVERTEMP.	La temperatura dell'ambiente del controller ha superato i 70 °C.	Attendere che la temperatura ritorni al di sotto della soglia.

<b>MESSAGGIO</b>	<b>DESCRIZIONE</b>	<b>AZIONE CORRETTIVA</b>
	Oppure La temperatura del radiatore del controller è superiore a 60 °C.	Premere due volte il pulsante START per riavviare la pompa.
TOO HIGH LOAD	La corrente assorbita dalla pompa è maggiore di quella programmata.	Verificare che il rotore della pompa abbia la possibilità di ruotare liberamente. Premere due volte il pulsante START per riavviare la pompa.
SHORT CIRCUIT	Durante il funzionamento normale (dopo la fase di avvio) la connessione di uscita è in corto circuito.	Verificare i collegamenti tra pompa e controller. Premere due volte il pulsante START per riavviare la pompa.
SYSTEM OVERRIDE	La pompa è stata fermata da un segnale di emergenza proveniente da un contatto remoto.	Staccare il cavo di alimentazione del controller e correggere la causa dell'emergenza. Ricollegare il cavo di alimentazione e premere due volte il pulsante START per riavviare la pompa.
OVERVOLTAGE	Si è verificato un guasto nella sezione di alimentazione del controller, o il controller ha ricevuto un segnale spurio.	Premere due volte il pulsante START per riavviare la pompa. Se il messaggio si ripresenta rivolgersi in Agilent per la manutenzione.
POWER FAIL	Guasto nella sezione di alimentazione della pompa, del controller.	Rivolgersi alla Agilent per Manutenzione.

**1 Istruzioni per l'uso**  
**Messaggi di errore**



## 2 Gebrauchsanleitung

Allgemeine Informationen	28
Lagerung	30
Installationsvorbereitung	30
Installation	31
Anwendung	33
Steuerungen, Anzeigen und Stecker	33
Beschreibung der Frontkonsole	34
Einschalten des Controllers	36
Pumpenstart	36
Pumpenstopp	36
Wartung	37
Entsorgung	37
Fehlermeldungen	38

Übersetzung der Originalanleitungen



## Allgemeine Informationen

Dieses Gerät ist zur Benutzung durch Fachleute gedacht. Vor seiner Anwendung sollte der Benutzer diese Gebrauchsanweisung und alles andere von Agilent gelieferten Informationsmaterial lesen. Agilent kann für nichts verantwortlich gemacht werden, was wegen auch nur teilweiser Nichtbeachtung dieser Anweisungen, falsche Benutzung durch nicht ausgebildetes Personal, nicht genehmigtes Eingreifen in das Gerät oder alle Aktionen im Widerspruch zu den spezifischen nationalen Normen verursacht wird.

Der Turbo-V Remote Panel wird von einem Mikroprozessor gesteuert und erlaubt den Betrieb einer Pumpe mit angebauter Steuerung über eine Frontkonsole:

- TV551 mit Navigator Controller
- TV701 mit Navigator Controller
- TV1001 mit Navigator Controller
- TV2k-G mit Controller on-board
- TV3k-T mit Controller on-board
- TV3k-G mit Controller on-board
- TV70 mit Navigator Controller
- TV81 mit Navigator Controller
- TV301 mit Navigator Controller
- TV303 mit Navigator Controller

Die folgenden Abschnitte enthalten alle notwendigen Informationen zur Garantie der Bedienersicherheit bei Benutzung des Geräts. Detaillierte Informationen finden sich im Anhang "Technical Information".

**Dies Handbuch benutzt die folgenden Standardkonventionen:**

**VORSICHT!**

Erscheint vor Prozeduren, bei deren Nichtbefolgen es zu schweren Schäden am Gerät kommen könnte.

---

**WARNUNG!**



Warnungen weisen den Bediener auf eine besondere Prozedur hin, bei deren Nichtbefolgen es zu schweren Verletzungen kommen könnte.

---

**HINWEIS**

Diese Hinweise enthalten wichtige Informationen aus dem Text.

---

## Lagerung

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

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

## Installationsvorbereitung

Der Turbo-V Remote Panel wird in einer speziellen Schutzverpackung geliefert. Wenn diese Schäden aufweist, die beim Transport entstanden sein können, das lokale Verkaufsbüro kontaktieren. Beim Auspacken darauf achten, dass das Gerät nicht herunterfällt oder gestoßen wird. Das Verpackungsmaterial vorschriftsmäßig entsorgen. Das Material kann zu 100 % recycelt werden und entspricht der EU-Richtlinie 85/399.

## Installation

Die Turbo-V Remote Panel muss mit dem Verbindungskabel (siehe die Tabelle der nachbestellbaren Teile) wie in der Abbildung weiter unten gezeigt an den Controller angeschlossen werden. Alle notwendigen Versorgungen und Signale liegen auf dem Kabel an. Daher ist keine externe Stromversorgung nötig.

Wenn der Anwender die Ventile zum Reinigen und Abblasen mit externen Befehlen betätigen will, die Kabelversion IP54 benutzen, den Controller in "valve setting" "on command" stellen und dann die Kabel wie in der Abbildung gezeigt anschließen. Die Ventile sind Öffner: sie öffnen sich bei Anlegen von 24 V zwischen den Vent Purge und Common Klemmen.

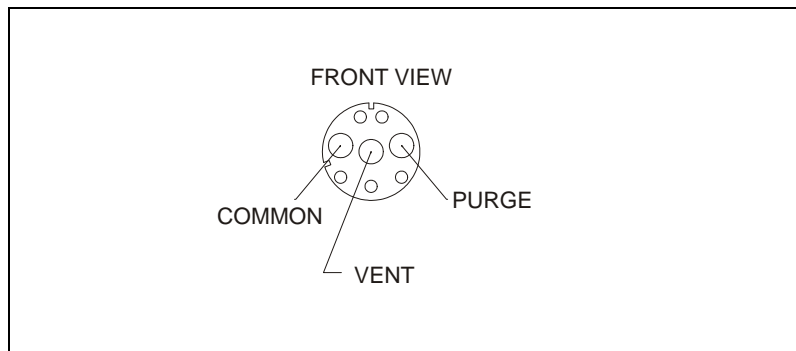
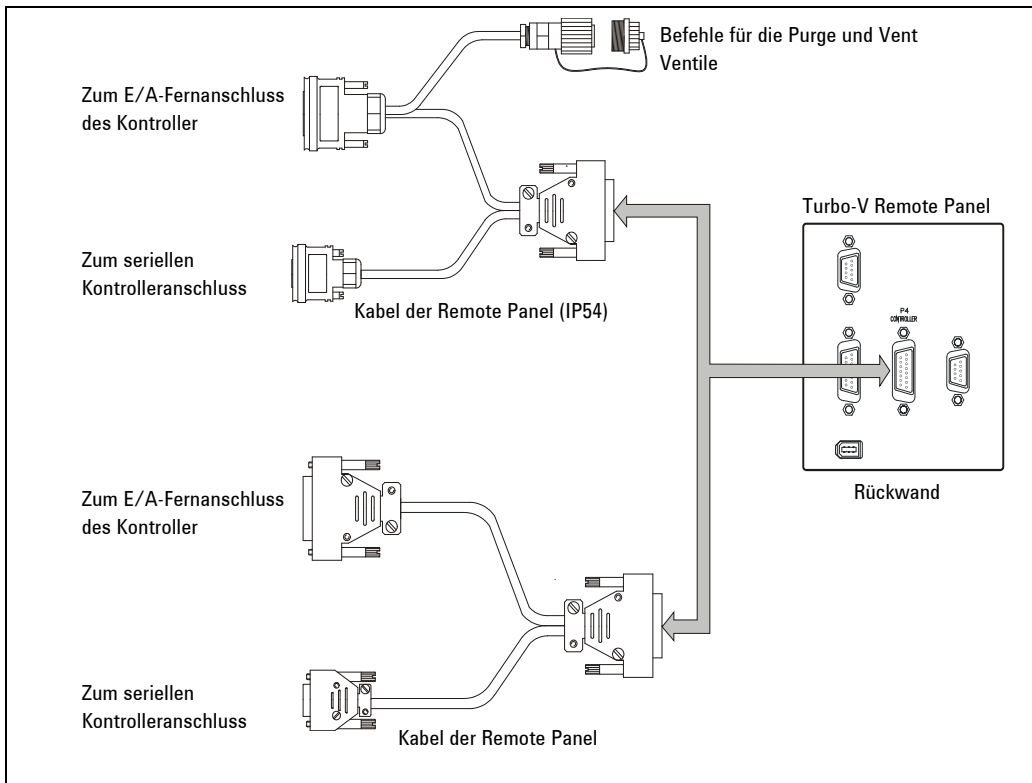


Abbildung 1

## 2 Gebrauchsanleitung

### Installation



**Abbildung 2** Anschlüsse

#### HINWEIS

Der Turbo-V Remote Panel kann als Tisch- oder Rackeinheit benutzt werden, aber sie muss so eingesetzt werden, dass die Luft frei durch die Löcher zirkulieren kann. Den Controller nicht in Umgebungen mit atmosphärischen Agenzien (Regen, Schnee, Eis), Staub, aggressiven Gasen oder explosions- und stark feuergefährdeten Umgebungen installieren oder benutzen.

Beim Betrieb müssen die folgenden Umgebungsbedingungen eingehalten werden:

- Temperatur: zwischen 5 °C und +45 °C
- rel. Luftfeuchtigkeit: 0 – 95 % (nicht kondensierend)

Siehe den Anhang " Technical Information" bezüglich den oben erwähnten und anderen Anschlüsse sowie der Installation von Optionen.

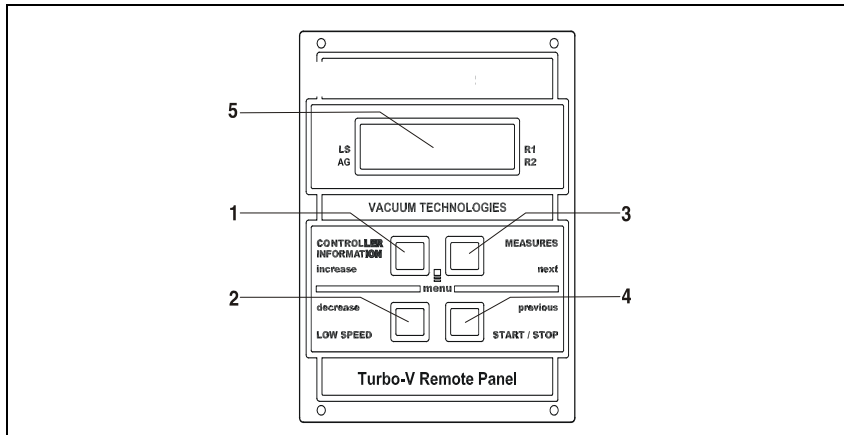
## **Anwendung**

Dieser Abschnitt beschreibt die Hauptbetriebsprozeduren. Im Einzelnen sind Informationen und Betriebsprozeduren für optionale Anschlüsse und Optionen im Abschnitt "Anwendungen" des Anhangs "Technical Information " beschrieben.

## **Steuerungen, Anzeigen und Stecker**

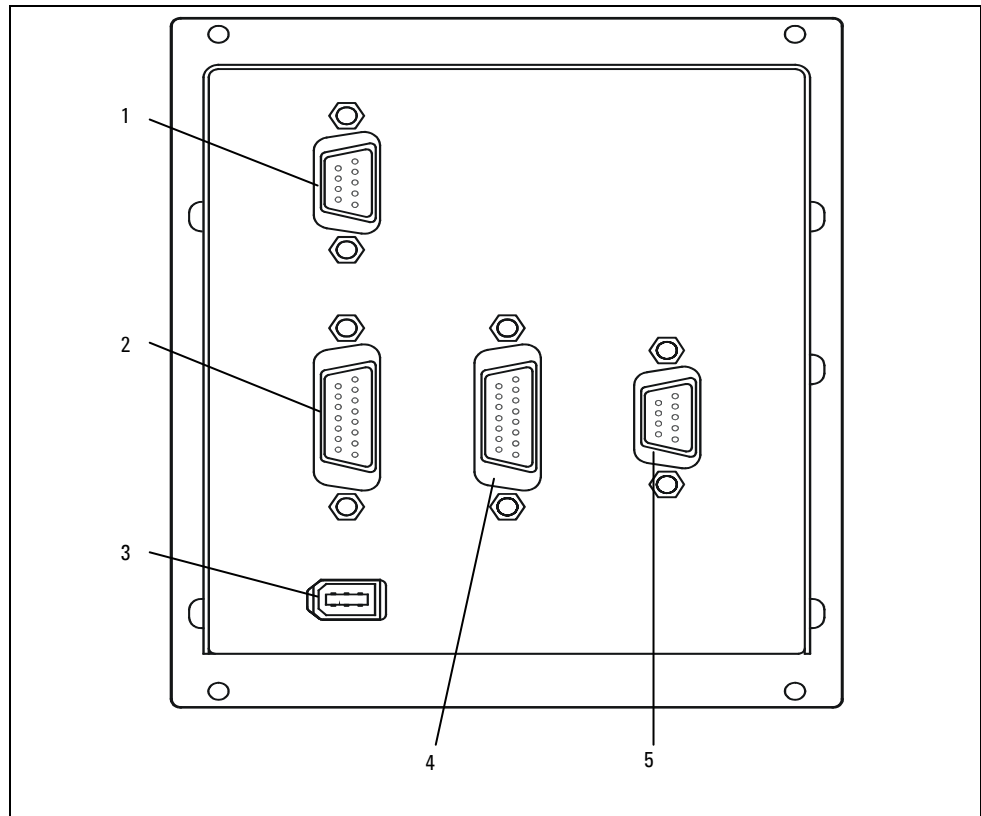
Der folgende Abschnitt beschreibt die Frontkonsole und die Rückwand des Kontroller. Weitere Einzelheiten finden sich im Anhang " Technical Information".

## Beschreibung der Frontkonsole



**Abbildung 3** Turbo-V Remote Panel

- 1 Taste zur Anzeige von Zyklusnummer, Zykluszeit und Pumpenlebensdauer auf dem Display
- 2 Taste zur Wahl des LANGSAM-MODUS. Sie ist nur bei Wahl der Frontkonsolenoperation aktiviert. Bei einmaligem Drücken läuft die Pumpe in Bereitschaftsgeschwindigkeit. Zum Abwählen des Modus die Taste wieder drücken.
- 3 Taste zur Anzeige von Pumpenstrom, Pumpentemperatur, Pumpleistung und Drehgeschwindigkeit auf dem Display. Sie ist immer unabhängig vom gewählten Betriebsmodus aktiviert. Bei gleichzeitigem Drücken der Tasten 3 und 4 für mehr als 2 Sekunden startet eine Kontrollerroutine, mit der einige Betriebsparameter programmiert werden können.
- 4 Taste zur Wahl der Modi START, STOP, RESET. Sie ist bei Wahl der Frontkonsolenoperation aktiviert. Bei einmaligem Drücken beginnt die Startphase, bei erneutem Drücken wird die Pumpe gestoppt. Wenn die Pumpe automatisch durch Fehler angehalten wurde, bewirkt einmaliges Drücken ein Reset des Controller und nochmaliges Drücken den Neustart der Pumpe.
- 5 Rückseitig beleuchtetes alphanumerisches LCD: 2 Zeilen mit je 16 Zeichen.



**Abbildung 4** Rückwand des Turbo-V Remote Panel

- 
- |   |                                                                                                                                  |
|---|----------------------------------------------------------------------------------------------------------------------------------|
| 1 | Eingangsanschlüsse der Folgesignale (der Verbindungsanschluss wird mit der dafür vorgesehenen Verschlussüberbrückung geliefert). |
|---|----------------------------------------------------------------------------------------------------------------------------------|
- 
- |   |                                                                                                              |
|---|--------------------------------------------------------------------------------------------------------------|
| 2 | Anschluss ausgehender Folgesignale und Überwachung der Pumpenfrequenz mit programmierbarem analogem Ausgang. |
|---|--------------------------------------------------------------------------------------------------------------|
- 
- |   |                                                         |
|---|---------------------------------------------------------|
| 3 | Anschluss des Versorgungsausgangs (24 Vcc pro Kaliber). |
|---|---------------------------------------------------------|
- 
- |   |                                                                                                                |
|---|----------------------------------------------------------------------------------------------------------------|
| 4 | Anschluss für Versorgungseingang vom Pumpenkontroller aus und serielle Kommunikation mit dem Pumpenkontroller. |
|---|----------------------------------------------------------------------------------------------------------------|
- 
- |   |                                        |
|---|----------------------------------------|
| 5 | Serieller Anschluss für Fernkontrolle. |
|---|----------------------------------------|
-

## **Einschalten des Controllers**

Zum Einschalten des Controllers genügt es, das Netzkabel an die Steckdose anzuschließen.

## **Pumpenstart**

Zum Starten der Pumpe den Verriegelungseingang bestätigen und die START Taste of der vorderen Konsole des Controllers drücken.

## **Pumpenstopp**

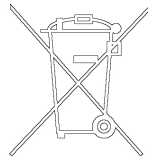
Zum Stoppen der Pumpe muß die STOPP-Taste an der Fronttafel gedrückt werden.

## Wartung

Die Controller der Serie Turbo-V 301 Rack sind wartungsfrei. Eventuell erforderliche Eingriffe müssen von dazu befugtem Fachpersonal ausgeführt werden. Bei einem Defekt kann der Agilent-Reparaturdienst bzw. der "Agilent advanced exchange service" in Anspruch genommen werden, der für die Erneuerung defekter Controller sorgt. Ein Controller muß unter Einhaltung der einschlägigen landesüblichen Vorschriften erfolgen.

## 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.



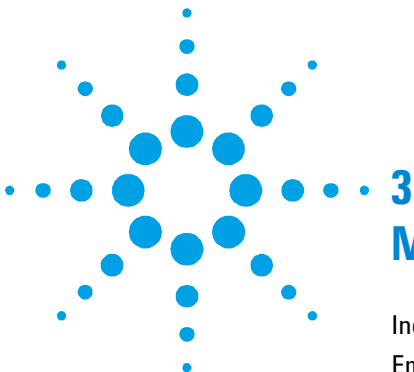
## Fehlermeldungen

In einigen Störungsfällen zeigt das Selbstdiagnosessystem des Controllers die in der nachstehenden Tabelle zusammengefaßten Meldungen an.

<b>MELDUNG</b>	<b>BESCHREIBUNG</b>	<b>BEHEBUNG</b>
COMMUNICATION ERROR	Das Remote Panel kann die Kommunikation nicht mit dem Controller starten.	Prüfen Sie den seriellen Anschluss zwischen dem Remote Panel und dem Controller.
COMMUNICATION BROKEN	Das Remote Panel kann die Kommunikation nicht mit dem Controller weiterführen.	Prüfen Sie den seriellen Anschluss zwischen dem Remote Panel und dem Controller.
XXXXXXX MODEL	Controller nicht unterstützt. Das Remote Panel funktioniert nur mit den folgenden Pumpen: TV551 mit Navigator Controller, TV701 mit Navigator Controller, TV1001 mit Navigator Controller, TV2k-G und TV3k-T.	Schließen Sie das Remote Panel an den unterstützten Controller an.
CHECK CONNECTION TO PUMP	Fehlfunktion der Pumpen-Controller Verbindung. oder Die Temperatur der Pumpe ist unter 0 °C.	Sicherstellen, daß das Verbindungskabel zwischen Pumpe und Controller an beiden Seiten korrekt befestigt ist und keine Unterbrechung vorliegt. Die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
WAITING INTERLOCK	Das Interlock-Signal auf dem Verbinder P1 ist wegen der Kurzschlußunterbrechung zwischen Pin 3 und Pin 8 des Verbinders K1 oder wegen der Öffnung des externen Interlock-Signals aktiv.	Den Kurzschluß zwischen Pin 3 und Pin 8 des Verbinders J1 rücksetzen oder das externe Interlock-Signal schließen.
PUMP OVERTEMP.	Die Temperatur des oberen Lagers bzw. der Pumpe hat 60 °C überschritten.	Warten bis die Temperatur unter den Schwellenwert gesunken ist.

<b>MELDUNG</b>	<b>BESCHREIBUNG</b>	<b>BEHEBUNG</b>
		Die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
CONTROLLER OVERTEMP.	Die Temperatur des Controller hat 70 °C überschritten. oder Die Temperatur des Controllerkühlkörpers ist über 60 °C.	Warten bis die Temperatur unter den Schwellenwert gesunken ist. Die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
TOO HIGH LOAD	Die Pumpe ist stromaufnahme größer als die vorgesehene.	Sicherstellen, daß der Pumpenrotor ungehindert drehen kann. Die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
SHORT CIRCUIT	Während des Normalbetriebs (nach der Startphase) erfolgt ein Kurzschluß der Ausgangsverbinding.	Die Verbindung zwischen Pumpe und Controller prüfen. Die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
SYSTEM OVERRIDE	Die Pumpe wurde durch ein von einem entfernten Kontakt kommendes Notsignal gestoppt.	Das Netzkabel des Controllers ausstecken und die Störungsursache beheben. Das Netzkabel wieder anschließen und die Pumpe durch zweimalige Betätigung der START-Taste neustarten.
OVERVOLTAGE	Defekt im Versorgungsbereich des Controllers bzw. der Controller hat ein falsches Signal erhalten	Die Pumpe durch zweimalige Betätigung der START-Taste neustarten. Erscheint die Meldung wieder sollte der Agilent-Wartungs-dienst gerufen werden.
POWER FAIL	Defekt im Pumpenstromkreis des Controllers.	Agilent für Wartung kontaktieren.

## **2** **Gebrauchsanleitung** **Fehlermeldungen**



### 3

## Mode d'emploi

Indications generales	42
Emmagasinage	44
Preparation pour l'installation	44
Installation	45
Utilisation	47
Commandes, indicateurs et connecteurs	47
Description du tableau avant	48
Allumage du Contrôleur	50
Mise en marche de la Pompe	50
Arrêt de la Pompe	50
Entretien	51
Mise au rebut	51
Messages d'erreur	52

Traduction de la mode d'emploi originale



## Indications générales

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'appareil. Agilent décline par conséquent 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.

Le Turbo-V Remote Panel est contrôlé par un microprocesseur et permet à l'utilisateur de commander une pompe du type "contrôleur embarqué" via un tableau avant :

TV551 avec Navigator controller  
TV701 avec Navigator controller  
TV1001 avec Navigator controller  
TV2k-G avec controller on-board  
TV3k-T avec controller on-board  
TV3k-G avec controller on-board  
TV70 avec Navigator controller  
TV81 avec Navigator controller  
TV301 avec Navigator controller  
TV303 avec Navigator controller

Les paragraphes suivants donnent toutes les indications nécessaires à garantir la sécurité de l'opérateur pendant l'utilisation de l'appareillage. Des renseignements plus détaillés se trouvent dans l'appendice "Technical Information".

Cette notice utilise les signes conventionnels suivants :

**ATTENTION!**

Les messages d'attention apparaissent avant certaines procédures qui, si elles ne sont pas observées, pourraient endommager sérieusement l'appareillage.

---

**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.

---

**NOTE**

Les notes contiennent des renseignements importants, isolés du texte.

---

## Emmagasinage

Pendant le transport et l'emmagasinage des contrôleurs, 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 Turbo-V Remote Panel est fourni dans un emballage de protection spécial; si l'on constate des marques de dommages pouvant s'être produits pendant le transport, contacter aussitôt le bureau de vente local. Pendant l'opération d'ouverture de l'emballage, veiller tout particulièrement à ne pas laisser tomber le contrôleur et à ne lui faire subir aucun choc. Ne pas jeter l'emballage dans la nature. Le matériel est entièrement recyclable et il est conforme aux directives CEE 83/399 en matière de protection de l'environnement.

## Installation

Le Turbo-V Remote Panel doit être connecté au contrôleur par le câble prévu à cet effet (voir tableau des pièces que l'on peut commander), comme illustré sur la figure suivante. La puissance/les signaux nécessaires étant fournis par ce câble, il n'y a besoin d'aucune autre alimentation extérieure.

Si l'opérateur souhaite commander les valves évent/purgeur via les commandes extérieures, utiliser la version avec câble IP54, régler le contrôleur sur "valve setting" "on command" et connecter les câbles comme illustré sur la figure. Les valves sont du type normalement fermées; si on fournit du 24 V entre les valves évent/purgeur et le terminal commun, les valves s'ouvriront.

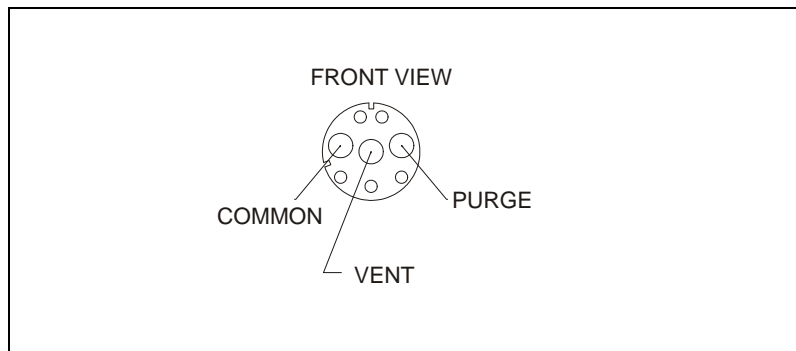
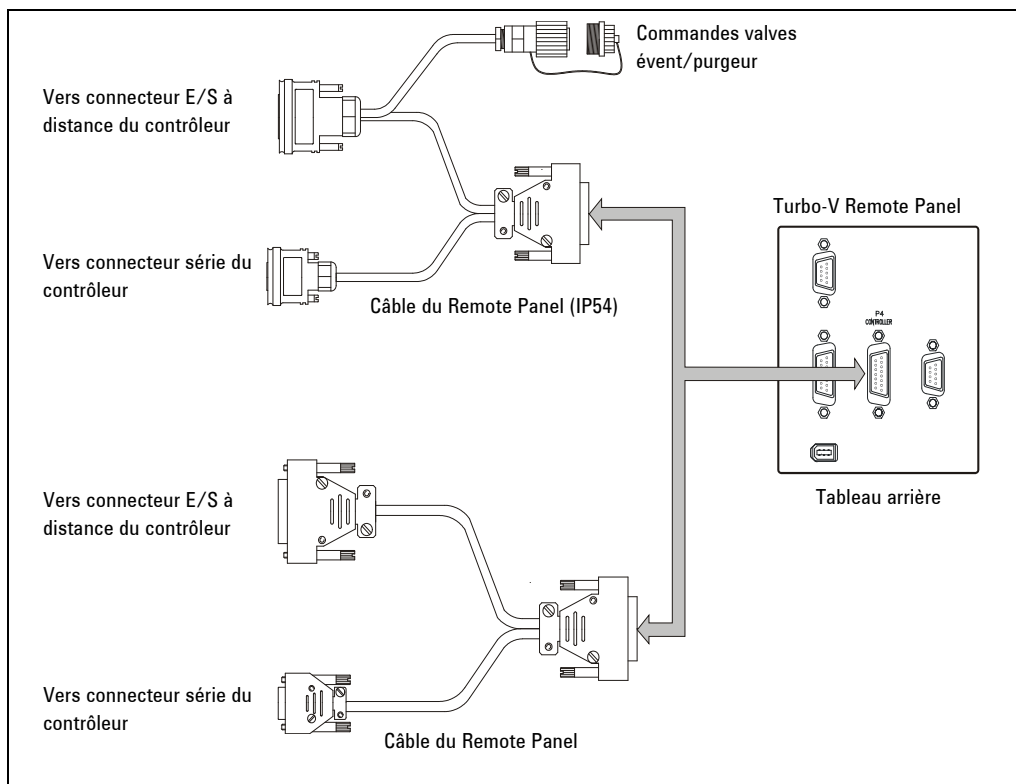


Figure 1

### 3 Mode d'emploi

#### Installation



**Figure 2** Connexions

#### NOTE

Le Turbo-V Remote Panel peut être installé sur un plateau d'appui ou à l'intérieur d'un rack prévu à cet effet. Il est en tout cas nécessaire que l'air de refroidissement puisse circuler librement à l'intérieur de l'appareil. Ne pas installer et/ou utiliser le contrôleur dans des milieux exposés à des agents atmosphériques (pluie, gel, neige), à des poussières, à des gaz de combat ainsi que dans des milieux explosifs ou à risque élevé d'incendie.

Pendant le fonctionnement, il est nécessaire de respecter les conditions environnementales suivantes :

- température : de 5 °C à +45 °C
- humidité relative : 0 – 95 % (non condensante).

Pour les autres connexions et pour l'installation des accessoires en option, voir la section "Technical Information".

## Utilisation

Dans ce paragraphe, on indique les principales procédures opérationnelles. Pour tous autres détails et pour les procédures concernant des connexions ou des éléments en option, se reporter au paragraphe "Use" de l'appendice "Technical Information".

## Commandes, indicateurs et connecteurs

On présente ci-dessous le tableau avant et le tableau arrière du contrôleur. Pour de plus amples détails, se reporter à la section "Technical Information".

## Description du tableau avant

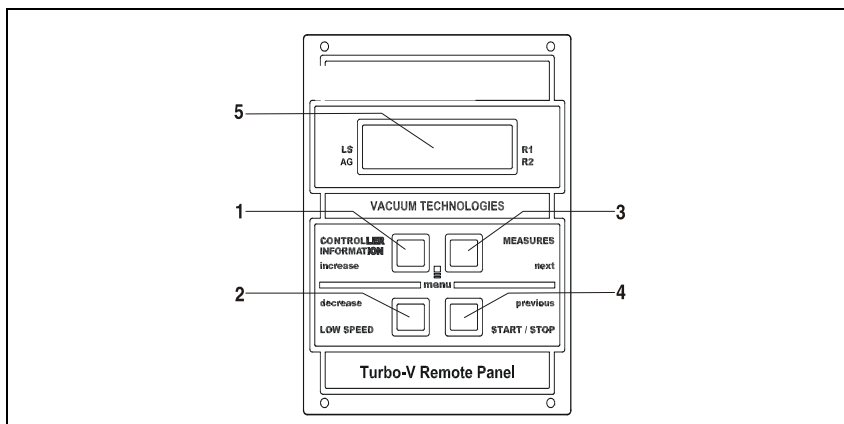
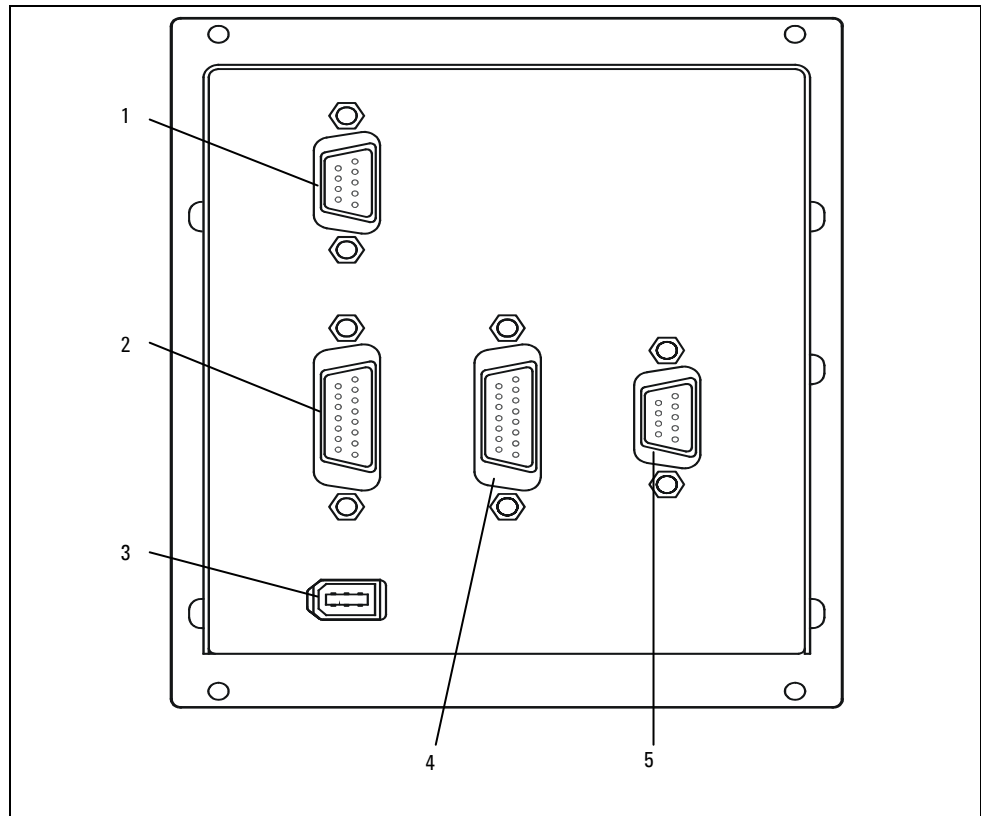


Figure 3 Turbo-V Remote Panel

- 1 Interrupteur rappelant sur l'afficheur les paramètres de cycle number, cycle time et pump life.
- 2 Interrupteur de sélection du mode LOW SPEED. Il n'est actif que lorsque le mode de commande est sélectionné depuis le tableau frontal. En le pressant une fois, la pompe tourne à vitesse "stand-by". En le pressant une deuxième fois, on désactive le mode LOW SPEED.
- 3 Interrupteur rappelant sur l'afficheur les paramètres de pump current, pump temperature, pump power et rotational speed. Il est toujours actif, indépendamment du mode de fonctionnement choisi. En pressant simultanément les interrupteurs 3 et 4 pendant 2 secondes au moins, on active un programme avec lequel il est possible de programmer certains paramètres opérationnels.
- 4 Interrupteur envoyant les commandes de START, STOP/RESET. Il n'est actif que lorsque le mode de commande est sélectionné depuis le tableau frontal. Une première pression de l'interrupteur active la phase de mise en marche; une deuxième pression provoque l'arrêt de la pompe. Si la pompe s'est arrêtée automatiquement à cause d'une panne, il faut presser cet interrupteur une première fois pour effectuer la mise à zéro du contrôleur et une deuxième fois pour remettre la pompe en marche.
- 5 Ecran alphanumérique à cristaux liquides: matrice de points, 2 lignes x 16 caractères.



**Figure 4** Panneau postérieur du Turbo-V Remote Panel

- |   |                                                                                                                      |
|---|----------------------------------------------------------------------------------------------------------------------|
| 1 | Connecteurs d'entrée des signaux logiques (le connecteur d'accouplement est fourni avec le pontet de fermeture).     |
| 2 | Connecteur signaux logiques en sortie et monitoring fréquence pompe plus sortie analogique programmable.             |
| 3 | Connecteur de sortie alimentation (24 Vcc par calibre).                                                              |
| 4 | Connecteur pour entrée alimentation du contrôleur de la pompe et communication série avec le contrôleur de la pompe. |
| 5 | Connecteur sériel pour contrôle à distance.                                                                          |

## **Allumage du Contrôleur**

Pour allumer le contrôleur, il suffit d'introduire le câble d'alimentation dans la prise du réseau.

## **Mise en marche de la Pompe**

Pour faire démarrer la pompe il faut habilitier l'entrée de interlock et appuyer sur la touche START du panneau avant.

## **Arrêt de la Pompe**

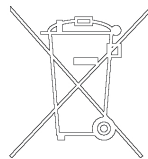
Pour arrêter la pompe, presser l'interrupteur STOP du tableau avant.

## Entretien

Les contrôleurs de la série Turbo-V 301 Rack n'exigent aucun entretien. Toute opération doit être effectuée par un personnel agréé. En cas de panne, il est possible de s'adresser au Service de réparation Agilent ou bien au "Agilent advance exchange service" qui permet d'obtenir un contrôleur régénéré à la place du contrôleur détraqué. En cas de mise au rebut du contrôleur, procéder à son élimination conformément aux réglementations nationales en la matière.

## Mise au rebut

**Signification du logo "WEEE" figurant 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

Dans certains cas de panne, l'ensemble de circuits d'autodiagnostic du contrôleur présente certains messages d'erreur indiqués dans le tableau ci-dessous.

MESSAGE	DESCRIPTION	INTERVENTION
COMMUNICATION ERROR	Le Remote Panel n'arrive pas à démarrer la communication avec le contrôleur.	Vérifier la connexion série entre le Remote Panel et le contrôleur.
COMMUNICATION BROKEN	Le Remote Panel n'arrive pas à continuer la communication avec le contrôleur.	Vérifier la connexion série entre le Remote Panel et le contrôleur.
XXXXXXX MODEL	Contrôleur sans support. le Remote Panel fonctionne uniquement avec les pompes suivantes: TV551 avec navigator controller, TV701 avec navigator controller, TV1001 avec navigator controller, TV2k-G et TV3k-T.	Connecter le Remote Panel au contrôleur de support .
CHECK CONNECTION TO PUMP	Dysfonctionnement de la connexion entre la pompe et le contrôleur. Ou bien La pompe a une température inférieure à 0 °C.	S'assurer que le câble de connexion entre la pompe et le contrôleur et le contrôleur est bien fixé aux deux extrémités et qu'il n'est pas coupé. Presser deux fois l'interrupteur START pour réactiver la pompe.
WAITING INTERLOCK	Le signal d'interlock situé sur le connecteur P1 est actif à cause de la coupure du court-circuit entre le pin 3 et le pin 8 du connecteur J1 ou à cause de l'ouverture du signal d'interlock extérieur.	Rétablir le court-circuit entre le pin 3 et le pin 8 du connecteur J1 ou fermer le signal d'interlock extérieur.
PUMP OVERTEMP.	La température du palier supérieur ou de la a dépassé 60 °C.	Attendre que la température retourne au-dessous du seuil. Presser deux fois l'interrupteur START pour remettre la pompe

MESSAGE	DESCRIPTION	INTERVENTION
CONTROLLER OVERTEMP.	La température du contrôleur a dépassé 70 °C. Ou bien La température du radiateur du contrôleur est supérieure à 60 °C.	en marche.  Attendre que la température retourne au-dessous du seuil. Presser deux fois l'interrupteur START pour remettre la pompe en marche.
TOO HIGH LOAD	Pendant le fonctionnement normal (après la phase de mise en marche), le courant absorbé par la pompe est plus grand que celui qui a été programmé.	S'assurer que le rotor de la pompe a la possibilité de tourner librement. Presser deux fois l'interrupteur START pour remettre la pompe en marche.
SHORT CIRCUIT	Pendant le fonctionnement normal (après la phase de mise en marche), la connexion de sortie est en court-circuit.	Vérifier les connexions entre la pompe et le contrôleur. Presser deux fois l'interrupteur START pour remettre la pompe en marche.
SYSTEM OVERRIDE	La pompe a été arrêtée par un signal d'alerte provenant d'un contact éloigné.	Débrancher le câble d'alimentation du contrôleur et corriger la cause de l'alerte. Reconnecter le câble d'alimentation et presser deux fois l'interrupteur START pour remettre la pompe en marche.
OVERVOLTAGE	Il s'est produit une panne de la section d'alimentation du contrôleur, ou bien le contrôleur a reçu un faux signal.	Presser deux fois l'interrupteur START pour remettre la pompe en marche.  Si le message se présente à nouveau, s'adresser à Agilent pour l'entretien.
POWER FAIL	Panne dans la section d'alimentation de la pompe, du contrôleur.	S'adresser à Agilent pour Maintenance.

### **3 Mode d'emploi**

#### **Messages d'erreur**



## 4 Manual de instrucciones

Información general	56
Almacenamiento	58
Preparación para la instalación	58
Instalación	59
Uso	61
Mandos, indicadores y conectores	61
Descripción del panel frontal	62
Encendido del controler	64
Puesta en marcha de la Bomba	64
Parada de la Bomba	64
Mantenimiento	65
Eliminación	65
Mensajes de error	66

Traducción de las instrucciones originales



## Información general

Este equipo se ha concebido para un uso profesional. El usuario deberá leer atentamente el presente manual de instrucciones y cualquier otra información suplementaria facilitada por Agilent antes de utilizar el equipo. Agilent se considera libre de cualquier responsabilidad debida al incumplimiento total o parcial de las instrucciones, al uso poco apropiado por parte de personal sin formación, a las operaciones no autorizadas o al uso que no cumpla con las normas nacionales específicas.

El Turbo-V Remote Panel está controlado por un microprocesador y permite al usuario dirigir una bomba de tipo “controller on-board” mediante un panel frontal:

TV551 con Navigator controller

TV701 con Navigator controller

TV1001 con Navigator controller

TV2k-G con controller on-board

TV3k-T con controller on-board

TV3k-G con controller on-board

TV70 con Navigator controller

TV81 con Navigator controller

TV301 con Navigator controller

TV303 con Navigator controller

En los apartados siguientes se ha incluido toda la información necesaria para garantizar la seguridad del operador durante la utilización del equipo. Si se desea información más detallada, se aconseja consultar el anexo "Technical Information".

**Este manual utiliza los siguientes símbolos convencionales:**

**¡ATENCIÓN!**

Los mensajes de atención se visualizan antes de procedimientos que, de no respetarse, podrían provocar daños al equipo.

---

**¡ADVERTENCIA!**



Los mensajes de advertencia atraen la atención del operador sobre un procedimiento o una operación específica que, de no realizarse correctamente, podría provocar graves lesiones personales.

---

**NOTA**

Las notas contienen información importante extraída del texto.

---

## **Almacenamiento**

Durante el transporte y el almacenamiento de los controlers se deberá cumplir con las condiciones ambientales siguientes:

- temperatura: de -20 °C a +70 °C
- humedad relativa: 0 – 95 % (no condensadora)

## **Preparación para la instalación**

El Turbo-V Remote Panel se suministra en un embalaje de protección especial; si se observan señales de daños, que podrían haberse producido durante el transporte, ponerse en contacto con la oficina de venta más cercana. Durante la operación de desembalaje, prestar una atención especial a no dejar caer el controler y evitarle golpes. No dispersar el embalaje en el medio ambiente. El material es totalmente reciclable y cumple con la directiva CEE 85/399 para la preservación del medio ambiente.

## Instalación

El Turbo-V Remote Panel debe conectarse al controler mediante el cable específico (ver tabla de las componentes que pueden ser ordenados), tal como se indica en la figura siguiente. Ya que la potencia y las señales requeridas son suministradas por este cable, no es necesaria ninguna otra alimentación externa.

Si el operador desea dirigir las válvulas de venteo y purga utilizando los mandos externos, será necesario utilizar la versión con cable IP54, programar el controler en "valve setting" "on command" y conectar los cables tal como indica la figura. Las válvulas son n.c. (normalmente cerradas); suministrando 24 V entre la válvula de venteo y purga y el terminal común, las válvulas se abrirán.

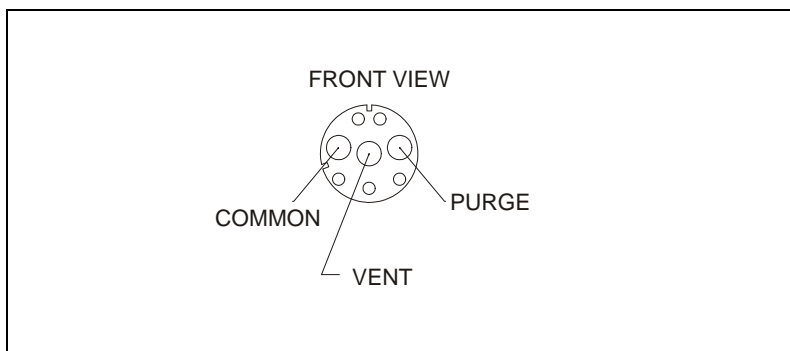
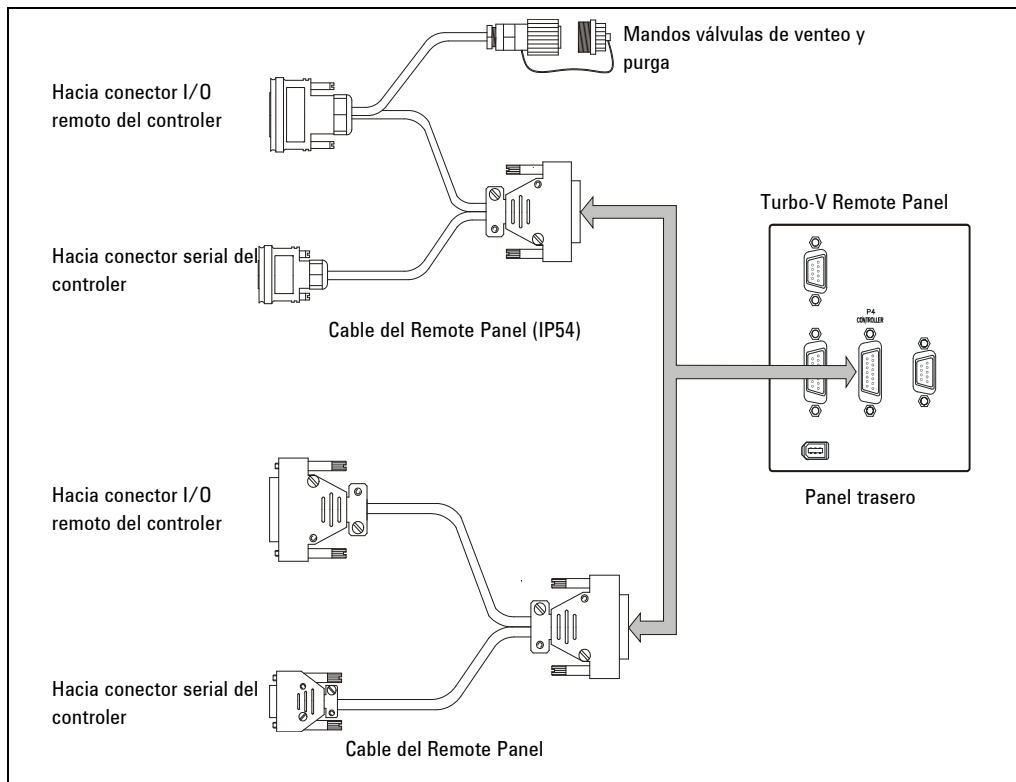


Figura 1

## 4 Manual de instrucciones

### Instalación



**Figura 2** Conexiones

#### NOTA

El Turbo-V Remote Panel puede ser instalado en una mesa o dentro de un Rack específico. En cualquier caso, es necesario que el aire de refrigeración pueda circular libremente alrededor del aparato. No instalar y/o utilizar el controler en ambientes expuestos a agentes atmosféricos (lluvia, hielo, nieve), polvo, gases agresivos, en ambientes explosivos o con alto riesgo de incendio .

Durante el funcionamiento es necesario que se respeten las siguientes condiciones ambientales:

- temperatura: de 5 °C a +45 °C
- humedad relativa: 0 – 95 % (no condensante).

Para otras conexiones y la instalación de los accesorios opcionales, véase la sección “Technical Information”.

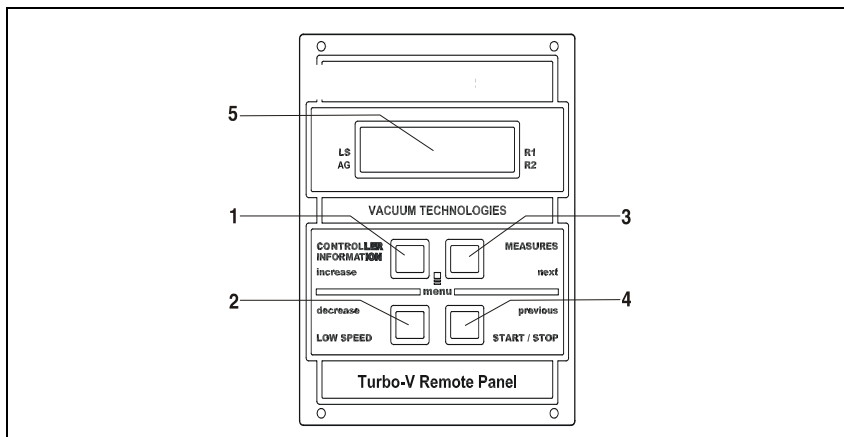
## Uso

En este apartado se citan los procedimientos operativos principales. Para más detalles y para procedimientos que impliquen conexiones u opcionales especiales, les remitimos al apartado “Use” del anexo “Technical Information”.

