

# Agilent TwisTorr 305 FS onboard Controller

## User Manual



# Notices

## Manual Part Number

87-901-060-01  
Edition A.00, June 2022

## Copyright

© Agilent Technologies, Inc. 2022

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.

Agilent Technologies Italia S.p.A.  
Vacuum Products Division  
Via F.lli Varian, 54  
10040 Leini (TO)  
ITALY

[www.agilent.com](http://www.agilent.com)

## Instrument Manufacturing

Manufactured by Agilent Technologies Italia S.p.A., Vacuum Products Division  
Via F.lli Varian, 54, 10040 Leini (TO),  
ITALY

Printed in Italy

## Operating Temperature

Operating Temperature: +5°C ÷ +45°C  
Storage Temperature: -20°C ÷ +70°C

## 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

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## 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.

# Contents

<b>1 Istruzioni per l'uso</b>	<b>11</b>
Informazioni su questo manuale	12
Validità	12
Definizioni e terminologia	13
Definizione di Attenzione, Avvertenza e Nota	13
Simboli di avvertenza	14
Sicurezza	16
Uso corretto	16
Uso improprio	17
Attrezzatura di protezione	18
Linee guida di sicurezza per i Controller di pompe turbomolecolari	18
Trasporto & Immagazzinamento	19
Descrizione del prodotto	19
Installazione	20
Preparazione per l'installazione	20
Set-up	21
Fissaggio del TwisTorr 305 FS onboard Controller	22
Avvio e funzionamento del Controller	
TwisTorr 305 FS onboard controller	23
Arresto del TwisTorr 305 FS onboard	23
Arresto di Emergenza	23
Manutenzione	24
Smaltimento	25
Servizio Post Vendita	26

## Contents

<b>2 Betriebsanleitung</b>	<b>27</b>
Informationen über diese Betriebsanleitung	28
Gültigkeit	28
Definitionen und Terminologie	29
Definition von Vorsicht, Warnung und Hinweis	29
Warnsymbole	30
Sicherheit	32
Sachgemäße Verwendung	32
Unsachgemäße Anwendung	33
Schutzausrüstung	34
Sicherheitsrichtlinien für Turbomolekularpumpen-Steuergeräte	34
Transport und Lagerung	35
Produktbeschreibung	35
Installation	36
Vorbereitung zum Einbau	36
Set-up	37
Montage von TwisTorr 305 FS onboard Controller	38
Inbetriebnahme und Betrieb des Controller TwisTorr 305 FS onboard	39
Abschalten des TwisTorr 305 FS onboard	39
Notabschaltung	39
Wartung	40
Entsorgung	41
Service	42

## Contents

<b>3</b>	<b>Mode d'emploi</b>	<b>43</b>
	À propos de ce manuel	44
	Validité	44
	Définitions et terminologie	45
	Signification des messages d'attention, d'avertissement et des notes	45
	Symboles d'avertissement	46
	Sécurité	48
	Utilisation appropriée	48
	Utilisation inappropriée	49
	Équipements de protection individuelle	50
	Consignes de sécurité relatives aux contrôleurs de pompes turbomoléculaires	50
	Transport et Emmagasinage	51
	Description du produit	51
	Installation	52
	Préparation pour l'installation	52
	Set-up	53
	Montage du Contrôleur TwisTorr 305 FS onboard	54
	Démarrage et utilisation du Contrôleur TwisTorr 305 FS onboard	55
	Arrêt du TwisTorr 305 FS onboard	55
	Arrêt d'urgence	55
	Entretien	56
	Élimination	57
	Service	58

## Contents

<b>4</b>	<b>Instrucciones de uso</b>	<b>59</b>
	Acerca del manual	60
	Validez	60
	Definiciones y terminología	61
	Definición de Precaución, Advertencia y Nota	61
	Símbolos de advertencia	62
	Seguridad	64
	Uso correcto	64
	Uso incorrecto	65
	Equipo de protección	66
	Pautas de seguridad para Controller de bombas turbomoleculares	66
	Transporte y almacenamiento	67
	Descripción del producto	67
	Instalación	68
	Preparación para instalación	68
	Puesta en marcha	69
	Montaje del TwisTorr 305 FS onboard Controller	70
	Puesta en marcha y funcionamiento del TwisTorr 305 FS onboard Controller	71
	Parada TwisTorr 305 FS onboard	71
	Parada de Emergencia	71
	Mantenimiento	72
	Eliminación	73
	Servicio	74