## Mandos, indicadores y conectores

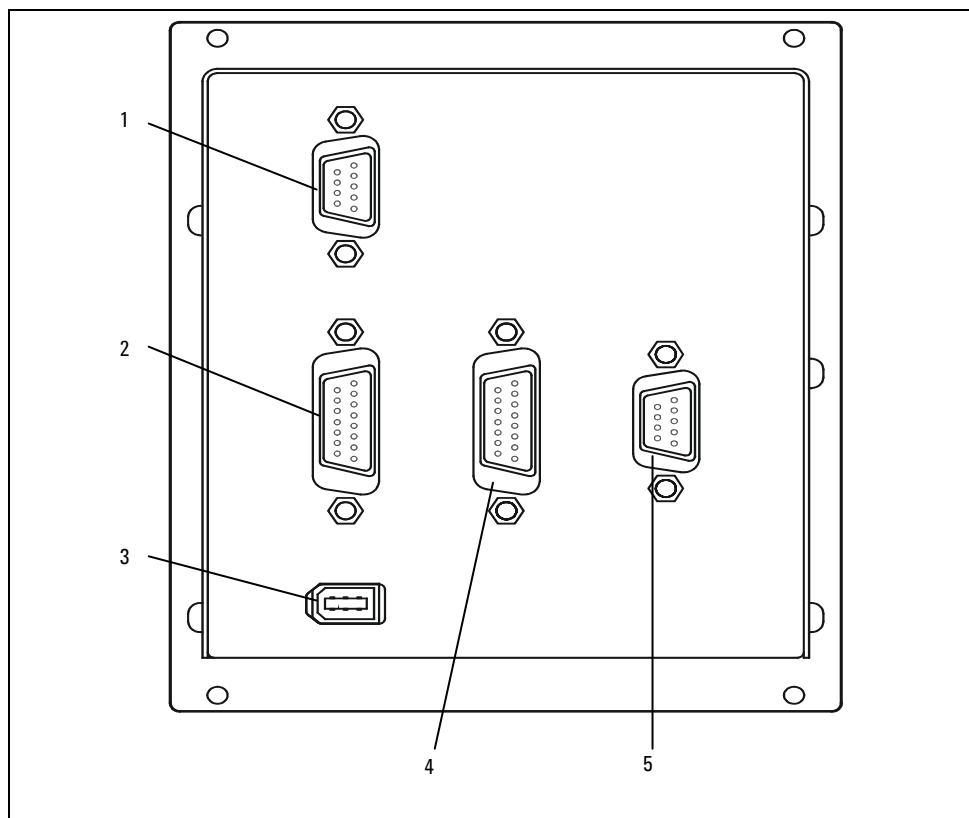
A continuación se ha ilustrado el panel frontal y el panel trasero del controler. Para más detalles consultar la sección “Technical Information”.

## Descripción del panel frontal



**Figura 3** Turbo-V Remote Panel

- 1 Pulsador para que aparezcan en el display los parámetros cycle number, cycle time y pump life.
- 2 Pulsador para la selección del modo LOW SPEED. Se activa sólo cuando se ha seleccionado el modo de mando del panel frontal. Apretándolo una vez, la bomba gira a velocidad "stand-by". Apretándolo otra vez se desactiva el modo LOW SPEED.
- 3 Pulsador para que aparezcan en el display los parámetros pump current, pump temperature, pump power y rotational speed. Está siempre activo independientemente del modo de funcionamiento seleccionado. Apretándolo junto con los pulsadores 3 y 4 durante 2 segundos, como mínimo, se activa un programa mediante el que es posible configurar algunos parámetros operativos.
- 4 Pulsador para enviar los mandos de START, STOP/RESET. Esta activo sólo cuando se ha seleccionado el modo de mando del panel frontal. Apretándolo una vez se activa la fase de puesta en marcha; apretándolo otra vez se para la bomba. Si la bomba se ha parado automáticamente a causa de una avería, será necesario apretar este pulsador primero una vez para efectuar el reset del controler y la segunda vez para volver a poner en marcha la bomba.
- 5 Display alfanumérico de cristales líquidos: matriz de puntos, 2 líneas x 16 caracteres.



**Figura 4** Panel posterior del Turbo-V Remote Panel

- 1 Conector de señales lógicas en entrada (el conector de acoplamiento se suministra con el correspondiente puente de cierre).
- 2 Conector de señales lógicas en salida y monitoreo de frecuencia bomba con salida analógica programable.
- 3 Conector de salida alimentación (24 Vcc para calibre).
- 4 Conector para entrada alimentación del controlador bomba y comunicación serial con controlador bomba.
- 5 Conector serial para control remoto.

## **Encendido del controler**

Para encender el controler es suficiente introducir el cable de alimentación en la toma de red.

## **Puesta en marcha de la Bomba**

Para activar la bomba es necesario habilitar la entrada de interlock y pulsar el botón START situado en el panel frontal.

## **Parada de la Bomba**

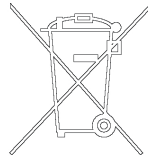
Para detener la bomba hay que apretar el pulsador STOP del panel frontal.

## Mantenimiento

Los controlers de la serie Turbo-V 301 Rack no necesitan ningún mantenimiento. Cualquier operación ha de ser efectuada por personal autorizado. En caso de avería es posible utilizar el servicio de reparación Agilent o del "Agilent advance exchange service", que permite obtener un controler regenerado en vez del averiado. En caso de que un controler se tenga que desguazar, efectuar su eliminación respetando las normas nacionales específicas.

## Eliminación

**Significado del logotipo "WEEE" presente en las etiquetas.** El símbolo que se indica a continuación, es aplicado en observancia de la directiva CE denominada "WEEE". Este símbolo (**válido sólo para los países miembros de la Comunidad Europea**) indica que el producto sobre el cual ha sido aplicado, NO debe ser eliminado junto con los residuos comunes sean éstos domésticos o industriales, y que, por el contrario, deberá ser sometido a un procedimiento de recogida diferenciada. Por lo tanto, se invita al usuario final, a ponerse en contacto con el proveedor del dispositivo, tanto si éste es la casa fabricante o un distribuidor, para poder proveer a la recogida y eliminación del producto, después de haber efectuado una verificación de los términos y condiciones contractuales de venta.



## Mensajes de error

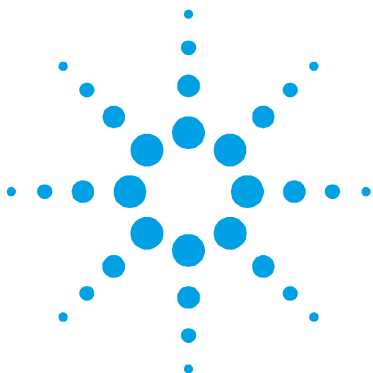
En algunos casos de avería los circuitos de autodiagnos del controler presenta algunos mensajes de error detallados en la tabla siguiente.

MENSAJE	DESCRIPCIÓN	ACCIÓN CORRECTIVA
COMMUNICATION ERROR	El Remote Panel no es capaz de establecer comunicación con el controlador.	Comprobar la conexión serial entre el Remote Panel y el controlador.
COMMUNICATION BROKEN	El Remote Panel no es capaz de continuar la comunicación con el controlador.	Comprobar la conexión serial entre el Remote Panel y el controlador.
XXXXXXX MODEL	Controlador no compatible. El Remote Panel funciona sólo con las siguientes bombas: TV551 con navigator controller, TV701 con navigator controller, TV1001 con navigator controller, TV2k-G y TV3k-T.	Conectar el Remote Panel al controlador compatible.
CHECK CONNECTION TO PUMP	Mal funcionamiento en la conexión entre la bomba y el Controler. O bien La bomba tiene una temperatura inferior a 0 °C.	Comprobar que el cable de conexión entra en la bomba y el controler está bien fijado por ambos extremos y no está interrumpido. Apretar dos veces el pulsador START para volver a poner en marcha la bomba.
WAITING INTERLOCK	Está activa la señal de interlock presente en el conector P1 a causa de la interrupción del cortocircuito entre el pin 3 y el pin 8 del conector J1, o a causa de la apertura de la señal de interlock externo.	Eliminar el cortocircuito entre el pin 3 y el pin 8 del conector J1, o cerrar la señal de interlock exterior.
PUMP OVERTEMP.	La temperatura del rodamiento superior o de la bomba ha superado los 60 °C.	Esperar a que la temperatura vuelva por debajo del umbral. Apretar dos veces el pulsador

MENSAJE	DESCRIPCIÓN	ACCIÓN CORRECTIVA
		START para volver a poner en marcha la bomba.
CONTROLLER OVERTEMP.	La temperatura del controler ha superado los 70 °C. O bien La temperatura del radiador del controler es superior a 60 °C.	Esperar a que la temperatura vuelva por debajo del umbral. Apertar dos veces el pulsador START para volver a poner en marcha la bomba.
TOO HIGH LOAD	La corriente absorbida por la bomba es superior a la programada.	Comprobar que el rotor de la bomba tiene la posibilidad de girar libremente. Apertar dos veces el pulsador START para volver a poner en marcha la bomba.
SHORT CIRCUIT	Durante el funcionamiento normal (tras la fase de puesta en marcha) la conexión de salida está en cortocircuito.	Comprobar las conexiones entre la bomba y el controler. Apertar dos veces el pulsador START para volver a poner en marcha la bomba.
SYSTEM OVERRIDE	La bomba ha sido parada por una señal de emergencia procedente de un contacto remoto.	Desenchufar el cable de alimentación del controler y corregir la causa de la emergencia. Volver a conectar el cable de alimentación y apertar dos veces el pulsador START para volver a poner en marcha la bomba.
OVERVOLTAGE	Se ha producido una avería en la sección de alimentación del controler o el controler ha recibido una señal espurio.	Apertar dos veces el pulsador START para volver a poner en marcha la bomba. Si el mensaje se vuelve a presentar dirigirse a Agilent para el mantenimiento.
POWER FAIL	Avería en la sección de alimentación de la bomba, del controler.	Ponerse en contacto con Agilent para el Mantenimiento.

## **4 Manual de instrucciones**

### **Mensajes de error**



## 5

# Manual de Instruções

Informações gerais	70
Armazenagem	72
Preparação para a instalação	72
Instalação	73
Utilização	75
Comandos, Indicadores e Conectores	75
Descrição do painel frontal	76
Acendimento do Controller	78
Activação da bomba	78
Paragem da bomba	78
Manutenção	79
Eliminação	79
Mensagens de erro	80

Tradução das instruções originais



## Informações gerais

Esta aparelhagem destina-se ao uso profissional. O utilizador deve ler atentamente o presente manual de instruções e todas as informações adicionais fornecidas pela Agilent antes de utilizar a aparelhagem. A Agilent não se responsabiliza pela inobservância total ou parcial das instruções, pelo uso indevido por parte de pessoas não treinadas, por operações não autorizadas ou pelo uso contrário às normas nacionais específicas.

O Turbo-V Remote Panel é controlado por um microprocessador e permite que o utilizador comande uma bomba de tipo “controller on-board” mediante um painel frontal:

- TV551 com Navigator controller
- TV701 com Navigator controller
- TV1001 com Navigator controller
- TV2k-G com controller on-board
- TV3k-T com controller on-board
- TV3k-G com controller on-board
- TV70 com Navigator controller
- TV81 com Navigator controller
- TV301 com Navigator controller
- TV303 com Navigator controller

Nos parágrafos seguintes, estão descritas todas as informações necessárias para garantir a segurança do operador durante o uso da aparelhagem. Informações pormenorizadas são fornecidas no apêndice "Technical Information".

**Este manual utiliza as seguintes convenções:**

**CUIDADO!**

As mensagens de cuidado são visualizadas antes de procedimentos que, se não observados, podem causar danos à aparelhagem.

---

**ATENÇÃO!**



As mensagens de atenção chamam a atenção do operador para um procedimento ou uma prática específica que, se não efectuada correctamente, pode provocar graves lesões pessoais.

---

**NOTA**

As notas contêm informações importantes destacadas do texto.

---

## **Armazenagem**

Durante o transporte e a armazenagem dos controllers, devem ser satisfeitas as seguintes condições ambientais:

- temperatura: da -20 °C a +70 °C
- humidade relativa: 0 – 95 % (não condensante)

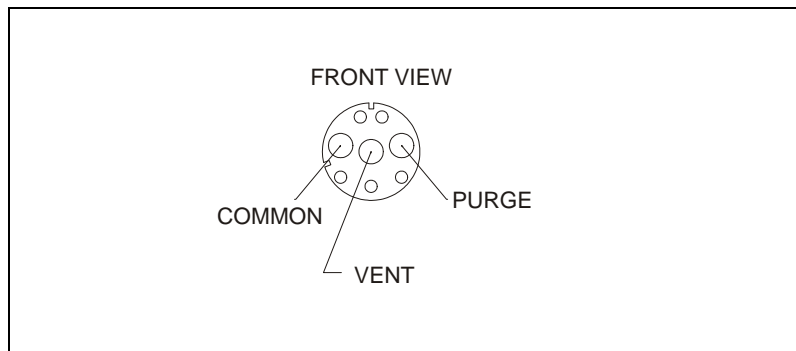
## **Preparação para a instalação**

O Turbo-V Remote Panel é fornecido numa embalagem protectora especial; se aparecerem sinais de danos, que podem ter ocorrido durante o transporte, entrar em contacto com o escritório de vendas local. Durante a retirada da embalagem, tomar muito cuidado para não deixar cair o controller e para não o submeter a choques. Não deitar a embalagem no meio ambiente. O material é completamente reciclável e corresponde à directriz CEE 85/399 para a protecção do meio ambiente.

## Instalação

O Turbo-V Remote Panel deve ser ligado ao controller mediante o respectivo cabo (ver tabela das peças a encomendar), como ilustrado na seguinte figura. Dado que a potência/os sinais necessários são fornecidos por este cabo, não é necessário nenhuma outra alimentação externa.

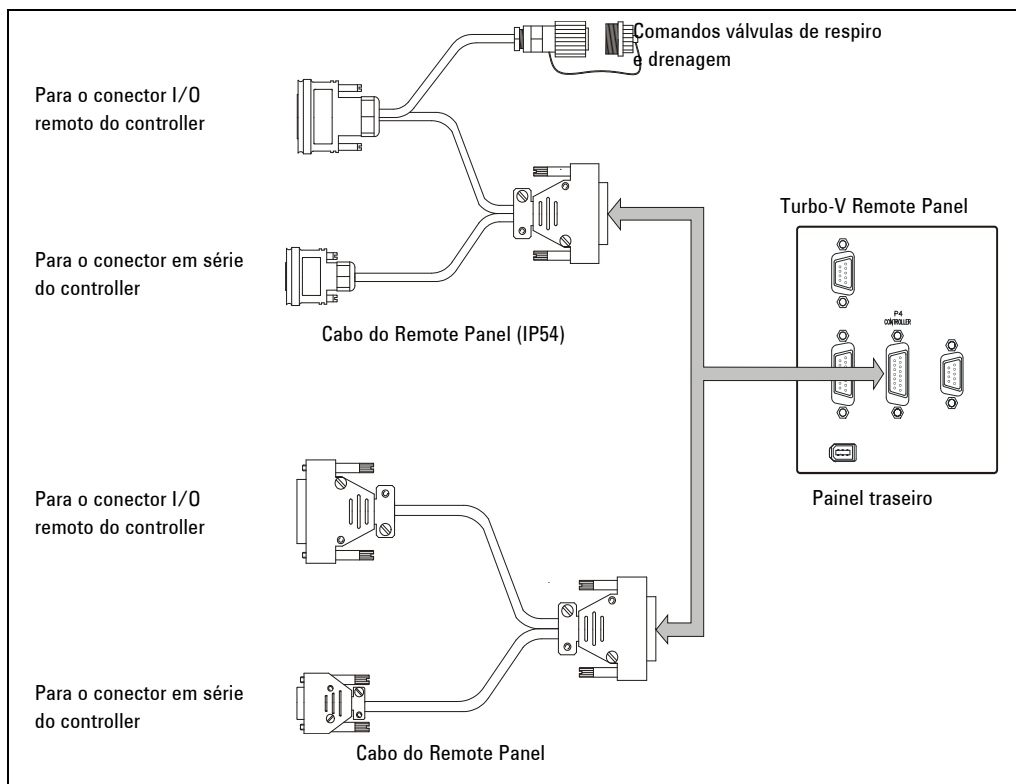
Se o operador desejar comandar as válvulas de respiro e drenagem utilizando os comandos externos, utilizar a versão com cabo IP54, definir o controller em "valve setting" "on command" e ligar os cabos como ilustrado na figura. As válvulas são n.c.; fornecendo 24 V entre as válvulas de respiro e drenagem e o terminal comum, as válvulas abrir-se-ão.



**Figura 1**

## 5 Manual de Instruções

### Instalação



**Figura 2** Ligações

#### NOTA

O Turbo-V Remote Panel pode ser instalado numa mesa ou no interior de um rack específico. Em todo o caso, é necessário que o ar de refrigeração possa circular livremente ao redor da aparelhagem. Não instalar e/ou utilizar o controller em ambientes expostos a agentes atmosféricos (chuva, gelo, neve), pós, gases agressivos ou em ambientes explosivos ou com elevado risco de incêndios.

Durante o funcionamento, é necessário que sejam respeitadas as seguintes condições ambientais:

- temperatura: da 5 °C a +45 °C
- humidade relativa: 0 – 95 % (não condensante).

Para as outras ligações e a instalação dos acessórios opcionais, ver a secção "Technical Information".

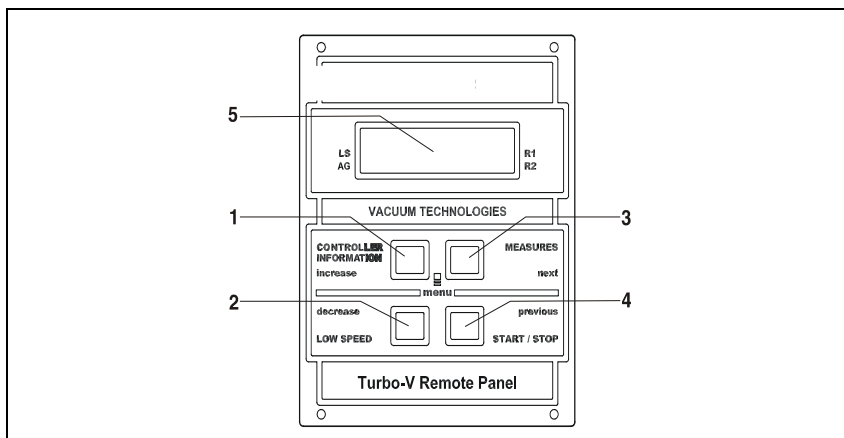
## Utilização

Neste parágrafo são descritos os principais procedimentos operativos. Para maiores detalhes e para procedimentos que envolvem ligações ou peças opcionais, consultar o parágrafo "Use" do apêndice "Technical Information".

## Comandos, Indicadores e Conectores

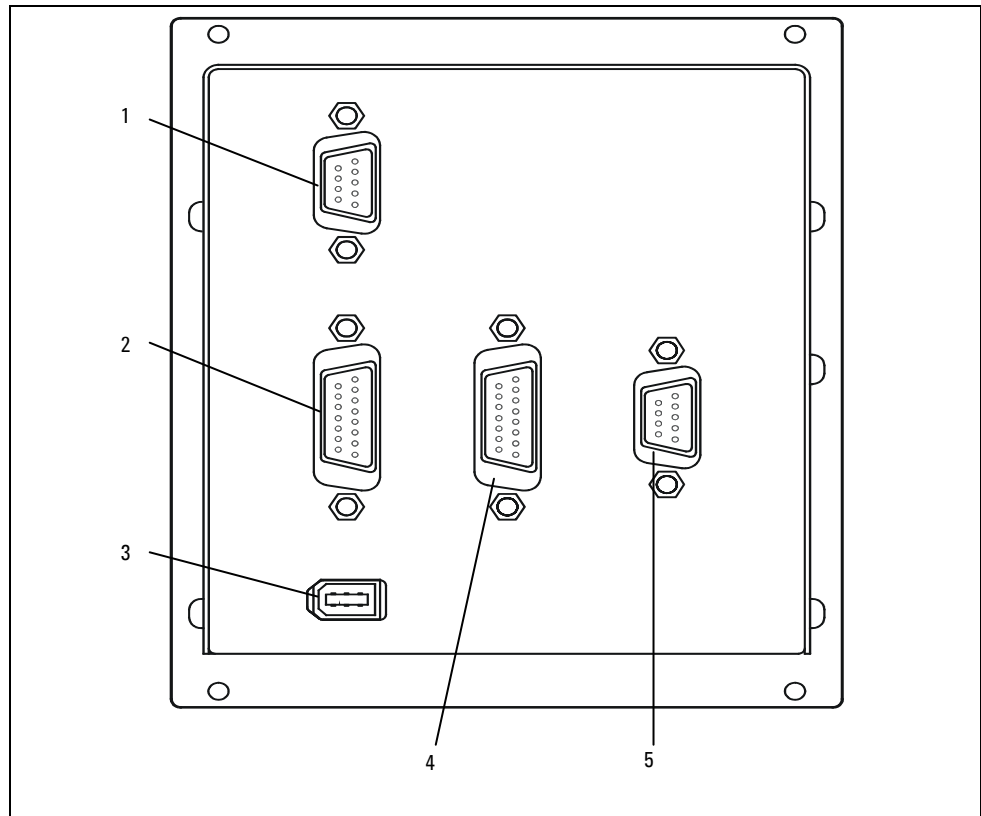
A seguir, são ilustrados os painéis frontal e posterior do controller. Para maiores detalhes, consultar a secção "Technical Information".

## Descrição do painel frontal



**Figura 3** Turbo-V Remote Panel

- 1 Botão para visualizar no visor os parâmetros cycle number, cycle time e pump life.
- 2 Botão para seleccionar o modo LOW SPEED. Está activo somente quando está seleccionado o modo de comando do painel frontal. Carregando nele uma vez, a bomba roda com velocidade "stand-by". Carregando nele novamente, desactiva-se o modo LOW SPEED.
- 3 Botão para visualizar no visor os parâmetros pump current, pump temperature, pump power e rotational speed. Está sempre activo independentemente do modo de funcionamento escolhido. Premindo os botões 3 e 4 em conjunto pelo menos 2 segundos, activa-se um programa com o qual é possível programar alguns parâmetros operativos.
- 4 Botão para enviar os comandos de START, STOP/RESET. Está activo somente quando estiver seleccionado o modo de comando do painel frontal. Carregando nele uma vez, activa-se a fase de arranque; carregando nele novamente, pára a bomba. Se a bomba tiver parado automaticamente por causa de uma avaria, é necessário carregar neste botão uma primeira vez para executar o reset do controller e uma segunda vez para reactivar a bomba.
- 5 Visor alfanumérico com cristais líquidos: matriz de pontos 2 linhas x 16 caracteres.



**Figura 4** Painel posterior do Turbo-V Remote Panel

- 1 Conectores de entrada dos sinais lógicos (o conector de acoplamento é dotado com a apropriada ponte de fechamento).
- 2 Conector de sinais lógicos de saída e monitoragem de frequência bomba mais saída analógica programável.
- 3 Conector de saída de alimentação (24 Vcc por calibragem).
- 4 Conector para entrada de alimentação do controller da bomba e comunicação serial com o controller da bomba.
- 5 Conector serial por controle remoto.

## **Acendimento do Controller**

Para ligar o controller é suficiente inserir o cabo de alimentação na tomada de rede.

## **Activação da bomba**

Para ligar a bomba ocorre habilitar a entrada de interlock e apertar o botão START do painel frontal.

## **Paragem da bomba**

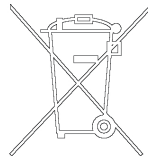
Para parar a bomba é necessário premer o botão STOP do painel frontal.

## Manutenção

Os controllers da série Turbo-V 301 Rack não requerem qualquer manutenção. Todas as operações devem ser efectuadas por pessoal autorizado. Em caso de defeito é possível utilizar o serviço de reparação Agilent ou o "Agilent advanced exchange service", que permite obter um controller regenerado que substitua o controller com defeito. Caso um controller deva ser destruído, proceder à sua eliminação respeitando as normas nacionais específicas.

## Eliminação

**Significado do logótipo "WEEE" presente nos rótulos.** O símbolo abaixo indicado é aplicado de acordo com a directiva CE denominada "WEEE". Este símbolo (**válido apenas para os países da Comunidade Europeia**) indica que o produto no qual está aplicado NÃO deve ser eliminado juntamente com os resíduos domésticos ou industriais comuns, mas deve ser dirigido a um sistema de recolha diferenciada. Portanto, convidamos o utilizador final a contactar o fornecedor do dispositivo, seja este o fabricante ou um revendedor, para encaminhar o processo de recolha e eliminação, após a oportuna verificação dos termos e condições do contrato de venda.



## Mensagens de erro

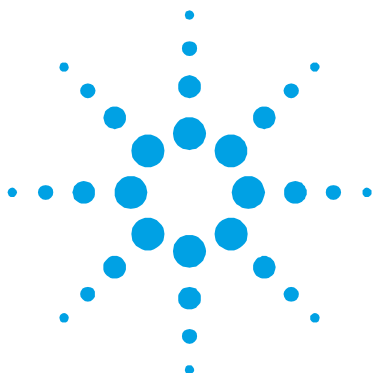
Em alguns casos de defeitos, os circuitos de autodiagnóstico do controller apresentam mensagens de erro relacionadas na tabela abaixo.

MENSAGEM	DESCRIÇÃO	ACÇÃO CORRECTIVA
COMMUNICATION ERROR	O Remote Panel não consegue arrancar a comunicação com o controller.	Verificar a conexão serial entre o Remote Panel e o controller.
COMMUNICATION BROKEN	O Remote Panel não consegue continuar a comunicação com o controller.	Verificar a conexão serial entre o Remote Panel e o controller.
XXXXXXX MODEL	Controller não suportado. O Remote Panel funciona somente com as seguintes bombas: TV551 com navigator controller, TV701 com navigator controller, TV1001 com navigator controller, TV2k-G e TV3k-T.	Ligar o Remote Panel ao controller suportado.
CHECK CONNECTION TO PUMP	Mau funcionamento na ligação entre a bomba e o controller. Ou A bomba possui temperatura inferior a 0 °C.	Verificar se o cabo de ligação entre a bomba e o controller está bem fixado em ambas as extremidades e não está interrompido. Premer duas vezes o botão START para reactivar a bomba.
WAITING INTERLOCK	É activado o sinal de interlock existente no conector P1 devido à interrupção do curto circuito entre o pin 3 e o pin 8 do conector J1 ou devido à abertura do sinal de interlock externo.	Restabelecer o curto circuito entre o pin 3 e o pin 8 do conector J1 ou fechar o sinal de interlock externo.
PUMP OVERTEMP.	A temperatura do rolamento superior ou da bomba superou os 60 °C.	Aguardar até que a temperatura volte ao limite estabelecido. Premer duas vezes o botão START para reactivar a bomba.
CONTROLLER OVERTEMP.	A temperatura do controller superou os 70 °C.	Aguardar até que a temperatura volte ao limite estabelecido.

<b>MENSAGEM</b>	<b>DESCRIÇÃO</b>	<b>ACÇÃO CORRECTIVA</b>
	<p>Ou</p> <p>A temperatura do radiador do controller é superior a 60 °C.</p>	<p>Premer duas vezes o botão START para reactivar a bomba.</p>
TOO HIGH LOAD	<p>Durante o funcionamento normal (após a fase de accionamento) a corrente absorvida pela bomba é maior do que a programada.</p>	<p>Verificar se o rotor da bomba pode rodar livremente.</p> <p>Premer duas vezes o botão START para reactivar a bomba.</p>
SHORT CIRCUIT	<p>Durante o funcionamento normal (após a fase de accionamento) a conexão de saída está em curto circuito.</p>	<p>Verificar as ligações entre a bomba e o controller.</p> <p>Premer duas vezes o botão START para reactivar a bomba.</p>
SYSTEM OVERRIDE	<p>A bomba parou por um sinal de emergência proveniente de um contacto remoto.</p>	<p>Remover o cabo de alimentação do controller e corrigir a causa da emergência. Ligar novamente o cabo de alimentação e premer duas vezes o botão START para reactivar a bomba.</p>
OVERVOLTAGE	<p>Verificou-se um defeito na secção de alimentação do controller, ou o controller recebeu um sinal falso.</p>	<p>Premer duas vezes o botão START para reactivar a bomba. Se a mensagem se reapresentar, dirigir-se à Agilent para a manutenção.</p>
POWER FAIL	<p>Avaria na seção de alimentação da bomba, do controller.</p>	<p>Dirigir-se a Agilent para Manutenção.</p>

## **5 Manual de Instruções**

### **Mensagens de erro**



## 6 Bedrijfshandleiding

Algemene Informatie	84
Opslag	86
Vorbereiding voor installatie	86
Installatie	87
Gebruik	89
Bedieningsorganen, controlelampjes en connectoren	89
Beschrijving frontpaneel	90
Inschakelen van de controller	92
Starten van de pomp	92
Stoppen van de pomp	92
Onderhoud	93
Afvalverwerking	93
Foutmeldingen	94

Vertaling van de originele instructies



## Algemene Informatie

Deze apparatuur is bestemd voor beroepsmatig gebruik. De gebruiker wordt verzocht aandachtig deze handleiding en alle overige door Agilent verstrekte informatie door te lezen alvorens het apparaat in gebruik te nemen. Agilent acht zich niet aansprakelijk voor de gevolgen van het niet of gedeeltelijk in acht nemen van de aanwijzingen, onoordeelkundig gebruik door niet hiervoor opgeleid personeel, reparaties waarvoor geen toestemming is verkregen of gebruik in strijd met de specifieke nationale wetgeving.

Het Turbo-V Remote Panel wordt gestuurd door een microprocessor en biedt de gebruiker de mogelijkheid een pomp van het type "controller on-board" te bedienen door middel van een frontpaneel:

TV551 met Navigator controller

TV701 met Navigator controller

TV1001 met Navigator controller

TV2k-G met controller on-board

TV3k-T met controller on-board

TV3k-G met controller on-board

TV70 met Navigator controller

TV81 met Navigator controller

TV301 met Navigator controller

TV303 met Navigator controller

In de volgende paragrafen wordt alle nodige informatie gegeven om de veiligheid van de operatie tijdens het gebruik van de apparatuur te verzekeren. Gedetailleerde informatie is te vinden in de bijlage "Technical Information".

Deze handleiding hanteert de volgende symbolen:

**VOORZICHTIG!**

Bij dit symbool staat tekst met procedures die, indien niet opgevolgd, schade aan apparatuur kunnen veroorzaken.

---

**WAARSCHUWING!**



Bij dit symbool staat tekst die de aandacht van de operator vestigt op een speciale procedure of methode die, indien niet correct uitgevoerd, ernstig lichamelijk letsel kan veroorzaken.

---

**OPMERKING**

De opmerkingen bevatten belangrijke informatie die uit de tekst is gelicht.

---

## Opslag

Tijdens het transport en de opslag van de controllers moeten de volgende omgevingscondities aanwezig zijn:

- Temperatuur: van -20 °C tot +70 °C
- Relatieve vochtigheid: 0 – 95 % (niet condenserend)

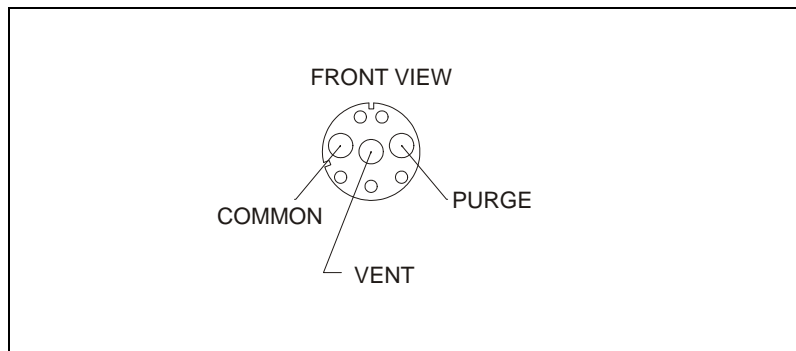
## Vorbereiding voor installatie

De Turbo-V Remote Panel wordt in een speciale beschermende verpakking geleverd; als er schade wordt geconstateerd die tijdens het transport veroorzaakt zou kunnen zijn, meteen contact opnemen met het plaatselijke verkoopkantoor. Zorg er bij het uitpakken voor dat de controller niet kan vallen of stoten te verduren krijgt. Laat de verpakking niet ergens buiten achter. Het verpakkingsmateriaal is volledig recyclebaar en voldoet aan de EEG-milieurichtlijn 85/399.

## Installatie

Het Turbo-V Remote Panel moet op de controller worden aangesloten met behulp van de speciale kabel (zie tabel bestelbare onderdelen), zoals getoond in onderstaande afbeelding. Aangezien het vermogen / de vereiste signalen door deze kabel worden geleverd, is geen andere externe voeding vereist.

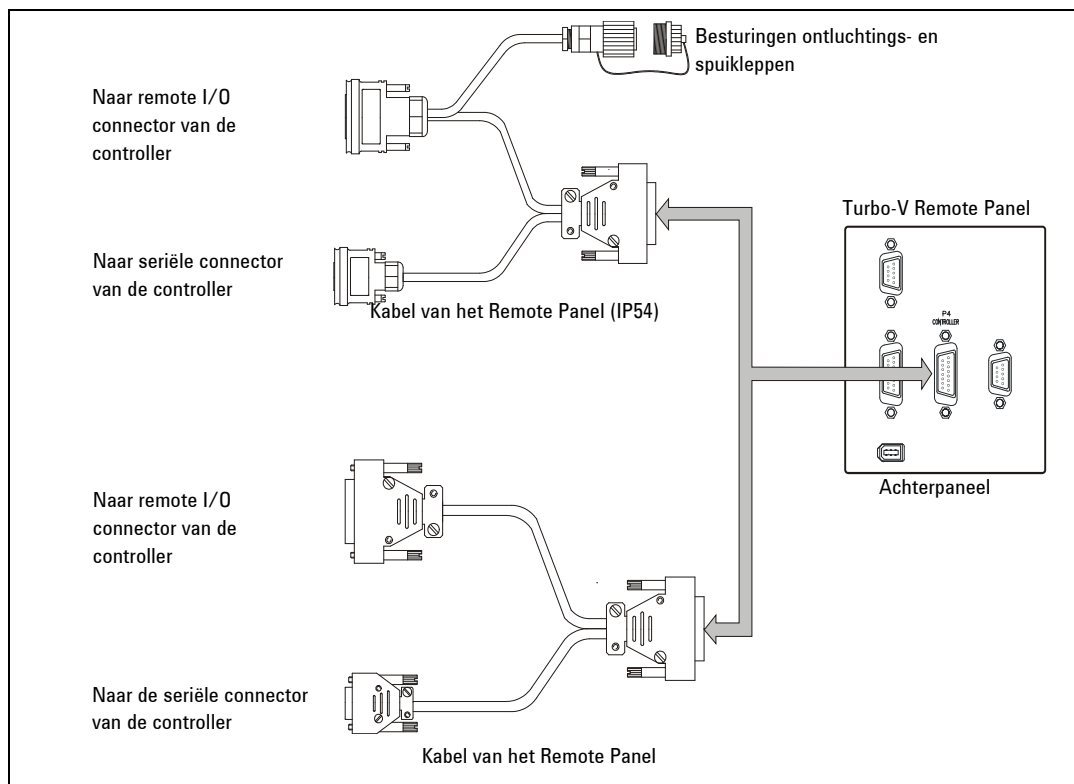
Indien de operator de ontluuchtings- en spuikleppen met gebruik van externe bedieningsorganen wenst te besturen, de uitvoering met IP54-kabewl gebruiken, de controller instellen op "valve setting" "on command" en de kabels aansluiten zoals in de afbeelding getoond. De kleppen zijn n.c.; bij leveren van 24 V tussen de ontluuchtings- en spuikleppen en het gezamenlijke eindcontact, gaan de kleppen open.



**Figuur 1**

## 6 Bedrijfshandleiding

### Installatie



**Figuur 2** Verbindingen

#### OPMERKING

Het Turbo-V Remote Panel kan op een tafel of in een speciaal rack worden geïnstalleerd. In ieder geval moet de koellucht vrij rondom het apparaat kunnen circuleren. De controller mag niet geïnstalleerd en/of gebruikt worden in ruimten die blootgesteld zijn aan weersinvloeden (regen, vorst, sneeuw), stof, agressieve gassen, of in ruimten met explosiegevaar of zeer hoog brandgevaar.

Tijdens de werking moeten de volgende omgevingscondities aanwezig zijn:

- temperatuur: van 5 °C tot +45 °C
- relatieve vochtigheid: 0 – 95 % (niet condenserend).

Voor de overige aansluitingen en de installaties van de optionele accessoires wordt verwezen naar het hoofdstuk "Technical Information".

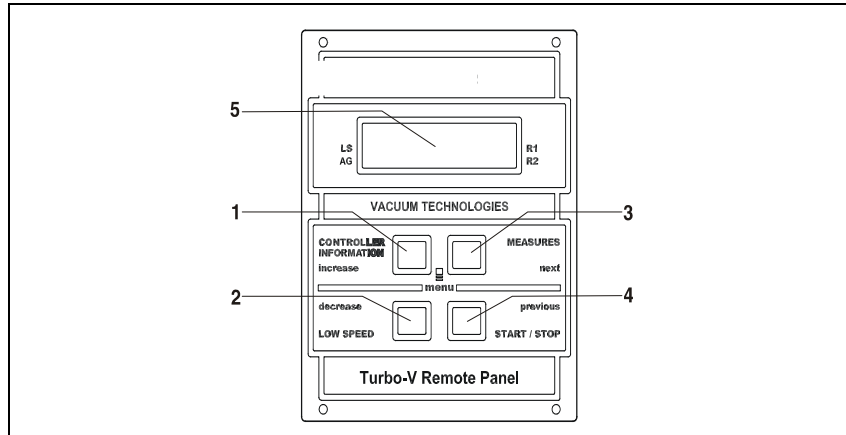
## Gebruik

In deze paragraaf worden de voornaamste bedieningswijzen uitgelegd. Voor meer informatie of procedures die aansluitingen of speciale opties betreffen, wordt verwezen naar de paragraaf "Use" van de bijlage "Technical Information".

## Bedieningsorganen, controlelampjes en connectoren

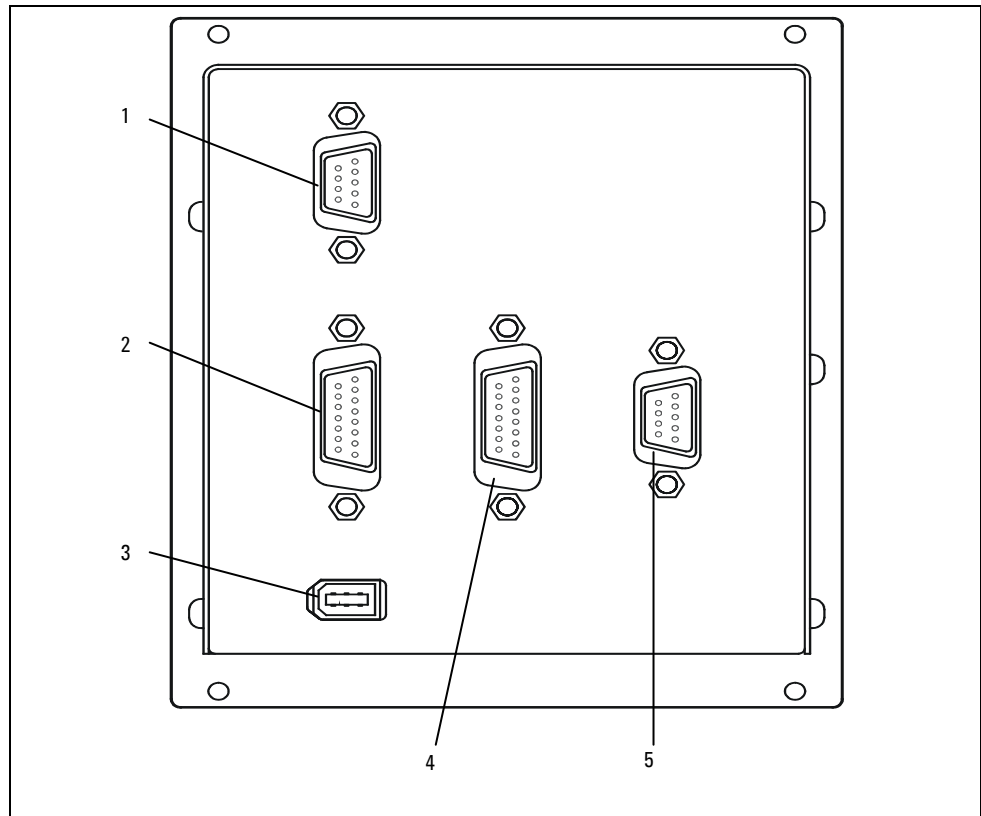
Hieronder worden het frontpaneel en het achterpaneel van de controller geïllustreerd. Voor meer informatie wordt verwezen naar het hoofdstuk "Technical Information".

## Beschrijving frontpaneel



**Figuur 3** Turbo-V Remote Panel

- 1 Drukknop voor weergave op het display van de parameters cycle number, cycle time en pump life.
- 2 Drukknop voor selectie van de modus LOW SPEED. Is alleen actief wanneer de bedieningsmodus vanaf het frontpaneel is geselecteerd. Door de knop eenmaal in te drukken, draait de pomp met "stand-by" snelheid. Door de knop opnieuw te bedienen wordt LOW SPEED geïnactiveerd.
- 3 Drukknop voor weergave op het display van de parameters pump current, pump temperature, pump power en rotational speed. Is altijd actief, onafhankelijk van de gekozen modus. Door de knoppen 3 en 4 ten minste 2 seconden gelijktijdig in te drukken, wordt een programma gestart waarmee enkele bedrijfsparameters kunnen worden geprogrammeerd.
- 4 Drukknop voor het zenden van de sturingen START, STOP/RESET. Is alleen actief wanneer de bedieningsmodus vanaf het frontpaneel is geselecteerd. Door de knop eenmaal in te drukken wordt de startfase geactiveerd; door de knop opnieuw te bedienen wordt de pomp gestopt. Als de pomp automatisch is gestopt als gevolg van een storing, moet deze knop een eerste maal worden bediend om de controller te resetten en een tweede maal om de pomp weer op te starten.
- 5 Alfanumeriek LCD-scherm: puntjespatroon, 2 regels x 16 tekens.



**Figuur 4** Achterpaneel van het Turbo-V Remote Panel

- 
- |   |                                                                                                               |
|---|---------------------------------------------------------------------------------------------------------------|
| 1 | Connectors voor ingang logische signalen (de te koppelen connector wordt met de speciale sluitklem geleverd). |
|---|---------------------------------------------------------------------------------------------------------------|
- 
- |   |                                                                                                           |
|---|-----------------------------------------------------------------------------------------------------------|
| 2 | Connector voor uitgang logische signalen en bewaking pompfrequentie plus programmeerbare analoge uitgang. |
|---|-----------------------------------------------------------------------------------------------------------|
- 
- |   |                                                |
|---|------------------------------------------------|
| 3 | Outputconnector voeding (24 Vdc voor kaliber). |
|---|------------------------------------------------|
- 
- |   |                                                                                         |
|---|-----------------------------------------------------------------------------------------|
| 4 | Inputconnector voeding vanaf pompcontroller en seriële communicatie met pompcontroller. |
|---|-----------------------------------------------------------------------------------------|
- 
- |   |                                           |
|---|-------------------------------------------|
| 5 | Seriële connector voor afstandsbediening. |
|---|-------------------------------------------|
-

## **Inschakelen van de controller**

Om de controller in te schakelen, de voedingskabel in de netcontactdoos inbrengen.

## **Starten van de pomp**

Om de pomp te starten, de interlock-ingang inschakelen en op de START-knop op het frontpaneel drukken.

## **Stoppen van de pomp**

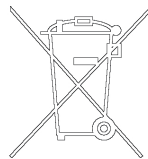
Voor het stoppen van de pomp de STOP-knop op het frontpaneel bedienen.

## Onderhoud

De controllers van de serie Turbo-V 301 zijn onderhoudsvrij. Eventuele werkzaamheden moeten door bevoegd personeel worden uitgevoerd. In geval van storing is het mogelijk om de reparatiedienst van Agilent of de "Agilent advanced exchange service" in te schakelen: zo krijgt men een ruilcontroller ter vervanging van de defecte controller. Mocht de controller gesloopt worden, ga dan overeenkomstig de specifieke nationale wetgeving te werk.

## Afvalverwerking

**Betekenis van het logo "WEEE" op de etiketten.** Het onderstaande symbool wordt aangebracht in overeenstemming met de EG-richtlijn "WEEE". Dit symbool (**alleen geldig voor de landen van de Europese Gemeenschap**) geeft aan dat het product waarop het is aangebracht, NIET mag worden afgevoerd samen met normaal huisvuil of industrieel afval, maar gescheiden moet worden ingezameld. De eindgebruiker wordt dus verzocht contact op te nemen met de leverancier van het apparaat, zij het de fabrikant of een wederverkoper, om het proces van gescheiden inzameling en verwerking in gang te zetten, na de van toepassing zijnde termen en voorwaarden van het verkoopcontract te hebben gecontroleerd.



## Foutmeldingen

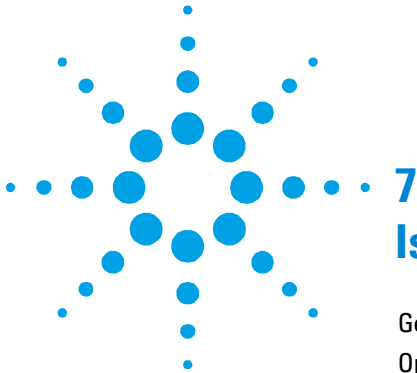
In geval van storingen wekt het zelfdiagnose-circuit van de controller enkele foutmeldingen op die in de volgende tabel zijn omschreven.

<b>BOODSCHAP</b>	<b>OMSCHRIJVING</b>	<b>REMEDIE</b>
COMMUNICATION ERROR	Het Remote Panel kan communicatie met de controller niet starten.	Controleer de seriële aansluiting tussen Remote Panel en controller.
COMMUNICATION BROKEN	Het Remote Panel kan communicatie met de controller niet standhouden.	Controleer de seriële aansluiting tussen Remote Panel en controller.
XXXXXXX MODEL	Controller niet ondersteund. Het Remote Panel werkt uitsluitend met de volgende pompen: TV551 met navigator controller, TV701 met navigator controller, TV1001 met navigator controller, TV2k-G en TV3k-T.	Sluit het Remote Panel aan op de ondersteunde controller.
CHECK CONNECTION TO PUMP	Foutieve verbinding tussen pomp en controller. Of De pomp heeft een lagere temperatuur dan 0 °C.	Controleren of de verbindingskabel tussen pomp en controller aan beide uiteinden goed bevestigd is en geen onderbrekingen vertoont. Bedien twee maal de START-knop om de pomp weer op te starten.
WAITING INTERLOCK	Het interlock-sigitaal op connector P1 is actief wegens onderbreking van de kortsluiting tussen pin 3 en pin 8 van connector J1 of wegens het openen van het externe interlock-sigitaal.	Herstel de kortsluiting tussen pin 3 en pin 8 van connector J1 of sluit het externe interlock-sigitaal.
PUMP OVERTEMP.	De temperatuur van het bovenste pomplager bedraagt meer dan 60 °C.	Wacht tot de temperatuur weer onder de drempelwaarde is gezakt. Bedien twee maal de START-knop om de pomp weer op te starten.
CONTROLLER OVERTEMP.	De temperatuur van de controller	Wacht tot de temperatuur weer

<b>BOODSCHAP</b>	<b>OMSCHRIJVING</b>	<b>REMEDIE</b>
	bedraagt meer dan 70 °C. Of De temperatuur van de radiator van de controller bedraagt meer dan 60 °C.	onder de drempelwaarde is gezakt. Bedien twee maal de START-knop om de pomp weer op te starten.
TOO HIGH LOAD	De stroomopname van de pomp is groter dan de geprogrammeerde waarde.	Controleer of de pomprotor vrij kan draaien. Bedien twee maal de START-knop om de pomp weer op te starten.
SHORT CIRCUIT	Tijdens normale werking (na startfase) is de uitgangsverbinding in kortsluiting.	Controleer de verbindingen tussen pomp en controller. Bedien twee maal de START-knop om de pomp weer op te starten.
SYSTEM OVERRIDE	De pomp is stilgelegd door een noodsignaal afkomstig van een afstandscontact.	Koppel de voedingskabel van de controller af en elimineer de oorzaak van de noodstop. Sluit de voedingskabel weer aan en bedien twee maal de START-knop om de pomp weer op te starten.
OVERVOLTAGE	Er is een storing opgetreden in de voedingssectie van de controller of de controller heeft een stoorsignaal ontvangen.	Bedien twee maal de START-knop om de pomp weer op te starten. Als de melding weer verschijnt zich voor onderhoud tot Agilent wenden.
POWER FAIL	Defect in het voedingsgedeelte van de pomp, van de controller.	Neem contact op met Agilent voor onderhoud.

## **6 Bedrijfshandleiding**

### **Foutmeldingen**



## 7 Istruktionsbog

Generel information	98
Opbevaring	100
Forberedelser før installation	100
Installation	101
Anvendelse	103
Betjeningsfunktioner, indikatorer og kontakter	103
Beskrivelse af frontpanel	104
Start af styreenheden	106
Start af pumpen	106
Stop af pumpen	106
Vedligeholdelse	107
Bortskaffelse	107
Fejlmeddelelser	108

Oversættelse af originalinstruktionerne



## Generel information

Dette materiel er beregnet til professionel anvendelse. Brugeren bør læse denne brugsanvisning og anden yderligere information fra Agilent, før udstyret anvendes. Agilent tager ikke ansvar for skader helt eller delvis som følge af tilsidesættelse af disse instruktioner, fejlagtig brug af personer uden tilstrækkelig kendskab, ukorrekt anvendelse af udstyret eller håndtering, der strider imod gældende lokale lovgivning.

Turbo-V Remote Panel er mikroprocessorstyret, så brugeren kan styre en pumpe af typen "on-board styreenhed " med et frontpanel:

- TV551 med Navigator styreenhed
- TV701 med Navigator styreenhed
- TV1001 med Navigator styreenhed
- TV2k-G med on-board styreenhed
- TV3k-T med on-board styreenhed
- TV3k-G med on-board styreenhed
- TV70 med Navigator styreenhed
- TV81 med Navigator styreenhed
- TV301 med Navigator styreenhed
- TV303 med Navigator styreenhed

De følgende afsnit indeholder al information der behøves, for at garantere operatørens sikkerhed under anvendelsen. Detaljeret information findes i bilaget "Technical Information".

I brugsanvisningen anvendes følgende standard-rubrikker:

**FORSIGTIG!**

Denne advarselsmeddelelse vises før procedurer, der skal følges nøje for at hindre maskinskader.

---

**ADVARSEL!**



Advarselsmeddelelserne informerer operatøren om, at en speciel procedure eller en vis type arbejde skal udføres præcist efter anvisningerne. I modsat fald er der risiko for alvorlige personskader.

---

**BEMÆRK**

Dette gør opmærksom på vigtig information i teksten

---

## Opbevaring

Følgende krav til omgivelserforholdene gælder ved transport og opbevaring af styreenhederne:

- temperatur: fra -20 °C til +70 °C
- relativ luftfugtighed: 0 – 95 % (ikke kondenserende)

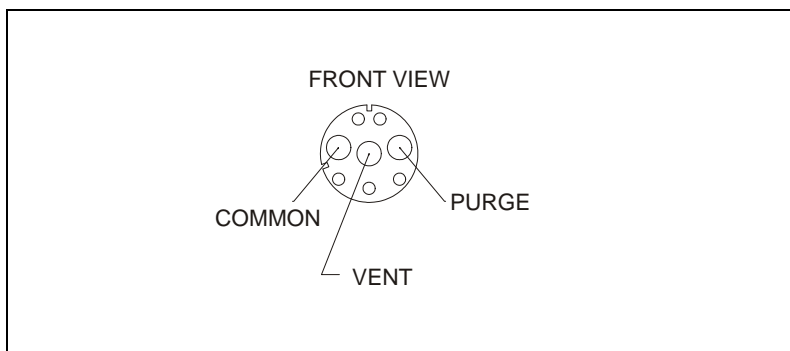
## Forberedelser før installation

Styreenheden leveres i en speciel beskyttende emballage. Kontakt den lokale forhandler, hvis emballagen viser tegn på skader, der kan være opstået under transporten. Sørg for at styreenheden ikke tabes eller udsættes for stød ved udpakningen. Smid ikke emballagen ud. Materialet kan genbruges 100 % og opfylder EU-direktiv 85/399 om miljøbeskyttelse.

## Installation

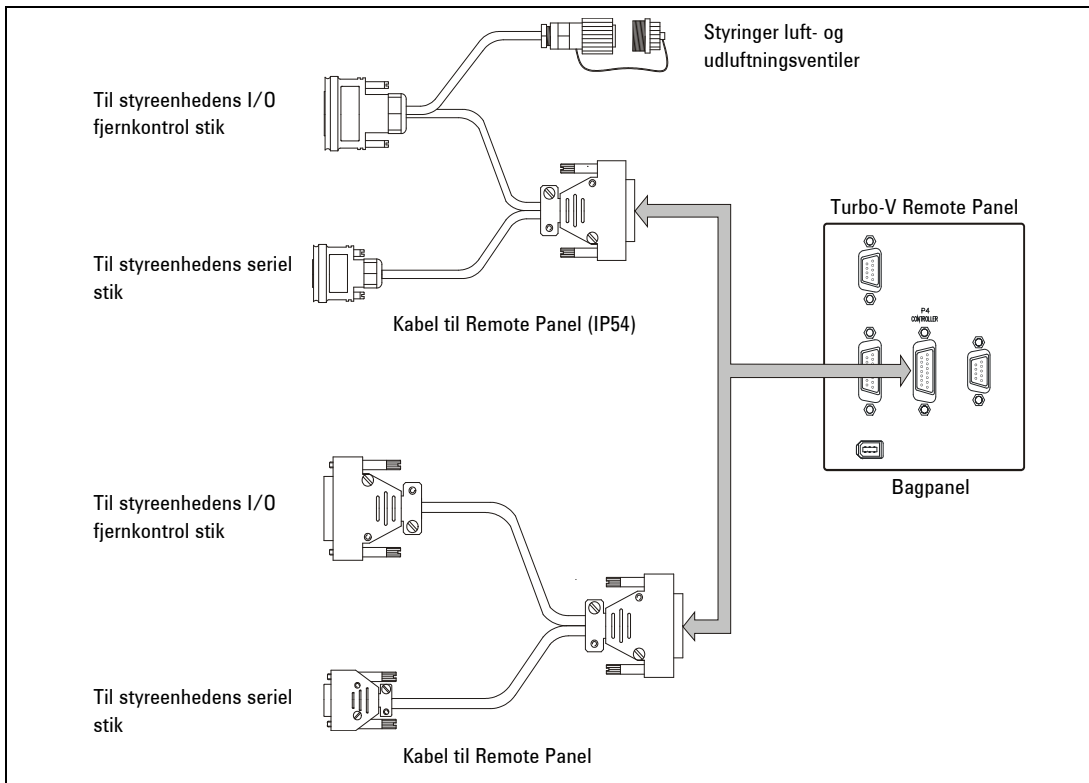
Turbo-V Remote Panel skal tilkobles styreenheden med det særlige kabel (se oversigt over reservedele, som kan bestilles), som vist på billedet. Da de nødvendige effektsignaler leveres af dette kabel, er anden ekstern strømtilførsel ikke nødvendig.

Hvis operatøren ønsker at styre luft- og udluftningsventilerne med de eksterne styringer, skal versionen med IP54 kabel anvendes. Indstil styreenheden på "valve setting" og tilkobl kablerne som vist på billedet. Ventilerne er n.l.; ved ydelse af 24 V mellem luft- og udluftningsventilen og den fælles klemme åbnes ventilerne.



Figur 1

## 7 Istruktionsbog Installation



**Figur 2** Tilslutninger

### BEMÆRK

Turbo-V Remote Panel kan installeres på et bord eller et velegnet stativ. I begge tilfælde skal der være plads nok til, at luft kan cirkulere frit omkring udstyret. Installér og anvend ikke styreenheden i miljøer, der udsættes for påvirkninger fra atmosfæren (regn, sne, is), damp, aggressive gasser, og ligeledes ikke i eksplosivt eller brandfarligt miljø.

Følgende krav til omgivelserforholdene gælder ved drift:

- temperatur: fra 5 °C til +45 °C
- relativ luftfugtighed: 0 – 95 % (ikke kondenserende).

For øvrige tilslutninger og installation af tilbehør henvises til afsnittet "Technical Information".

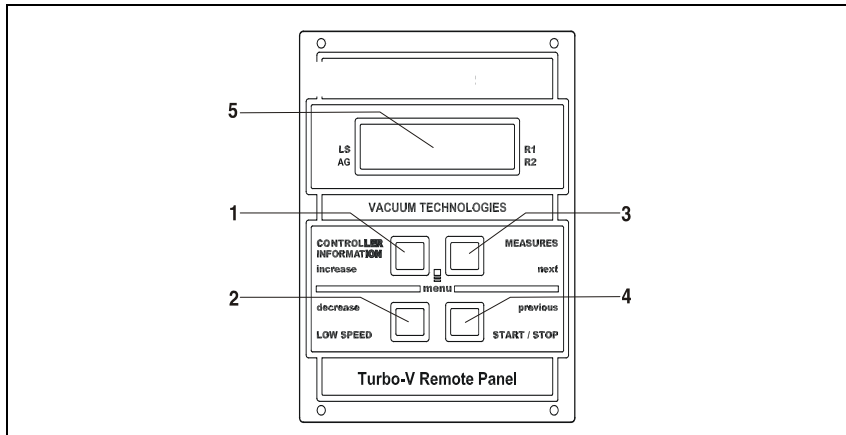
## Anvendelse

Dette afsnit beskriver de vigtigste driftsprocedurer. For en detaljeret beskrivelse samt procedurer, der involverer tilslutninger eller tilbehør, henvises til afsnittet "Use" i bilag "Technical Information".

## Betjeningsfunktioner, indikatorer og kontakter

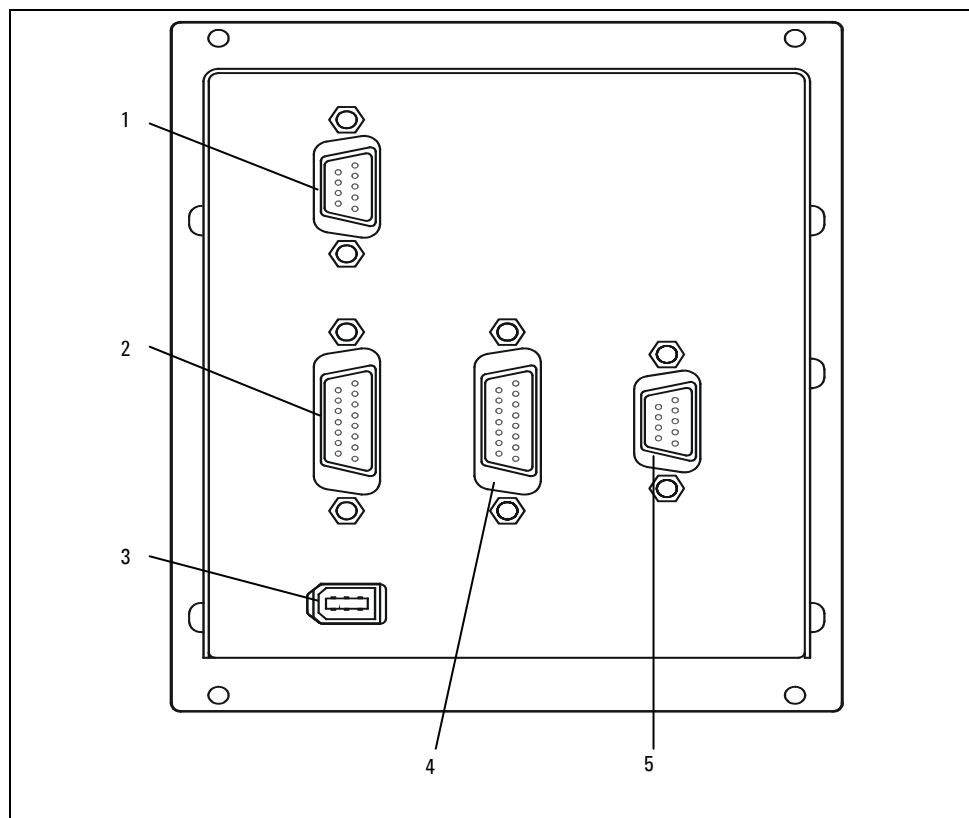
Nedenfor vises styreenhedens front- og bagpanel. For yderligere information henvises til bilag "Technical Information".

## Beskrivelse af frontpanel



**Figur 3** Turbo-V Remote Panel

- 1 Trykknop til at hente cycle number, cycle time og pump life frem på displayet.
- 2 Trykknop til installering af LOW SPEED driftsvalg. Trykknappen fungerer kun, når der vælges fra frontpanelet. Tryk på knappen en gang for at få pumpen til at rotere med "standby" hastighed. Tryk på knappen igen for at gå ud af LOW SPEED driften.
- 3 Trykknop til at hente pump current, pump temperature, pump power og rotational speed frem på displayet. Trykknappen fungerer altid, uanset den valgte styretype. Når trykknapperne 3 og 1 holdes trykket ind samtidigt i mindst 2 sekunder, starter et program, der tillader, at programmere visse driftsparametre.
- 4 Trykknop til overføring af betjeningsfunktionerne START, STOP/RESET. Trykknappen fungerer kun, når der vælges fra frontpanelet. Tryk på knappen en gang for at starte pumpen, tryk på knappen igen for at standse pumpen. Hvis pumpen standser automatisk ved en fejl, tryk på denne knap en gang for at nulstille styreenheden, og yderligere en gang for at starte pumpen igen.
- 5 LCD-display med tal og bogstaver: punktmatrice, 2 linier med 16 tegn.



**Figur 4** Turbo-V Remote Panel bagpanelet

- 1 Til de logiske signalers indgangsstik (koblingsstikket leveres med den pågældende lukningsbro).
- 2 Stik til logiske signaler i udgang og overvågning af pumpefrekvens samt programmerbar analogisk udgang.
- 3 Udgangsstik til strømforsyningen (24 Vdc pr. kaliber).
- 4 Stik til strømforsyningens indgang fra pumpens styreenhed og serial kommunikation med pumpens styreenhed.
- 5 Serial stik til fjernkontrollen.

## **Start af styreenheden**

Styreenheden startes ved at sætte strømkablet i vægudtaget.

## **Start af pumpen**

Pumpen startes ved at aktivere interlock indgangen og trykke på START-knappen på frontpanelet.

## **Stop af pumpen**

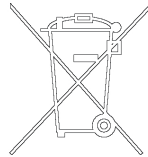
Pumpen stopper ved at trykke på STOP-trykknappen på frontpanelet.

## Vedligeholdelse

Styreenhederne i Turbo-V 301 Rack serien behøver ikke nogen vedligeholdelse. Ethvert indgreb på pumpen skal foretages af autoriseret personale. Hvis styreenheden går i stykker, kan man benytte sig af Agilents reparations-service eller Agilent udvekslingsservice, hvor man kan få en repareret pumpe i bytte for den, der er gået i stykker. Skrotning af styreenheden skal foregå i overensstemmelse med det pågældende lands særlige love.

## Bortskaffelse

**Betydningen af "WEEE" logoet på mærkaterne.** Nedenstående symbol anvendes i overensstemmelse med det såkaldte EU-direktiv "WEEE". Symbolet (**kun gældende for EU-landene**) viser, at produktet, som det sidder på IKKE må bortskaffes sammen med affald fra private husholdninger eller industriel affald men skal indleveres på en godkendt affaldsstation. Vi opfordrer derfor slutbrugeren til at kontakte leverandøren af anordningen, enten fabrikken eller en forhandler, for igangsættelse af afhentnings- og bortskaffelsesprocessen efter nøje at have kontrolleret betingelserne i salgskontrakten.



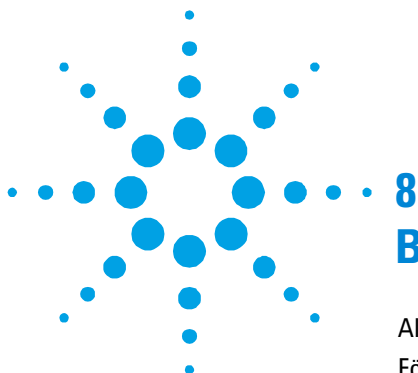
## Fejlmeddelelser

Når visse fejl opstår, viser styreenheden ved selvdiagnose aktuelle fejl på displayet. De mulige meddelelser listes i følgende tabel.

<b>MEDDELELSE</b>	<b>BESKRIVELSE</b>	<b>KONTROL</b>
COMMUNICATION ERROR	Remote Panel kan ikke starte kommunikationen med styreenheden.	Kontrollér den serielle forbindelse mellem Remote Panel og styreenheden.
COMMUNICATION BROKEN	Remote Panel kan ikke fortsætte kommunikationen med styreenheden.	Kontrollér den serielle forbindelse mellem Remote Panel og styreenheden.
XXXXXXX MODEL	Styreenhed ikke understøttet. Remote Panel virker kun med følgende pumper: TV551 med navigator styreenhed, TV701 med navigator styreenhed, TV1001 med navigator styreenhed, TV2k-G og TV3k-T.	Forbind Remote Panel med den understøttede styreenhed.
CHECK CONNECTION TO PUMP	Tilslutning mellem pumpe og styreenhed er defekt. eller Temperaturen på pumpen ligger under 0 °C.	Kontrollér at tilslutningskabel mellem pumpe og styreenhed er korrekt monteret samt at ingen afbrydninger forekommer. Tryk to gange på START for at starte pumpen igen.
WAITING INTERLOCK	Interlocksignalet findes på kontakt P1 p.g.a. kortslutning mellem stift 3 og stift 8 i kontakten J1 eller p.g.a. at det eksterne interlocksignal er åbent.	Tilbagestil kortslutningen mellem stift 3 og stift 8 på kontakt J1 eller sluk for det eksterne interlocksignal.
PUMP OVERTEMP.	Temperaturen på det øverste leje eller pumpen overskrider 60 °C.	Vent på at temperaturen falder til under tærskelværdi. Tryk to gange på START for at starte pumpen igen.
CONTROLLER OVERTEMP.	Temperaturen på styreenhedens overskrider 70 °C. eller Temperaturen på styreenhedens køler overskrider 60 °C.	Vent på at temperaturen falder til under tærskelværdi. Tryk to gange på START for at starte pumpen igen.

<b>MEDDELELSE</b>	<b>BESKRIVELSE</b>	<b>KONTROL</b>
TOO HIGH LOAD	Under normal drift (efter startfasen) forbruger pumpen en større effekt end den programmerede værdi.	Kontrollér om pumpens rotor kan rotere frit. Tryk to gange på START for at starte pumpen igen.
SHORT CIRCUIT	Under normal drift (efter startfasen) er udgangseffekten kortsluttet. Strømmen, som optages af pumpen overskrider den programmerede værdi.	Kontrollér forbindelserne mellem pumpe og styreenhed. Tryk to gange på START for at starte pumpen igen.
SYSTEM OVERRIDE	Pumpen er blevet standset af et nødstopsignal fra en fjernkontakt.	Frakobl styreenhedens strømkabel og kontrollér årsagen til nødstopet. Sæt derefter strømkablet i. Tryk to gange på START, for at starte pumpen igen.
OVERVOLTAGE	Der er opstået fejl i styreenhedens fødespænding eller styreenheden har fået et falsk signal.	Tryk to gange på START-knappen for at starte pumpen igen. Hvis meddelelsen kommer igen tag kontakt med Agilent for nødvendig vedligeholdelse.
POWER FAIL	Der er opstået fejl i sektionen for strømtilførsel til pumpen, styreenheden.	Tag kontakt med Agilent for Vedligeholdelse.

## **7** **Istruktionsbog** **Fejlmeddelelser**



## 8

# Bruksanvisning

Allmän information	112
Förvaring	114
Förberedelser för installation	114
Installation	115
Användning	117
Styrningar, indikatorer och kontakter	117
Beskrivning av frontpanel	118
Start av styrenheten	120
Start av pumpen	120
Stopp av pumpen	120
Underhåll	121
Bortskaffning	121
Felmeddelanden	122

Översättning av originalinstruktionerna



## Allmän information

Utrustningen är avsedd för yrkesmässig användning. Användaren bör noga läsa igenom denna bruksanvisning, samt övrig dokumentation från Agilent före användning av utrustningen. Agilent tar inget ansvar för skador som helt eller delvis orsakats av åsidosättande av instruktionerna, olämplig användning av person utan tillräcklig kunskap, obehörigt bruk av utrustningen eller hantering som strider mot gällande lokala föreskrifter.

Turbo-V Remote Panel styrs via en mikroprocessor och möjliggör styrning av en pump av typ "styrenhet ombord" genom en frontpanel:

TV551 med Navigator controller

TV701 med Navigator controller

TV1001 med Navigator controller

TV2k-G med styrenhet ombord

TV3k-T med styrenhet ombord

TV3k-G med styrenhet ombord

TV70 med Navigator controller

TV81 med Navigator controller

TV301 med Navigator controller

TV303 med Navigator controller

I avsnitten nedan ges all information som erfordras för att garantera operatörens säkerhet under användningen av utrustningen. Detaljerad information tillhandahålls i bilagan "Technical Information".

I bruksanvisningen används följande standard-rubriker:

**OBSERVER!**

Detta varningsmeddelande visas framför procedurer som måste följas exakt för att undvika skador på utrustningen.

---

**VARNING!**



Varningsmeddelandena informerar operatören om att en speciell procedur eller en viss typ av arbete måste utföras exakt enligt anvisningarna. I annat fall finns risk för svåra personskador.

---

**OBSERVERA**

Detta visar på viktig information i texten.

---

## Förvaring

Följande krav på omgivningsförhållanden gäller vid transport och förvaring av styrenheten:

- temperatur: från -20 °C till +70 °C
- relativ fuktighet: 0 – 95 % (utan kondens)

## Förberedelser för installation

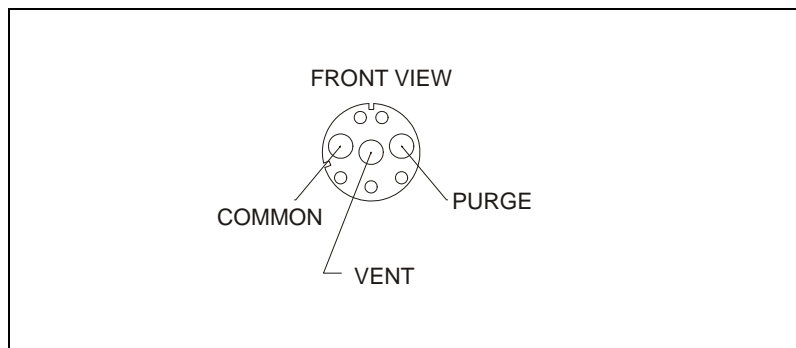
Styrenheten levereras i ett särskilt skyddande emballage. Kontakta det lokala försäljningskontoret om emballaget visar tecken på skador som kan ha uppstått under transporten. Se till att styrenheten inte tappas eller utsätts för stötar vid uppackningen. Kasta inte packmaterialet i soporna. Materialet är återvinningsbart till 100 % och uppfyller EU-direktiv 85/399 om miljöskydd.

## Installation

Turbo-V Remote Panel måste anslutas till styrenheten genom den särskilt avsedda kabeln (se tabellen som visar de delar som kan beställas), enligt bilden nedan. Eftersom den här kabeln leder både ström och signaler kommer ingen annan extern strömförsörjning att behövas.

I det fall operatören önskar reglera luft- och tömningsventilerna med hjälp av de externa styrningarna, bör versionen med IP54 kabeln användas, ställ in styrenheten på "valve setting" "on command" och anslut kablarna enligt vad som visas på bilden.

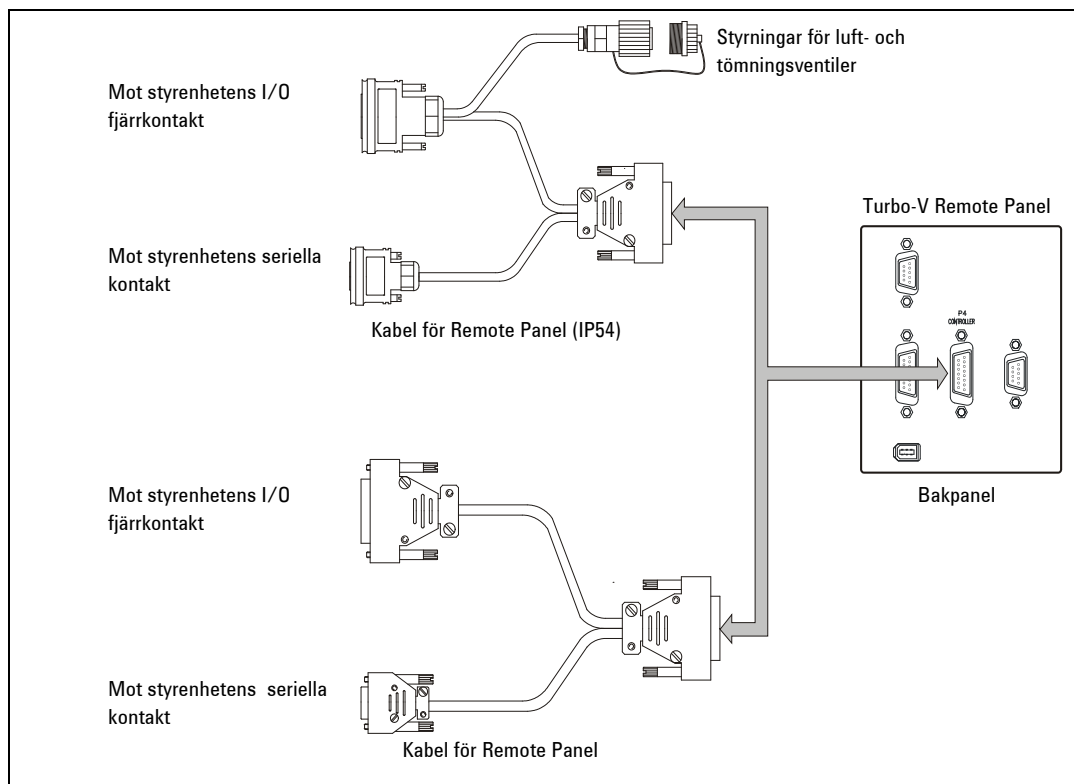
Ventilerna är normalt stängda (nc); genom tillförsel av 24 V mellan luft- och tömningsventilen och det gemensamma uttaget, kommer ventilerna att öppnas.



Figur 1

## 8 Bruksanvisning

### Installation



**Figur 2** Anslutningar

#### OBSERVERA

Turbo-V Remote Panel kan installeras på ett bord eller inuti ett därtill avsett rack. I samtliga fall måste dock kyl luften kunna cirkulera fritt kring apparaten. Installera och/eller använd inte styrenheten i miljöer som utsätts för påverkan från atmosfären (regn, snö, is), damm, aggressiva gaser, och inte heller i explosiv eller brandfarlig miljö.

Följande krav på omgivningsförhållanden gäller vid drift:

- temperatur: från 5 °C till +45 °C
- relativ fuktighet: 0 – 95 % (utan kondens).

Beträffande övriga anslutningar och installation av tillbehör hänvisas till avsnittet "Technical Information".

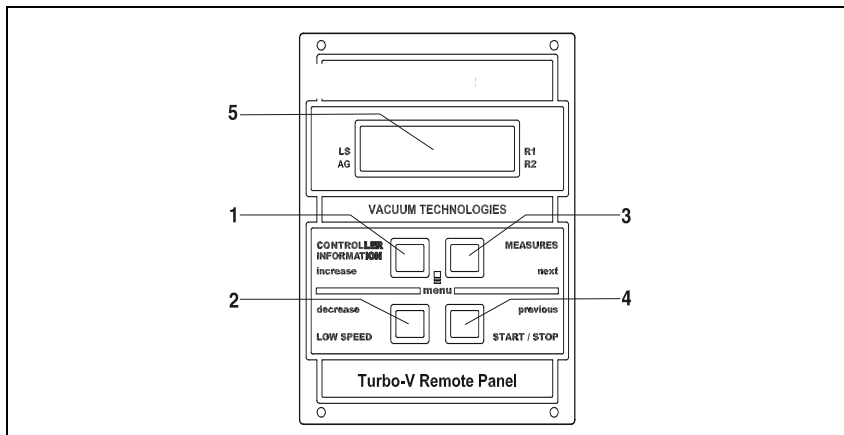
## Användning

Detta avsnitt beskriver de viktigaste driftmomenten. För en detaljerad beskrivning samt procedurer som involverar anslutningar eller tillbehör hänvisas till avsnittet "Use" i bilagan "Technical Information".

## Styrningar, indikatorer och kontakter

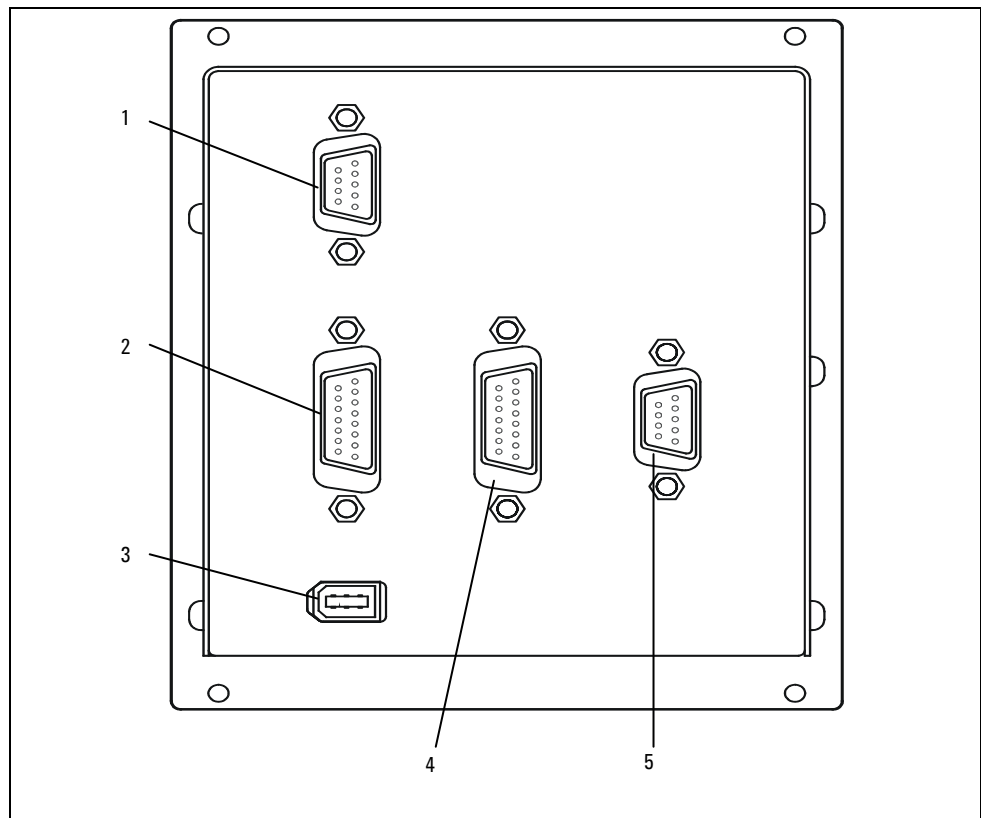
Nedan visas styrenhetens front- och bakpanel. För ytterligare information hänvisas till bilagan "Technical Information".

## Beskrivning av frontpanel



**Figur 3** Turbo-V Remote Panel

- 1 Tryckknapp för att hämta parametrarna cycle number, cycle time och pump life på displayen.
- 2 Tryckknapp för inställning av driftsättet LOW SPEED. Tryckknappen fungerar endast när styrsätt från frontpanelen har valts. Genom att trycka på knappen en gång kommer pumpen att rotera vid "stand-by" hastighet. Tryck på knappen igen för att gå ur driftsättet LOW SPEED.
- 3 Tryckknapp för att hämta parametrarna pump current, pump temperature, pump power och rotational speed. Tryckknappen fungerar alltid oavsett det valda styrsättet. När tryckknapparna 3 och 4 hålls samtidigt intryckta i minst 2 sekunder, startas ett programmera vissa driftparametrar.
- 4 Tryckknappen för överföring av START, STOP/RESET. Tryckknappen fungerar endast när styrsätt från frontpanelen har valts. Tryck på knappen en gång för att starta pumpen, tryck på knappen igen för att stanna pumpen. Om pumpen stannat automatiskt i samband med ett fel, tryck på denna knapp en gång för att återställa styrenheten, och ytterligare en gång för att starta om pumpen.
- 5 LCD Display med siffror och bokstäver: Punktmatris, 2 rader x 16 tecken.



**Figur 4** Bakpanel för Turbo-V Remote Panel

- |   |                                                                                                               |
|---|---------------------------------------------------------------------------------------------------------------|
| 1 | Kontakter för ingående logiksignaler (kopplingskontakten levereras med avsedd stängningsbrygga).              |
| 2 | Kontakt för utgående logiksignaler och kontroll av pumpfrekvens plus programmerbar analogisk utgång.          |
| 3 | Kontakt för utgående strömförsörjning (24 Vdc för mätinstrument).                                             |
| 4 | Kontakt för ingående strömförsörjning från pumpens styrenhet och seriell kommunikation med pumpens styrenhet. |
| 5 | Seriell kontakt för fjärrstyrning.                                                                            |

## **Start av styrenheten**

Styrenheten startas enkelt genom att sätta strömkabeln i vägguttaget.

## **Start av pumpen**

Vid start av pumpen måste ingången till interlock aktiveras och START-knappen tryckas in på frontpanelen.

## **Stopp av pumpen**

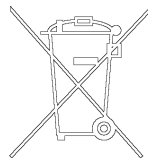
Pumpen stoppas genom att trycka på tryckknappen STOPP på frontpanelen.

## Underhåll

Styrenheterna i Turbo-V 301 Rack serien är under-hållsfria. Allt servicearbete måste utföras av auktoriserad personal. Om styrenheten havererar, kontakta Agilent reparationsverkstad eller Agilent utbytesservice, som kan ersätta styrenheten med en renoverad styrenhet. Skrotning av styrenheten skall ske enligt gällande lagstiftning.

## Bortskaffning

**Betydelse av logotypen "WEEE" på etiketterna.** Symbolen som visas nedan har tillämpats i enlighet med CD-direktivet som har betecknats som "WEEE". Den här symbolen (**gäller endast i de länder som tillhör den Europeiska Unionen**) indikerar att produkten på vilken symbolen har applicerats INTE får skaffas bort tillsammans med vanliga hushålls- eller industriavfall, men att däremot ett differentierat uppsamlingsystem måste upprättas. Vi ekommanderar därför att slutanvändaren tar kontakt med leverantören av anordningen, oberoende om det handlar om moderföretaget eller återförsäljaren, för att kunna starta uppsamlings- och bortskaffningsprocessen, detta efter lämplig kontroll av kontraktsevenliga tidsgränser och försäljningsvillkor.



## Felmeddelanden

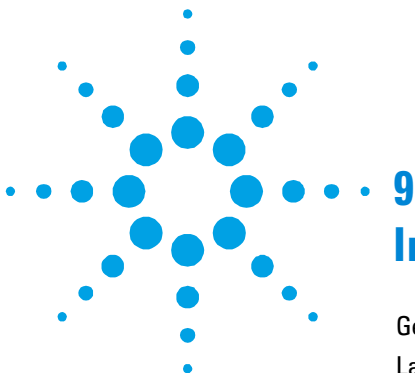
När vissa fel uppstår visar styrenhetens självdiagnoskrets aktuellt felmeddelande på displayen. De möjliga meddelandena listas i följande tabell.

MEDELLENDE	BESKRIVNING	ÅTGÄRD
COMMUNICATION ERROR	Remote Panel kan inte starta kommunikationen med styrenheten.	Kontrollera den seriella anslutningen mellan Remote Panel och styrenheten.
COMMUNICATION BROKEN	Remote Panel kan inte fortsätta kommunikationen med styrenheten.	Kontrollera den seriella anslutningen mellan Remote Panel och styrenheten.
XXXXXXX MODEL	Inget stöd till styrenheten. Remote Panel fungerar endast med följande pumpar: TV551 med navigator controller, TV701 med navigator controller, TV1001 med navigator controller, TV2k-G och TV3k-T.	Anslut Remote Panel till styrenheten som ger stöd.
CHECK CONNECTION TO PUMP	Anslutningen mellan pump och styrenhet är defekt. eller också Pumpens temperatur ligger under 0 °C.	Kontrollera att anslutningskabeln mellan pump och styrenhet är ordentligt monterad samt att inget avbrott förekommer. Tryck två gånger på knappen START för att starta om pumpen.
WAITING INTERLOCK	Interlock-signalen finns på kontakt P1 på grund av kortslutning mellan stift 3 och stift 8 i kontakten J1, eller på grund av att den externa interlock-signalen är öppen.	Åtgärda kortslutningen mellan stift 3 och stift 8 på kontakt J1 eller stäng den externa interlock-signalen.
PUMP OVERTEMP.	Temperaturen på det övre lagret eller pumpen överskrider 60 °C.	Vänta tills temperaturen sjunker under tröskelvärdet. Tryck två gånger på knappen START för att starta om pumpen.
CONTROLLER OVERTEMP.	Temperaturen på styrenhetens överskrider 70 °C. eller också	Vänta tills temperaturen sjunker under tröskelvärdet. Tryck två gånger på knappen

MEDELANDE	BESKRIVNING	ÅTGÄRD
	Temperaturen på styrenhetens kylare överskrider 60 °C.	START för att starta om pumpen.
TOO HIGH LOAD	Strömmen som absorberas av pumpen överstiger det värde som programmerats.	Kontrollera att pumpens rotor kan rotera fritt. Tryck två gånger på knappen START för att starta om pumpen.
SHORT CIRCUIT	Under normal drift (efter startmomentet) har utgången kortslutits.	Kontrollera anslutningarna mellan pump och styrenhet. Tryck två gånger på knappen START för att starta om pumpen.
SYSTEM OVERRIDE	Pumpen har stannats av en nödstoppsignal från en fjärrkontakt.	Frånkoppla styrenhetens strömkabel och kontrollera nödstoppets orsak. Sätt därefter i strömkabeln. Tryck två gånger på knappen START för att starta om pumpen.
OVERVOLTAGE	Ett fel i styrenhetens matningsdel har uppstått eller styrenheten har fått en falsk signal.	Tryck två gånger på knappen START för att starta om pumpen. Om meddelandet visas igen, ta kontakt med Agilent för nödvändigt underhåll.
POWER FAIL	Fel i sektionen för strömtillförsel till pumpen, styrenheten.	Ta kontakt med Agilent för Underhåll.

## **8 Bruksanvisning**

### **Felmeddelanden**



## 9

# Instruksjon Manual

Generell informasjon	126
Lagring	128
Forberede installasjonen	128
Installasjon	129
Bruk	131
Styringer, indikatorer og kontakter på styreenheten	131
Beskrivelse av frontpanelet	132
Starte styreenheten	134
Starte pumpen	134
Stoppe pumpen	134
Vedlikehold	135
Eliminering	135
Feilmeldinger	136

Oversetting av den opprinnelige samsvarserklæringen



## Generell informasjon

Dette apparatet er beregnet til bruk av profesjonelle brukere. Brukeren må lese denne brukerveiledningen og all annen ekstra informasjon fra Agilent før apparatet tas i bruk. Agilent fratar ansvar for hendelser som skjer på grunn av manglende oppfølging, selv delvis, av disse instruksjonene, feilaktig bruk av utrenet personell, av inngrep på utstyret som ikke er autorisert eller handlinger som på noen måte er i strid med nasjonale bestemmelser.

Turbo-V Remote Panel er styrt av en mikroprosessor som gjør det mulig for brukeren å styre en pumpe av type "controller on-board" fra et frontpanelet:

- TV551 med Navigator controller
- TV701 med Navigator controller
- TV1001 med Navigator controller
- TV2k-G med controller on-board
- TV3k-T med controller on-board
- TV3k-G med controller on-board
- TV70 med Navigator controller
- TV81 med Navigator controller
- TV301 med Navigator controller
- TV303 med Navigator controller

De følgende avsnittene inneholder all informasjon som er nødvendig for å sikre brukeren når utstyret er i bruk. For mer detaljert bruk vises det til tillegget "Technical Information".

**Denne veiledningen bruker følgende standard-protokoll:**

**FORSIKTIG!**

Denne advarselen vises foran fremgangsmåter som, dersom de ikke følges, kan føre til at utstyret skades.

---

**ADVARSEL!**



Disse meldingene skal tiltrekke seg brukerens oppmerksomhet til en spesiell fremgangsmåte eller praksis som, hvis den ikke følges, kan medføre alvorlige skader.

---

**MERK**

Merknadene inneholder viktig informasjon som er hentet fra teksten.

---

## Lagring

Når styreenhetene transporteres eller lagres, må følgende værforhold være oppfylt:

- temperatur: fra - 20 °C til +70 °C
- relativ fuktighet: 0 – 95 % (uten kondens)

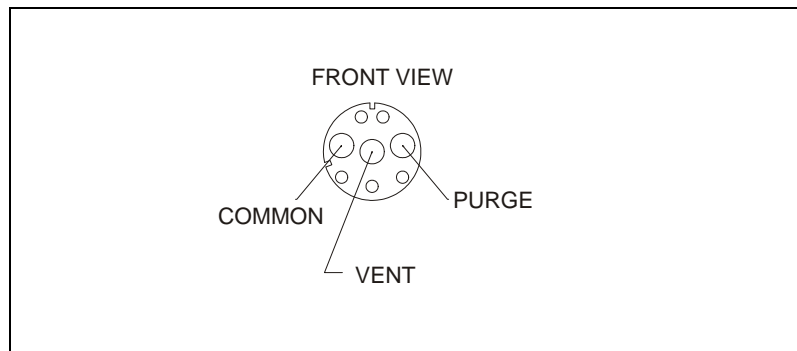
## Forberede installasjonen

Styreenheten leveres i en spesiell beskyttelses-emballasje. Viser denne tegn på skader som kan ha oppstått under transporten, må du ta kontakt med det lokale salgskontoret. Når styreenheten pakkes ut, må du passe på at den ikke slippes ned eller utsettes for støter. Emballasjen må ikke kastes ut i omgivelsen. Alle materialer er fullstendig resirkulerbare og er i samsvar med EUDirektiv 85/399 om miljøbeskyttelse.

## Installasjon

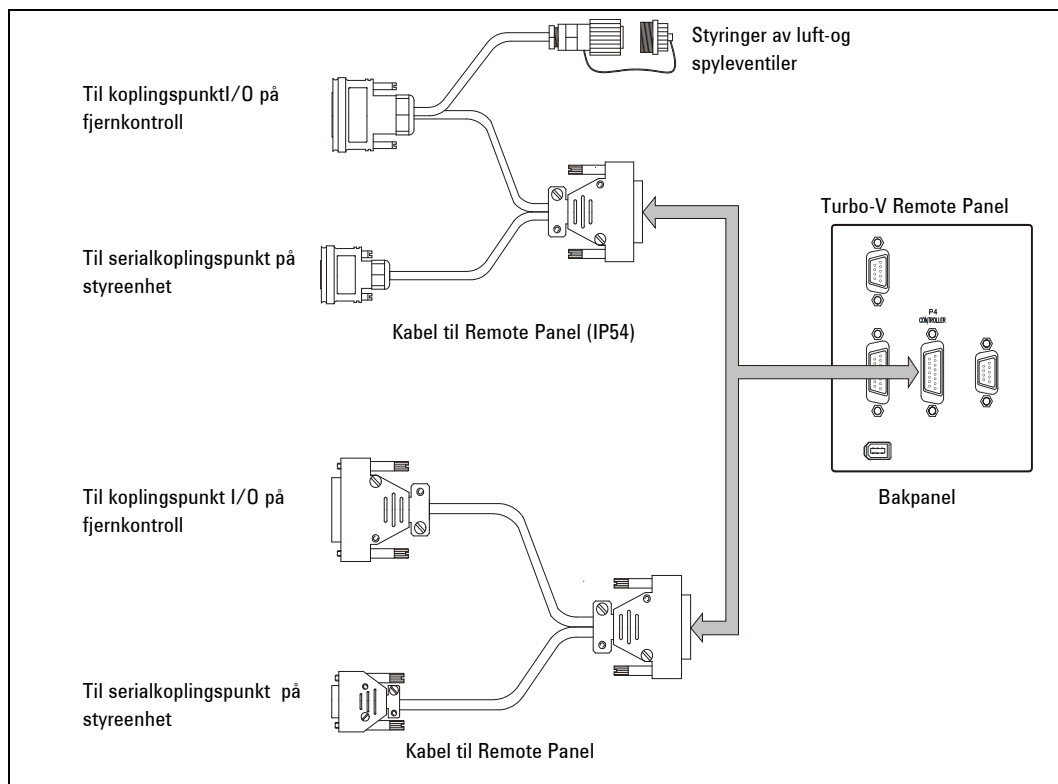
Turbo-V Remote Panel må koples til styreenheten med en spesiell kabel (se tabell over deler som kan bestilles), som illustrert på figuren som følger. Siden kraft/signaler som kreves, blir gitt av denne kablet, er det ikke nødvendig med en ekstern mating.

Hvis operatøren vil styre luft-og spyleventilene ved å bruke de eksterne styringene, må versjon med IP54-kabel brukes. Innstill styreenheten på "valve setting" "on command" og velge kablene som illustrert på figuren. Ventilene er n.c.: ved tilføsel av 24 volt mellom luft- og spyleventilen og terminalen, åpnes ventilene.



Figur 1

## 9 Instruksjon Manual Installasjon



**Figur 2** Koblere

### MERK

Turbo-V Remote Panel kan installeres på et bord eller inne i et passende stativ. Uansett så må luft for avkjøling kunne sirkulere fritt rundt apparatet. Ikke installer eller bruk styreenheten i miljøer som utsettes for regn, snø eller is, støv, aggressive gasser, eksplosjonsfarlige miljøer eller i miljøer med stor brannfare.

Under bruk må følgende værforhold respekteres:

- temperatur: fra 5 °C til +45 °C
- relativ fuktighet: 0 – 95 % (uten kondens)

Når det gjelder andre tilkoblinger og installasjon av valgfritt tilbehør vises det til avsnittet "Technical Information".

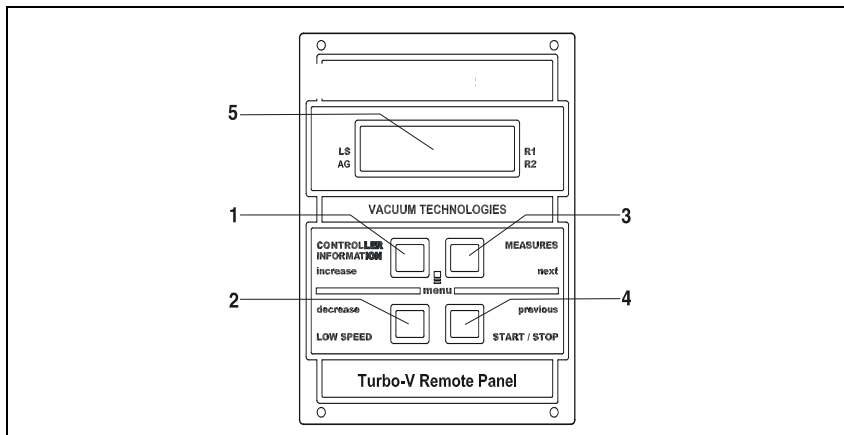
## Bruk

Dette avsnittet beskriver de viktigste operative prosedyrene. For en detaljert beskrivelse og prosedyrer som omfatter tilkoblinger eller særskilte valgfrie utstyr vises det til avsnittet "Use" i vedlegget "Technical Information".

## Styringer, indikatorer og kontakter på styreenheten

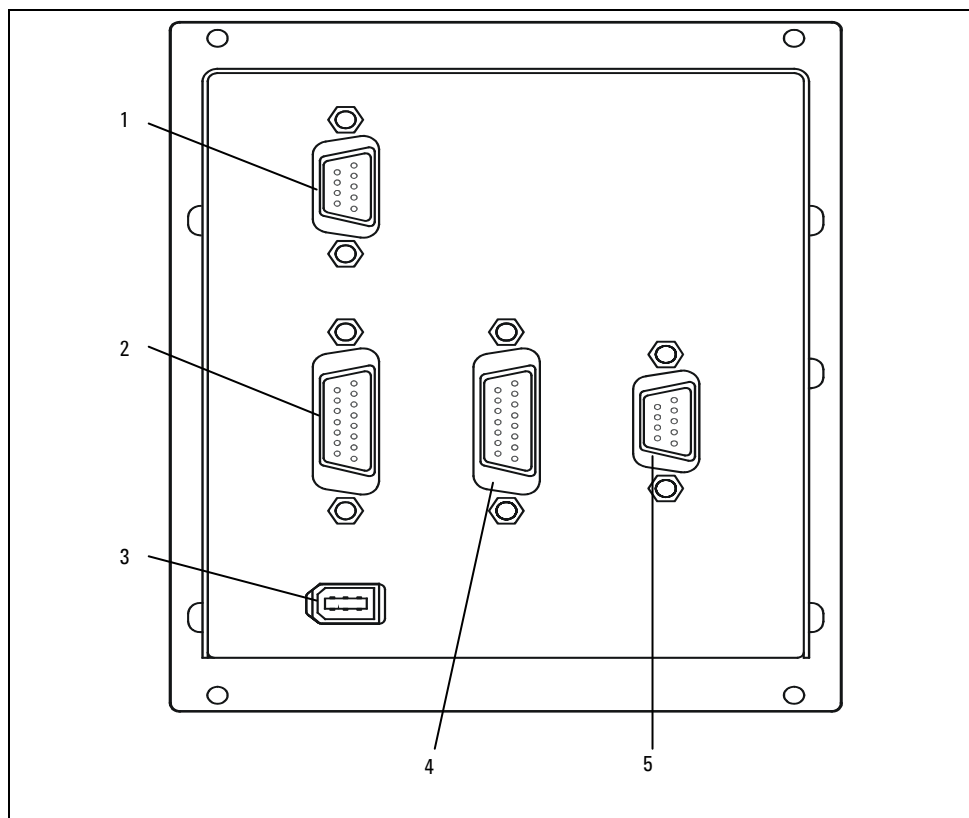
Nedenunder er frontpanelet og bakpanelet på styreenheten illustrert. For ytterligere informasjon vises det til avsnittet "Technical Information".

## Beskrivelse av frontpanelet



**Figur 3** Turbo-V Remote Panel

- 1 Trykknapp for å vise parametrene cycle number, cycle time og pump life på displayet.
- 2 Trykknapp for innstilling av LOW SPEED drift. Trykknappen fungerer kun når det er valgt styremåte fra frontpanelet. Ved å trykke på knappen én gang, roterer pumpen i "stand-by" hastighet. Trykk på knappen igjen for å gå ut av LOW SPEED drift.
- 3 Trykknapp for å vise parametrene pump current, pump temperature, pump power og rotational speed på displayet. Trykknappen fungerer alltid, uavhengig av styremåten. Når trykknappene 3 og 4 trykkes ned samtidig i minst to sekunder, startes et program som tillater programmering av visse driftsparametre.
- 4 Trykknapp for overføring av START, STOP/RESET. Trykknappen fungerer kun når det er valgt styremåte fra frontpanelet. Trykk på knappen en gang for å starte pumpen, trykk på knappen igjen for å stanse pumpen. Har pumpen stanset automatisk i forbindelse med en feil, trykk en gang på denne knappen for å tilbakestille styreenheten, en gang til for å starte pumpen.
- 5 LCD-display med siffer og bokstaver: punkt-matrise, 2 rader med 16 tegn.



**Figur 4** Turbo-V Remote Panels bakpanel

- 
- |   |                                                                                               |
|---|-----------------------------------------------------------------------------------------------|
| 1 | Inngangskoblingsstykker for logiske signaler (koblingsstykket leveres med passende lukkebro). |
|---|-----------------------------------------------------------------------------------------------|
- 
- |   |                                                                                                                      |
|---|----------------------------------------------------------------------------------------------------------------------|
| 2 | Koblingsstykke ofr logiske signaler i utgang og overvåking av pumpens frekvens pluss programmerbar analogisk utgang. |
|---|----------------------------------------------------------------------------------------------------------------------|
- 
- |   |                                                                |
|---|----------------------------------------------------------------|
| 3 | Utgangskoblingsstykke for strømtilførsel (24 V dc pr kaliber). |
|---|----------------------------------------------------------------|
- 
- |   |                                                                                                                      |
|---|----------------------------------------------------------------------------------------------------------------------|
| 4 | Koblingsstykke for inngang av strømtilførsel fra pumpens controller og seriell kommunikasjon med pumpens controller. |
|---|----------------------------------------------------------------------------------------------------------------------|
- 
- |   |                                          |
|---|------------------------------------------|
| 5 | Seriellkoblingsstykke for fjernkontroll. |
|---|------------------------------------------|
-

## **Starte styreenheten**

Styreenheten startes ved å sette strømkabelen i veggkontakten.

## **Starte pumpen**

For å starte pumpen må man preparere interlockinngangen og trykke på knappen START på frontpanelet.

## **Stoppe pumpen**

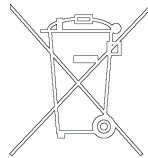
Pumpen stoppes ved å trykke på knappen STOPP på frontpanelet.

## Vedlikehold

Turbo-V 301 Rack seriens styreenheter er vedlikeholdsfrie. Alt arbeid på styreenheten må kun utføres av autorisert personell. Dersom styreenheten stanser, må du ta kontakt med Agilents reparasjonsservice eller med Agilents avanserte bytteservice, som kan tilby overholte styreenheter til erstatning for den ødelagte styreenheten. Dersom en styreenhet skal kasseres, må dette skje i henhold til nasjonale bestemmelser.

## Eliminering

**Betydelsen av symbolet på logo "WEEE" på etikettene.** Symbolet nedenunder som finnes, er anvendt i henhold til EC-direktiv kalt "WEEE". Dette symbolet (**som bare gjelder for land i Det europeiske fellesskap**), viser at produktet som det sitter på, **IKKE** må behandles som vanlig hus-industriavfall, men må legges i kildesortering. Derfor oppfordrer man den sluttelige brukeren av anordningen å henvende seg til leverandøren av anordningen, som kan være et firma eller en forhandler, som sørger for oppsamling og eliminering etter å ha kontrollert avtal og betingelser i kjøpekontrakten.



## Feilmeldinger

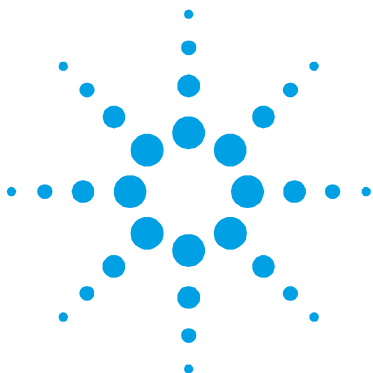
Når det oppstår visse feil viser selvdiagnosekretsen i styreenheten den aktuelle feilmeldingen i displayet. De aktuelle feilmeldingen fremgår av tabellen nedenfor.

MELDING	BESKRIVELSE	FORHOLDSREGEL
COMMUNICATION ERROR	Remote Panel klarer ikke å starte opp kommunikasjonen med controller.	Kontrollere den serielle kommunikasjonen mellom Remote Panel og controller.
COMMUNICATION BROKEN	Remote Panel klarer ikke å fortsette kommunikasjonen med controller.	Kontrollere den serielle kommunikasjonen mellom Remote Panel og controller.
XXXXXXX MODEL	Controller støttes ikke. Remote Panel fungerer bare med følgende pumper: TV551 med navigator controller, TV701 med navigator controller, TV1001 med navigator controller, TV2k-G og TV3k-T.	Koble Remote Panel til controller som støttes.
CHECK CONNECTION TO PUMP	Defekt kopling mellom pumpe og styreenhet. Eller Pumpen har en temperatur under 0 °C.	Kontroller at tilkoplingskabelen mellom pumpe og styreenhet er skikkelig montert samt at kabelen ikke er skadet. Trykk to ganger på knappen START for å starte pumpen.
WAITING INTERLOCK	Låsesignalet for kontakt P1 skyldes en kortslutning mellom stift 3 og stift 8 i kontakten J1 eller fordi det eksterne låsesignalet er åpent.	Tilbakestill kortslutningen mellom stift 3 og stift 8 på kontakt J1 eller steng det eksterne låsesignalet.
PUMP OVERTEMP.	Temperaturen på det øvre lageret eller pumpen er over 60 °C.	Vent til temperaturen synker under terskelverdien. Trykk to ganger på knappen START for å starte pumpen.
CONTROLLER OVERTEMP.	Temperaturen på styreenhetens er over 70 °C. Eller Temperaturen på radiatoren på styreenheten er over 60 °C.	Vent til temperaturen synker under terskelverdien. Trykk to ganger på knappen START for å starte pumpen.

MELDING	BESKRIVELSE	FORHOLDSREGEL
TOO HIGH LOAD	Den absorberte strømmen av pumpen er større enn den som er programmert	Kontroller om pumpens rotor kan rotere fritt. Trykk to ganger på knappen START for å starte pumpen.
SHORT CIRCUIT	Ved normal drift (etter startmomentet) er utgangen kortsluttet.	Kontroller tilkoplingene mellom pumpe og styreenhet. Trykk to ganger på knappen START for å starte pumpen.
SYSTEM OVERRIDE	Pumpen har stanset av et nødstoppsignal en fra fjernkontakt.	Kople fra styreenhetens strømkabel og finn frem til årsaken til nødstoppen. Kople deretter maskinen til strømmettet igjen. Trykk to ganger på knappen START for å starte pumpen.
OVERVOLTAGE	Det har oppstått en feil i styreenhetens matedel, eller så har styreenheten fått et falskt signal.	Trykk to ganger på knappen START for å starte pumpen igjen. Vises feilmeldingen om igjen, må du ta kontakt med Agilent for nødvendig vedlikehold.
POWER FAIL	Feil på matesektoren på pumpen, på styreenheten.	Henvend deg til Agilent for vedlikehold.

## **9 Instruksjon Manual**

### **Feilmeldinger**



## 10 Ohjekäsikirja

Yleistä	140
Säilytys	142
Asennuksen valmistelu	142
Asennus	143
Käyttö	145
Ohjaimet, ilmaisimet ja liittimet	145
Beskrivelse av frontpanelet	146
Valvojan päälle pano	148
Pumpun käynnistys	148
Pumpun pysäyttäminen	148
Huolto	149
Hävittäminen	149
Vianetsintä	150

Alkuperäisen ohjeiden käännös



## Yleistä

Tämä laite on tarkoitettu ammattimaiseen käyttöön. Ennen laitteen käyttöä käyttäjän tulee lukea huolellisesti tämä käyttöopas ja kaikki muut Agilentin toimittamat tiedot. Agilent ei vastaa seurauksista, jotka ovat johtuneet laitteen käyttöohjeiden täydellisestä tai osittaisesta laiminlyönnistä, laitteen virheellisestä käytöstä tehtävään perehdyttämättömien henkilöiden taholta, laitteeseen luvattomasti tehdyistä toimenpiteistä tai erillisten maakohtaisten säädösten vastaisesta käytöstä.

Turbo-V Remote Panel toimii mikroprosessorin avulla. Käyttäjä voi taulun ansiosta ohjata "controller on board" tyyppistä pumppua etupaneelin välityksellä:

TV551 varustettu Navigator controller-säätimellä

TV701 varustettu Navigator controller-säätimellä

TV1001 varustettu Navigator controller-säätimellä

TV2k-G varustettu controller on board-säätimellä

TV3k-T varustettu controller on board-säätimellä

TV3k-G varustettu controller on board-säätimellä

TV70 varustettu Navigator controller-säätimellä

TV81 varustettu Navigator controller-säätimellä

TV301 varustettu Navigator controller-säätimellä

TV303 varustettu Navigator controller-säätimellä

Seuraavissa kappaleissa annetaan kaikki tarvittava tieto käyttäjän turvallisuuden takaamiseksi laitteen käytön aikana. Yksityiskohtaisemmat tiedot löytyvät liitteestä "Technical Information".

Tässä ohjeessa on käytetty seuraavia merkkejä:

**HUOMIO!**

Huomioviestejä on merkitty sellaisten toimenpiteiden edelle, jotka voivat virheellisesti suoritettuina aiheuttaa vaurioita laitteistolle.

---

**VAROITUS!**



Vaaraviestien tarkoituksena on kiinnittää käyttäjän huomio tiettyyn toimenpiteeseen tai käytäntöön, joka voi virheellisesti suoritettuna johtaa vakaviin henkilövahinkoihin.

---

**HUOM**

Huomautuksissa on tekstistä poimittua tärkeää tietoa

---

## Säilytys

Säädinten kuljetuksen ja varastoinnin aikana on huolehdittava siitä, että ympäristöolosuhteet ovat seuraavien kriteerien mukaiset:

- lämpötila: -20 °C – +70 °C
- suhteellinen kosteus: 0–95 % (ei kondensoituva)

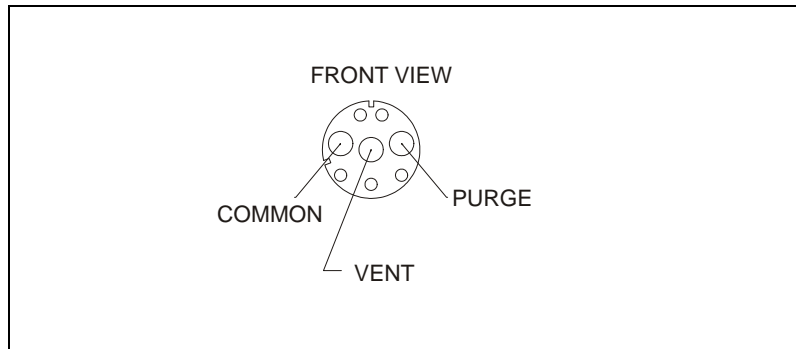
## Asennuksen valmistelu

Säädin toimitetaan erityisessä suojapakkauksessa. Jos siinä näkyy merkkejä vaurioista, jotka ovat saattaneet syntyä kuljetuksen yhteydessä, ota yhteyttä paikalliseen jälleenmyyjään. Kun purat säädintä pakkauksesta, varo ettei se pääse putoamaan tai osumaan mihinkään. Älä heitä pakkausta luontoon. Pakkausmateriaali on täysin kierrätyskelpoista ja ympäristönsuojelua koskevan direktiivin 85/399/ETY vaatimusten mukaista.

## Asennus

Turbo-V Remote Panel on liitettävä säätimeen tähän tarkoitettulla kaapelilla (ks. tilaamalla saatavien osien luettelo) seuraavassa kuvassa esitetyllä tavalla. Teho ja tarvittavat signaalit siirtyvät tätä kaapelia pitkin, joten muuta ulkoista virransyöttöä ei tarvita.

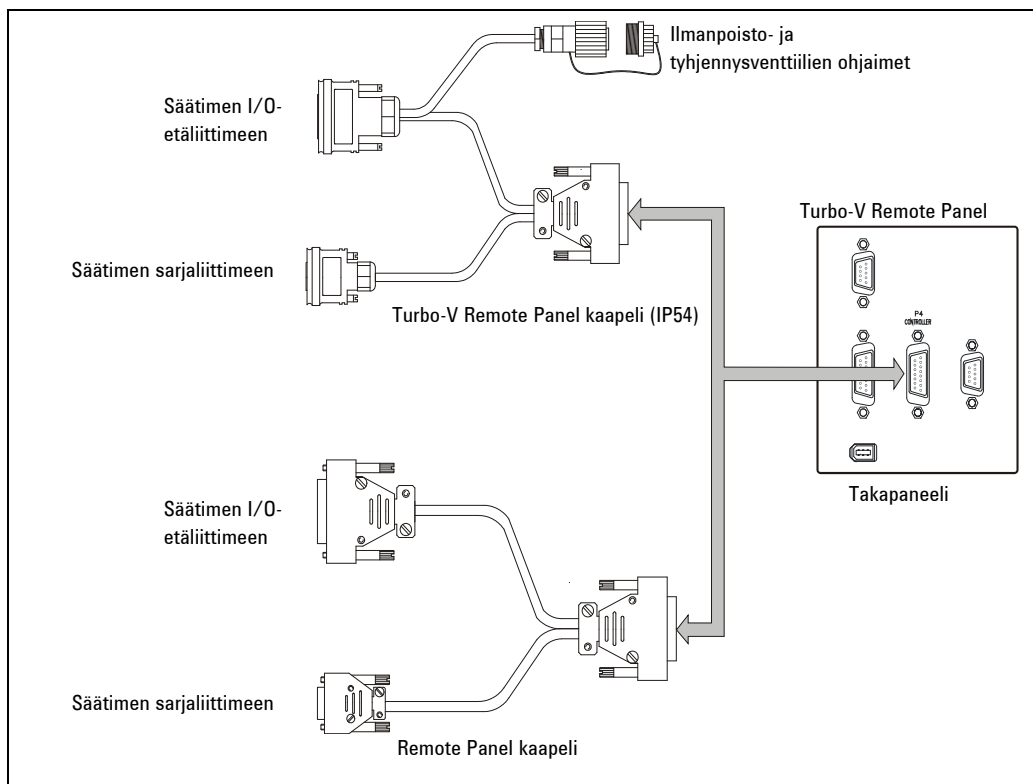
Jos käyttäjä haluaa ohjata ilmanpoisto- ja tyhjennysventtiileitä ulkoisia ohjaimia käyttämällä, on käytettävä versiota, jossa on IP54-kaapeli, asetettava säädin asentoon "valve setting" "on command" ja liitettävä kaapelit kuvassa esitetyllä tavalla. Venttiilit ovat tyypiltään normaalisti suljettuja: kun ilmanpoisto- ja tyhjennysventtiilin ja yhteisen päätteen välille syötetään 24 voltin virtaa, venttiilit avautuvat.



Kuva 1

## 10 Ohjekäsikirja

### Asennus



**Kuva 2** Liitännät

#### HUOM

Turbo-V Remote Panel voidaan asentaa pöydälle tai erilliseen telineeseen. Jäähdytysilman täytyy joka tapauksessa päästä kiertämään vapaasti laitteen sisällä. Älä asenna ja/tai käytä säädintä paikassa, jossa se altistuu ankarille sääolosuhteille (sade, pakkanen, lumi), pölylle tai haitallisille kaasuille, älä myöskään räjähdysalttiissa tai erityisen palovaarallisissa paikoissa.

Toiminnan aikana on ehdottomasti noudatettava seuraavia ympäristöolosuhteita:

- lämpötila: 5 °C – 45 °C
- suhteellinen kosteus: 0–95 % (ei kondensoituva).

Tietoa muista liitännöistä ja lisävarusteiden asentamisesta löydät osiosta ”Technical Information”.

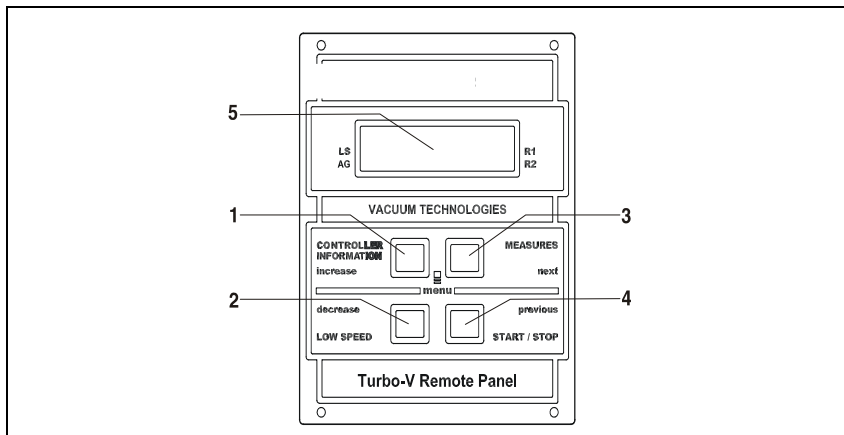
## Käyttö

Tässä kappaleessa esitellään keskeiset toiminnot. Tarkempia tietoja ja tietoa liitännöihin tai lisävarusteisiin liittyvistä toimenpiteistä löydät liitteen ”Technical Information” kappaleesta ”Use”.

## Ohjaimet, ilmaisimet ja liittimet

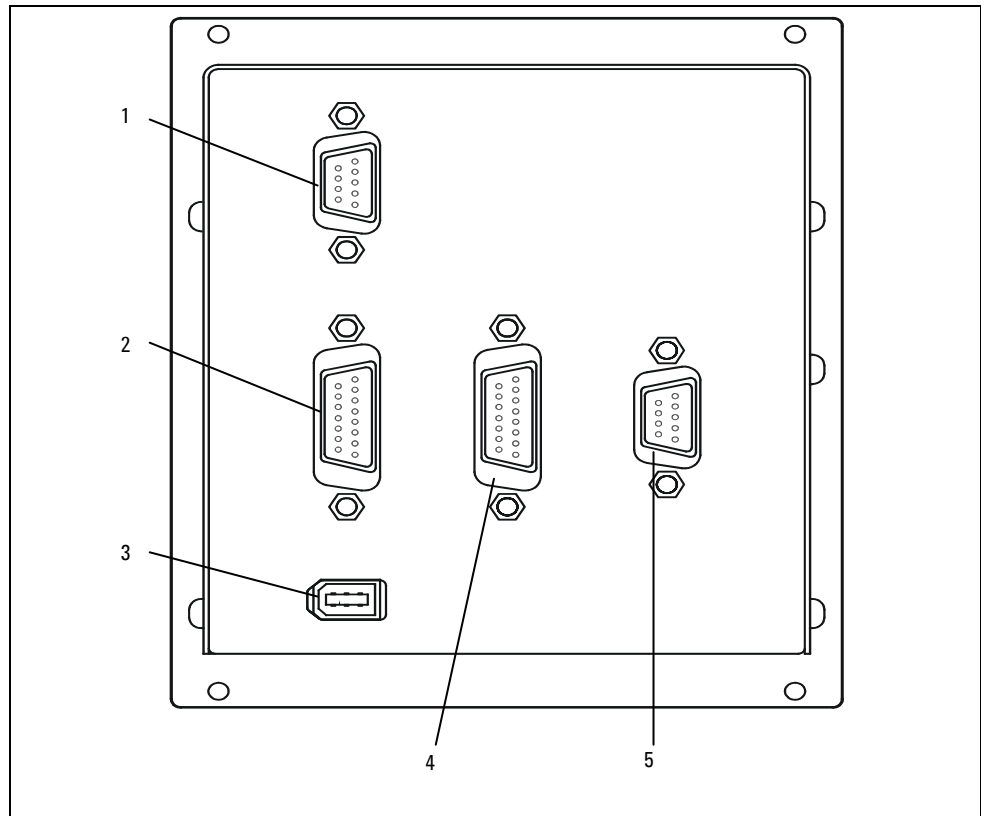
Seuraavassa on kuvattu säätimen etu- ja takapaneeli. Lisätietoa löydät osiosta ”Technical Information”.

## Beskrivelse av frontpanelet



**Kuva 3** Turbo-V Remote Panel

- 1 Painike, jolla näytölle haetaan parametrit cycle number, cycle time ja pump life.
- 2 Painike, jolla valitaan toimintatapa LOW SPEED. Käytössä ainoastaan silloin, kun ohjaustapa on valittu etupaneelistä. Jos painiketta painetaan kerran, pumppu pyörii standby-nopeudella. Jos painiketta painetaan uudelleen, LOW SPEED -toimintatapa poistuu käytöstä.
- 3 Painike, jolla näytölle haetaan parametrit pump current, pump temperature, pump power ja rotational speed. Se on aina käytössä riippumatta valitusta toimintatavasta. Kun painikkeita 3 ja 4 painetaan yhtä aikaa vähintään 2 sekunnin ajan, aktivoituu ohjelma, jolla voidaan ohjelmoida muutamia toimintaparametrejä.
- 4 Painike START- ja STOP/RESET-komentojen lähettämiseksi. Käytössä ainoastaan silloin, kun ohjaustapa on valittu etupaneelistä. Kun painiketta painetaan kerran, käynnistysvaihe aktivoituu; kun painiketta painetaan uudestaan, pumppu pysähtyy. Jos pumppu on pysähtynyt automaattisesti vian vuoksi, tätä painiketta on painettava ensin kerran säätimen nollaamiseksi ja sitten toisen kerran pumpun käynnistämiseksi uudelleen.
- 5 Aakkosnumeerinen nestekidenäyttö: pistematriisi, 2 riviä x 16 merkkiä.



**Kuva 4** Turbo-V Remote Panel takalevy

- 1 Loogisten signaalien sisääntuloliitin (kytkentäliitin toimitetaan asianmukaisella sulkukappaleella varustettuna).
- 2 Loogisten signaalien ulostuloliitin ja pumpun taajuusvalvonta sekä ohjelmoitava analoginen ulostulo.
- 3 Virransyötön ulostuloliitin (24 V tasavirta/anturi)
- 4 Virransyötön sisääntuloliitin pumpun säätimestä ja pumpun säätimen sarjaliitin.
- 5 Etäohjauksen sarjaliitin.

## **Valvojan päälle pano**

Valvoja käynnistyy asettamalla virtakaapeli pistorasiaan.

## **Pumpun käynnistys**

Käynnistääksenne pumpun pitää ottaa käyttöön interlock-sisääntulo ja painaa etupaneelin painiketta START.

## **Pumpun pysäyttäminen**

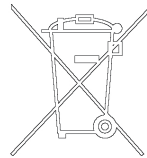
Pumppu pysähtyy painamalla etupanelissa olevaa STOP painonappia.

## Huolto

Turbo-V 301 Rack sarjan valvojat eivät kaipaa minkäänlaista huoltoa. Mahdolliset valvojaan tehtävät toimenpiteet tulee jättää aina valtuutetun henkilön tehtäviksi. Toimintahäiriön sattuessa on mahdollista käyttää Agilentin korjauspalvelua tai "Agilent advance exchange service" -palvelua, jolloin on mahdollista vaihtaa rikkiönyt valvoja ladattuun valvojaan. Mikäli valvoja täytyy romuttaa, toimikaa sen hävittämisessä kansallisten säädösten ja normien määrävällä tavalla.

## Hävittäminen

**Pakkausmerkinnöissä olevan WEEE-logon merkitys.** Alla näkyvä merkki on lisätty pakkaukseen EY:n ns. WEEE-direktiivin mukaisesti. Merkki (**koskee ainoastaan Euroopan Unionin jäsenmaita**) tarkoittaa, että tuotetta EI saa hävittää tavallisen kotitalous- tai teollisuusjätteen mukana, vaan se on toimitettava erilliseen keräyspisteeseen. Loppukäyttäjää kehoitetaan sen vuoksi ottamaan keräys- ja hävittämisprosessia varten yhteyttä laitteen toimittajaan, olipa se sitten laitteen valmistaja tai jälleenmyyjä, tarkastettuaan ensin kaupan sopimusehdot.



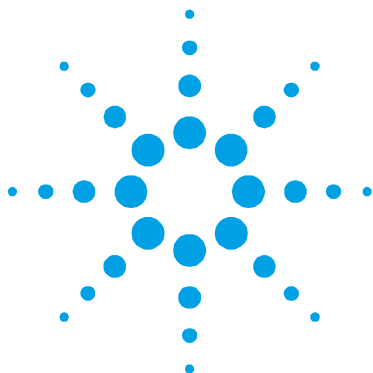
## Vianetsintä

Joidenkin toimintahäiriöiden yhteydessä valvojan itsemäärittelypiiri analysoi virheen, joka näkyy viesteinä, jotka on kuvailtu seuraavassa taulukossa.

VIESTI	VIKA	KORJAUSTOIMENPITEET
COMMUNICATION ERROR	Remote Panel ei kykene käynnistämään yhteyttä säätimen kanssa.	Tarkista Remote Panelin ja säätimen välinen sarjaliitäntä.
COMMUNICATION BROKEN	Remote Panel ei kykene jatkamaan yhteyttä säätimen kanssa.	Tarkista Remote Panelin ja säätimen välinen sarjaliitäntä.
XXXXXXX MODEL	Järjestelmä ei tue säädintä. Remote Panel toimii ainoastaan seuraavien pumppumallien kanssa: TV551, TV701 ja TV1001, joissa navigator controller sekä pumput TV2k-G ja TV3k-T.	Kytke Remote Panel järjestelmän tukemaan säätimeen.
CHECK CONNECTION TO PUMP	Toimintahäiriö pumpun ja valvojan liitännässä. Tai Pumpun lämpötila on alempi kuin 0 °C.	Tarkistakaa että pumpun ja valvojan välinen yhteyskaapeli on hyvin kiinnitetty päistään eikä sen varrella ole esteitä. Painakaa kaksi kertaa painonappia START jolloin pumppu käynnistyy.
WAITING INTERLOCK	Lukitussignaali (interlock) liittimessä P1 on aktiivinen johtuen liittimen J1 neulojen 3 ja 8 välillä tapahtuneen oikosulun keskeytyksestä tai ulkoisen lukitussignaalin (interlock) avautumisesta.	Palauttakaa liittimen J1 neulojen 3 ja 8 välinen oikosulku tai sulkekaa ulkopuolinen lukitussignaali (interlock).
PUMP OVERTEMP.	Ylemmän laakerin tai pumpun lämpötila on ylittänyt 60 °C.	Odottakaa että lämpötila putoaa kynnysarvon alapuolelle. Painakaa painonappia START kaksi kertaa jolloin pumppu käynnistyy.
CONTROLLER OVERTEMP.	Valvojan muuntajan lämpötila on 70 °C. Tai Valvojan lämmittimen lämpötila on korkeampi kuin 60 °C.	Odottakaa, että lämpötila putoaa kynnysarvon alapuolelle. Painakaa painonappia START kaksi kertaa jolloin pumppu

VIESTI	VIKA	KORJAUSTOIMENPITEET
		käynnistyy.
TOO HIGH LOAD	Pumpun absorptiovirta ylittää ohjelmoidun arvon.	Tarkistakaa, että pumpun roottori pyörii vapaasti. Painakaa painonappia START kaksi kertaa jolloin pumppu käynnistyy.
SHORT CIRCUIT	Normaalityöskentelyn aikana (käynnistyksen jälkeen) poistoliitin on oikosulussa.	Tarkistakaa pumpun ja valvojan välinen liitos. Painakaa painonappia START kaksi kertaa jolloin pumppu käynnistyy.
SYSTEM OVERRIDE	Kaukokytimestä tuleva hälytysignaali on pysäyttänyt pumpun.	Irroitakaa valvojan sähkökaapeli verkkovirrasta ja korjatkaa hälytyksen aiheuttaja. Kytkekää sähkökaapeli uudelleen ja painakaa painonappia START kaksi kertaa jolloin pumppu käynnistyy.
OVERVOLTAGE	Valvojan sähkövirran syötössä on ilmennyt ongelma tai valvojan saama signaali on väärä.	Painakaa painonappia START kaksi kertaa, jolloin pumppu käynnistyy. Mikäli viesti näkyy uudelleen kääntykää Agilent huoltopalvelun puoleen.
POWER FAIL	Vika pumpun, valvojan virransyöttöosassa.	Kääntykää Agilent-yhtiön puoleen Huoltoa tarvittaessa.

## **10 Ohjekäsikirja** **Vianetsintä**



## 11 Felhasználói Kézikönyv

Általános információk	154
Tárolás	156
A telepítésre való előkészítés	156
Telepítés	157
Használat	159
Vezérlések, kijelzők és csatlakozók	159
Az előlap leírása	160
A vezérlő bekapcsolása	162
A szivattyú beindítása	162
A szivattyú leállítása	162
Karbantartás	163
Megsemmisítés	163
Hibaüzenetek	164

Az eredeti utasítás fordítása



## Általános információk

Ez a berendezés hivatásos használatra tervezett. A felhasználónak, a berendezés használatát megelőzően, gondosan el kell olvasnia a felhasználói kézikönyvet és minden egyéb, a Agilent által szolgáltatott információt. A Agilent elhárít magától minden esetleges felelősséget, mely a felhasználói utasítások teljes vagy részleges figyelmen kívül hagyásához, a nem kiképzett személyzet által történő nem megfelelő használathoz, meghatalmazás nélküli beavatkozásokhoz, vagy a specifikus nemzeti irányelvekkel ellentétes használathoz kötődik.

A mikroprocesszor által vezérelt Turbo-V Remote Panel segítségével a felhasználó előlapról vezérelhet egy "controller on-board" típusú szivattyút:

- TV551 Navigator vezérlővel
- TV701 Navigator vezérlővel
- TV1001 Navigator vezérlővel
- TV2k-G on-board vezérlővel
- TV3k-T on-board vezérlővel
- TV3k-G on-board vezérlővel
- TV70 Navigator vezérlővel
- TV81 Navigator vezérlővel
- TV301 Navigator vezérlővel
- TV303 Navigator vezérlővel

Az elkövetkező szakaszok tartalmazzák mindazokat az ismereteket, melyek a kezelő biztonságát hivatottak garantálni a berendezés használata során. Részletes információk találhatóak a "Technical Information" című mellékletben.

**E kézikönyv az alábbi megegyezésszerű jelöléseket alkalmazza:**

**FIGYELEM!**

A figyelmeztető üzenetek azon eljárások előtt kerülnek megjelenítésre, melyek figyelmen kívül hagyása a berendezés károsodásához vezethet.

---

**VESZÉLY!**



A veszélyt jelző üzenetek olyan eljárásra vagy specifikus tevékenységre hívják fel a kezelő figyelmét, melyek, ha nem megfelelő módon kerülnek végrehajtásra, súlyos személyi sérüléseket okozhatnak.

---

**MEGJEGYZÉS**

A megjegyzések fontos, a szövegből kivonatolt információkat tartalmaznak.

---

## Tárolás

A vezérlők szállítása és tárolása során az alábbi környezeti feltételeket kell biztosítani:

- hőmérséklet: -20 °C és +70 °C között
- relatív nedvességtartalom: 0 – 95 % (nem lecsapódó)

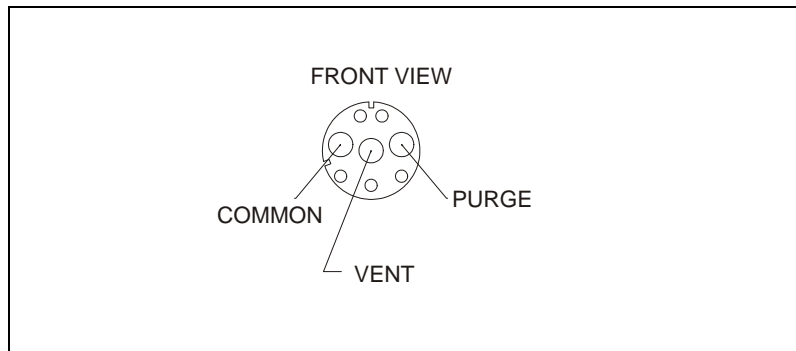
## A telepítésre való előkészítés

A vezérlő speciális védőcsomagolásban kerül leszállításra; ha ez károsodás jeleit mutatja, melyek a szállítás során keletkezettek, kapcsolatba kell lépni a helyi forgalmazó irodával. A kicsomagolás során különleges figyelmet kell fordítani arra, hogy a vezérlő le ne essék, és ne szenvedjen ütődéseket. A csomagolóanyagot nem szabad szétszórni a környezetben. Az anyag teljes mértékben újrafelhasználható, s megfelel a környezetvédelemmel kapcsolatos EGK 85/399-es irányelvnek.

## Telepítés

A Turbo-V Remote Panel a vezérlővel az e célra szolgáló kábellel kell összekötni (lásd a rendelhető alkatrészek táblázatát), az alábbi ábrán látható módon. Minthogy a szükséges feszültséget / jeleket e kábel biztosítja, nincs szükség semmiféle külső táplálásra.

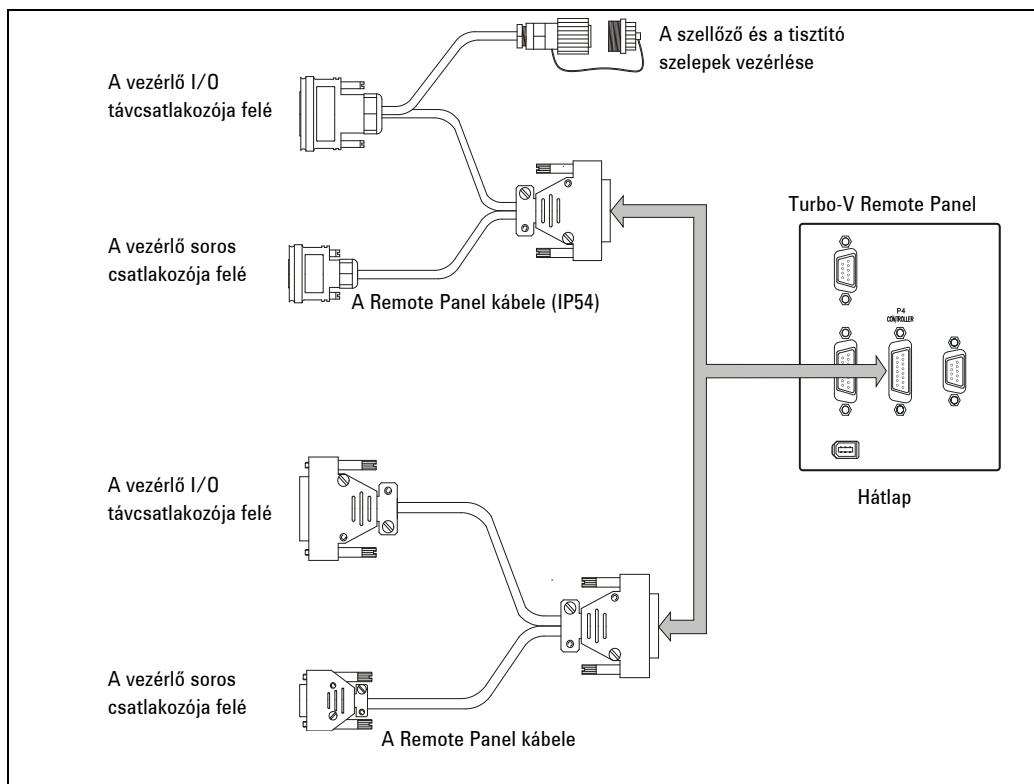
Ha a kezelő a szelőző és tisztító szelepeket külső vezérléssel kívánja irányítani, az IP54 kábellel ellátott változatot kell használni, a vezérlőt a "valve setting" "on command" állásba kell helyezni és a kábeleket az ábrán látható módon bekötni. A szelepek alaphelyzetben zártak (n.c.); 24V-ot adva a szellőző és a tisztító szelepek valamint a közös lezárás közé, a szelepek nyitott állapotba kerülnek.



Ábra 1

## 11 Felhasználói Kézikönyv

### Telepítés



Ábra 2 Bekötések

#### MEGJEGYZÉS

A vezérlő telepítésre kerülhet egy asztalon, vagy egy megfelelő rack belsejében. Minden esetben szükséges azonban, hogy a hűtőlevegő szabadon áramolhasson a berendezés belsejében. A vezérlőt nem szabad légköri hatásoknak (eső, fagy, hó) kitett, poros, maró hatású gázt tartalmazó, robbanás- vagy fokozottan tűzveszélyes környezetben telepíteni és/vagy használni.

Üzem közben az alábbi környezeti feltételeket kell biztosítani:

- hőmérséklet: 5 °C és +45 °C között
- relatív nedvességtartalom: 0 – 95 % (nem lecsapódó).

A többi összekötéssel és az opcionális kiegészítő felszerelések telepítésével kapcsolatosan lásd a “Technical Information” című alfejezetet.

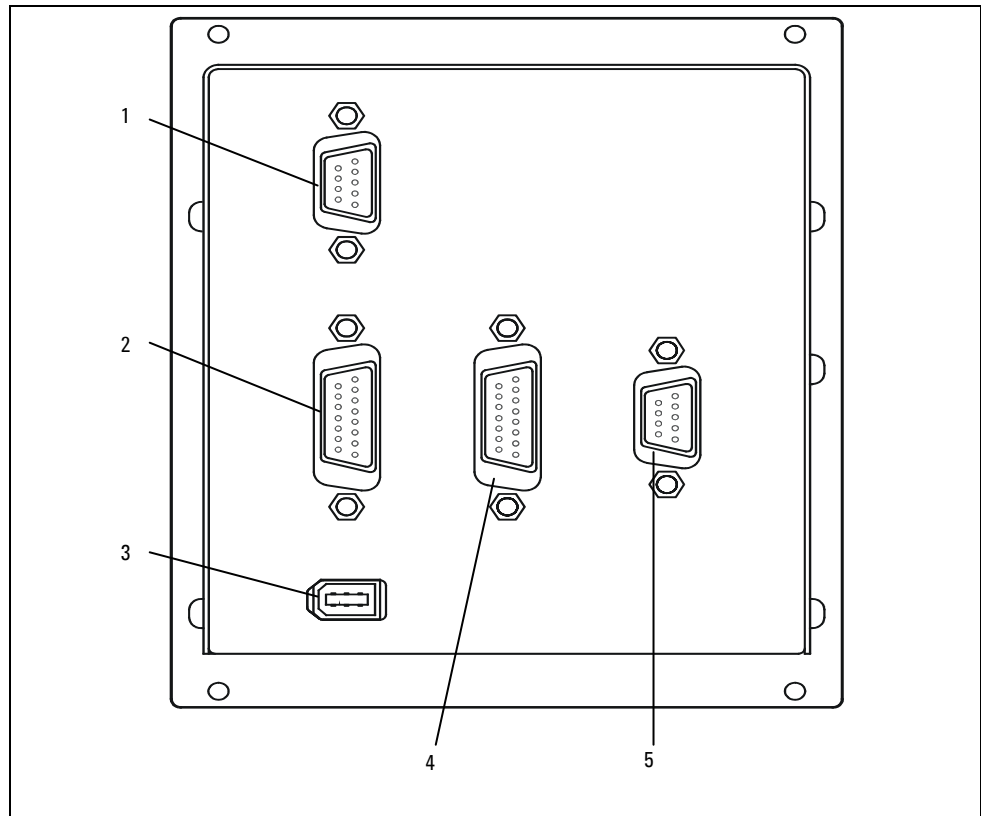
## Használat

Ebben a szakaszban a fontosabb működési eljárások kerülnek ismertetésre. További részletek és az összekötéseket vagy az opcionális elemeket érintő eljárások találhatóak a “Technical Information” melléklet “Use” című szakaszában.

## Vezérlések, kijelzők és csatlakozók

Az alábbiakban bemutatásra kerül a vezérlő elő- és hátlapja. További részletek találhatóak a “Technical Information” című szakaszban.





**Ábra 4** A Turbo-V Remote Panel hátlapja

- 1 Logikai jelek bemeneti csatlakozói (az összekötő csatlakozó megfelelő lezáró átkötéssel van ellátva).
- 2 Csatlakozó a kimeneti logikai jelek, a szivattyúfrekvencia ellenőrzése, valamint programozható analóg kimenet számára.
- 3 Kimeneti tápcsatlakozó (24 V egyenfeszültség a vakuum-mérőhöz (gauge)).
- 4 A szivattyú vezérlőjének bemeneti tápcsatlakozója, illetve csatlakozó a szivattyú vezérlőjével való soros kommunikációhoz.
- 5 Soros csatlakozó távvezérléshez.

## A vezérlő bekapcsolása

A vezérlő bekapcsolásához elegendő a hálózati kábelt a hálózati aljzatba illeszteni.

## A szivattyú beindítása

A szivattyú beindításához engedélyezni kell az interlock bemenetet és megnyomni az előlapon a START nyomógombot.

## A szivattyú leállítása

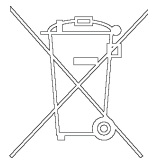
A szivattyú leállításához meg kell nyomni az előlapon található STOP nyomógombot.

## Karbantartás

A Turbo-V 301 Rack sorozatú vezérlők semmiféle karbantartást sem igényelnek. Valamennyi beavatkozást meghatalmazott személyzet hajthat csak végre. Meghibásodás esetén igénybe vehető a Agilent javítási szolgáltatása, vagy a "Agilent advanced exchange service", mely a meghibásodott helyett, egy javított vezérlőt biztosít. Amikor egy vezérlőt le kell selejtezni, megsemmisítésekor a tárgykörbe tartozó nemzeti normáknak megfelelően kell eljárni.

## Megsemmisítés

**A címkén jelenlévő "WEEE" logo jelentése.** Az alább látható szimbólum az EK "WEEE" elnevezésű irányelvvel összhangban kerül alkalmazásra. Ez a szimbólum (**mely csak az Európai Közösség országaiban érvényes**), azt jelzi, hogy a termék, melyen megtalálható, NEM kerülhet közönséges háztartási vagy ipari hulladékkal együtt megsemmisítésre, hanem azt egy szelektív hulladékgyűjtő rendszerbe kell továbbítani. A végfelhasználónak, a gyűjtési és megsemmisítési eljárás beindítása céljából, az adásvételi szerződés határidőinek és feltételeinek alapos áttanulmányozása után, fel kell vennie a kapcsolatot a berendezés szállítójával, legyen ez a gyártó vagy egy viszonteladó.



## Hibaüzenetek

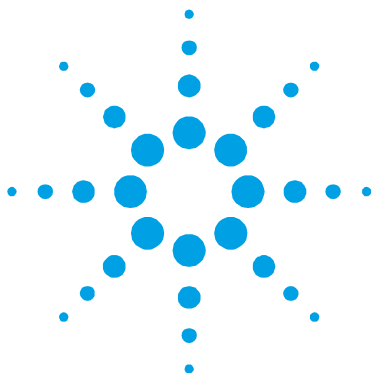
Néhány meghibásodási esetben a vezérlő öndiagnosztikai áramkörei hibaüzenetet jelenítenek meg (lásd az alábbi táblázatot).

ÜZENET	LEÍRÁS	JAVÍTÁSI TEVÉKENYSÉG
COMMUNICATION ERROR	A Remote Panel nem tud kommunikációba lépni a vezérlővel.	Ellenőrizze a Remote Panel és a vezérlő között a soros kapcsolat meglétét.
COMMUNICATION BROKEN	A Remote Panel nem tudja folytatni a kommunikációt a vezérlővel.	Ellenőrizze a Remote Panel és a vezérlő között a soros kapcsolat meglétét.
XXXXXXX MODEL	Inkompatibilis vezérlő. A Remote Panel csak az alábbi szivattyúkkal működik: TV551 navigator vezérlővel, TV701 navigator vezérlővel, TV1001 navigator vezérlővel, TV2k-G e TV3k-T.	A Remote Panelt csatlakoztassa kompatibilis vezérlőhöz.
CHECK CONNECTION TO PUMP	Összeköttetési hiba a szivattyú és a vezérlő között, vagy a szivattyú hőmérséklete 0 °C alatt van.	Ellenőrizni, hogy a szivattyú és a vezérlő közti összekötő kábel mindkét végén jól van-e rögzítve és nincs e megszakadva. A szivattyú újraindításához kétszer megnyomni a START nyomógombot.
WAITING INTERLOCK	A P1 csatlakozó reteszelés jelen jele aktív a J1 csatlakozó 3. és 8. pin-je közti rövidzár megszakadása, vagy a külső reteszelő jel nyitott állapota következtében.	Visszaállítani a J1 csatlakozó 3. és 8. pin-je közti rövidzárát, vagy zárni a külső reteszelő jelet.
PUMP OVERTEMP.	A felső csapágy vagy a szivattyú hőmérséklete meghaladta a 60 °C-t.	Várakozni, míg a hőmérséklet a küszöbérték alá esik. A szivattyú újraindításához kétszer megnyomni a START nyomógombot.
CONTROLLER OVERTEMP.	A vezérlő ának hőmérséklete túllépte a 70 °C-t, vagy	Várakozni, míg a hőmérséklet a küszöbérték alá esik. A szivattyú újraindításához kétszer megnyomni a START

ÜZENET	LEÍRÁS	JAVÍTÁSI TEVÉKENYSÉG
	a vezérlő hűtésének hőmérséklete túllépte a 60 °C-t.	nyomógombot.
TOO HIGH LOAD	A szivattyú által felvett áram meghaladja a programozott értéket.	Ellenőrizni, hogy a szivattyú forgórésze szabadon foroghat-e. A szivattyú újraindításához kétszer megnyomni a START nyomógombot.
SHORT CIRCUIT	Normál üzem közben (az indítási fázist követően) a kimeneti összeköttetés zártos.	Ellenőrizni az összeköttetéseket a szivattyú és a vezérlő között. A szivattyú újraindításához kétszer megnyomni a START nyomógombot.
SYSTEM OVERRIDE	A szivattyú leállt egy távérintkezőtől származó vészjel hatására.	Kihúzni a vezérlő hálózati csatlakozóját és megszüntetni a vészjelzés okát. Újra csatlakoztatni a hálózati kábelt, és kétszer megnyomni a START nyomógombot a szivattyú újraindításához.
OVERVOLTAGE	Meghibásodás történt a vezérlő tápellátási részegységében, vagy a vezérlő egy hamis jelet kapott.	A szivattyú újraindításához kétszer megnyomni a START nyomógombot. Ha az üzenet ismétlődik, a Agilent-től karbantartási beavatkozást kérni.
POWER FAIL	Meghibásodás a szivattyú vagy a vezérlő tápellátási részegységében.	A Agilent-től karbantartási beavatkozást kérni.

## **11 Felhasználói Kézikönyv**

### **Hibaüzenetek**



## 12 Podrecznik Instrukcji

Informacje ogólne	168
Magazynowanie	170
Przygotowanie do instalacji	170
Instalacja	171
Użytkowanie	173
Sterowniki, Wskazniki i Łączniki	173
Opis pulpitu frontального	174
Zaswiecenie kontrolera	176
Uruchomienie pompy	176
Zatrzymanie pompy	176
Konserwacja	177
Przetworstwo odpadów	177
Błędne informacje	178

Tłumaczenie instrukcji oryginalnej



## Informacje ogólne

Urządzenie to jest przeznaczone do użytku zawodowego. Przed przystąpieniem do korzystania użytkownik powinien wnikliwie przesledzić zarówno ten podrecznik zawierający instrukcje jak również każda inna dostarczona przez Agilent informację. Agilent uchyła się od jakiegokolwiek odpowiedzialności w przypadku całkowitego lub częściowego nie przestrzegania swoich instrukcji, przeznaczenia urządzenia niezgodnie z jego przeznaczeniem i jego obsługi przez personel uprzednio nie przeszkolony, naprawy urządzenia bez odpowiedniej autoryzacji lub przeznaczenia niezgodnego z obowiązującymi, specyficznymi normami krajowymi. Turbo-V Remote Panel jest kontrolowany przy pomocy mikroprocesora i pozwala użytkownikowi sterować pompą typu “controller on-board” poprzez pulpit frontalny:

TV551 z kontrolerem sterowanym

TV701 z kontrolerem sterowanym

TV1001 z kontrolerem sterowanym

TV2k-G z controller on-board

TV3k-T z controller on-board

TV3k-G z controller on-board

TV70 z kontrolerem sterowanym

TV81 z kontrolerem sterowanym

TV301 z kontrolerem sterowanym

TV303 z kontrolerem sterowanym

Następne paragrafy zawierają wszystkie niezbędne informacje dotyczące gwarancji bezpieczeństwa personelu obsługującego urządzenie. Bardziej szczegółowe informacje są zawarte w dodatku do ”Technical Information”.

**Podrecznik ten opiera się na następujących konwencjach:**

**UWAGA!**

Sygnaly wzrokowe ukazujące się przed procedurą jeżeli zostaną zlekceważone mogą spowodować uszkodzenia urządzenia.

---

**ZAGROZENIE!**



Sygnaly ostrzegające o zagrożeniu powinny zwrócić uwagę operatora na odpowiednią specyficzną procedurę którą jeżeli nie jest przestrzegana w sposób jednoznaczny może przyczynić się do poważnych obrażeń osoby obsługującej.

---

**PRZYPIS**

Przypisy zawierają najważniejsze informacje wyciągnięte z tekstu.

---

## Magazynowanie

Kontrolery podczas transportu i przechowywania w magazynie powinny mieć zapewnione następujące warunki otoczenia:

- temperatura: od -20 °C do +70 °C
- wilgotność względna: od 0-95 %(nie kondensująca)

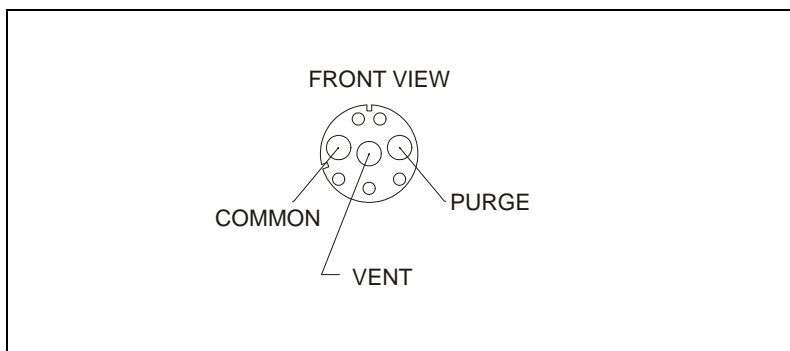
## Przygotowanie do instalacji

Turbo-V Remote Panel zostaje dostarczony w specjalnym, ochronnym opakowaniu; jeżeli opakowanie prezentuje jakiegokolwiek uszkodzenia które mogłyby powstać podczas transportu wówczas należy niezwłocznie powiadomić najbliższy, lokalny punkt sprzedaży. Podczas operacji rozpakowywania należy zachować specjalną ostrożność żeby nie spowodować jego upadku lub nie narazić kontrolera na uderzenia. Nie wyrzucać opakowania ponieważ jest ono całkowicie recyklowalne i odpowiada normom CEE 85/399 dotyczącym ochrony środowiska

## Instalacja

Turbo-V Remote Panel należy podłączyć do kontrolera przy pomocy odpowiedniego kabla ( patrz tabela części dostarczanych na zamówienie), w sposób zilustrowany na poniższym rysunku. Ponieważ moc/i i wymagane sygnały są dostarczane przez ten kabel zbytecznym jest jakiekolwiek dodatkowe zasilanie zewnętrzne.

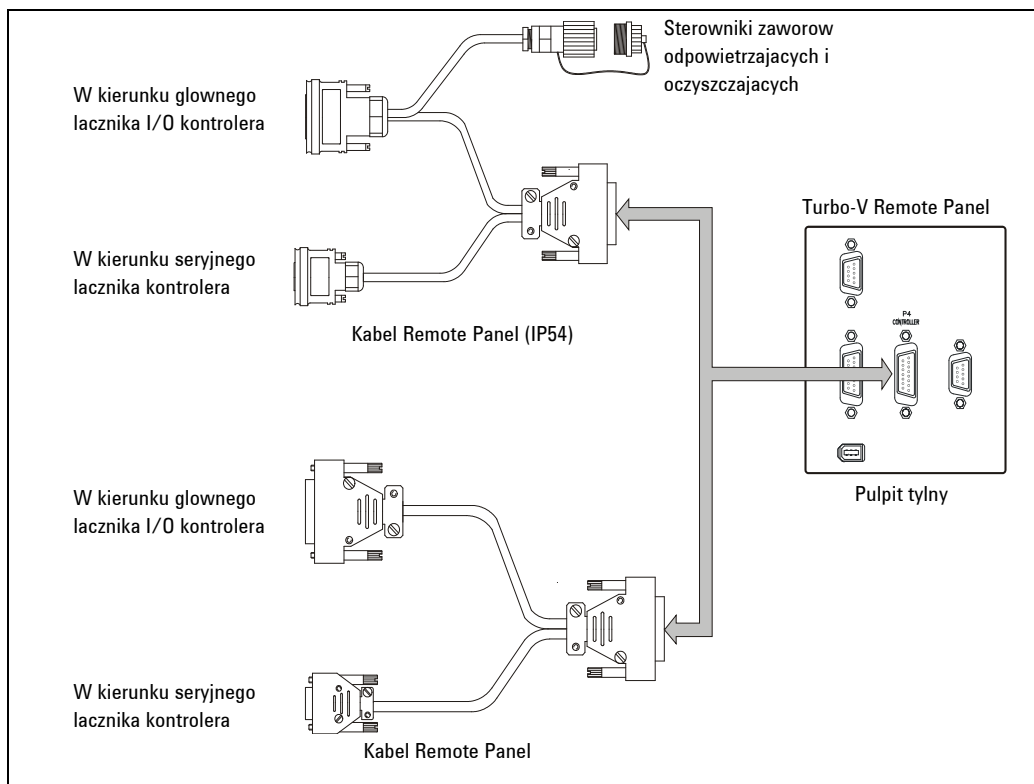
Jeżeli operator zamierza sterować zaworem odpowietrzającym i zaworem oczyszczającym złogi powinien posłużyć się do tej operacji kablem IP54, ustawiając kontroler na “valve setting” “on command” w sposób przedstawiony na szkicu. Zawory są nie podłączone; dostarczając napięcie 24V pomiędzy zawór odpowietrzający, zawór oczyszczający i wspólnym zacisk zawory się otworzą.



Posta 1

## 12 Podrecznik Instrukcji

### Instalacja



#### Posta 2 Connections

##### PRZYPIS

Turbo-V Remote Panel moze byc zainstalowany na stole lub wewnatrz odpowiedniego Rack. W obydwu przypadkach nalezy pamietac aby wokol kontrolera zapewnic swobodny przeplyw powietrza dla prawidlowego chlodzenia. Nie instalowac i/lub uzytkowac kontrolerow w miejscach narazonych na dzialanie warunkow atmosferycznych (jak deszcz, mroz, snieg), pylow, gazow agresywnych w srodowisku wybuchowym lub o duzym stopniu zagrozenia pozarowego.

Podczas pracy niezbedne jest zachowanie nastepujacych warunkow otoczenia:

- temperatura: od 5 °C do +45 °C;
- wilgotnosc wzgledna: 0 – 95 % (nie kondensujaca).

Sposoby innego rodzaju polaczenia oraz instalowania akcesoriow dodatkowych opisane sa w jednym z rozdzialow "Technical Information".

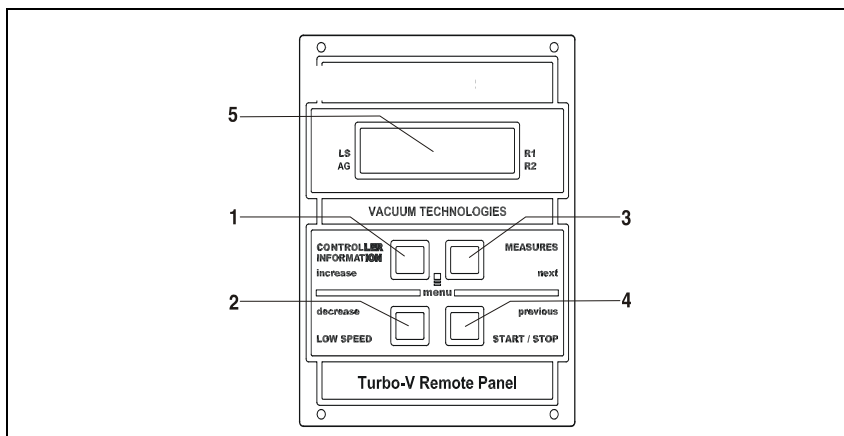
## Uzytkowanie

Paragraf ten zawiera podstawowe procedury operatywne. Dla uzyskania dodatkowych bardziej szczegolowych informacji lub dostepu do procedur dotyczacych polaczen lub dodatkowych akcesoriow konsultowac paragraf „Use“ znajdujacy sie w dodatku do „Technical Information“.

## Sterowniki, Wskazniki I Laczniki

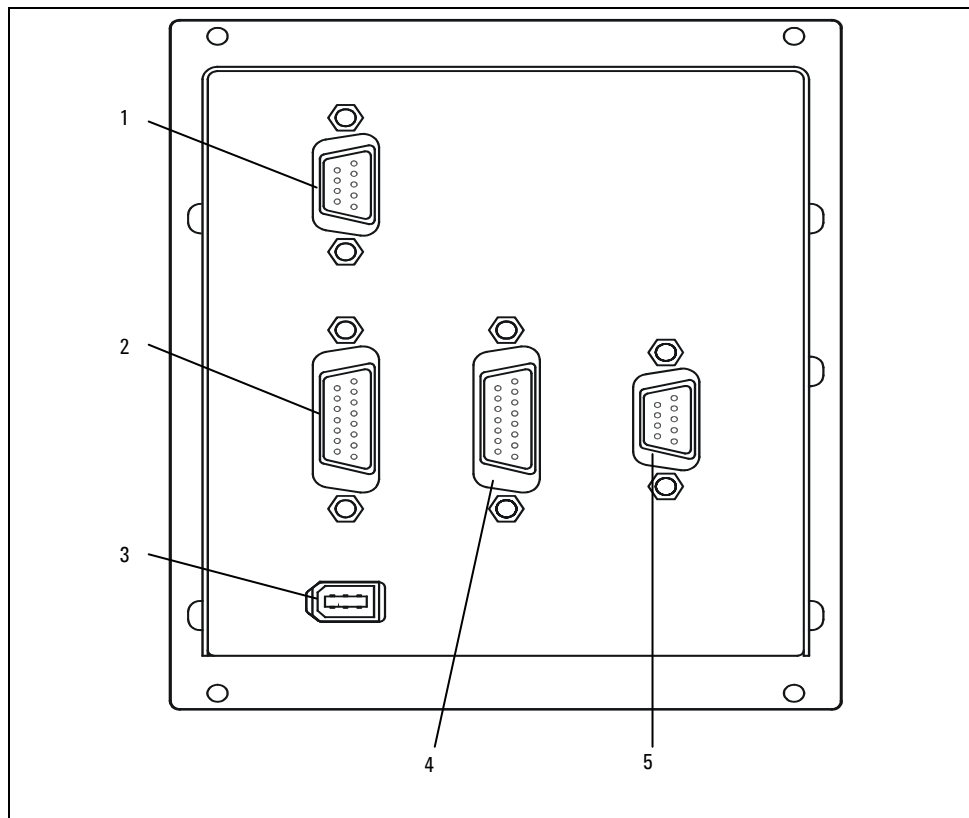
Ponizej jest przedstawiony pulpit sterowniczy przedni i tylny kontrolera. Dla uzyskania dokladniejszych szczegolow skonsultowac odpowiedni rozdzial "Technical Information".

## Opis pulpitu frontalnego



### Posta 3 Turbo-V Remote Panel

- 1 Przycisk sluzacy do przywołania na display parametrow cycle number, cycle time i pump life.
- 2 Przycisk do wybierania sposobu LOW SPEED. Dziala tylko wtedy kiedy sposob sterowania jest wybierany na pulpicie frontalnym. Naciskajac jeden raz pompa obraca sie z szybkościa "stand-by". Naciskajac ponownie następuje wylaczenie sposobu LOW SPEED.
- 3 Przycisk sluzacy do przywołania na display parametrow pump current, pump temperature, pump power i rotational speed. Jest zawsze aktywny niezaleznie od wybranego sposobu dzialania. Naciskajac jednoczesnie przyciski 3 i 4 przez conajmniej 2 sekundy uruchamia sie program przy pomocy ktorego jest mozliwe programowanie niektorych parametrow operatywnych.
- 4 Przycisk do przesyłania sygnalu START, STOP lub RESET. Dziala tylko wtedy kiedy sposob sterowania jest wybierany na pulpicie frontalnym. Naciskajac jeden raz uaktywnia sie faza uruchomienia; naciskajac ponownie zatrzymuje sie pompe. Jezeli pompa zatrzyma sie automatycznie z powodu uszkodzenia nalezy nacisnac przycisk jeden raz zeby zresetowac kontrolera i poraz drugi zeby uruchomic pompe.
- 5 Display alfanumeryczny na cieklych kryształach: matryca punktowana, 2 linie x 16 czcionek.



**Posta 4** Pulpit tylni Turbo-V Remote Panel

- 1 Łącznik wejściowy sygnałów logicznych (łącznik bliźniaczy jest dostarczony ze specjalnym mostkiem zamykającym).
- 2 Łącznik wyjściowy sygnałów logicznych I nadzór nad częstotliwością na wyjściu podczas zaprogramowanej wielokrotnej pracy pompy.
- 3 Łącznik zasilania na wyjściu (kalibru 24 Vcc ).
- 4 Łącznik wejściowy zasilania kontrolera pompy oraz łącznik seryjny wraz z kontrolem pompy.
- 5 Łącznik seryjny zdalnie sterowany.

## **Zaswiecenie kontrolera**

Do zaswiecenia kontrolera wystarczy wprowadzić kabel zasilający do gniazdka sieci.

## **Uruchomienie pompy**

Zeby uruchomic pompe nalezy odblokowac wejscie interlock i nacisnac znajdujacy sie na pulpicie frontalnym przycisk START .

## **Zatrzymanie pompy**

Do zatrzymania pompy wystarczy nacisnac przycisk STOP znajdujacy sie na pulpicie frontalnym.

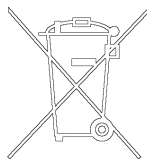
## Konservacja

Kontroler z serii Turbo-V 301 Rack nie wymaga zadnej konserwacji. Jakakolwiek interwencja moze byc dokonywana wytlacalnie przez osoby autoryzowane. W przypadku uszkodzenia mozliwe jest korzystanie z serwisow naprawczych Agilent lub "Agilent advance exchange service", w ktorym mozna otrzymac kontroler zregenerowany zastepujac go tym uszkodzonym. W przypadku decyzji zlomowania nalezy przeprowadzic operacje respektujac krajowe specyficzne normy.

## Przetworstwo odpadow

### Objasnienie znajdujacego sie na etykiecie znaczenia logo "WEEE".

Uzyty ponizej symbol jest zgodny z wymogiem zarzadzenia "WEEE" Unii Europejskiej. Symbol ten (**prawomocny tylko w krajach Unii Europejskiej**) oznacza, ze wyrob ktory nim zostal oznaczony NIE moze byc przetworzony jako odpad razem z innymi domowymi lub przemyslowymi natomiast musi byc skladowany w miejscu przeznaczonym dla odpadow do przerobki zroznicowanej . Dlatego tez poleca sie uzytkownikowi, , po uprzedniej weryfikacji terminu i warunkow zawartych w kontrakcie sprzedazy, nawiazac kontakt z dostawca lub sprzedawca urzadzenia w celu uruchomienia procesu zbiorke i przerobu.



## Bledne informacje

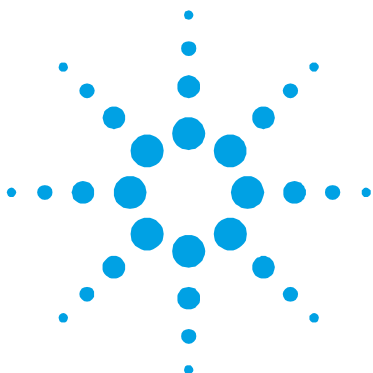
W niektórych przypadkach uszkodzen obwod do autodiagnozy kontrolera pokazuje niektore blendne informacje wymienione w ponizszej tabeli.

<b>INFORMACJA</b>	<b>OPIS</b>	<b>AKCJA KORYGUJACA</b>
COMMUNICATION ERROR	Il Remote Panel nie jest w stanie nawiązać łączności z kontrolerem	Sprawdzić łącze seryjne pomiędzy Remote Panel i kontrolerem
COMMUNICATION BROKEN	Il Remote Panel nie jest w stanie kontynuować łączności z kontrolerem	Sprawdzić łącze seryjne pomiędzy Remote Panel i kontrolerem
XXXXXXX MODEL	Kontroler bez wspomagania. Remote Panel działa jedynie w połączeniu z następującymi pompami: TV551 z kontrolerem sterowanym, TV701 z kontrolerem sterowanym, TV1001 z kontrolerem sterowanym, TV2k-G e TV3k-T.	Podłączyć Remote Panel do kontrolera z zasilaniem
CHECK CONNECTION TO PUMP	Zakłócenia na połączeniu między pompa i kontrolerem. Lub Temperatura pompy jest poniżej 0 °C.	Zweryfikować czy kabel łączący pompe z kontrolerem jest dobrze zamocowany na obydwu końcach I czy nie jest przerwany. Nacisnąć dwa razy przycisk START dla ponownego uruchomienia pompy.
WAITING INTERLOCK	Jest aktywny sygnał Interlock na łączniku P1 spowodowany zwarcie między pin 3 i pin 8 łącznika J1, Lub na skutek otwarcia sygnału zewnętrznego interlock.	Dokonać naprawy obwodu elektrycznego pomiędzy pin 3 i pin 8 łącznika J 1 lub zamknąć sygnał interlock zewnętrzny.
PUMP OVERTEMP.	Temperatura górnego łożyska lub pompy przekroczyła 60 °C.	Odczekać aż temperatura się obniży poniżej ustalonego poziomu. Nacisnąć dwa razy przycisk START dla ponownego uruchomienia pompy.
CONTROLLER OVERTEMP.	Temperatura w kontrolerze	Odczekać aż temperatura się

INFORMACJA	OPIS	AKCJA KORYGUJACA
	przekroczyła 70 °C. Lub Temperatura grzejnika nie przekracza 60 °C.	obnizy ponizej ustalonego poziomu. Nacisnac dwa razy przycisk START dla ponownego uruchomienia pompy.
TOO HIGH LOAD	Absorpcja pradu przez pompe jest wyzsza od zaprogramowanej.	Sprawdzic czy wirnik nie ma przeszkod w swobodnym obrocie. Nacisnac dwa razy przycisk START dla ponownego uruchomienia pompy.
SHORT CIRCUIT	Podczas normalnej pracy (po fazie uruchomienia) na wyjsciu sa zwarcia.	Sprawdzic polaczenie pompy z kontrolerem Nacisnac dwa razy przycisk START dla ponownego uruchomienia pompy.
SYSTEM OVERRIDE	Pompa sie zatrzymala na skutek sygnalu zagrozenia pochodzacego od pierwotnego kontaktu.	Odlaczyc kabel zasilania kontrolera i sprawdzic przyczynę powstania sygnalu zagrozenia. Podlaczyc ponownie kabel zasilajacy i nacisnac dwa razy przycisk START dla ponownego uruchomienia pompy.
OVERVOLTAGE	Nastapilo uszkodzenie w sekcji zasilania kontrolera lub kontroler otrzymal sygnal nieprawidlowy.	Nacisnac dwa razy przycisk START dla ponownego uruchomienia pompy. Jezeli sygnal sie powtorzy zwrocic sie do Agilent o przeprowadzenie konserwacji.
POWER FAIL	Uszkodzenie w sekcji zasilania pompy kontrolera.	Zwrocic sie do Agilent w celu naprawy.

## **12 Podrecznik Instrukcji**

### **Bledne informacje**



## 13 Návod k Použití

Všeobecné informace	182
Uskladnění	184
Příprava k instalaci	184
Instalace	185
Použití	187
Ovládání, indikátory a konektory	187
Popis čelního panelu	188
Zapalování Kontroléru	190
Spouštění čerpadla	190
Zastavení čerpadla	190
Údržba	191
Likvidace	191
Chybné zprávy	192

Překlad originálního návodu



## Všeobecné informace

Tento přístroj je určen pro profesionální potřebu. Uživatel si musí tuto příručku před použitím velmi pozorně přečíst, ale i každou další aktualizaci, kterou podnik Agilent bude předávat. V důsledku nedostatečného i částečně nesprávného použití, nebo v případě chyb nevyškoleného personálu či neoprávněných zásahů a neodpovídajících specifickým státním normám, Agilent nepřijímá a nenese žádnou zodpovědnost.

Turbo-V Remote Panel je kontrolovaný mikro-procesorem a uživatel může na čelním panelu řídit čerpadlo typu "controller on-board":

TV551 s Navigator controller

TV701 s Navigator controller

TV1001 s Navigator controller

TV2k-G s controller on-board

TV3k-T s controller on-board

TV3k-G s controller on-board

TV70 s Navigator controller

TV81 s Navigator controller

TV301 s Navigator controller

TV303 s Navigator controller

V následujících odstavcích jsou uvedeny veškeré nutné informace, týkající se ochrany operátora a bezpečnostních podmínek během provozu přístroje. Detailní informace se nacházejí v dodatku "Technical Information".

Tato příručka používá následující běžná označení:

**POZOR!**

Tyto zprávy se objevují před jakoukoliv procedurou, je třeba se jimi řídit, aby nedošlo k poškození přístroje.

---

**NEBEZPEČÍ!**



Odkazy, které upozorňují na nebezpečí a donutí operátora, aby se řídil dle procedury nebo specifického způsobu, což nebude-li řádně vykonáno, může ohrozit vážným způsobem jeho zdraví.

---

**POZNÁMKA**

Poznámky, vyjmuté z textu obsahují velmi důležité informace.

---

## Uskladnění

Během dopravy a skladování kontrolérů, se musí dbát na následující podmínky prostředí:

- teplota: od -20 °C do +70 °C
- relativní vlhkost: 0 – 95 % (nekondenzující)

## Příprava k instalaci

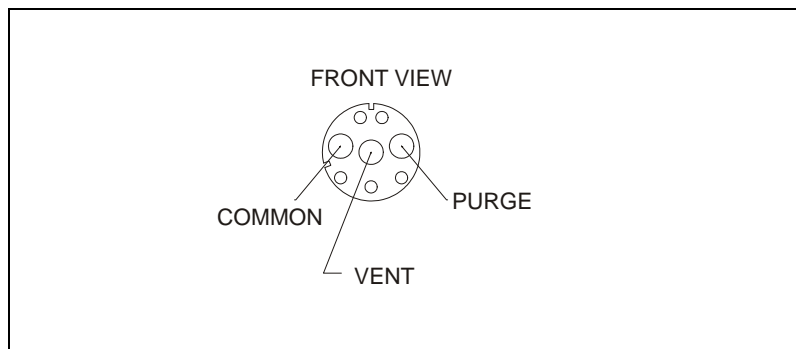
Turbo-V Remote Panel je dodáván ve speciálním ochranném balení; objeví-li se následkem dopravy nějaké poškození, je třeba se okamžitě obrátit na místní prodejní servis. Během rozbalování se obzvláště musí dávat pozor, aby kontrolér nespadol nebo aby neutrpěl nárazy. Balicí materiál se nesmí odkládat v přírodě. Tento materiál se může kompletně recyklovat, neboť se shoduje se směrnicemi EU, CEE 85/399, které se týkají životního prostředí.

## Instalace

Turbo-V Remote Panel musí být napojen na kontrolér vhodným kabelem (viz tabulku součástí, které lze objednat), viz následující obrázek. Vzhledem k tomu, že požadovaný výkon/signály jsou poskytovány uvedeným kabelem, není zapotřebí vnějšího napájení.

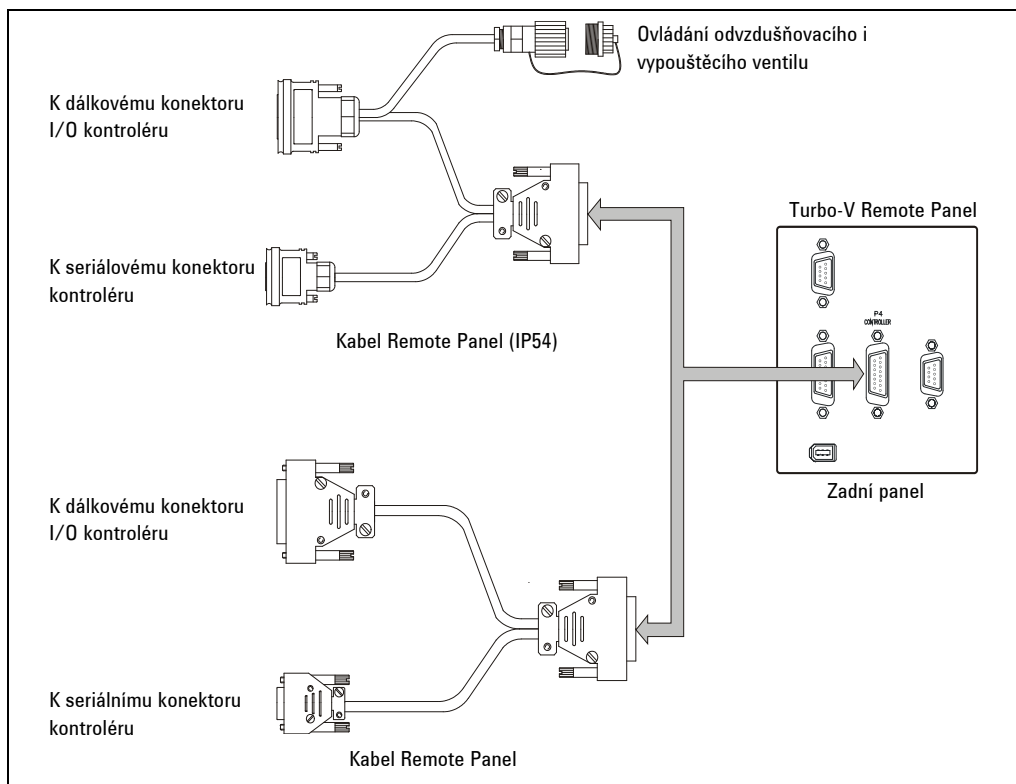
V případě, že operátor si zvolí řízení odvěšovací a vypouštěcí ventilů používajíc vnější ovládání, je třeba použít verzi s kabelem IP54, nařídit kontrolér na "valve setting" "on command" a kabely spojit dle obrázku.

Ventily jsou n. z. ; ventily se otevřou přivedí-li se 24 V do odvěšovacího i vypouštěcího ventilu a společné koncovky.



Postava 1

## 13 Návod k Použití Instalace



**Postava 2** Spojení

### POZNÁMKA

Turbo-V Remote Panel může být nainstalován na stole nebo uvnitř příslušného rack. V každém případě je třeba, aby okolo přístroje vzduch proudil volně. Kontrolér se nesmí instalovat v místech vystavených atmosferickým vlivům (déšť, mráz, sníh), prachu, agresivním plynům, na výbušných místech nebo na místech kde existuje velké riziko požárů.

Během provozu je třeba, aby se respektovaly následující podmínky, týkající se operativního prostředí:

- teplota: od 5 °C do +45 °C;
- relativní vlhkost: 0 – 95 % (nesmí kondenzovat).

Pro další spojení a instalaci opčného příslušenství je třeba se podívat do sekce "Technical Information".

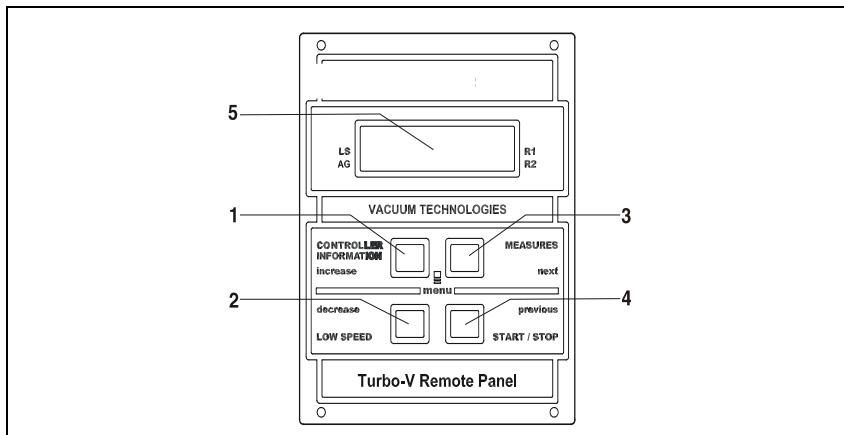
## Použití

V tomto odstavci jsou uvedeny hlavní operativní postupy. Pro další detaily i pro procedury týkající se spojení nebo dalších opčných detailů, je třeba se obracet na odstavec "Use" v příloze "Technical Information".

## Ovládání, indikátory a konektory

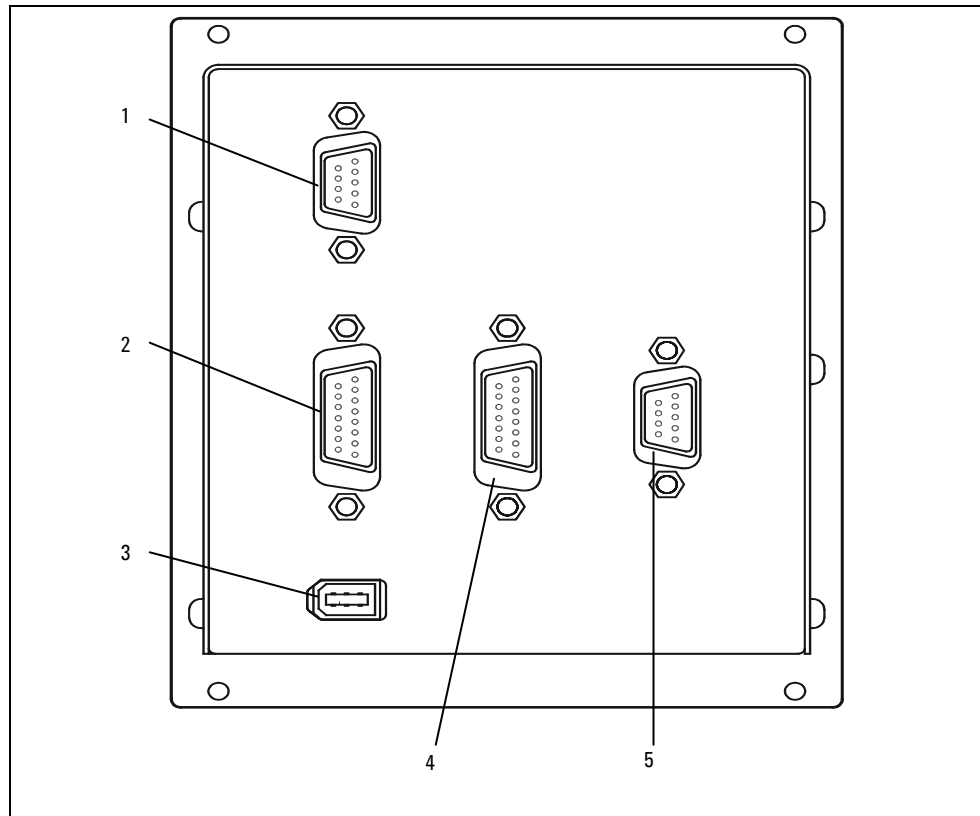
V následujícím je ukázán čelní a zadní panel kontroléru. Pro další detaily je nutno se odkazovat na sekci "Technical Information".

## Popis čelního panelu



**Postava 3** Turbo-V Remote Panel

- 1 Tlačítko na vyvolání parametrů na displeji: cycle number, cycle time a pump life.
- 2 Tlačítko na volbu způsobem LOW SPEED. Je aktivní jen při zvoleném způsobu ovládání z čelního panelu. Stiskne-li se jen jednou, čerpadlo pracuje rychlostí "stand-by". A stiskne-li se ještě jednou, deaktivuje se způsob LOW SPEED.
- 3 Tlačítko na přivolání parametrů na displeji: pump current, pump power, rotational speed. Nezávisle od zvoleného provozního způsobu, je neustále aktivní. Jestliže se tlačítka 3 a 1 stisknou společně alespoň na 2 vteřiny, bude aktivován program, dle kterého se mohou naprogramovat některé operativní parametry.
- 4 Tlačítko ovládání START, STOP/RESET. Je aktivním jen je-li zvolen způsob ovládání na čelním panelu. Stiskne-li se jen jednou, spouštěcí fáze bude aktivována; jestliže se znovu stiskne, čerpadlo se zastaví. Jestliže se čerpadlo automaticky zastaví z důvodu nějaké poruchy, uvedené tlačítko se musí napoprvé stisknout, aby se uskutečnil reset kontroléru a po druhé aby se čerpadlo znovu spustilo.
- 5 Alfanumerický displej na tekuté krystaly: matice bodů, 2 linky x 16 charakterů.



**Postava 4** Zadní panel Turbo-V Remote Panel

- 
- |   |                                                                                                        |
|---|--------------------------------------------------------------------------------------------------------|
| 1 | Konektory vstupních logických signálů (spojovací konektor je dodáván s příslušným uzavíracím můstkem). |
|---|--------------------------------------------------------------------------------------------------------|
- 
- |   |                                                                                                              |
|---|--------------------------------------------------------------------------------------------------------------|
| 2 | Konektor výstupních logických signálů a monitorování frekvence čerpadla a analogický programovatelný výstup. |
|---|--------------------------------------------------------------------------------------------------------------|
- 
- |   |                                                 |
|---|-------------------------------------------------|
| 3 | Konektor výstupního přívodu (24 Vcc na kalibr). |
|---|-------------------------------------------------|
- 
- |   |                                                                                               |
|---|-----------------------------------------------------------------------------------------------|
| 4 | Konektor vstupního přívodu z kontroléru čerpadla a sériové komunikace s kontrolérem čerpadla. |
|---|-----------------------------------------------------------------------------------------------|
- 
- |   |                                        |
|---|----------------------------------------|
| 5 | Sériový konektor na dálkovou kontrolu. |
|---|----------------------------------------|
-

## **Zapalování Kontroléru**

Kontrolér se zapne vložením kabele do zásuvky na síti.

## **Spouštění čerpadla**

Ke spuštění čerpadla, je třeba oprávnit vstup interlocku a stisknout tlačítko START na čelním panelu.

## **Zastavení čerpadla**

Čerpadlo se zastaví stisknutím tlačítka STOP na čelním panelu.

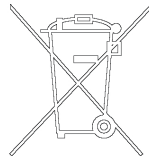
## Údržba

Kontrolér série Turbo-V 301 Rack nevyžadují žádnou údržbu. Jakýkoliv zásah musí být vykonáván jen oprávněným personálem. V případě poruchy se může využít opravářského servisu Agilent nebo "Agilent advanced exchange service", kde poškozený kontrolér bude vyměněn za nový. Bude-li zapotřebí kontrolér likvidovat, musí se postupovat dle platných specifických státních norem

## Likvidace

**Význam loga "WEEE" nacházejícího se na štítku.** Níže uvedený symbol odpovídá směrnici CE pojmenovaným "WEEE". Tento symbol (**platný jen pro státy Evropské Unie**) určuje, že výrobek, který je takto označen NESMÍ být likvidován společně s ostatními domácími nebo průmyslovými odpady, ale je nutno ho předat do patřičných provozních sběrů, kde musí být tříděn a likvidován odděleně od městského odpadu.

Uživateli, který míní výrobek likvidovat doporučujeme, aby se nakontaktoval přímo s výrobcem nebo prodejcem, který se po patřičném prověření termínů a smluvních podmínek postará o kompletní likvidaci uvedeného výrobku.



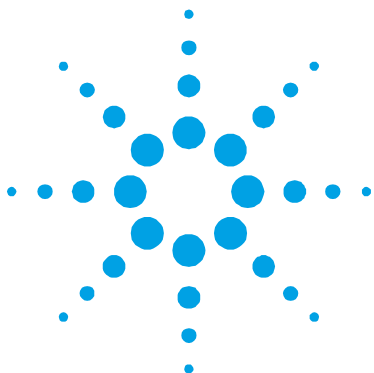
## Chybné zprávy

V některých případných poruchách na autodiagnostických obvodech, kontrolér předává několik chybných odkazů, které jsou vyjmenovány v tabulce, která se nachází na následující stránce.

ZPRÁVA	POPIS	KOREKCE
COMMUNICATION ERROR	Remote Panel není schopen zahájit komunikaci s kontrolérem.	Prověřit sériové spojení mezi Remote Panelem a kontrolérem.
COMMUNICATION BROKEN	Remote Panel není schopen pokračovat v komunikaci s kontrolérem.	Prověřit sériové spojení mezi Remote Panelem a kontrolérem.
XXXXXXX MODEL:	Kontrolér bez podpory. Remote Panel funguje jen s následujícími čerpadly: TV551 s navigator controller, TV701 s navigator controller, TV1001 s navigator controller, TV2k-G a TV3k-T.	Spojit Remote Panel a kontrolér s podporou.
CHECK CONNECTION TO PUMP	Chybné fungování spojení mezi čerpadlem a kontrolérem (P31). Nebo Čerpadlo má nižší teplotu než je 0 °C.	Prověřit zda kabelové spojení mezi pumpou a kontrolérem je správně upevněné na obou dvou koncích a že nedošlo k přerušení. Stisknout tlačítko START dvakrát pro opětné spuštění čerpadla.
WAITING INTERLOCK	Aktivní signál interlock se nachází na konektoru P1, neboť došlo k přerušení zkratu mezi pin 3 a pin 8 del konektoru J1, nebo kvůli otevření signálu vnějšího interlock .	Znovu obnovit spojení nakrátko mezi pin 3 a pin 8 konektoru J1, nebo uzavřít signál vnějšího interlock.
PUMP OVERTEMP.	Teplota horního ložiska nebo čerpadla převýšila 60 °C.	Počkat až se teplota opět vrátí pod mezní hodnotu. Stisknout dvakrát tlačítko START pro opětné spuštění čerpadla.
CONTROLLER OVERTEMP.	Teplota kontroléru převýšila 70 °C. Nebo Teplota radiátoru kontroléru je vyšší než 60 °C.	Vyčkat až se teplota opět vrátí pod mezní hodnotu. Stisknout dvakrát tlačítko START pro opětné spuštění čerpadla.
TOO HIGH LOAD	Absorbovaný proud čerpadla je vyšší	Prověřit, zda rotor čerpadla se

ZPRÁVA	POPIS	KOREKCE
	než ten naprogramovaný.	může volně otáčet. Stisknout dvakrát tlačítko START pro opětné spuštění čerpadla.
SHORT CIRCUIT	Během normálního provozu (po spouštěcí fázi) výstupní spojení je ve zkratu.	Prověřit spojení mezi čerpadlem a kontrolérem. Stisknout dvakrát tlačítko START pro opětné spuštění čerpadla.
SYSTEM OVERRIDE	Pumpa byla zastavena nouzovým signálem prostřednictvím dálkového kontaktu.	Odpojit napájecí kabel kontroléru a napravit příčinu poplachu. Znovu připojit napájecí kabel a stisknout dvakrát tlačítko START pro opětné spuštění čerpadla.
OVERVOLTAGE	Vznikla porucha v napájecí sekci kontroléru, nebo kontrolér dostal falešný signál.	Stisknout dvakrát tlačítko START pro opětné spuštění čerpadla. Po opětné zprávě ohledně údržby je nutno se obrátit na Agilent .
POWER FAIL	Porucha v napájecí sekci čerpadla, kontroléru.	Ohledně údržby je třeba se obrátit na Agilent.

## **13** **Návod k Použití** Chybné zprávy



## 14 Návod na Obsluhu

Všeobecné informácie	196
Uskladňovanie	198
Príprava pre inštaláciu	198
Inštalácia	199
Použitie	201
Povely, ukazovatele a prípojky	201
Popis čelného panelu	202
Zapalování Kontroléru	204
Naštartovanie čerpadla	204
Zastavenie čerpadla	204
Údržba	205
Likvidácia	205
Oznamy väd	206

Preklad originálneho návodu



## Všeobecné informácie

Tento prístroj slúži k profesionálnemu použitiu. Užívateľ si pred použitím tohto prístroja musí pozorne prečítať tento návod k použitiu a jeho ďalšie doplnky, dodané podnikom Agilent. Podnik Agilent nepreberá na seba žiadnu zodpovednosť za úplné, alebo čiastočné nedodržanie návodu, za nesprávne používanie prístroja nekompetentnými osobami, za nepovolené zásahy a za používanie, ktoré nie je v zhode s príslušnými národnými normami.

Turbo-V Remote Panel je kontrolovaný mikro-procesorom a čelným panelom a umožní užívateľovi riadiť čerpadlo typu "controller on-board":

TV551 s Navigator controller

TV701 s Navigator controller

TV1001 s Navigator controller

TV2k-G s controller on-board

TV3k-T s controller on-board

TV3k-G s controller on-board

TV70 s Navigator controller

TV81 s Navigator controller

TV301 s Navigator controller

TV303 s Navigator controller

Ďalej sú popísané všetky potrebné informácie, ktorých úlohou je zabezpečiť bezpečnosť pracovníka počas používania prístroja. Podrobné informácie sú uvedené v dodatku "Technical Information".

**Tento návod používa nasledujúce signály:**

**POZOR!**

Signály “pozor” sú pred tými postupmi, ktoré, v prípade ich nedodržania, môžu poškodiť prístroj.

---

**NEBEZPEČÍ!**



Signál nebezpečie upozorňuje pracovníka na určitý pracovný postup, ktorý, keď nie je prevedený správne, môže zapríčiniť vážne osobné poranenia.

---

**POZNÁMKA**

Poznámky zahrňujú dôležité informácie, vytiahnuté z textu.

---

## Uskladňovanie

Počas prepravy a uskladňovania “controllers” musia byť dodržané tieto podmienky prostredia:

- teplota: od -20 °C do +70 °C
- relatívna vlhkosť: 0 – 95 % (nekondenzujúca)

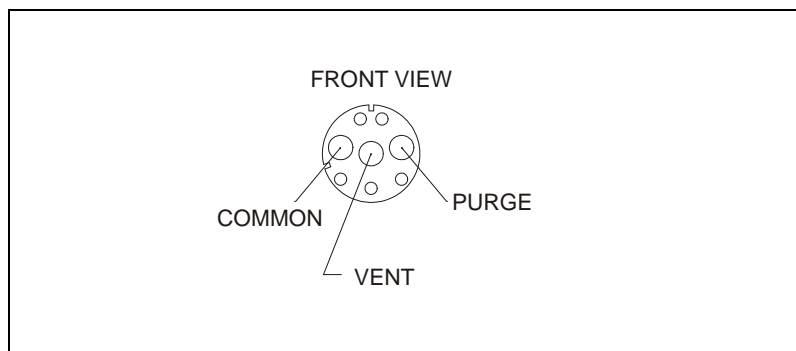
## Príprava pre inštaláciu

Turbo-V Remote Panel je dodaný v špeciálnom ochrannom obale; keby ste objavili škodu, spôsobenú prepravou, skontaktujte sa s kanceláriou miestneho predaja. Pri rozbaľovaní dávajte pozor, aby Vám “controller” nespadol, alebo niekde nenarazil. Obal nezahadzujte. Je vyrobený z recyklovateľného materiálu a zodpovedá smernici CEE 85/399 pre ochranu životného prostredia.

## Inštalácia

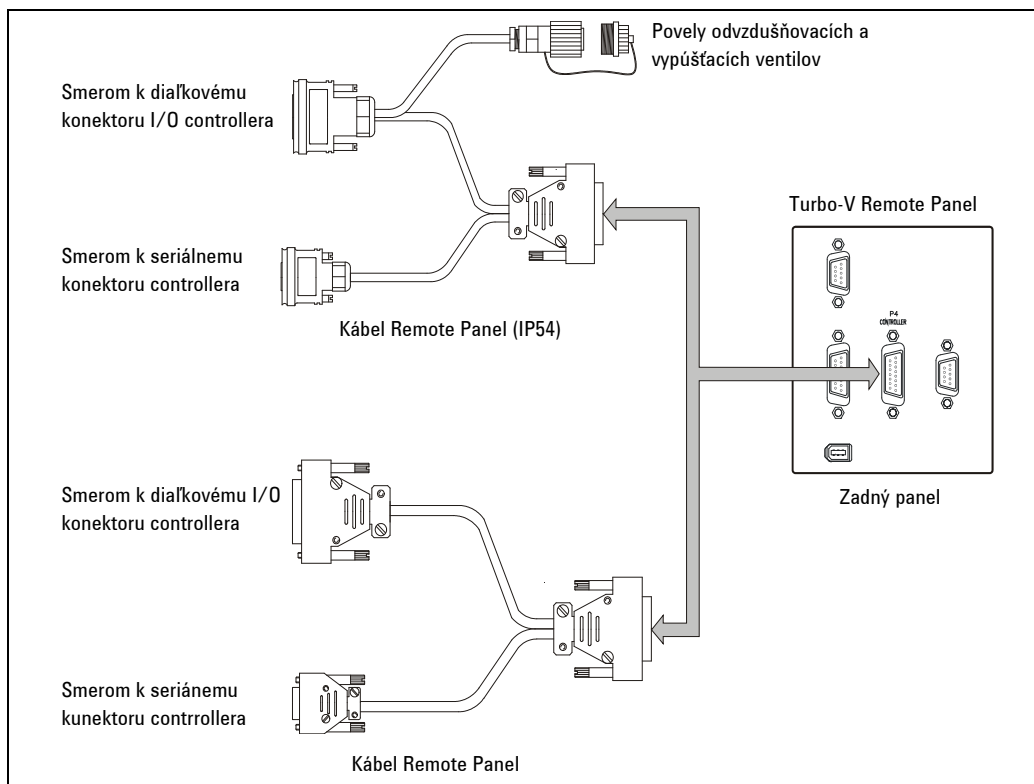
Turbo-V Remote Panel musí byť napojený na controller príslušným káblom (viď tabuľku objednávajúcich súčiastok), tak, ako je to znázornené na nasledujúcom obrázku. Pretože výkon a požadované signály dodáva tento kábel, nie je potrebné ďalšie vonkajšie napojenie.

Keď chce pracovník riadiť odvzdušňovacie a vypúšťacie ventily pomocou vonkajších povelov, použije verziu s káblom IP54, controller nastaví na "valve setting" "on command" a káble spojí tak, ako je to znázornené na obrázku. Ventily sú n.c.; dodaním 24 V medzi odvzdušňovací a vypúšťací ventil a spoločnú koncovku, sa ventily otvoria.



Postava 1

## 14 Návod na Obsluhu Inštalácia



**Postava 2** Napojenia

### POZNÁMKA

Turbo-V Remote Panel môže byť inštalovaný na pracovnom stole alebo vo vnútri rack. V každom prípade je potrebné, aby chladiaci vzduch mohol voľne krúžiť okolo prístroja. Controller nepoužívajte a neinštalujte v prostredí, vystavenom atmosferickým vplyvom (dážď, mráz, sneh), prachu, agresívnym plynom a v prostredí s vysokým rizikom požiaru.

Počas chodu je dôležité, aby boli rešpektované tieto podmienky prostredia:

- teplota: od 5 °C do +45 °C
- relatívna vlhkosť: 0 – 95 % (nekondenzujúca).

Pre ostatné napojenia a inštaláciu zvoleného príslušenstva vid' časť "Technical Information".

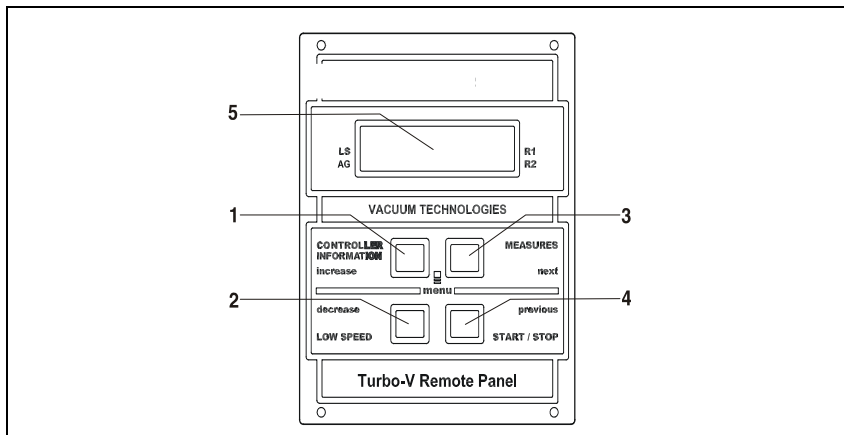
## Použitie

V tomto odstavci sú uvedené hlavné operatívne postupy. Pre ostatné detaily a pre postupy, ktoré sa vzťahujú na napojenia alebo na zvolené príslušenstvo, postupujte podľa časti "Use" v dodatku "Technical Information".

## Povely, ukazovatele a prípojky

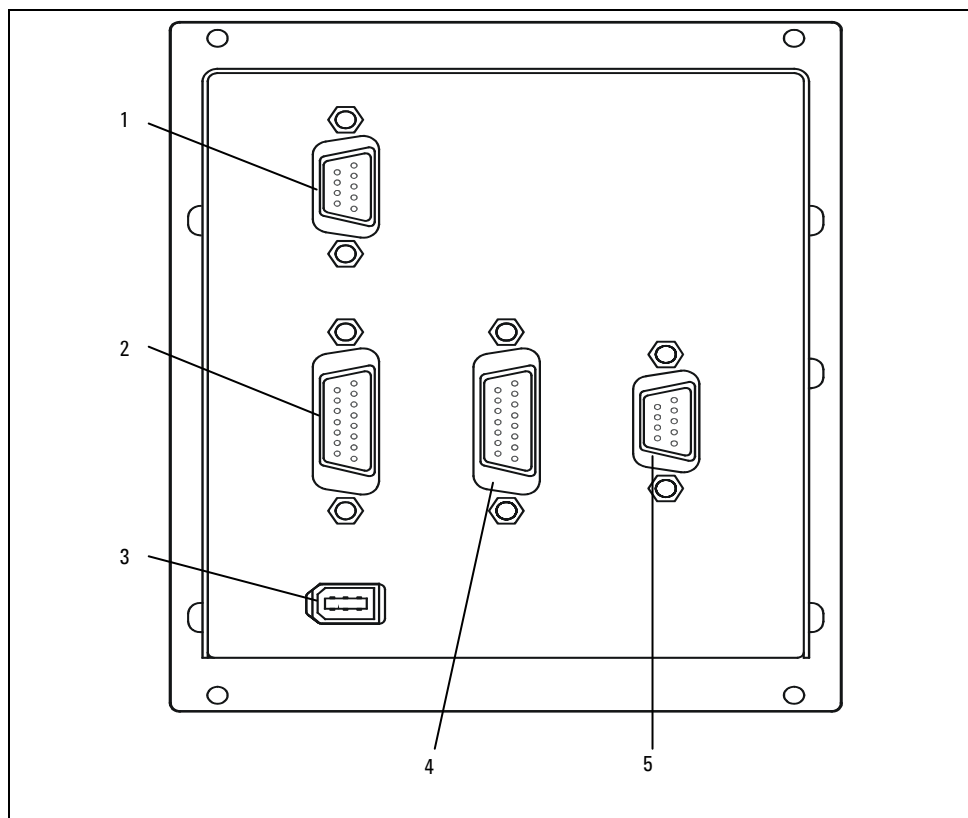
Ďalej je zobrazený čelný a zadný panel controllera. Pre ďalšie detaily konzultujte časť "Technical Information".

## Popis čelného panelu



**Postava 3** Turbo-V Remote Panel

- 1 Tlačidlo pre vyvolanie parametrov cycle number, cycle time a pump life.
- 2 Tlačidlo pre voľbu LOW SPEED. Je aktívne iba vtedy, keď povel vychádza z čelného panelu. Keď sa stlačí raz, čerpadlo sa točí rýchlosťou "stand-by". Stlačiac ho druhý raz, sa vypne LOW SPEED.
- 3 Tlačidlo pre vyvolanie parametrov pump current, pump temperature, pump power a rotational speed. Je stále aktívne, nezávisle od voľby funkcie. Stlačiac spolu tlačidlá 3 a 4 na aspoň 2 sekundy sa spustí program, s ktorým je možné zvoliť niektoré operatívne parametre.
- 4 Tlačidlo povelov START, STOP/RESET. Je aktívne iba vtedy, keď povel vychádza z čelného panelu. Keď sa stlačí raz, spustí sa fáza naštartovania; stlačiac ho znovu sa zastaví čerpadlo. Keď sa čerpadlo v prípade poruchy zastaví automaticky, treba stlačiť tlačidlo prvý krát pre resetovanie controllera a druhý krát pre znovunaštartovanie čerpadla.
- 5 Alfanumerický displej z tekutých kryštáľkov: bodová matica, 2 línie x 16 písmem. i.



**Postava 4** Zadný panel Turbo-V Remote Panel

- 1 Konektory výstupných logických signálov (spojovací konektor je dodaný s príslušným uzáverovým mostikom).
- 2 Konektor logických výstupových signálov a monitorovanie kmitočtu čerpadla a programovateľný analogový výstup.
- 3 Výstupný napájací konektor (24 Vcc na kaliber).
- 4 Vstupný napájací konektor čerpadlového controllera a seriálna komunikácia s čerpadlovým controllerom.
- 5 Seriálny konektor pre diaľkovú kontrolu.

## Zapnutie “controller”

Pre zapnutie “controller” stačí zapojiť kábel na elektrickú sieť a vypínač siete dať do polohy I.

## Naštartovanie čerpadla

Pre naštartovanie čerpadla je potrebné sprístupniť vstup Interlock a stlačiť tlačidlo START na čelnom paneli.

## Zastavenie čerpadla

Pre zastavenie čerpadla treba stlačiť tlačidlo STOP na čelnom paneli.

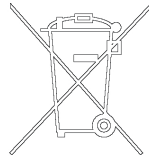
## Údržba

“Controllers” série Turbo-V 301 Rack nevyžadujú žiadnu údržbu. Akýkoľvek zásah musí previesť poverený personál. V prípade vady je možné využiť opravnú službu Agilent alebo "Agilent advanced exchange service", ktorá vymení vadný “controller”. V prípade, že “controller” musí ísť do zberu, riadte sa platnými príslušnými národnými normami.

## Likvidácia

**Význam loga "WEEE" nachádzajúceho sa na štítkoch.** Aplikovanie doluo značeného symbolu dodržiava smernicu EÚ s názvom "WEEE".

Tento symbol (**platný iba pre štáty Európskej Únie**) znamená, že výrobok s týmto štítkom NEMIE byť odstránený spolu s bežným domácim alebo priemyselným odpadom, ale sa musí odstrániť ako delený odpad. Vyzývame preto konečného užívateľa, aby sa skontaktoval s dodávateľom prístroja, či už je to výrobca alebo predajca za účelom jeho likvidácie podľa zmluvných podmienok predaja.



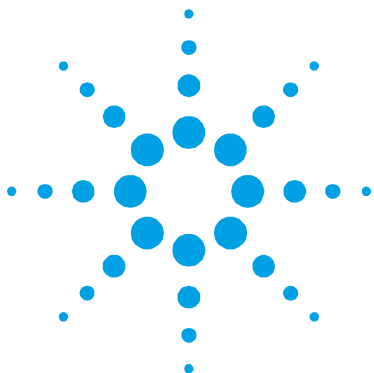
## Oznamy vąd

V niektorých prípadoch, týkajúcich sa vąd, obvod samodiagnostiky “controller” signalizuje niekoľko oznamov vąd, uvedených na ďalšej strane.

OZNAM	POPIS	ZÁSAH OPRAVY
COMMUNICATION ERROR	Remote Panel nie je schopný naštartovať komunikáciu s controller.	Skontrolujte seriálne spojenie medzi Remote Panel a controller.
COMMUNICATION BROKEN	Remote Panel nie je schopný pokračovať v komunikácii s controller.	Skontrolujte seriálne spojenie medzi Remote Panel a controller.
XXXXXXX MODEL	Controller nie je podporený. Remote Panel funguje iba s týmito čerpadlami: TV551 s navigator controller, TV701 s navigator controller, TV1001 s Navigator Controller, TV2k-G a TV3k-T.	Napojte Remote Panel na podporený controller.
CHECK CONNECTION TO PUMP	Nefunguje napojenie medzi čerpadlom a “controller”. Alebo Čerpadlo má nižšiu teplotu ako 0 °C.	Skontrolujte, či je spojovací kábel medzi čerpadlom a “controller” správne zasunutý na oboch koncoch a či nie je prerušený. Stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla.
WAITING INTERLOCK	je aktívny signál interlock na spojke P1 z dôvodu prerušenia skratu medzi pin 3 a pin 8 spojky J1, alebo z dôvodu otvorenia signálu vonkajšieho interlock.	Obnoviť skrat medzi pin 3 a pin 8 spojky J1, alebo zavrieť signál vonkajšieho interlock.
PUMP OVERTEMP.	Teplota horného ložiska alebo čerpadla prekročila 60 °C.	Vyčkať, až sa teplota opäť vráti pod hranicu. Stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla.
CONTROLLER OVERTEMP.	Teplota “controller” prekročila 70 °C. Alebo Teplota radiátora controllera presahuje 60 °C.	Počkať, kým teplota neklesne pod hranicu. Stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla
TOO HIGH LOAD	Počas bežného chodu (po naštartovaní) čerpadlom absorbovaný prúd prevyšuje ten naprogramovaný.	Skontrolujte, či má rotor čerpadla možnosť točiť sa voľne. Stlačte dvakrát tlačidlo START pre

OZNAM	POPIS	ZÁSAH OPRAVY
SHORT CIRCUIT	Počas bežného chodu (po naštartovaní) východzie napojenie je v skrate.	znovunaštartovanie čerpadla Skontrolujte napojenia medzi čerpadlom a "controller". Stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla
SYSTEM OVERRIDE	Núdzový signál, vychádzajúci z diaľkového kontaktu zastavil čerpadlo.	Vytiahnite šnúru z elektrickej siete "controller" a ostráňte dôvod alarmu. Zasuňte napájaciu šnúru a stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla.
OVERVOLTAGE	Objavila sa vada v sekcii napájania "controller", alebo "controller" dostal falošný signál.	Stlačte dvakrát tlačidlo START pre znovunaštartovanie čerpadla. Keď sa tento oznam objaví znovu, obráťte sa na firmu Agilent pre údržbu.
POWER FAIL	Vada v napájacom systéme čerpadla, controllera.	Pre údržbu sa obráťte na Agilent.

## **14** **Návod na Obsluhu** Oznamy vád



## 15 Priločnik za Navodila

Splošne informacije	210
Skladiščenje	212
Priprava za montažo	212
Montaža	213
Uporaba	215
Komande, indikatorji in konektorji	215
Opis prednjega odbora	216
Vklonitev Controllera	218
Zagon črpalke	218
Zaustavitev črpalke	218
Vzdrževanje	219
Odlaganje odpadkov	219
Obvestilo o napaki	220

Prevod navodil v izvorniku



## Splošne informacije

Naprava je namenjena samo za strokovno rabo. Uporabnik mora temeljito prebrati priročnik in pri tem upoštevati vsako dodatno informacijo, ki jo nudi Agilent. Agilent ne odgovarja v slučaju celotnega ali delnega nespoštovanja navodil, pri nepravilni uporabi strokovnega osebja, v slučaju nedovoljenih posegov ali pri zanemarjenju specifičnih nacionalnih meril.

Turbo-V Remote Panel upravlja mikrokrmilnik in dovoli uporabniku, da upravlja črpalko kot je "controller on board" s pomočjo prednjega odbora:

- TV551 z Navigator controller
- TV701 z Navigator controller
- TV1001 z Navigator controller
- TV2k-G s controller on board
- TV3k-T s controller on board
- TV3k-G s controller on board
- TV70 z Navigator controller
- TV81 z Navigator controller
- TV301 z Navigator controller
- TV303 z Navigator controller

V naslednjih odstavkih so navedene vse informacije potrebne pri zagotavljanju sigurnosti uporabnika med rokovanjem s to napravo. Podrobnejše informacije dobite v priponki "Technical Information".

V tem priročniku so varnostne informacije razvrščene v dva razredai:

**POZOR!**

Znak 'pozor' se pojavi pred postopke, ki če jih uporabnik zanemari, lahko napravo poškoduje.

---

**SVARILO!**



Znak svarila opozarja upravnika, da mora pri določenem postopku ali pri posebnem delovanju paziti. Kršitev svarila lahko privede do lažjih ali hudih telesnih poškodb.

---

**OPOMBA**

Opombe vsebujejo vse najbolj pomembne informacije iz besedila.

---

## Skladiščenje

Med transportom in skladiščenjem controllerjev morate zagotoviti naslednje pogoje okolice:

- temperatura: od -20 °C do +70 °C
- Relativna vlaga: 0 – 95 % (ne-kondenzna)

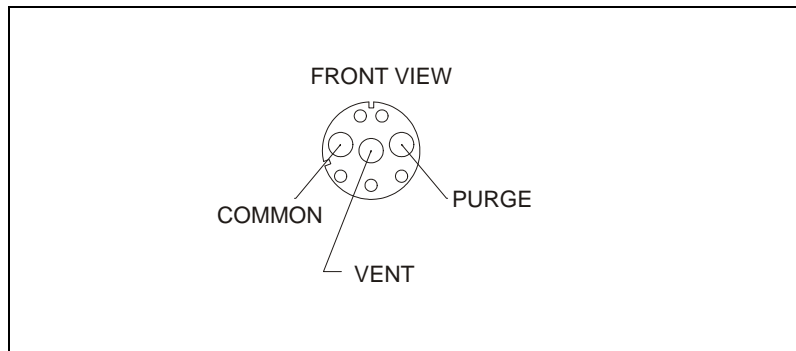
## Priprava za montažo

Turbo-V Remote Panel dobite v posebni embalaži; če so na embalaži znaki poškodbe, ki naj bi nastale med transportom obrnite se krajevnemu prodajalnemu biroju. Pri odstranitvi embalaže morate še posebno paziti, da controller ne pade na tla ali da ne podleigne udarcem. Ne zavržite embalažo v okolico. Material lahko v celoti recikirate in ustreza direktivi CEE 85/399 za zaščito okolja.

## Montaža

Turbo-V Remote Panel morate priključiti h controllerju s pomočjo ustrezne žice (oglejte si tudi preglednico naročljivih delov), kot prikazano na spodnji sliki. Potrebno moč/potrebne znake dobite dobavljene s to žico, zaradi tega ni potrebno nobeno drugo zunanje napajanje.

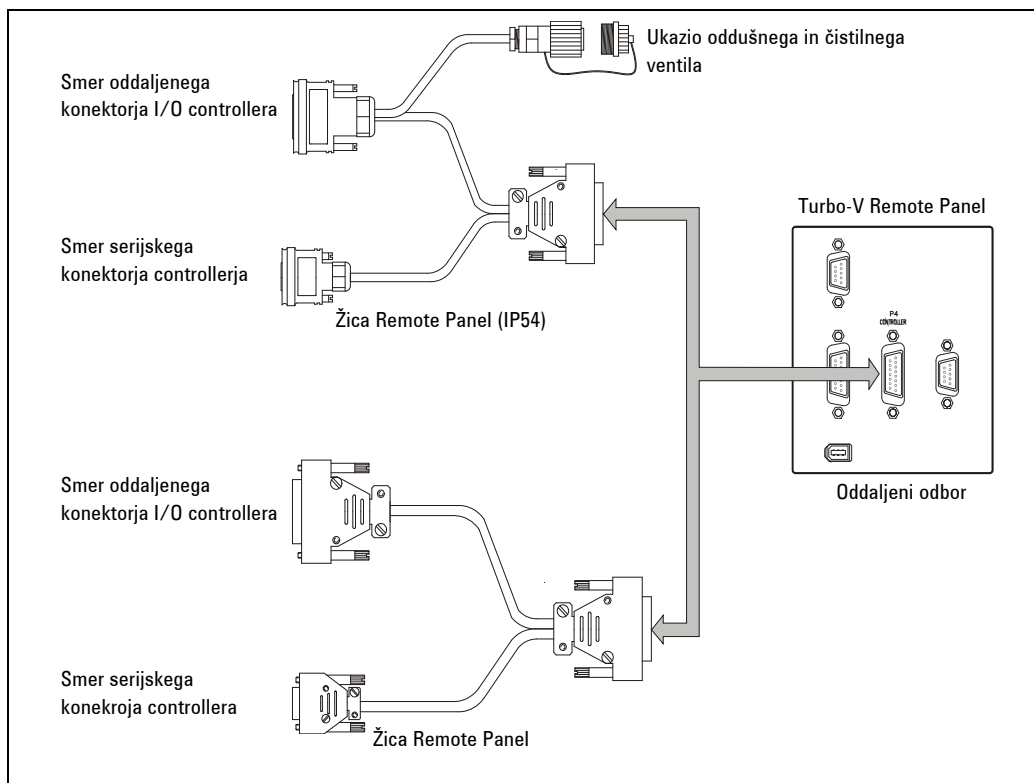
Če želi uporabnik upravljati oddušne in čistilne ventile s pomočjo zunanjih ukazov, mora koristiti verzijo z žico IP54, controller namestiti na "valve setting" "on command" in priključiti žice kot je prikazano na sliki. Ventili so navadno zaprti; če priključimo na 24 V oddušni in čistilni ventil ter skupni terminal, se bodo ventili odprli.



Slika 1

## 15 Priročnik za Navodila

### Montaža



**Slika 2** Priključki

#### OPOMBA

Turbo-V Remote Panel lahko vgradite na mizo ali znotraj ustreznega rack-a. V vsakem slučaju morate poskrbeti, da bo zrak za hlajenje prosto krožil skozi napravo. Controllera ne smete montirati ali uporabiti v okolici, ki je izpostavljena klimatskim dejavnikom (dežju, ledu, snegu), prahu, agresivnim plinom, v okolici, kjer obstaja nevarnost vnetja ali požara.

Med delovanjem morajo biti zagotovljeni naslednji pogoji okolice:

- temperatura: od 5 °C do +45 °C
- relativna vlaga: 0 – 95 % (ne-kondenzna).

Kar se tiče ostalih priključkov in montaže opcionalnih delov, si lahko ogledate priponko "Technical Information".

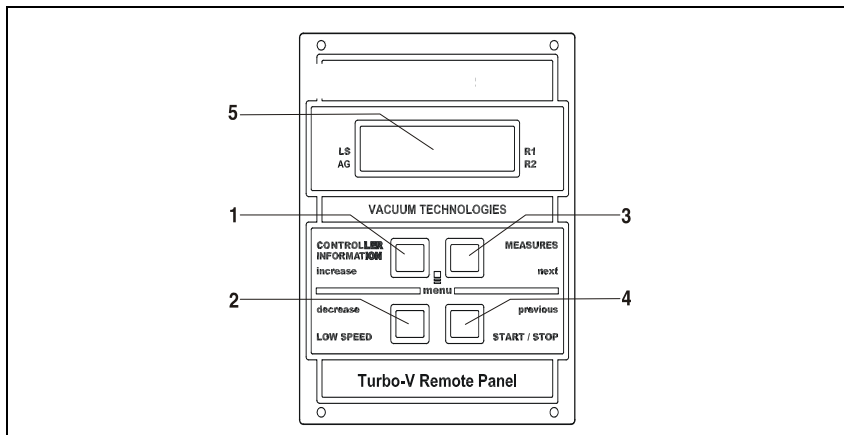
## Uporaba

V tem odstavku so navedeni vsi glavni operativni postopki. Za dodatne detalje in postopke, ki zadevajo priključitve ali posamezne opcije, si lahko ogledate odstavek "Use" v priponki "Technical Information".

## Komande, indikatorji in konektorji

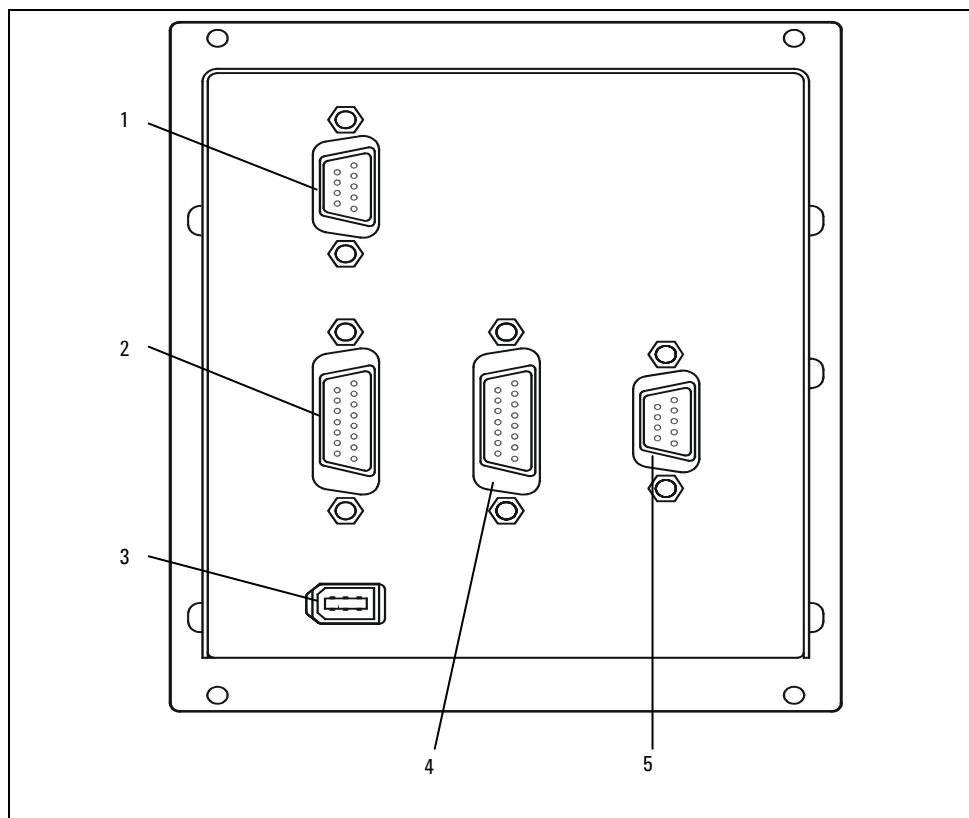
Sledi prikaz prednje in ozadnjega odbora controllera. Za podrobnejše informacije si ogledajte paragraf "Technical Information".

## Opis prednjega odbora



**Slika 3** Turbo-V Remote Panel

- 1 Tipka za preklic na zaslonu parametrov cycle number, cycle time in pump life.
- 2 Tipka za LOW SPEED modus, je aktivna samo nato, ko ste nastavili ta ukaz na sprednjem odboru. Če gumb pritisnete enkrat, se črpalka vrti s hitrostjo "stand-by". Če še enkrat pritisnete gumb, izključite LOW SPEED modus.
- 3 Tipka za preklic na zaslonu parametrov pump current, pump temperature, pump power in rotational speed. Ta tipka je vedno aktivna neodvisno od nastavljenega modusa delovanja. Če skupaj pritisnete gumb 3 in 4 za vsaj 2 sekundi boste aktivirali program, s katerim je mogoče nastaviti nekatere operativne parametre.
- 4 Tipka za zagon komand START, STOP/RESET. Aktivna je samo nato, ko ste to komando nastavili na prednjem odboru. Če gumb enkrat pritisnete boste pognali črpalko; če ga še enkrat pritisnete, boste črpalko zaustavili. Če se je črpalka zaradi kvara sama zaustavila, morate to tipko še enkrat pritisniti, da bi lahko controller resetirali in še drugič za ponoven zagon črpalke.
- 5 Alfaštevilčni zaslon iz tekočih kristalov: prebijalo bodov, 2 linije x 16 znakov.



**Slika 4** Zadnji pult Turbo-V Remote Panel

- 1 Vhodni konektor za logične signale (konektor za zduževanje je opremljen z ustreznim zapiralnim mostičkom).
- 2 Konektor logičnih signalov na izhodu in monitoriranje frekvence črpalke, in nastavljivi analogični izhod.
- 3 Izhodni konektor za napajanje (24 Vcc za kaliber).
- 4 Vhodni konektor za napajanje iz controller-a črpalke in serijska komunikacija s controller-jem črpalke.
- 5 Serijski konektor za daljinski nadzor.

## **Vklopitev Controllera**

Za vklopitev controllera, priključite napajalni kabel v omrežno vtičnico.

## **Zagon črpalke**

Da bi zagnali črpalko morate usposobiti vhod interlock-a in pitisniti tipko START na prednji plošči.

## **Zaustavitev črpalke**

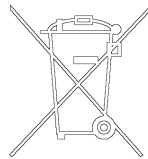
Zaustavite črpalko s tipko STOP na prednji plošči.

## Vzdrževanje

Controllerji serije Turbo-V 301 Rack ne potrebujejo vzdrževanja. Samo pooblaščen osebje lahko rokuje s to napravo. V slučaju okvare se lahko obrnete servisu Agilent ali "Agilent advanced exchange service", ki vam stavlja na razpolago obnavljeni controller-a Controller zavržite kot industrijski odpadek v skladu z državnimi predpisi.

## Odlaganje odpadkov

**Pomen znamke "WEEE" na etiketah.** Spodaj navedeni simbol je v skladu z direktivo ES znano pod imenom "WEEE". Ta simbol (**ki velja samo v državah Evropske Skupnosti**) pomeni, da izdelek NE SMETE ODSTRANITI skupaj z ostalimi komunalnimi ali pa industrijskimi odpadki, temveč morate poskrbeti za njihovo primerno ločevanje. Zato pozivamo uporabnike, da se ali pri prodajalnem centru ali pa pri prodajalcu seznanijo o postopku ločevanja in odstranitve odpadkov, šele nato, ko se je seznanil s pogoji in z merili kupopordajne pogodbe.



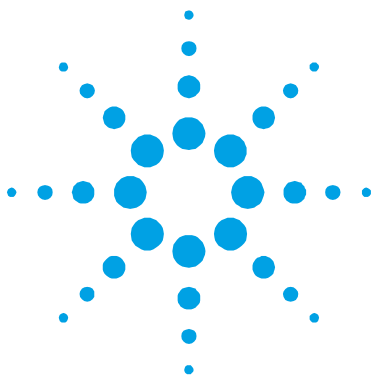
## Obvestilo o napaki

V nekaterih slučajih kvara pokaže električno vezje avtodiagnoze controller-a nekaj obvestil o napakah kot so opisane v spodnji tabeli.

OBVESTILO	OPIS	POPRAVA
COMMUNICATION ERROR	Remote Panel ne zaganja komunikacijo s controller-jem.	Preverite serijski priključek med Remote Panel in controller-jem.
COMMUNICATION BROKEN	Remote Panel ne nadaljuje komunikacijo s controller-jem.	Preverite serijski priključek med Remote Panel in controller-jem.
XXXXXXX MODEL	Controller ni med opcijami. Remote Panel deluje le z naslednjimi črpalkami: TV551 z navigator controller, TV701 z navigator controller, TV1001 z navigator controller, TV2k-G in TV3k-T	Priključite Remote Panel k vgrajenemu controllerju.
CHECK CONNECTION TO PUMP	Napaka v priključitvi črpalke in controller-a. Ali Črpalka ima nižjo temperaturo od 0 °C.	Preverite, da je priključni kabel med črpalko in controller-em na obeh koncih trdno pritrjen in da ni prekinjen. Dvakrat pritisnite tipko START za ponoven zagon črpalke.
WAITING INTERLOCK	Signal interlocka prisoten na konektorju P1 zaradi prekinitve kratkega stika med pinom 3 in pinom 8 konektorja J1, ali ker se je odprl signal zunanjega interlocka.	Ponovno nastavite kratki stik med pinom 3 in pinom 8 konektorja J1, ali zaprite signal zunanjega interlocka.
PUMP OVERTEMP.	Temperatura gornjega ležišča ali črpalke je preseгла i 60 °C.	Počakajte, da se temperatura vrne pod normalnim vrednostnim pragom. Pritisnite dvakrat tipko START za ponoven zagon črpalke.
CONTROLLER OVERTEMP.	Temperatura controllera je preseгла 70 °C. Ali Temperatura controllerskega radiatorja je višja od 60 °C.	Počakajte, da se temperatura vrne pod normalnim vrednostnim pragom. Dvakrat pritisnite tipko START za ponoven zagon črpalke.
TOO HIGH LOAD	Tok, ki ga črpalka prejme je večji od programiranega.	Preverite, da lahko rotor črpalke prosto vrti. Dvakrat pritisnite tipko START za ponoven zagon črpalke.

<b>OBVESTILO</b>	<b>OPIS</b>	<b>POPRAVA</b>
SHORT CIRCUIT	Med navadnim delovanjem (po zagonu) je priklopitev na izhodu v kratkem stiku.	Preverite priključke črpalke in controllera. Dvakrat pritisnite tipko START za ponoven zagon črpalke.
SYSTEM OVERRIDE	Delovanje črpalke je zaustavil signal zasilnosti iz oddaljenega kontakta.	Izklopite napajalni kabel controllera in popravite napako, ki se je pojavila s signalom zasilnosti. Ponovno priključite napajalni kabel in pritisnite dvakrat tipko START za ponoven zagon črpalke.
OVERVOLTAGE	Kvar v oddelku za napajanje controllera, ali je controller dobil nepristen znak.	Dvakrat pritisnite tipko START za ponoven zagon črpalke. Če se to obvestilo spet pojavi, obrnite se Agilent za pravilno vzdrževanje.
POWER FAIL	Kvar v oddelku napajanja črpalke, controllera.	Obrnite se Agilent za pravilno vzdrževanje.

**15** Priročnik za Navodila  
Obvestilo o napaki



## 16 Instructions for Use

General Information	224
Storage	226
Preparation for Installation	226
Installation	227
Use	229
Controls, Indicators and Connectors	229
Front Panel Description	230
Switch-on the Remote Panel	232
Starting the Pump	232
Pump Shutdown	232
Maintenance	233
Disposal	233
Error Messages	234

Original Instructions



## General Information

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

The Turbo-V Remote Panel is micro-processor-controlled, and allows the user to operate one of the following pumps using a front panel:

TV551 with Navigator controller

TV701 with Navigator controller

TV1001 with Navigator controller

TV2k-G with on board controller

TV3k-T with on board controller

TV3k-G with on board controller

TV70 with Navigator controller

TV81 with Navigator controller

TV301 with Navigator controller

TV303 with Navigator controller

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

**This manual uses the following standard protocol:**

**CAUTION!**

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

---

**WARNING!**



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

---

**NOTE**

The notes contain important information taken from the text.

---

## Storage

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

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

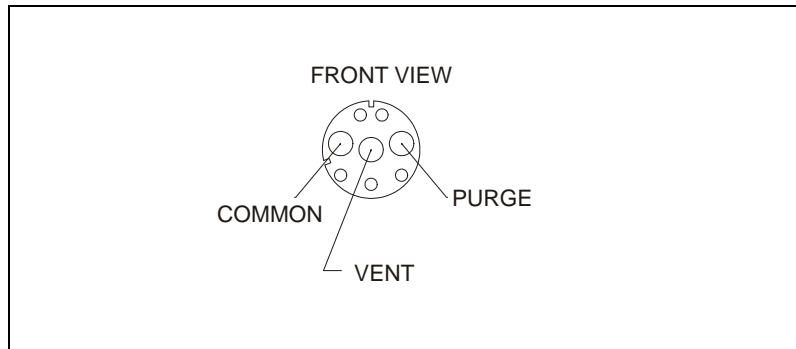
## Preparation for Installation

The Turbo-V Remote Panel 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, ensure that it is not dropped or subjected to any form of impact. Do not dispose of the packing materials in an unauthorized manner. The material is 100 % recyclable and complies with EEC Directive 85/399.

## Installation

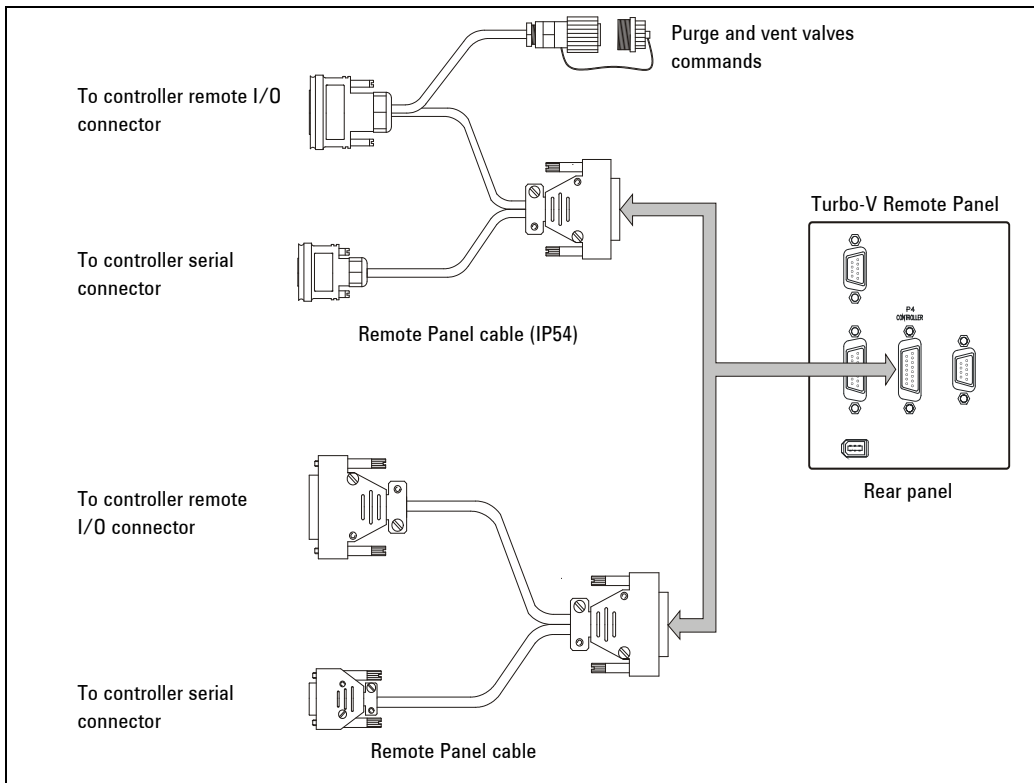
The Turbo-V Remote Panel must be connected to the controller by connection cable (see orderable parts table), as shown in the picture below. All the necessary power/signals are present on this cable, so any external power supply is not necessary.

In order to drive the purge and vent valves using external commands, use the IP54 cable version, enter the configuration menu, set the “valve setting” on the value “on command”, and connect the cables as shown in the figure. The valves are n.c., providing 24V between vent purge and common terminal the valves will open.



**Figure 1**

## 16 Instructions for Use Installation



**Figure 2** Connections

### NOTE

The Turbo-V Remote Panel can be used as a bench unit or a rack module, but it must be positioned so that free air can flow through the holes. Do not install or use the controller in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

During operation, the following environmental conditions must be respected:

- temperature: from 5 °C to +45 °C
- relative humidity: 0 – 95 % (without condensation)

See the appendix "Technical Information" for detailed information about the above mentioned and the other connections, and about the options installation.

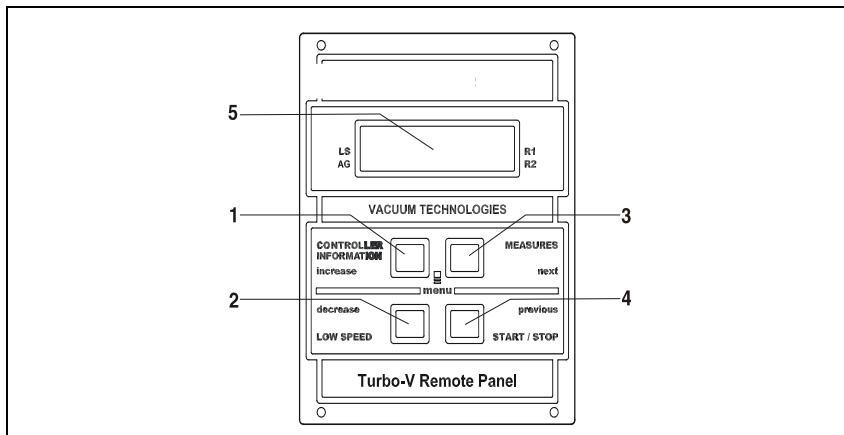
## Use

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".

## Controls, Indicators and Connectors

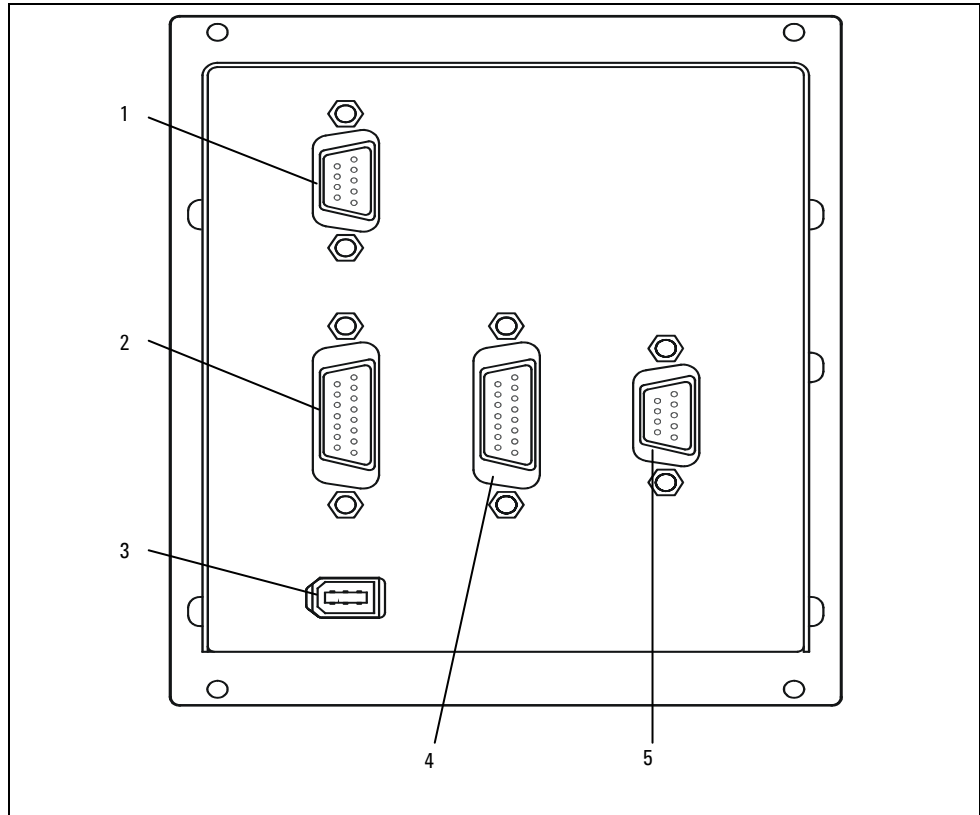
The following paragraph illustrates the Controller front panel and rear panel. More details are contained in the appendix "Technical Information".

## Front Panel Description



**Figure 3** Turbo-V Remote Panel

- 1 Keyboard push-button to recall on the display the cycle number, cycle time and pump life.
- 2 Keyboard push-button for LOW SPEED mode selection. It is active only when the front panel operation has been selected. Pressed once, the pump runs at “stand-by” speed. To unselect the mode, press the push-button again.
- 3 Keyboard push-button to recall on the display the pump current, pump temperature, pump power and rotational speed. It is always active regardless of the operating mode selected. Push-buttons 3 and 4, if pressed together for at least 2 seconds, put the controller in a routine where it is possible to program some operation parameters.
- 4 Keyboard push-button for START, STOP, RESET mode selection. It is active only when the front panel operation has been selected. By pressing once the starting phase begins; if pressed again it stops the pump. If the pump has been stopped automatically by a fault, this push-button must be pressed once to reset the controller and a second time to restart the pump.
- 5 LCD back-lighted alphanumeric display: dot matrix 2 lines x 16 characters.



**Figure 4** Turbo-V Remote Panel Rear panel

- 
- |   |                                                                       |
|---|-----------------------------------------------------------------------|
| 1 | Logic input signals connectors (mating connector supplied with link). |
|---|-----------------------------------------------------------------------|
- 
- |   |                                                                                            |
|---|--------------------------------------------------------------------------------------------|
| 2 | Logic output signals connector and pump frequency monitor plus programmable analog output. |
|---|--------------------------------------------------------------------------------------------|
- 
- |   |                                            |
|---|--------------------------------------------|
| 3 | Power output connector (24 Vdc for Gauge). |
|---|--------------------------------------------|
- 
- |   |                                                                                                       |
|---|-------------------------------------------------------------------------------------------------------|
| 4 | Connector for power input from the pump controller and serial communication with the pump controller. |
|---|-------------------------------------------------------------------------------------------------------|
- 
- |   |                                  |
|---|----------------------------------|
| 5 | Remote control serial connector. |
|---|----------------------------------|
-

## **Switch-on the Remote Panel**

To switch-on the Turbo-V Remote Panel plug the connection cable to the controller on both sides and then power on the controller.

## **Starting the Pump**

To start the pump, simply press the START pushbutton on the Turbo-V Remote Panel.

## **Pump Shutdown**

To shutdown the pump press the STOP pushbutton on the Turbo-V Remote Panel.

## Maintenance

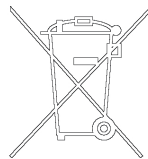
The Turbo-V Remote Panel does not require any maintenance. Any work performed on the controller must be carried out by authorized personnel.

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.

If a controller is to be scrapped, it must be disposed of in accordance with the specific national standards.

## 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.



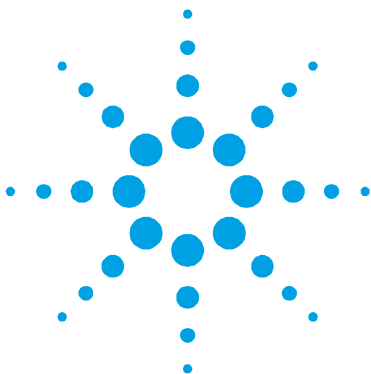
## Error Messages

For a certain type of failure, the Turbo-V Remote Panel will show a list of error message described in the following table.

MESSAGE	DESCRIPTION	REPAIR ACTION
COMMUNICATION ERROR	The Remote Panel is unable to start a communication with the controller.	Check serial connection between the Remote Panel and the controller.
COMMUNICATION BROKEN	The Remote Panel is unable to continue the communication with the controller.	Check serial connection between the Remote Panel and the controller.
XXXXXXX MODEL	Controller not supported. The Remote Panel works only with the following pumps: TV551 with navigator controller, TV701 with navigator controller, TV1001 with navigator controller, TV2k-G and TV3k-T.	Connect the Remote Panel to a supported controller.
CHECK CONNECTION TO PUMP	Wrong connection between the pump and the controller. Or The pump's temperature is below 0 °C.	Check connection between controller and pump. Press the START push-button twice to start the pump.
WAITING INTERLOCK	The interlock signal of P1 connector is activated by an interruption of the link between pin 3 and 8 of J1 connector, or because the external interlock signal is open.	Reset the short circuit between pin 3 and pin 8 of J1 connector, or close the external interlock signal.
PUMP OVERTEMP.	The upper bearing/pump temperature exceeds 60 °C.	Wait until the temperature decrease below threshold value. Press the START push-button twice to start the pump.
CONTROLLER OVERTEMP.	The controller environment temperature exceeds 70 °C. Or The controller's radiator temperature is above 60 °C.	Wait until the temperature decrease below threshold value. Press the START push-button twice to start the pump.

<b>MESSAGE</b>	<b>DESCRIPTION</b>	<b>REPAIR ACTION</b>
TOO HIGH LOAD	The current drawn by the pump is higher than programmed.	Check that the pump rotor is free to rotate. Press the START push-button twice to start the pump.
SHORT CIRCUIT	After the starting phase the output connection is shorted.	Check connections and shortages between pump and controller. Press the START push-button twice to start the pump.
SYSTEM OVERRIDE	The pump is stopped by an emergency stop signal provided via a remote contact.	Check the emergency condition which is causing the override signal on the P1 connector. Remove the override signal on P1 connector, then press the START push-button twice to start the pump.
OVERVOLTAGE	Controller power supply circuitry is faulty, or the Controller received a spike.	Press the START push-button twice to start the pump. Should the message still be present, call the Agilent service.
POWER FAIL	Failure in the controller's pump power supply section.	Contact Agilent for Maintenance.

## **16** Instructions for Use Error Messages



## 17 Technical Information

Specifications	239
Controller Outline	240
Controller Connection	242
P1 – Input	242
J1 – Output	244
J5 – Gauge Connector	251
J2 – Serial Connector	252
Serial Communication Descriptions	254
Window Protocol Description	255
Description	255
Communication Format	255
Communication Protocol	256
Window Meanings	262
How to use by Front Panel	266
Use	272
General	272
Startup	273

Original Instructions



## 17 Technical Information

Programming	278
Configuration Menu	278
Mode Menu	280
Pump Setting Menu for TV2k-G, TV3k-T	281
Pump Setting Menu for TV551, TV701 and TV1001	283
INPUT/OUTPUT Menu	285
Setpoint R1 and Setpoint R2 Submenus	288
Gauge Menu	290
Serial Menu	291
How to use by Remote I/O	292
How to use in Serial Mode	292
Error Messages	294
Orderable Parts	297

## Specifications

Input:	
Voltage	24 Vdc
Power	2 W
Environment storage	-20 °C to +70 °C 0-95 %
Operating	+5 - +45 °C 0 – 90 %
Applied norms.	EN61010-1 EN61000-4-2 EN61000-4-3 EN61000-4-4
Weight:	0.8 Kg
Pollution Degree	2
Installation Category	II

## Controller Outline

The outline dimensions for the Turbo-V Remote Panel and cables are shown in the following figures.

**NOTE**

The Turbo-V Remote Panel 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.

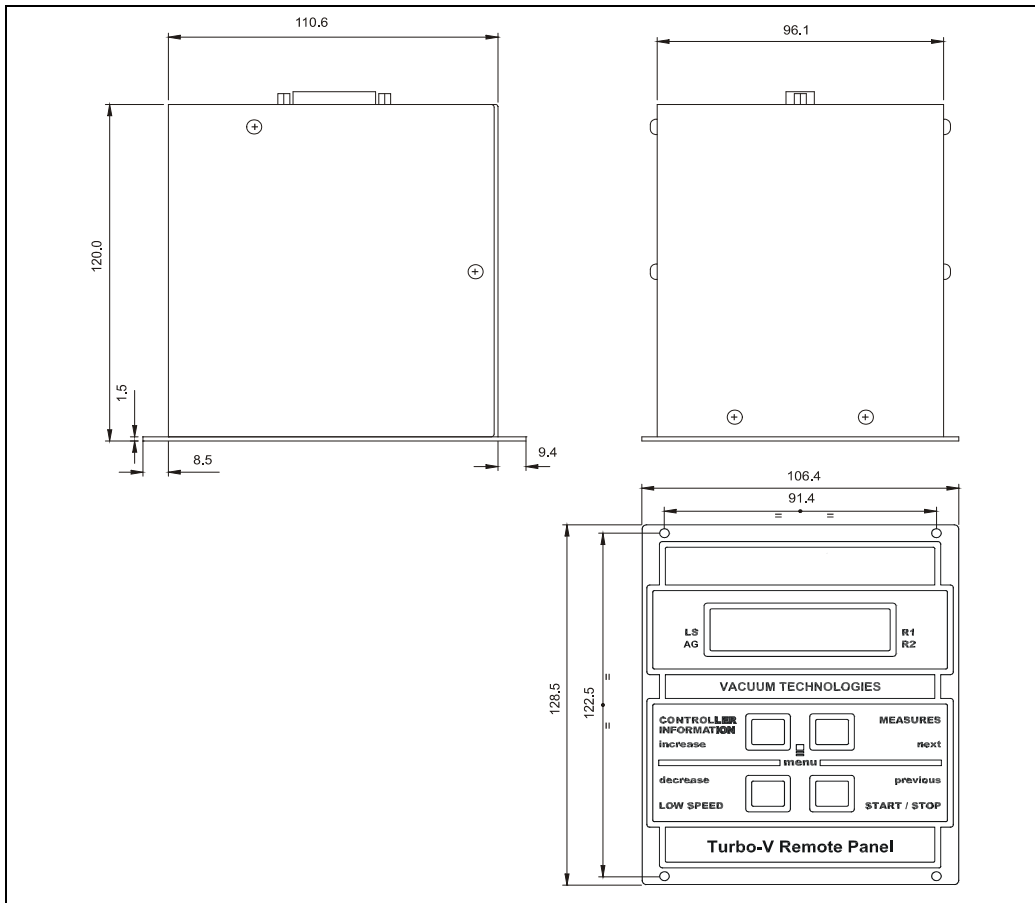
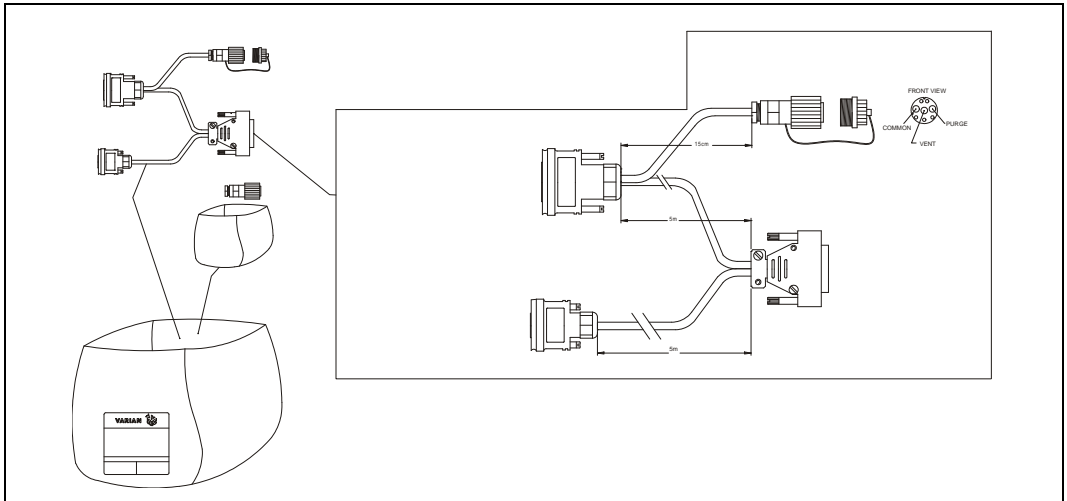
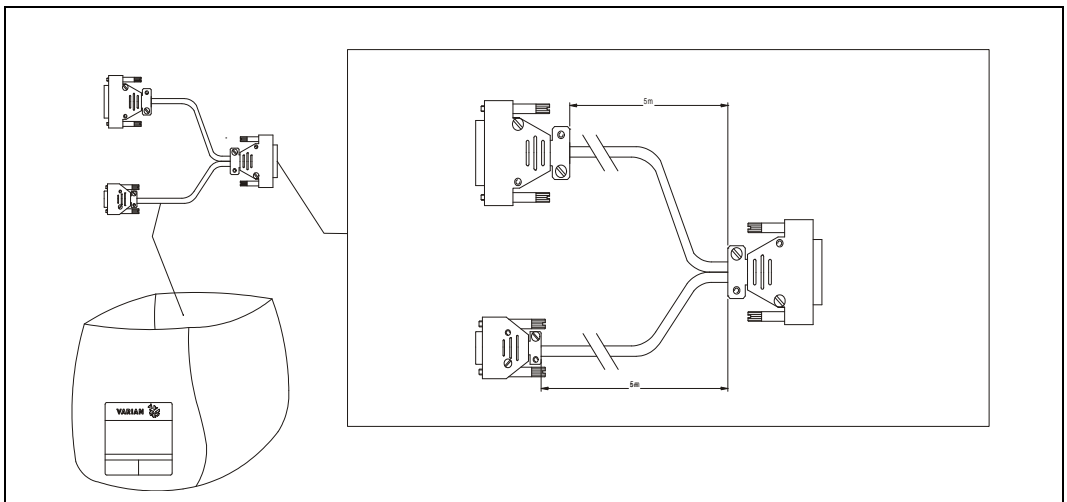


Figure 1 Outline



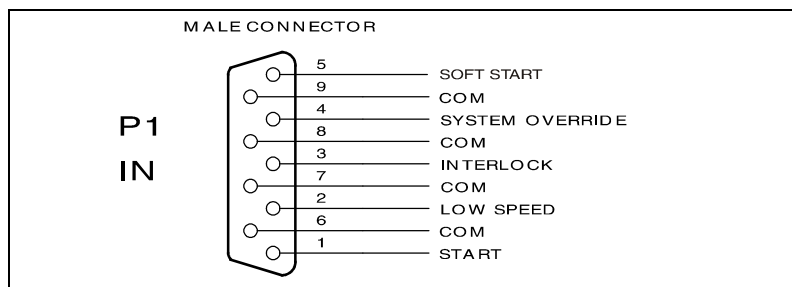
**Figure 2** Remote panel IP54 Cable – 969-9945



**Figure 3** Remote panel Cable – 969-9946

## Controller Connection

### P1 – Input



**Figure 4** P1 Input Connector

All the logic input to the controller **must** be connected at J1 mating connector.

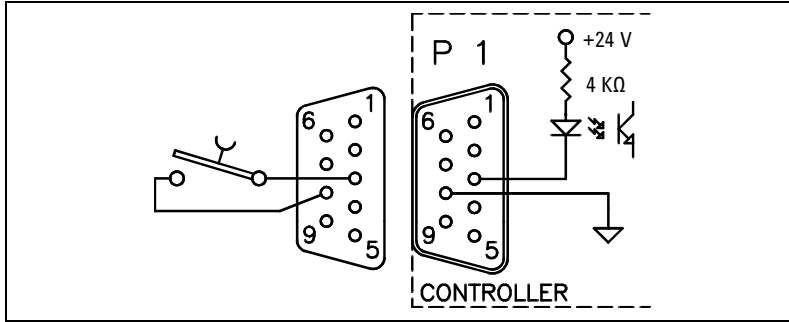
With the provided J1 mating connector (shipped with pin 3 and pin 8 shorted) make the connections with AWG 24, (0.24 mm<sup>2</sup>) or smaller wire to the pins indicated in the figure to obtain the desired capability.

The following table describes the signals available on the connector.

Pins	Name	Description	Electrical	Note
1-com	Start/Stop	Contact closed = Start Contact open = Stop	Optically isolated Vlow, max = 10V Imax = 5mA Imin (on) = 2mA	Active only in Remote Mode operation
2-com	Low Speed	Contact closed = Low Speed Contact open = High Speed	Optically isolated Vlow, max = 10v Imax = 5mA	Active only in Remote Mode operation
3-com	Interlock	Contact closed = Active Contact open = Inactive If active allows the pump running, if inactive stops the pump.	Optically isolated Vlow, max = 10V Imax = 5mA Imin (on) = 2mA	Active in all modes
4-com	System Override	Contact closed = Active Contact open = Inactive If inactive allows the pump running, if active stops the pump and all the others connected devices (Fan, Vent Valve...).	Optically isolated Vlow, max = 10V Imax = 5mA Imin (on) = 2mA	Active in all modes
5-com	Soft Start	Contact closed= Active Contact open= Inactive If active before the start command activates the soft start procedure. To be used only at first pump run-up.	Optically isolated Vlow, max = 10V Imax = 5mA Imin (on) = 2mA	Active only in Remote Mode operation, and in STOP status.

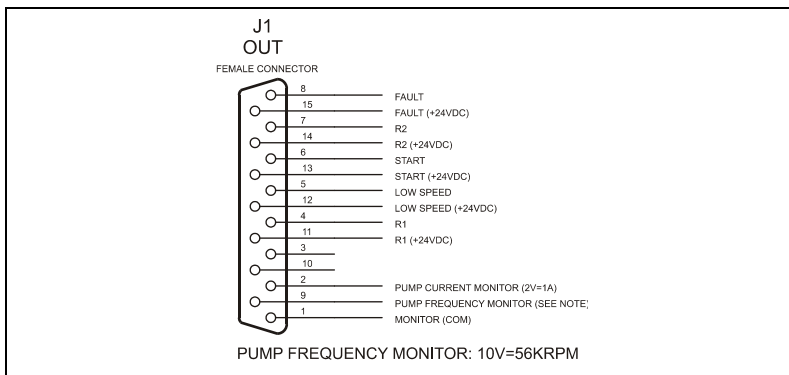
**17 Technical Information**  
**Controller Connection**

The following figure shows a typical contact logic input connection and the related simplified circuit of the controller.



**Figure 5** Typical Logic Input Connection

## J1 – Output



**Figure 6** Logic Output Connector

All the logic output from the controller must be connected at P2 mating connector. The following table describes the signals available on the connector.

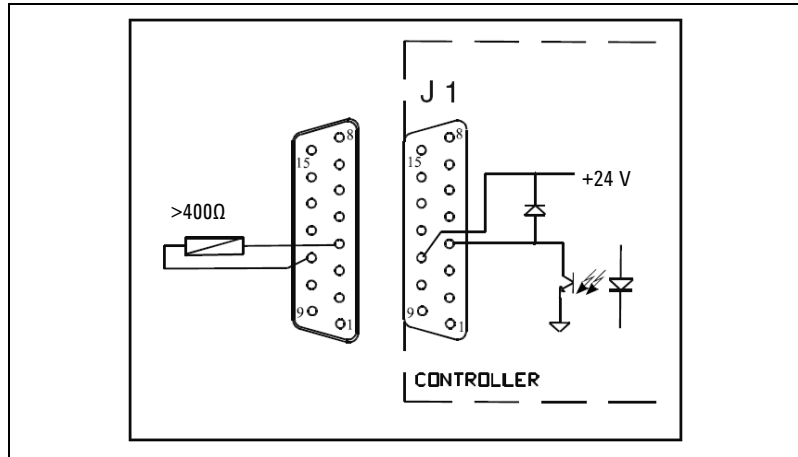
Pins	Name	Description	Electrical	Note
15-8	Fault	0V = no fail 24V = fail	Optically isolated Vmax (=open) = 24v Vlow, max = 1,5v Imax = 60mA Pin-8 negative Pin-15 positive	Connect a 10kohm resistor between the 2 pins before measuring voltage A STOP command (by front Panel or serial or Start/stop input) reset the FAIL status.
13-6	Starting / Running	Configurable output. Default setting (win.177 = 0): 24V pump in RAMP state, 0V = other state. If win.177 = 1: 0V = pump in stop or FAIL status, 24V = other state	Optically isolated Vmax (=open) = 24v Vlow, max = 1,5v Imax = 60mA Pin-6 negative Pin-13 positive	Connect a 10kohm resistor between the 2 pins before measuring voltage
12-5	Low Speed	24V = low speed active 0V = low speed not active	Optically isolated Vmax (=open) = 24v Vlow, max = 1,5v Imax = 60mA Pin-5 negative Pin-12 positive	Connect a 10kohm resistor between the 2 pins before measuring voltage
11-4	R1	Programmable Set Point Out Can be related to: Frequency, Power, Time Status or Pressure. See following figures for more details. All the setting can be done by: front panel or serial line (win.101 to 105 and 162).	Optically isolated Vmax (=open) = 24v Vlow, max = 1,5v Imax = 60mA Pin-5 negative Pin-12 positive	Connect a 10kohm resistor between the 2 pins before measuring voltage

## 17 Technical Information

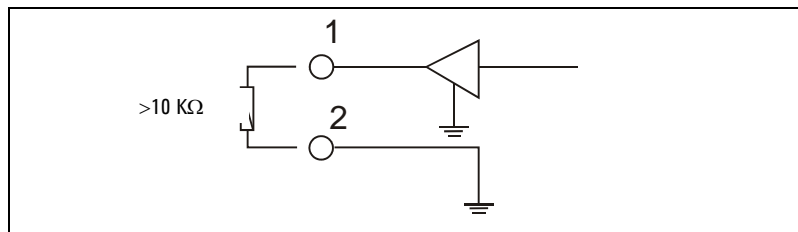
### Controller Connection

Pins	Name	Description	Electrical	Note
14-7	R2	<p>Programmable Set Point Out</p> <p>Can be related to: Frequency, Power, Time Status or Pressure.</p> <p>See following figures for more details. All the setting can be done by: front panel or serial line (win.171 to 176).</p>	<p>Optically isolated</p> <p><math>V_{max} (=open) = 24v</math></p> <p><math>V_{low, max} = 1,5v</math></p> <p><math>I_{max} = 60mA</math></p> <p>Pin-7 negative</p> <p>Pin-14 positive</p>	<p>Connect a 10kohm resistor between the 2 pins before measuring voltage</p>
2-1	Programmable Analog Output	<p>Programmable analog output.</p> <p>Can be related to : frequency, power, pump temperature (linear) or pressure (logarithmic).</p> <p>The setting can be done by: front panel or serial line (win.111).</p> <p>0V = 0Hz, 0W, 0 °C</p> <p>10V = 963Hz, 150W, 100 °C</p> <p>See following chart for more details</p>	<p>Optically isolated</p> <p><math>R_{load, min} = 10k\Omega</math></p> <p>Ripple <math>V_{pp,max} =</math> 200mv</p> <p>Precision = 5 %</p> <p>Resolution = 40mv</p> <p>Pin-1 negative</p> <p>Pin-2 positive</p>	
9-1	Frequency Analog Output	<p>Analog output proportional to frequency (linear).</p> <p>0V = 0Hz</p> <p>10V = 963Hz</p>	<p>Optically isolated</p> <p><math>R_{load, min} = 10k\Omega</math></p> <p>Ripple <math>V_{pp,max} =</math> 200mv</p> <p>Precision = 1 %</p> <p>Resolution = 40mv</p> <p>Pin-1 negative</p> <p>Pin-9 positive</p>	

The following figure shows a typical logic output connection (relay coil) but any other device may be connected e.g. a LED, a computer, etc., and the related simplified circuit of the controller.

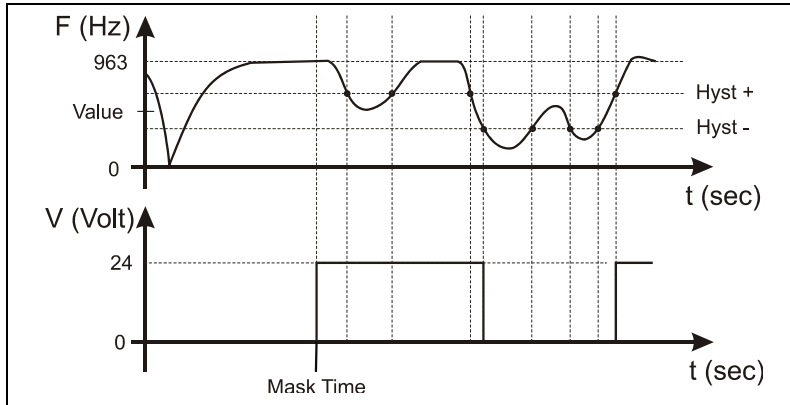


**Figure 7** Digital Output Connection Example

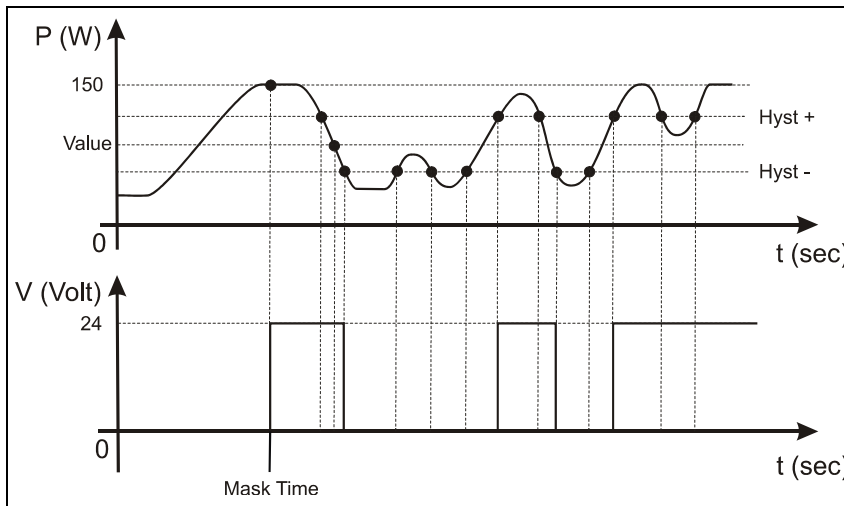


**Figure 8** Analog Output Connector Example

**R1 – R2 Output Diagrams**



**Figure 9** R1 or R2 related to frequency



**Figure 10** R1 or R2 related to power

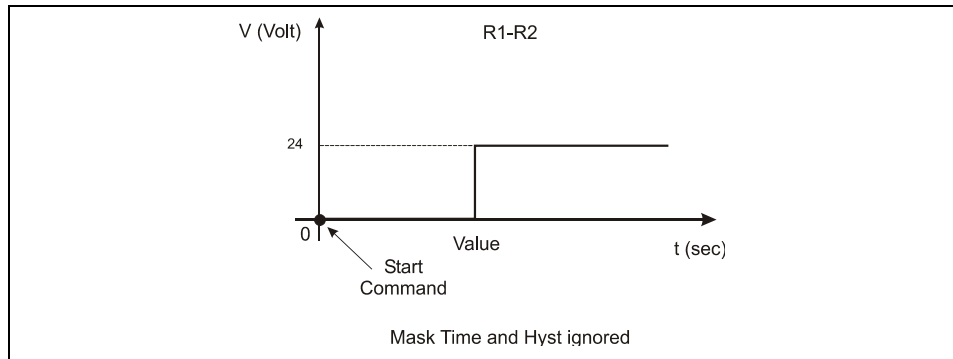


Figure 11 R1 or R2 related to time

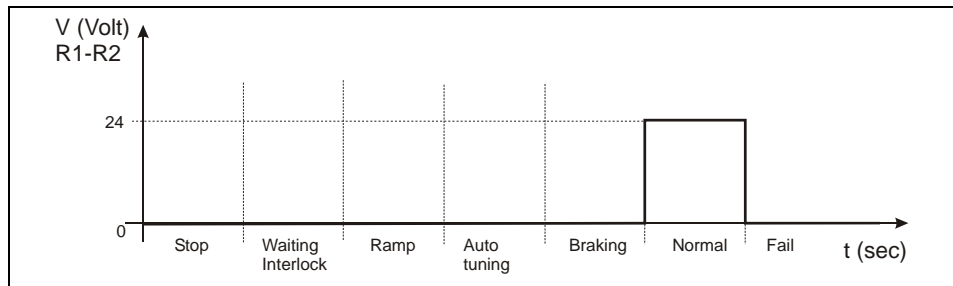


Figure 12 R1 or R2 related to Normal status

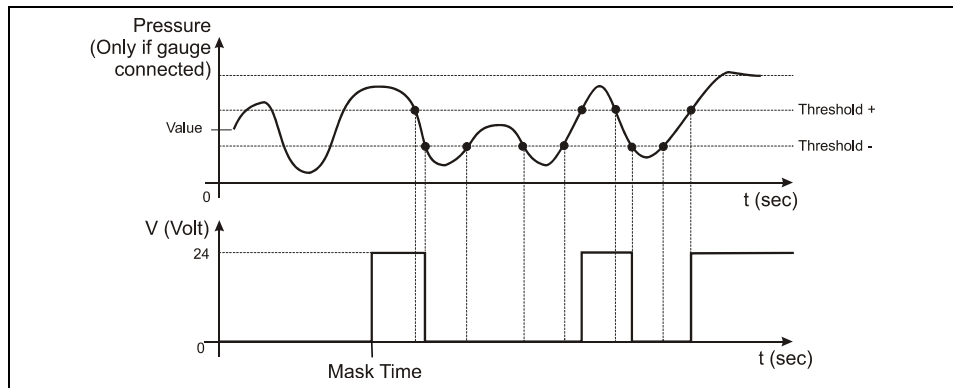


Figure 13 R1 or R2 related to pressure reading (if gauge connected)

### Programmable Analog Output Diagrams

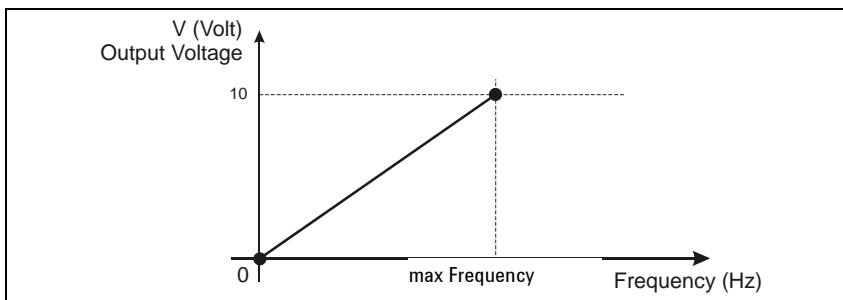


Figure 14 Programmable analog output related to frequency

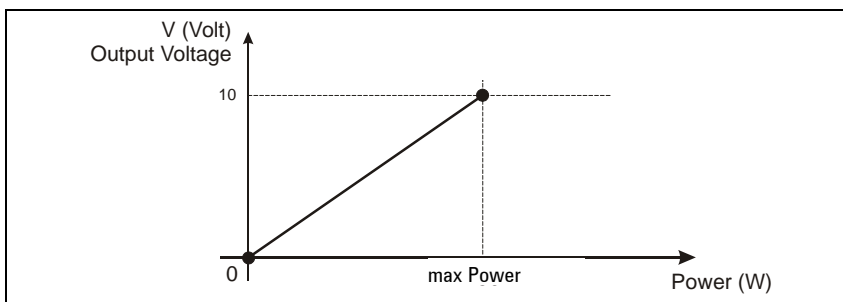


Figure 15 Programmable analog output related to power

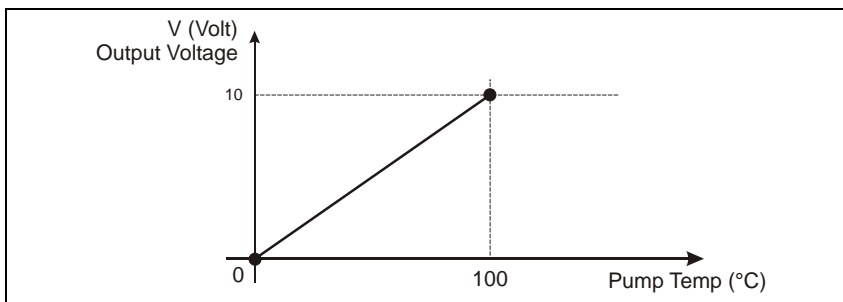
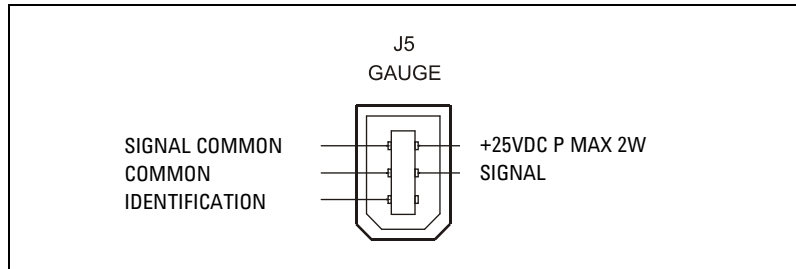


Figure 16 Programmable analog output related to pump temperature

## J5 – Gauge Connector

This connector is available to connect “EyeSys Mini-IMG” gauge and EyeSys ConvecTor® gauge with their cables (see ordering parts table).

The pressure can be read by display or by serial line (win. 224) or by REMOTE I/O on Programmable Analog Out (in this case the Remote Panel replicates the signal coming from the gauge).



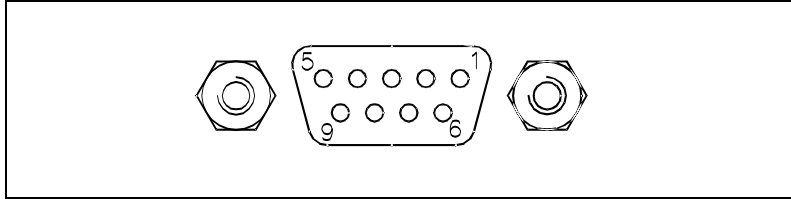
**Figure 17** Gauge Connector

**NOTE**

The Mini-IMG gauge is intended for high vacuum use, so with pressure above 50 mTorr the signal is misinterpreted and a “false” pressure is displayed. Use a transition gauge to understand if the reading is really under 50 mTorr or not.

## J2 – Serial Connector

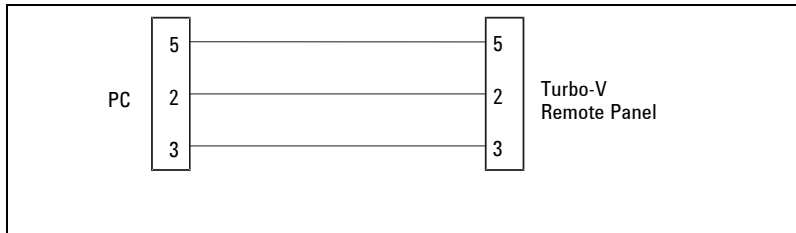
This connector provides the connection for RS – 232 and RS – 485 serial line.



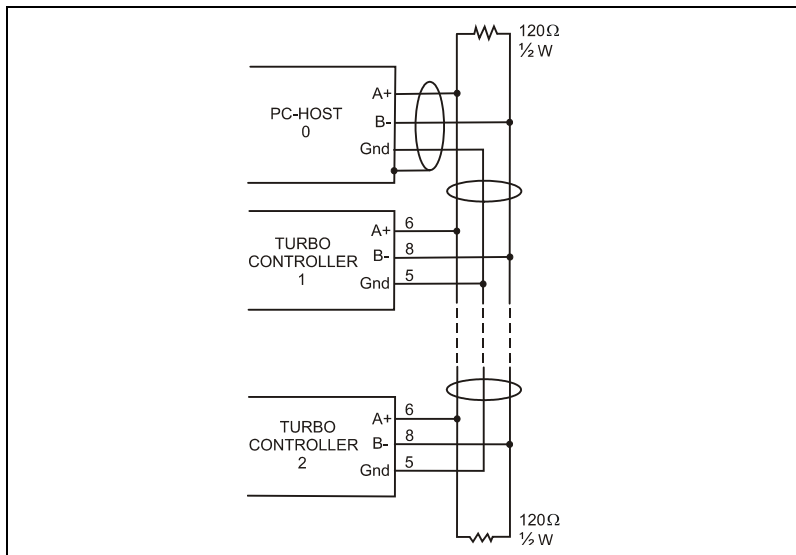
**Figure 18** J2 Pin-out

PIN N.	SIGNAL NAME
1	+5 V (OUT) (Reserved)
2	TX (RS232)
3	RX (RS232)
4	Not Connected
5	GND
6	A + (RS485)
7	Not Connected
8	B – (RS485)
9	Reserved

**Connector Examples**



**Figure 19** RS – 232 Connection



**Figure 20** RS – 485 Connection

The communication port mating connector is supplied with the RS 232 PCB (AMP/Cannon or equivalent 19-pin "D" type male connector).

## 17 Technical Information

### J2 – Serial Connector

For example, the Transmit data signal from controller (pin 2) must be connected to the host computer's receive data line (pin 2) and vice versa. Consult the host computer's instruction manual for its serial port connections.

#### NOTE

Agilent cannot guarantee compliance with FCC regulations for radiated emissions unless all external wiring is shielded, with the shield being terminated to the metal shroud on the 0-subconnector. The cable should be secured to the connector with screws.

---

## Serial Communication Descriptions

This unit can communicate by the “Window” protocol. This protocol can be used as well with 232 or 485 media.

# Window Protocol Description

## Description

The remote panel acts as a gateway between the host and the controller. Some command are managed directly by the Remote Panel, which sends the appropriate commands to the controller. Other commands are just passed to the controller and then the answer is sent back to the host. This section of the manual describes the former commands, while for the latter commands refer to the manual of the controller the remote panel is connected to. In case of duplicated commands the first types ones have always the priority.

### NOTE

The Window 008 of the controller to which the remote panel is attached to is always fixed to the value 0 (Serial), in order to allow the Remote Panel to write data on the controller. Trying to set this window to any other value has no effect (the controller will turn it back to 0).

---

## Communication Format

- 8 data bit
- no parity
- 1 stop bit
- baud rate: 600/1200/2400/4800/9600 programmable

## Communication Protocol

The communication protocol is a MASTER/SLAVE type where:

- Host = MASTER
- Controller = SLAVE

The communication is performed in the following way:

1. the host (MASTER) send a MESSAGE + CRC to the controller (SLAVE);
2. the controller answer with an ANSWER + CRC to the host.

The MESSAGE is a string with the following format:

<STX>+<ADDR>+<WIN>+<COM>+<DATA>+<ETX>+<CRC>

where:

### NOTE

When a data is indicated between two quotes ('...') it means that the indicated data is the corresponding ASCII character.

- <STX> (Start of transmission) = 0x02
- <ADDR> (Unit address) = 0x80 (for RS 232)
- <ADDR> (Unit address) = 0x80 + device number (0 to 31) (for RS 485)
- <WIN> (Window) = a string of 3 numeric character indicating the window number (from '000' to '999'); for the meaning of each window see the relevant paragraph.
- <COM> (Command) = 0x30 to read the window, 0x31 to write into the window
- <DATA> = an alphanumeric ASCII string with the data to be written into the window. In case of a reading command this field is not present. The field length is variable according to the data type as per the following table:

Data Type	Field Length	Valid Characters
Logic (L)	1	'0' = OFF '1' = ON
Numeric (N)	6	'-', '.', '0' ... '9' right ustified with '0'
Alphanumeric (A)	10	from blank to '_' (ASCII)

- <ETX> (End of transmission) = 0x03
- <CRC> = XOR of all characters subsequent to <STX> and including the <ETX> terminator. The value is hexadecimal coded and indicated by two ASCII character.

The addressed SLAVE will respond with an ANSWER whose structure depends from the MESSAGE type. When the MESSAGE is a reading command, the SLAVE will respond transmitting a string with the same structure of the MESSAGE.

**NOTE**

Using the RS 485 interface, the message structure remains identical to the one used for the RS 232 interface, the only difference being that the value assigned to the ADDRESS <ADDR>.

## 17 Technical Information

### Window Protocol Description

The controller can answers with the following response types:

Response Type	Response Length	Response Value	Description
Logic	1 byte	-	after a read instruction of a logic window
Numeric	6 bytes	-	after a read instruction of a numeric window
Alphanumeric	10 bytes	-	after a read instruction of an alphanumeric window
ACK	1 byte	(0x6)	the command execution has been successfully completed
NACK	1 byte	(0x15)	the command execution has been failed
Unknown Window	1 byte	(0x32)	the specified window in the command is not a valid window
Data Type Error	1 byte	(0x33)	the data type specified in the command (Logic, Numeric or Alphanumeric) is not accorded with the specified Window
Out of Range	1 byte	(0x34)	the value expressed during a write command is out of the range value of the specified window
Win Disabled	1 byte	(0x35)	the specified window is Read Only or temporarily disabled (for example you can't write the Soft Start when the Pump is running)

**Examples:**

**Command: START**

Source: PC

Destination: Controller

02	80	30	30	30	31	31	03	42	33
STX	ADDR	WINDOW			WR	ON	ETX	CRC	

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

**Command: STOP**

Source: PC

Destination: Controller

02	80	30	30	30	31	30	03	42	32
STX	ADDR	WINDOW			WR	OFF	ETX	CRC	

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

**17 Technical Information**  
**Window Protocol Description**

**Command: LOW SPEED ACTIVATION (YES)**

Source: PC

0280 3030 3131 3003 4233

Destination: Controller

02	80	30	30	31	31	31	03	42	32
STX	ADDR	WINDOW			WR	YES	ETX	CRC	

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

**Command: LOW SPEED ACTIVATION (NO)**

Source: PC

Destination: Controller

02	80	30	30	31	31	31	03	42	33
STX	ADDR	WINDOW			WR	NO	ETX	CRC	

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

**Command: READ PUMP STATUS**

Source: PC

Destination: Controller(with address = 3)

02	83	32	30	35	30	03	38	37
STX	ADDR	WINDOW			RD	ETX	CRC	

Source: Controller(with address = 3 in stop status)

Destination: PC

02	83	32	30	35	30	30	30	30	30	03	38	37
STX	ADDR	WINDOW			DATA (STATUS)					ETX	CRC	

**Command: READ SERIAL TYPE**

Source: PC

Destination: Controller(with address = 3 in 485 mode)

02	83	35	30	34	30	03	38	37
STX	ADDR	WINDOW			RD	ETX	CRC	

Source: Controller

Destination: PC

02	83	35	30	34	30	31	03	42	30
STX	ADDR	WINDOW			RD	DATA	ETX	CRC	

## Window Meanings

<b>N.</b>	<b>Read/ Write</b>	<b>Data Type</b>	<b>Description</b>	<b>Admitted Values</b>
000	R/W	L	Start/Stop (in remote/ Front mode the window is a read only)	Start = 1 Stop = 0
001	R/W	L	Low Speed Activation	No = 0 Yes = 1 (default = 0)
107	R/W	N	Front (default)/ Remote/ Serial configuration	Front = 0 Remote = 1 Serial = 2 (default = 0)
101	R/W	N	R1 Set Point type	0 = Frequency 1 = Power 2 = Time 3 = Normal 4 = Pressure (default = 3)
103	R/W	N	R1 Set Point value (expressed in Hz, W or s)	
104	R/W	N	Set Point delay: time between the pump start and the set point check (seconds)	0 to 99999 (default = 0)
105	R/W	N	Set point hysteresis (in % of value)	0 to 100 (default = 2)
108	R/W	N	Baud rate	600 = 0 1200 = 1 2400 = 2 4800 = 3 9600 = 4 (default = 4)

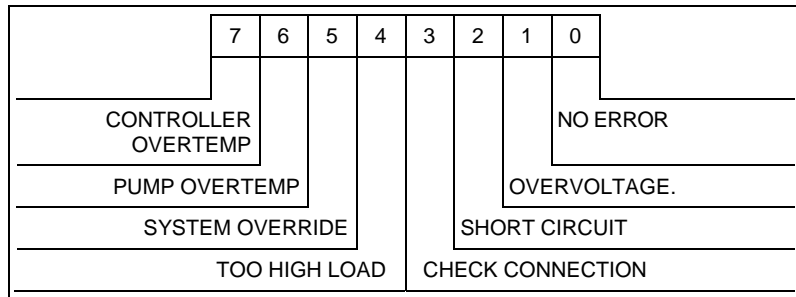
<b>N.</b>	<b>Read/ Write</b>	<b>Data Type</b>	<b>Description</b>	<b>Admitted Values</b>
110	R/W	L	Interlock type (default = 1)	Impulse = 0 Continuous = 1
111	R/W	L	Analog output type: output voltage signal	0 = frequency 1 = power 2 = pump temperature 3 = pressure (default = 1)
161	R/W	N	Pressure Reading Correction Factor See "GAUGE CONNECTOR" for more details	0 to 10 0 = 0 10 = 1 = N <sub>2</sub>
162	R/W	A	R1 Set Point Pressure Value Valid if win. 101 = 4 Format X.X E. XX Where X = 0 to 9 s = + or -	
163	R/W	N	Pressure unit of measure	0 = mBar 1 = Pa 2 = Torr
171	R/W	N	R2 Set Point Type	0 = Freq 1 = Power 2 = Time 3 = Normal 4 = Pressure
172	R/W	N	R2 Set Point Value(Hz, W, s)	
173	R/W	N	Set Point delay: time between the pump start and the set point check (seconds)	
174	R/W	L	R2 Set Point Signal Activation Type	0 = high level active 1 = low level active
175	R/W	N	R2 Set front Hysteresis (in % of R2 Valve)	

## 17 Technical Information

### Window Protocol Description

N.	Read/ Write	Data Type	Description	Admitted Values
176	R	A	R2 Set Point Pressure Value Valid in win 171 = 4 Format X.X E X.X Where: X= 0 to 9 s = + or -	
200	R	N	Pump current in mA dc	
203	R	N	Driving frequency in Hz	
205	R	N	Pump status	Stop = 0 Waiting intlk = 1 Starting = 2 Normal = 3 High Load = 4 Fail = 5 Approaching low speed = 6
206	R	N	Error code	Bit description: see the following figure
206	R	N	Controller Heatsink Temperature ( °C)	
207	R	L	R1 Set Point Status	Off: 0 On: 1
208	R	L	R2 Set Point Status	Off: 0 On: 1
221	R	L	R1 Set Point Status	Off: 0 On: 1
224	R	A	Pressure reading Format = X.X E XX	
225	R	L	R2 Set Point Status	Off: 0 On: 1
408	R	A	Firmware revision	

N.	Read/ Write	Data Type	Description	Admitted Values
503	R/W	N	RS 485 address	0 to 31 (default = 0)
504	R/W	L	Serial type select	0 = RS 232 1 = RS 485 (default = 0)



**Figure 21** Window N. 206 Bit Description

## How to use by Front Panel

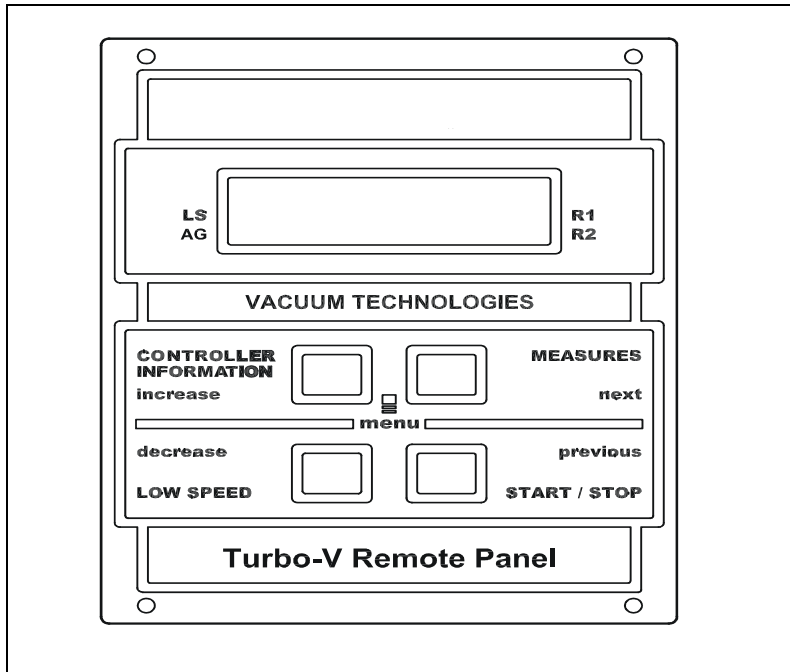


Figure 22

### The main Panel

After startup the following panel is displayed.

X	X	X	X	X	X	X	X			X	X	X	X	H	z
					X	X	X	X	X	X					

Where:

**XXXXXX:** indicates the status of the pump

**XXXHz:** is the current frequency of rotation of the pump. This value can be modified by the user according to needs, proceeding as described in the “PROGRAMMING” paragraph. The unit of measurement (HZ / Krpm, default = Hz) is selected by the user from the INPUT/OUTPUT->SPEED menu.

The operating mode of the controller is shown at the bottom in the center. This may be:

**FRONT:** front mode (the controller accepts START/STOP and LOW SPEED commands from the front panel)

**REMOTE:** remote mode (the controller accepts START/STOP and LOW SPEED commands via the rear connector)

**RS232:** the controller accepts commands from the serial port according to the RS232 standard.

Black rectangles as shown below may be displayed at any time in the first and last box of the bottom line of the display (opposite the LS / AG / R1 / R2 captions). The meaning of these rectangles is as follows:

**LS** → the controller is at LOW SPEED

**AG** → the Gauge is connected to the controller

**R1** → relay (set point ) R1 is active

**R2** → relay (set point ) R2 is active

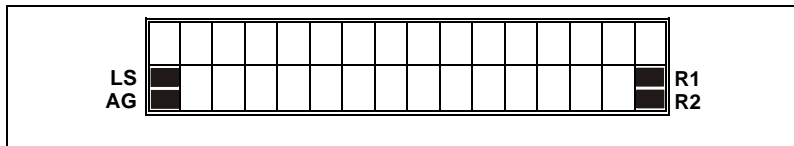


Figure 23

## 17 Technical Information

### How to use by Front Panel

#### Start Stop and Reset Fail button

Press once for the START command and again for the STOP command. In the case of controller's failure, press once to reset the fail condition.

#### NOTE

The front panel START/STOP function is available only if the controller is set to "FRONT mode". If it is set to "REMOTE mode", when the key is pressed an error message is shown indicating "COMMAND LOCKED IN REMOTE", if it is set to "SERIAL" mode, the "COMMAND LOCKED IN SERIAL MODE" error message is displayed.

The "Reset Fail" function is always available regardless of controller mode.

---

#### Low Speed

Press once to activate the LOW SPEED command; press again to de-activate.

#### NOTE

The front panel LOW SPEED function is available only if the controller is set to "FRONT"; otherwise an error message is returned.

If it is set to "REMOTE" mode, the COMMAND LOCKED IN REMOTE error message is displayed while if it is in "SERIAL" mode the "COMMAND LOCKED IN SERIAL MODE" error message is returned.

---

### Controller Information

Pressing the first time, the following information is displayed:

		<b>C</b>	<b>Y</b>	<b>C</b>	<b>L</b>	<b>E</b>		<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>		
	<b>T</b>	<b>I</b>	<b>M</b>	<b>E</b>				<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>m</b>

Where

**CYCLE**    XXXX    Number of start/stop cycles

**TIME**    XXXXXm    Time of last cycle (minutes)

Pressing a second time, the following information is displayed:

		<b>P</b>	<b>U</b>	<b>M</b>	<b>P</b>		<b>L</b>	<b>I</b>	<b>F</b>	<b>E</b>			
		<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>h</b>				

#### **PUMP LIFE**

**XXXXXX**    total operating time (hours) of the pump

Pressing a third time, the following message is displayed:

	<b>M</b>	<b>N</b>	<b>:</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>
	<b>S</b>	<b>N</b>	<b>:</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>	<b>Y</b>

**MN**    Model Number of the controller

**SN**    Serial Number of the controller

On pressing the key again, the main panel is redisplayed.

**17 Technical Information**  
 How to use by Front Panel

**Measures**

Pressing once, the following is obtained:

<b>P</b>	<b>=</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>W</b>					<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>H</b>	<b>z</b>
					<b>T</b>	<b>=</b>	<b>X</b>	<b>X</b>	<b>°</b>	<b>C</b>					

Where:

**P= XXX W** is the current power supplied to the pump

**T=XX °C** is the current temperature of the pump

**XXXX Hz** is the current frequency of rotation of the pump. This value can be modified by the user according to needs, proceeding as described in the “PROGRAMMING” paragraph. The unit of measurement (HZ / Krpm, default = Hz) is selected by the user from the INPUT/OUTPUT->SPEED menu.

Pressing twice, the following is obtained:

<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>					<b>X</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>H</b>	<b>z</b>
	<b>I</b>	<b>=</b>	<b>X</b>	<b>.</b>	<b>X</b>	<b>X</b>	<b>A</b>		<b>V</b>	<b>=</b>	<b>X</b>	<b>X</b>	<b>X</b>	<b>V</b>	

Where:

**XXXXXX**: indicates the status of the controller

**XXXHz**: see description “Pressing once” **Measures** key

**I=X.XXA**: indicates the current absorbed by the pump

**V= XXX V**: indicates the pump driving voltage

Pressing 3 times, the following is obtained:

P	=	X	X	X	W					X	X	X	X	H	z
		X	.	X	X	E	-	Y	Y	m	B	a	r		

Where:

**P= XXXW:** indicates the current power supplied to the pump

**XXXHz:** see description “Pressing once” Measures key

**x.xxE-yy mbar:** indicates the pressure currently measured by the gauge.

Other possible indications:

P	=	X	X	X	W					X	X	X	X	H	z
		G	A	U	G	E		E	R	R	O	R			

Where:

**GAUGE\_ERROR:** gauge not connected or broken

P	=	X	X	X	W					X	X	X	X	H	z
		P	R	E	S	S	U	R	E	U	N	D	E	R	

Where:

**PRESS UNDER:** indicates that the pressure read is below the minimum measurement threshold set (5 E-9 mBar)

P	=	X	X	X	W					X	X	X	X	H	z
		P	R	E	S	S	U	R	E	O	V	E	R		

Where:

**PRESS OVER:** indicates that the pressure measured is above the maximum measurement threshold set (1000 mBar)

The unit of measurement of the pressure can be selected (GAUGE menu→Unit) from “mBar”, “Pa”, “torr”.

Pressing 4 times the main panel is displayed again.

# Use

## General

Make all vacuum manifold and electrical connections and refer to Turbo-V pump instruction manual before to operating the Turbo-V Remote Panel.

---

### **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 Turbo-V pump if the pump inlet is not connected to the system or blanked off.**

---

### **NOTE**

The input signal P1 connector should be left in position including the shipping links if no external connections are made. The forepump and the Turbo-V pump can be switched on at the same time.

---

## Startup

- Plug the controller power cable into a suitable power source.
- The display lights up for approx. 2 seconds and shows:

				A	U	T	O	T	E	S	T				
							O	K							

### NOTE

If the Remote Panel is not able to communicate with the controller after the displaying for 2 seconds

				A	U	T	O	T	E	S	T				

the following message is displayed:

	C	O	M	M	U	N	I	C	A	T	I	O	N		
					E	R	R	O	R						

This can be due to a missing or bad RS232 connection. From now on if the Remote Panel cannot any more communicate with the controller the following message is displayed:

*	C	O	M	M	U	N	I	C	A	T	I	O	N	*	
*					B	R	O	K	E	N				*	

## 17 Technical Information

### Use

If the remote Panel is connected to an unsupported controller the following message is displayed:

				A	U	T	O	T	E	S	T				
X	X	X	X	X	X	X		M	O	D	E	L	:		

Where:

**XXXXXXXX**: is the model number of the controller connected to the Remote Panel

#### NOTE

If the interlock on the P1 connector is not closed (pin 3 and 8 shorted) the following message is displayed:

				W	A	I	T	I	N	G					
			I	N	T	E	R	L	O	C	K				

Use the mating connector provided with the Remote Panel or make a short circuit on the pin 3 and 8 of the P1 connector.

#### NOTE

At startup the Remote Panel auto detects the baud rate of the controller in order to have a proper communication.

- After 2 seconds, the display shows the following screen page:

S	T	O	P						.	.	.	.	H	z
				F	R	O	N	T						

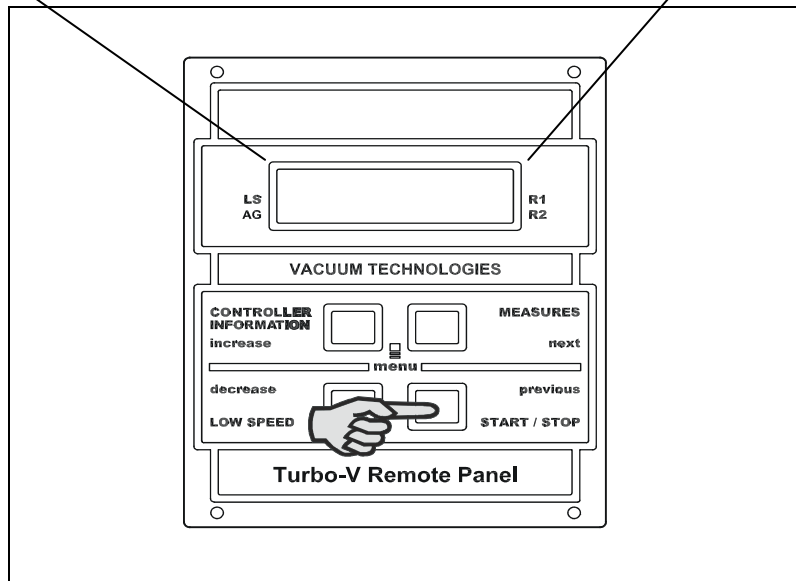
**NOTE**

If the pump is not connected, the display will be as follows:

C	H	E	C	K	C	O	N	N	E	C	T	I	O	N
				T	O	P	U	M	P					

- Pressing the START button, the pump is started and the display shows the following screen page (if the interlock is closed).

S	T	A	R	T	I	N	G			X	X	X	X	H	z
						F	R	O	N	T					



**Figure 24**

This screen page will be visible throughout the acceleration phase of the pump. The increasing value of the rotational frequency will be displayed in the boxes marked X; the user may decide to express this value in Hz (default) or in KRMP.

**17 Technical Information**  
Use

Once operating speed has been reached, the following screen page is displayed:

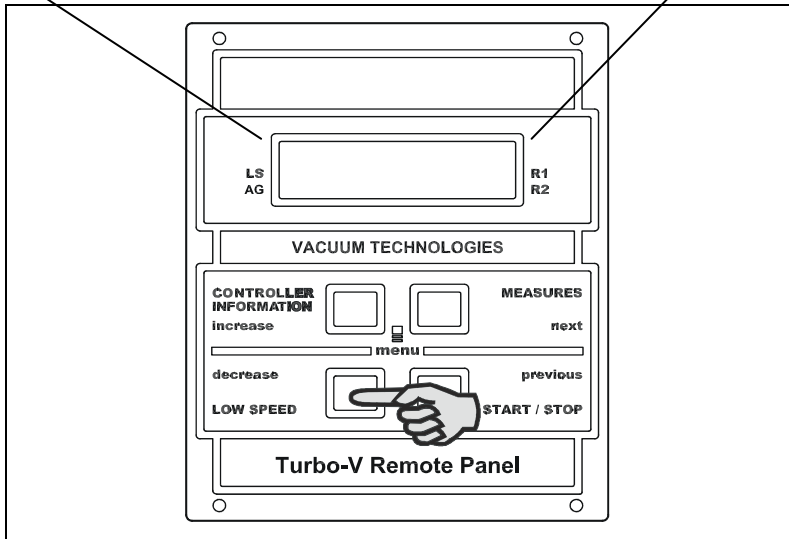
N	O	R	M	A	L					X	X	X	H	z
					F	R	O	N	T					

Where:

**XXXHz** is the current frequency of rotation of the pump.

If the LOW SPEED key is pressed and the pump connected is either a TV2k-G or TV3k-T, the above screen page changes to:

B	R	A	K	I	N	G				X	X	X	X	H	z
■					F	R	O	N	T						



**Figure 25**

The following screen page is displayed on reaching the pre-established speed:

N	O	R	M	A	L					X	X	X	H	z
■					F	R	O	N	T					

Where:

**XXXHz** is the current frequency of rotation of the pump.

The rotation value set in both NORMAL and LOW SPEED mode can be modified by the user as described in the Programming section of this manual.

## Programming

### Configuration Menu

To access the configuration menu, press the COUNTERS + MEASURES buttons at the same time for at least 2 sec.

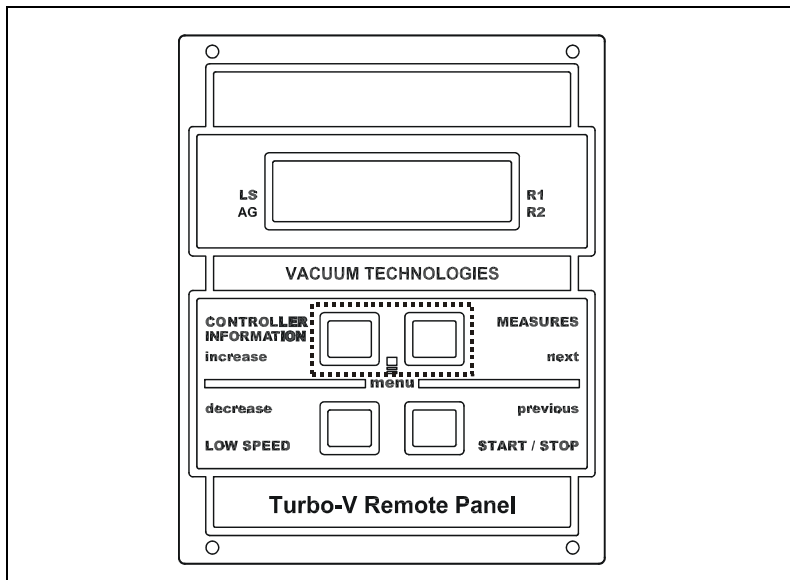


Figure 26

In programming mode, the meaning of the 4 buttons changes as follows:

- **COUNTERS/LOW SPEED** speed become “**INCREASE** and **DECREASE**” and make it possible to select the next or previous value of a parameter (in the case of compulsory parameters) or to increase/decrease the individual digit of a parameter (in the case of numeric parameters)
- The “**MEASURES** and **START/STOP**” buttons become “**NEXT** and **PREVIOUS**” and make it possible to select the next parameter (or next digit of a numeric parameter) or previous parameter.

Switching from one parameter to the next (using the “**NEXT**” key) indicates confirmation of any changes made to the specific parameter. If the changes cannot be accepted by the controller, the “**Data out of limits**” error message is displayed, the changes are cancelled and the value of the parameter prior to the modification is reposed.

Switching from a parameter to the previous parameter (using the “**PREVIOUS**” key) indicates the intention to abort any changes made to the specific parameter. Abort is indicated by the caption “**Changes aborted**” that is displayed for approx. 1 second before moving to the previous parameter.

The configuration environment has a tree-like structure and permits “circular” navigation; on reaching the end of a branch, the user is returned automatically to the start of this. Navigation can proceed in a forwards (**NEXT** button) or backwards (**PREVIOUS** button) direction

It is possible to enter a sub-branch by pressing the “**INCREASE**” button.

If no buttons are pressed for more than 1 minute, the controller automatically quits the programming environment canceling any modifications made to the current parameter.

## 17 Technical Information

### Programming

The figure below shows the first level of the configuration menu and the links to the respective second level menus (indicated below).

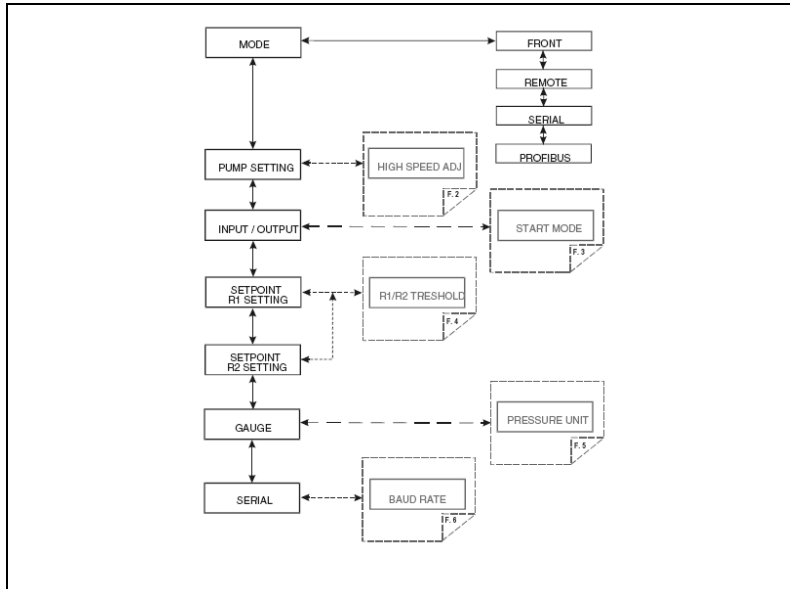


Figure 27 Flow – 1/6

## Mode Menu

- **MODE:** permits selection (see Flow 1/6) of the operating mode of the controller from those indicated below:
  - **FRONT:** accepts commands from the front panel
  - **REMOTE:** accepts commands from the remote input connector
  - **SERIAL:** accepts commands from the serial interface

## Pump Setting Menu for TV2k-G, TV3k-T

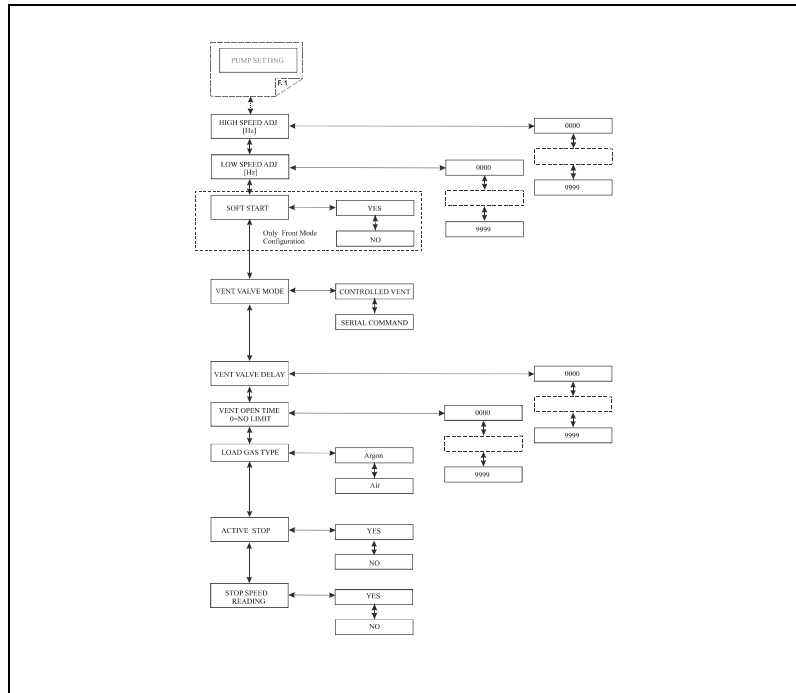


Figure 28 Flow - 2/6

This menu contains all the parameters that have a direct effect on driving of the pump.

- **HIGH SPEED:** Sets (in Hz) the speed of rotation of the pump when the low-speed function is not active.
- **LOW SPEED:** Sets (in Hz) the speed of rotation of the pump when the low-speed function is active.
- **SOFT START:** enables (YES) or disables (NO) the soft start function. This parameter is visible only if the controller is configured in FRONT mode.
- **VENT VALVE MODE:** makes it possible to select driving of the vent valve

## 17 Technical Information

### Programming

- **CONTROLLER VENT:** the valve is controlled by the controller
- **SERIAL COMMAND:** the valve is controlled by the serial port.
- **VENT VALVE DELAY:** Sets the delay, expressed in seconds, between stopping of the pump and opening of the vent valve.
- **VENT VALVE OPEN TIME:** sets the time, expressed in seconds, of opening of the vent valve. Enter 0 to be keep the valve opened indefinitely (until the next start).
- **LOAD GAS TYPE:** selects the type of gas used in the process. Therefore, the controller limits the power supplied to the pump according to the setting of this parameter.
- **ACTIVE STOP:** enables (YES) or disables (NO) active braking.
- **STOP SPEED READING:** enables (YES) or disables (NO) the pump rotational frequency reading function also during the slowing phase that follows a stop command

## Pump Setting Menu for TV551, TV701 and TV1001

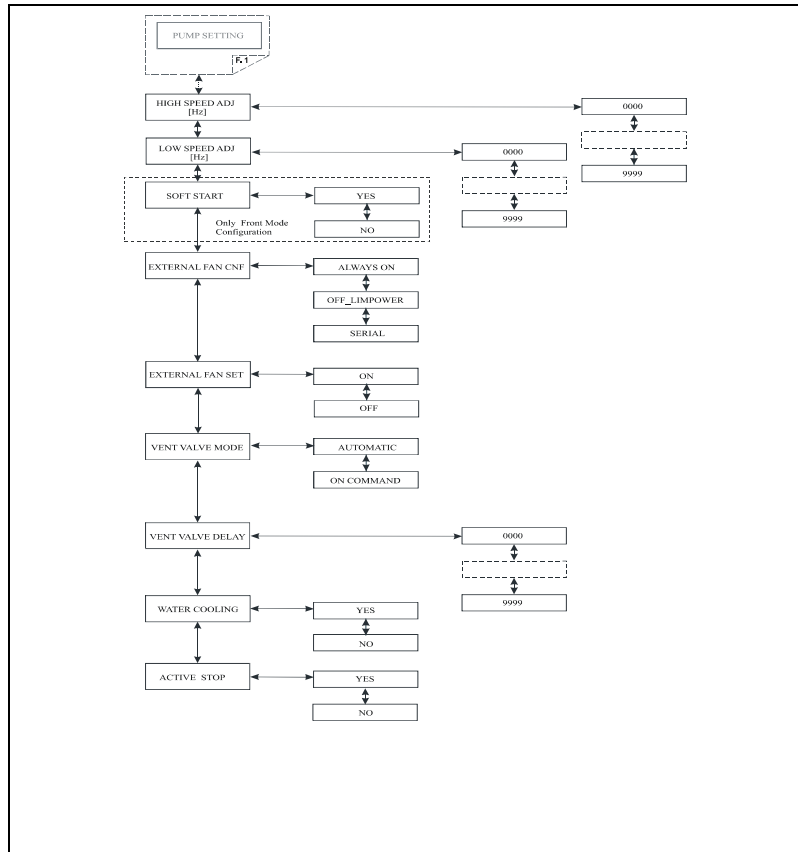


Figure 29 Flow – 2/6

## 17 Technical Information

### Programming

This menu contains all the parameters that have a direct effect on driving of the pump.

- **HIGH SPEED:** Sets (in Hz) the speed of rotation of the pump when the low-speed function is not active.
- **LOW SPEED:** Sets (in Hz) the speed of rotation of the pump when the low-speed function is active.
- **SOFT START:** enables (YES) or disables (NO) the soft start function. This parameter is visible only if the controller is configured in FRONT mode.
- **EXTERNAL FAN CNF:** Makes it possible to select driving mode of the external cooling fan. Possible options:
  - ALWAYS ON: the cooling fan is always ON
  - OFF\_LIMPOWER: 10 seconds after the pump has reached NORMAL status, the fan is switched off and the driving power of the pump is reduced to 50 W
  - SERIAL: the cooling fan is controlled directly via the serial port.
- **EXTERNAL FAN SET:** enables/disables the possibility to drive the external cooling fan. The function EXTERNAL FAN CFG is useless when EXTERNAL FAN SET is set to OFF.
- **VENT VALVE MODE:** makes it possible to select driving of the vent valve
  - AUTOMATIC: the valve is controlled by the controller
  - ON COMMAND: the valve is controlled by the serial port.
- **VENT VALVE DELAY:** Sets the delay, expressed in seconds, between stopping of the pump and opening of the vent valve.
- **WATER COOLING:** indicates whether the pump is cooled with water or not. The controller limits the power supplied to the pump according to this setting.
- **ACTIVE STOP:** enables (YES) or disables (NO) active braking.

## INPUT/OUTPUT Menu

This menu contains parameters that determine the operating mode of the inputs and outputs of the controller.

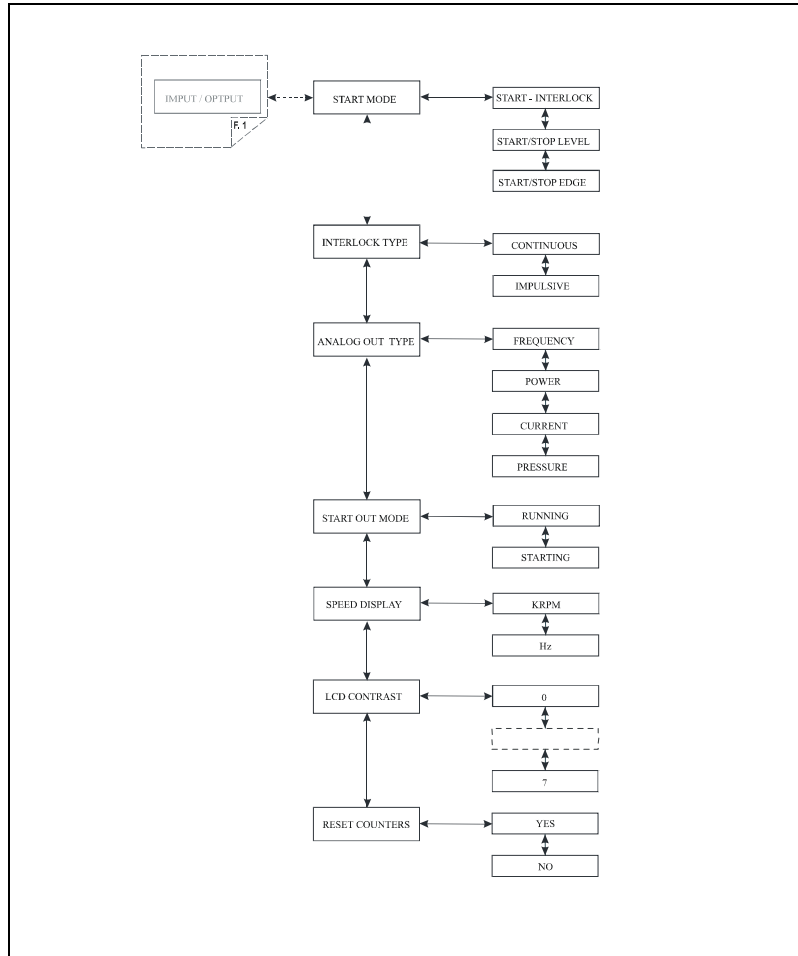


Figure 30 Flow – 3/6

- **START MODE:** Permits selection of one of the three types of functioning of the START/STOP and INTERLOCK inputs:
  - START/INTERLOCK: The pump is started if the START/STOP input is activated and is stopped if the START/STOP input is de-activated. The interlock input must be constantly active (INTERLOCK TYPE parameter = continuous) or may be active only at the time of starting (INTERLOCK TYPE parameter = impulsive)
  - START/STOP LEVEL The START/STOP input acts as only START command (active on the level) The INTERLOCK input acts as STOP command (active on the level) In the case of concurrent activation of the two inputs, the STOP command takes priority
  - START/STOP EDGE The START/STOP input acts as START command (active on the edge). The INTERLOCK input acts as STOP command (active on the edge).
- **INTERLOCK TYPE:** This parameter is displayed only when the previous parameter (START MODE) is equal to START/INTERLOCK and permits selection of one of the following two operating modes of the INTERLOCK input:
  - CONTINUOUS: the INTERLOCK input must be active continuously so that the pump can be driven by the controller. If, with the controller in start, the input INTERLOCK is de-activated, the controller switches to WAITING INTERLOCK status and interrupts driving of the pump
  - IMPULSIVE: it is sufficient for the input to be active when the START command is sent to the pump. Subsequently, the input can also be de-activated. If the interlock is not active when the pump is started, the controller switches too WAITING INTERLOCK status.

- **ANALOG OUT TYPE:** permits selection of the type of signal to be sent to the programmable analogue output. One of the following signals can be selected:
  - **FREQUENCY:** the drive frequency of the pump is shown
  - **POWER:** the power supplied to the pump is shown
  - **TEMPERATURE:** the temperature of the pump is shown, this selection is valid only for Tv2k-G and TV3k-T pumps
  - **PRESSURE:** the pressure read by the gauge is shown.
- **START OUT MODE:** Selects the operating mode of the START output. Possible options:
  - **RUNNING:** the output is active when the pump is driven
  - **STARTING:** the output is active only during acceleration ramp-up.
- **SPEED DISPLAY (KRPM/Hz):** permits selection of the unit of measurement of the speed of rotation shown on the display of the controller. This setting acts only on the measurement shown on the display. Possible options:
  - KRPM
  - Hz
- **LCD CONTRAST:** sets the LCD contrast.
- **RESET COUNTERS:** resets all the counters of the pump (cycles, time and pump life).

## Setpoint R1 and Setpoint R2 Submenus

This submenu contains the settings relating to the output with relay R1 (R2).

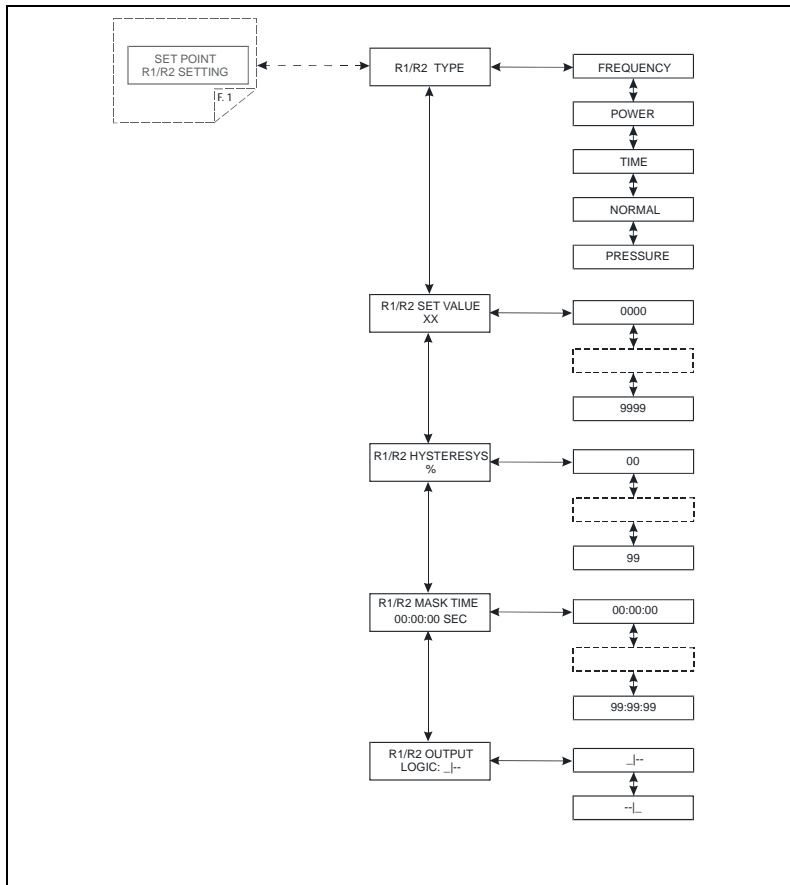


Figure 31 Flow – 4/6

- **R1 (R2) TYPE:** selects the type of reference measurement to drive the relay. Possible types:
    - Frequency: driving frequency of the pump
    - Power: power supplied to the pump
    - Time: time from start
    - Normal: achievement of normal condition
    - Pressure: pressure read by the gauge, if present
  - **R1 (R2) SET VALUE:** Sets the reference value at which the relay R1 (R2) is activated. This parameter is not available if SETPOINT TYPE is set to NORMAL.
  - **R1 (R2) HISTERESYS:** Sets, in percentage, the hysteresis of the relay. This parameter is not available if SETPOINT TYPE is set to TIME or NORMAL.
  - **R1 (R2) MASK TIME:** sets the masked time (in hh:mm:ss), measured from start, within which the SETPOINT R1 (R2) is not driven even if the threshold value has been reached. This parameter is not available if SETPOINT TYPE is set to TIME.
  - **R1 (R2) OUTPUT LOGIC:** Selects the control logic of the SETPOINT output.
    - \_|-- positive logic (the relay is energized when the measurement exceeds the reference)
    - |\_ negative logic (the relay is energized when the measurement is below the reference)
- This parameter is not available if SETPOINT TYPE is set to TIME or NORMAL.

## Gauge Menu

This menu contains the parameters relating to functioning of the active gauge.

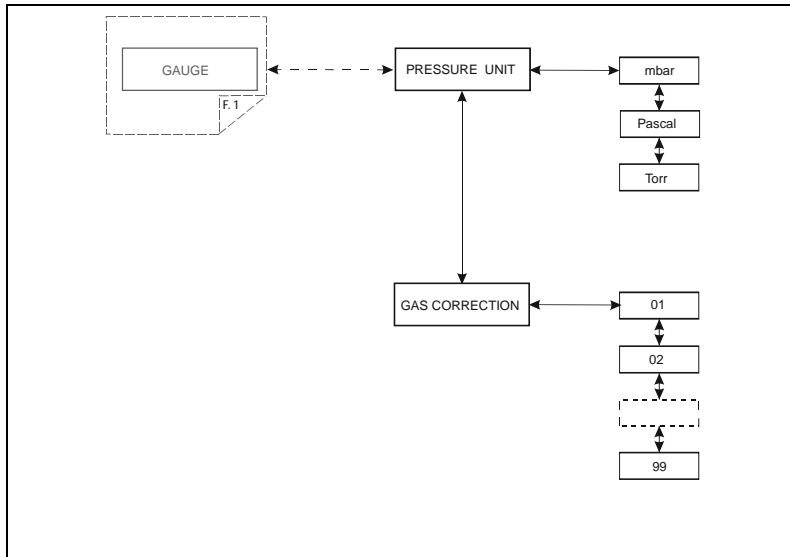


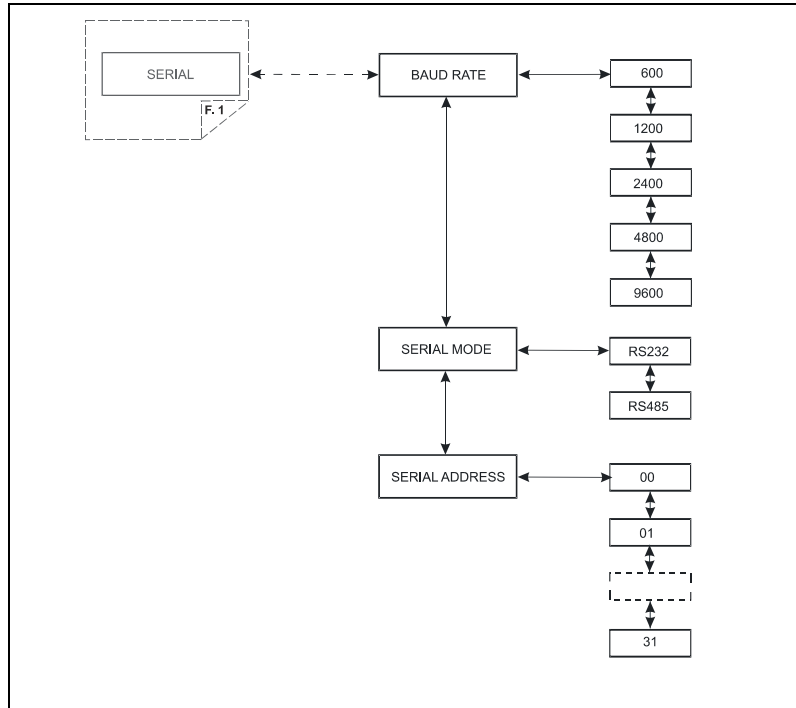
Figure 32 Flow – 5/6

- **PRESSURE UNIT:** selects the unit of measurement of the pressure used to show this on the display, for display on the serial line and for setting of the setpoint and autopressure thresholds. Possible options:
  - mbar
  - Pascal
  - Torr
- **GAS CORRECTION:** sets the conversion factor of the pressure measurement in the case in which the process gas is not nitrogen. The pressure indicated by the instrument is equal to:

$$P \text{ indicated} = P \text{ measured} \times \text{Gas correction} / 10$$

## Serial Menu

This menu contains the settings for management of the RS232/485 serial line.



**Figure 33** Flow - /6

- **BAUD RATE:** permits selection of the communication speed between the host (PC) and the remote panel; the possible values are: 600, 1200, 4800, 9600 baud
- **SERIAL MODE:** permits setting of the type of interface to RS232 (point to point) or RS485 (multidrop)
- **SERIAL ADDRESS:** sets the address of the node in the case of the RS485 interface. This parameter is not visible if the RS232 is selected.

## How to use by Remote I/O

1. Power ON the controller
2. Press the “COUNTERS” and “MEASURES” button together.
3. Select MODE = REMOTE from the “INPUT/OUTPUT” menu
4. Exit from configuration menu
5. Connection example:

## How to use in Serial Mode

1. Power on the controller
2. Connect the serial cable
3. Select by front panel if you want to work in 232 or 485, and if necessary, the serial address.
4. Select SERIAL from the INPUT/OUTPUT menu
5. Open the T+ software (see orderable parts table)
6. Ready to operate by SERIAL connections.

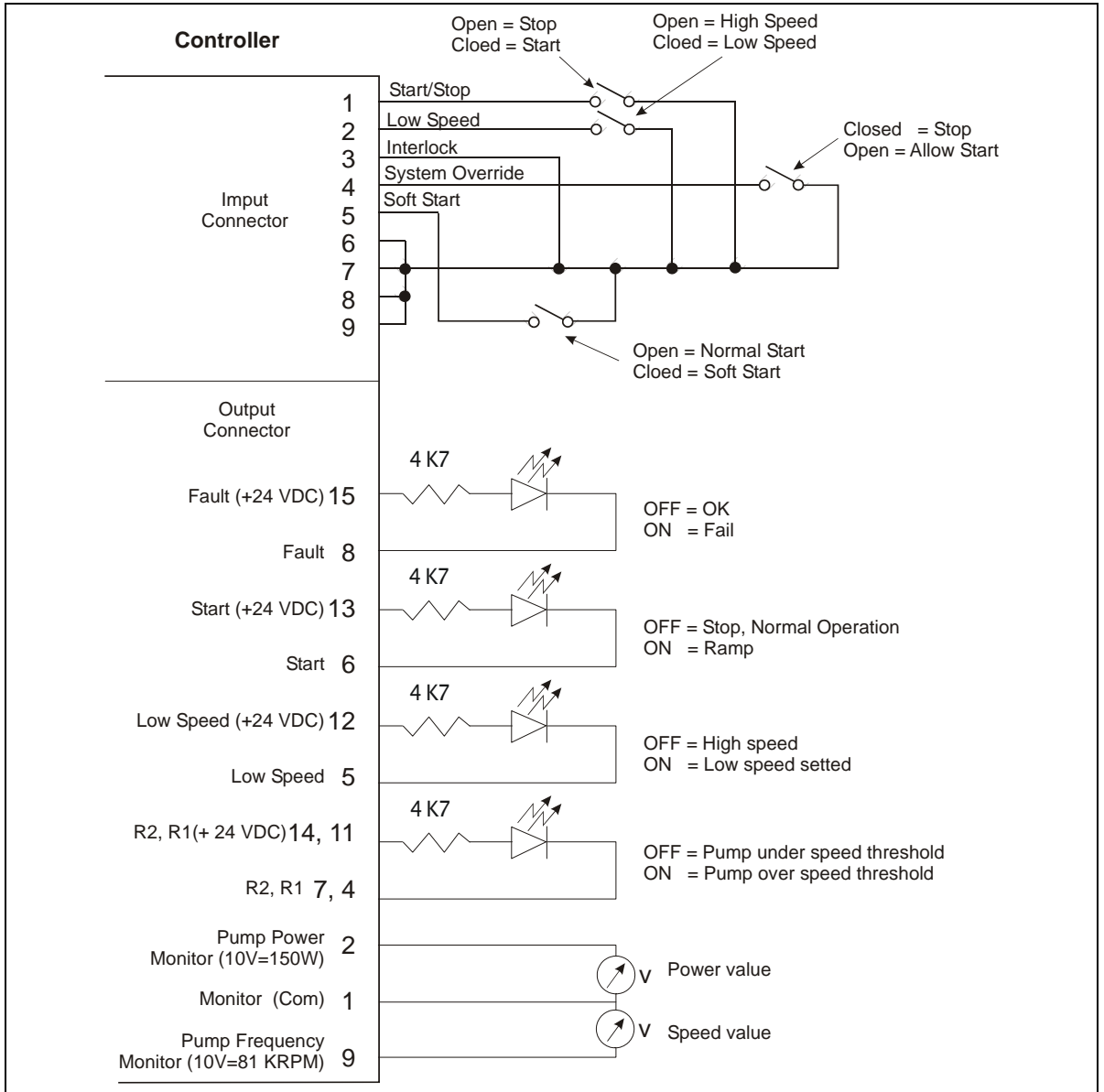


Figure 34

## Error Messages

For a certain type of failure, the controller will self-diagnose the error and the following messages will be displayed.

	C	O	M	M	U	N	I	C	A	T	I	O	N		
					E	R	R	O	R						

At startup there's a missing or faulty connection between the serial port of the Remote Panel and the serial port of the controller.

*	C	O	M	M	U	N	I	C	A	T	I	O	N	*
*					B	R	O	K	E	N				*

After a first successful communication the Remote Panel isn't any more able to communicate with the controller.

				A	U	T	O	T	E	S	T				
X	X	X	X	X	X	X		M	O	D	E	L	:		

The Remote Panel is connected to a not supported controller. Remote Panel can be connected only to the following pumps: TV551 with navigator controller, TV701 with navigator controller, TV1001 with navigator controller, TV2k-G with on-board controller and TV3k-T with on-board controller.

If the pump is not connected, the display will be as shown in the following figure.

C	H	E	C	K		C	O	N	N	E	C	T	I	O	N
			T	O		P	U	M	P						

Check connection between controller and pump, then press STOP push-button twice to start the pump.

If the P1 input connector is not in position with the link or the external interlock connections are open, when the START pushbutton is pressed the display will be as shown in the following figure.

					W	A	I	T	I	N	G			
					I	N	T	E	R	L	O	C	K	

Disconnect connector P1 and check the link or the external interlock, then install the connector to start the pump.

If within each step of the soft start mode the rotational speed of the pump do not reach the planned value within 15 minutes

**OR**

if in normal operation (after the starting phase) the current drawn by the pump is higher than programmed (1.8 A), the pump and the inter-connected devices are switched off and the display will be as shown in the following figure.

	T	O	O		H	I	G	H		L	O	A	D	

Verify that systems has no leaks than press STOP RESET pushbutton twice to start the pump.

Check that pump rotor is free to rotate then press the STOP RESET pushbutton twice to start the pump.

If the upper bearing/pump temperature exceeds 60 °C, the pump is shut off, and the display will be as shown in the following figure.

						P	U	M	P					
						O	V	E	R	T	E	M	P	

## 17 Technical Information

### Error Messages

The message will stay on until the temperature decreases below threshold value. Press the STOP RESET pushbutton twice to start the pump.

If the controller transformer temperature exceeds 90 °C, the pump is shut off, and the display will be as shown in the following figure.

			C	O	N	T	R	O	L	L	E	R		
			O	V	E	R	T	E	M	P				

The message will stay on until the temperature decreases below threshold value. Press the STOP RESET pushbutton twice to start the pump.

			P	O	W	E	R		F	A	I	L		

Indicates a fault in the circuit of the controller used to power the pump.

If the pump is stopped by an emergency stop signal provided via a remote contact, the display will be as shown in the following figure.

			S	Y	S	T	E	M						
			O	V	E	R	R	I	D	E				

Remove the controller power cable and check the emergency condition.

Then reconnect the power cable and press the START pushbutton to start the pump.

			O	V	E	R	V	O	L	T	G	E		

Controller power supply circuitry is faulty, or the Controller received a spike.

Press the START push-button twice to start the pump. Should the message still be present, call the Agilent service.

	S	H	O	R	T		C	I	R	C	U	I	T		

Check connections and shortages between pump and controller, then press the STOP RESET pushbutton twice to start the pump.

## Orderable Parts

DESCRIPTION	PART NUMBER
Turbo-V Remote Panel	969-8850
Remote Panel Connection Cable with valves remote inputs and IP-54 protection grade (Use with: TV2KG and TV3KT with on-board controller)	969-9945
Remote Panel Connection Cable (Use with: TV551, TV701 and TV1001 with navigator controller)	969-9946
EyeSys Mini-IMG Gauge cable (1 m)	969-9943
EyeSys Mini-IMG Gauge cable (special)	R1400301
EyeSys ConvecTor® Gauge cable (1 m)	969-9947
Gauge extention cable (5 m)	969-9944
Serial cable and Navigator Software	969-9883

**17 Technical Information**  
**Orderable Parts**



**Agilent Technologies**

***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,*

**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 Leinì (TO) – Italy

E-MAIL: [vpd-qualityassurance\\_pdl-ext@agilent.com](mailto: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.

- 1) Complete the attached Request for Return form and send it to Agilent Technologies (see below), taking particular care to identify all products that have pumped or been exposed to any toxic or hazardous materials.
- 2) After evaluating the information, Agilent Technologies will provide you with a Return Authorization (RA) number via email or fax, as requested.

**Note:** Depending on the type of return, a Purchase Order may be required at the time the Request for Return is submitted. We will quote any necessary services (evaluation, repair, special cleaning, eg).

**3) Important steps for the shipment of returning product:**

- Remove all accessories from the core product (e.g. inlet screens, vent valves).
- Prior to shipment, drain any oils or other liquids, purge or flush all gasses, and wipe off any excess residue.
- If ordering an Advance Exchange product, please use the packaging from the Advance Exchange to return the defective product.
- Seal the product in a plastic bag, and package product carefully to avoid damage in transit. You are responsible for loss or damage in transit.
- Agilent Technologies is not responsible for returning customer provided packaging or containers.
- **Clearly label package with RA number.** Using the shipping label provided will ensure the proper address and RA number are on the package. Packages shipped to Agilent without a RA clearly written on the outside cannot be accepted and will be returned.

- 4) Return only products for which the RA was issued.
- 5) **Product being returned under a RA must be received within 15 business days.**
- 6) **Ship to the location specified on the printable label, which will be sent, along with the RA number, as soon as we have received all of the required information.** Customer is responsible for freight charges on returning product.
- 7) Return shipments must comply with all applicable **Shipping Regulations** (IATA, DOT, etc.) and carrier requirements.

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

<b>EUROPE:</b>	<b>NORTH AMERICA:</b>	<b>PACIFIC RIM:</b>
Fax: 00 39 011 9979 330	Fax: 1 781 860 9252	please visit our website for individual
Fax Free: 00 800 345 345 00	Toll Free: 800 882 7426, Option 3	office information
Toll Free: 00 800 234 234 00	<a href="mailto:vpl-ra@agilent.com">vpl-ra@agilent.com</a>	<a href="http://www.agilent.com">http://www.agilent.com</a>
<a href="mailto:vpt-customer@agilent.com">vpt-customer@agilent.com</a>		



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

1) CUSTOMER INFORMATION

Form with fields: Company Name, Contact Name, Tel, Email, Fax, Customer Ship To, Customer Bill To, Europe only: VAT reg. Number, USA/Canada only: Taxable, Non-taxable

2) PRODUCT IDENTIFICATION

Table with 4 columns: Product Description, Agilent P/N, Agilent S/N, Original Purchasing Reference

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

- 3A. Non-Billable, Billable, New PO #
3B. Exchange, Repair, Upgrade, Consignment/Demo, Calibration, Evaluation, Return for Credit

4) HEALTH and SAFETY CERTIFICATION

AGILENT TECHNOLOGIES CANNOT ACCEPT ANY PRODUCTS CONTAMINATED WITH BIOLOGICAL OR EXPLOSIVE HAZARDS, RADIOACTIVE MATERIAL, OR MERCURY AT ITS FACILITY. Call Agilent Technologies to discuss alternatives if this requirement presents a problem. The equipment listed above (check one): HAS NOT pumped or been exposed to any toxic or hazardous materials. OR HAS pumped or been exposed to the following toxic or hazardous materials. If this box is checked, the following information must also be filled out. Check boxes for all materials to which product(s) pumped or was exposed: Toxic, Corrosive, Reactive, Flammable, Explosive, Biological, Radioactive. List all toxic/hazardous materials. Include product name, chemical name, and chemical symbol or formula: NOTE: If a product is received at Agilent which is contaminated with a toxic or hazardous material that was not disclosed, the customer will be held responsible for all costs incurred to ensure the safe handling of the product, and is liable for any harm or injury to Agilent employees as well as to any third party occurring as a result of exposure to toxic or hazardous materials present in the product. Print Name: Authorized Signature: Date:

5) FAILURE INFORMATION:

Failure Mode (REQUIRED FIELD. See next page for suggestions of failure terms): Detailed Description of Malfunction: (Please provide the error message) Application (system and model):

I understand and agree to the terms of Section 6, Page 3/3. Print Name: Authorized Signature: Date:



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

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

TURBO PUMPS and TURBO CONTROLLERS

Table with 3 columns: APPARENT DEFECT/MALFUNCTION, POSITION, and PARAMETERS. Lists various failure modes like 'Does not start', 'Noise', 'Vertical', 'Horizontal', etc.

ION PUMPS/CONTROLLERS

Table listing failure modes for Ion Pumps/Controllers: Bad feedthrough, Vacuum leak, Error code on display, Poor vacuum, High voltage problem, Other.

VALVES/COMPONENTS

Table listing failure modes for Valves/Components: Main seal leak, Solenoid failure, Damaged sealing area, Bellows leak, Damaged flange, Other.

LEAK DETECTORS

Table listing failure modes for Leak Detectors: Cannot calibrate, Vacuum system unstable, Failed to start, No zero/high background, Cannot reach test mode, Other.

INSTRUMENTS

Table listing failure modes for Instruments: Gauge tube not working, Communication failure, Error code on display, Display problem, Degas not working, Other.

SCROLL AND ROTARY VANE PUMPS

Table listing failure modes for Scroll and Rotary Vane Pumps: Pump doesn't start, Doesn't reach vacuum, Pump seized, Noisy pump (describe), Over temperature, Other.

DIFFUSION PUMPS

Table listing failure modes for Diffusion Pumps: Heater failure, Doesn't reach vacuum, Vacuum leak, Electrical problem, Cooling coil damage, Other.

Section 6) ADDITIONAL TERMS

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

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

## Sales and Service Offices

### United States

**Agilent Technologies  
Vacuum Products Division**  
121 Hartwell Avenue  
Lexington, MA 02421 - USA  
Tel.: +1 781 861 7200  
Fax: +1 781 860 5437  
Toll-Free: +1 800 882 7426

### Benelux

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

### Canada

**Central coordination through: Agilent Technologies  
Vacuum Products Division**  
121 Hartwell Avenue  
Lexington, MA 02421 - USA  
Tel.: +1 781 861 7200  
Fax: +1 781 860 5437  
Toll-Free: +1 800 882 7426

### China

**Agilent Technologies (China) Co. Ltd  
Vacuum Products Division**  
No.3, Wang Jing Bei Lu,  
Chao Yang District,  
Beijing, 100102  
China  
Tel.: +86 (10) 6439 7718  
Toll-Free: 800 820 6556

### France

**Agilent Technologies France  
Vacuum Products Division**  
7 Avenue des Tropiques  
Z.A. de Courtaboeuf - B.P. 12  
91941 Les Ulis cedex - France  
Tel.: +33 (0) 1 69 86 38 84  
Fax: +33 (0) 1 69 86 29 88  
Toll free: 00 800 234 234 00

### Germany and Austria

**Agilent Technologies  
Vacuum Products Division**  
Alsfelder Strasse 6 Postfach 11  
14 35  
64289 Darmstadt – Germany  
Tel.: +49 (0) 6151 703 353  
Fax: +49 (0) 6151 703 302  
Toll free: 00 800 234 234 00

### India

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

### Italy

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

### Japan

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

### Korea

**Agilent Technologies Korea, Ltd.  
Vacuum Products Division**  
Shinsa 2nd Bldg. 2F 966-5 Daechi-dong  
Kangnam-gu, Seoul  
Korea 135-280  
Tel.: +82 2 3452 2452  
Fax: +82 2 3452 2451  
Toll-Free: 080 222 2452

### Mexico

**Agilent Technologies  
Vacuum Products Division**  
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,  
Vacuum Products Division**  
Agilent Technologies Building,  
1 Yishun Avenue 7,  
Singapore 768923  
Tel : (65) 6215 8045  
Fax : (65) 6754 0574

### Southeast Asia

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

### Taiwan

**Agilent Technologies Taiwan Limited  
Vacuum Products Division (3F)**  
20 Kao-Shuang Rd.,  
Pin-Chen City, 324  
Taoyuan Hsien , Taiwan, R.O.C.  
Tel. +886 34959281  
Toll Free: 0800 051 342

### UK and Ireland

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

### Other Countries

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

### Customer Support & Service

#### **NORTH AMERICA:**

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

#### **EUROPE:**

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

#### **PACIFIC RIM:**

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

#### **Worldwide Web Site, Catalog and Order**

##### **On-line:**

[www.agilent.com](http://www.agilent.com)  
Representative in most countries  
12/10

© Agilent Technologies, Inc. 2011

Printed in ITALY

11/2011

Publication Number: 87-900-987-01 (E)



# Agilent Technologies