## Contents

<b>5</b>	<b>使用说明</b>	<b>75</b>
	关于本手册	76
	有效性	76
	<b>定义与术语</b>	77
	小心、警告和注意的定义	77
	<b>警告符号</b>	78
	<b>安全</b>	80
	正确使用	80
	使用不当	81
	<b>防护装备</b>	82
	<b>涡轮分子泵控制器安全指南</b>	82
	<b>运输及存储</b>	83
	<b>般说明</b>	83
	<b>安装</b>	84
	<b>安装准备</b>	84
	<b>设置</b>	85
	TwisTorr 305 FS 遥控器安装	86
	TwisTorr 305 FS onboard Controller 的启动和运行	87
	TwisTorr 305 FS <b>板载控制器的关闭</b>	87
	<b>紧急停止</b>	87
	<b>维护</b>	88
	<b>处置</b>	89
	<b>服务</b>	90

## Contents

<b>6</b>	<b>取扱説明書</b>	<b>91</b>
	この説明書について	92
	有効期	92
	定義と用語	93
	警告・注意・注の定義	93
	警告の記号	94
	安全性	96
	適切な使用	96
	不適切な使用	97
	保護装具	98
	運輸及存储	99
	概要情報	99
	取り付け	100
	取り付けの準備	100
	取り付け	101
	TwisTorr 305 FS の取り付け	102
	TwisTorr 305 FS onboard Controller コンパクトの起動と操作	103
	TwisTorr 305 FS オンボードのスイッチを OFF にします	103
	緊急停止	103
	メンテナンス	104
	処分	105
	サービス	106

## Contents

<b>7</b>	<b>Instructions for Use</b>	<b>107</b>
	About this manual	108
	Validity	108
	Definitions and terminology	109
	Definition of Caution, Warning and Note	109
	Warning Symbols	110
	Safety	112
	Proper use	112
	Improper use	113
	Protective equipment	114
	Safety guideline for Turbomolecular Pump Controllers	114
	Transport & Storage	115
	Product description	115
	Installation	116
	Preparation for installation	116
	Set-up	117
	TwisTorr 305 FS onboard Controller Mounting	118
	Startup and Operation of the TwisTorr 305 FS onboard Controller	119
	Switching off the TwisTorr 305 FS onboard	119
	Emergency Stop	119
	Maintenance	120
	Disposal	121
	Service	122

## Contents

<b>8</b>	<b>Technical Information</b>	<b>123</b>
	Description of the TwisTorr 305 FS onboard Controller	124
	Technical specification	125
	Controller Outline	127
	TwisTorr 305 onboard Controller Installation	129
	Bottom Mounting	129
	Side Mounting	130
	Interconnections	131
	P3 – Vent	131
	P5 – External Fan	132
	J5 – In-Out	132
	Signal Description	134
	How to Connect the Open Collector Inputs of the Controller	137
	How to Connect the Outputs of the Controller	141
	J6 – Serial	141
	Input power connection on the model X3507-64130	143
	RS 232/RS 485 communication description	143
	Communication Format	143
	Communication Protocol	144
	Examples	146
	Window Meanings	148
	Air Cooling Kit Installation	152
	Use in the Presence of Magnetic Fields	155
	Accessories and Spare Parts	155

# 1

## Istruzioni per l'uso

Informazioni su questo manuale	12
Validità	12
Definizioni e terminologia	13
Definizione di Attenzione, Avvertenza e Nota	13
Simboli di avvertenza	14
Sicurezza	16
Uso corretto	16
Uso improprio	17
Attrezzatura di protezione	18
Linee guida di sicurezza per i Controller di pompe turbomolecolari	18
Trasporto & Immagazzinamento	19
Descrizione del prodotto	19
Installazione	20
Preparazione per l'installazione	20
Set-up	21
Fissaggio del TwisTorr 305 FS onboard Controller	22
Avvio e funzionamento del Controller	
TwisTorr 305 FS onboard controller	23
Arresto del TwisTorr 305 FS onboard	23
Arresto di Emergenza	23
Manutenzione	24
Smaltimento	25
Servizio Post Vendita	26

## Informazioni su questo manuale

### Validità

Questo manuale elenca le istruzioni per gli utenti del TwisTorr 305 FS onboard Controller, con particolare riferimento alle nozioni relative a sicurezza, funzionamento e manutenzione di primo livello, limitatamente alle operazioni di manutenzione di cui l'utente è responsabile.

Le operazioni di manutenzione, illustrate nelle sezioni specifiche, con disposizioni specifiche relative al livello più elevato di manutenzione (personale appositamente addestrato per le operazioni di manutenzione) non devono essere eseguite dall'utente.

Per una corretta installazione e avvio/arresto, consultare il paragrafo "Installazione". Per un uso tecnico più dettagliato, consultare il capitolo "Technical Information".

#### NOTA

- 1** Questo manuale contiene informazioni utili affinché tutto il personale che utilizza il TwisTorr 305 FS onboard Controller possa utilizzarla in sicurezza e garantire la perfetta efficienza, per tutta la sua durata.
  - 2** Conservare questo manuale, insieme a tutte le pubblicazioni ad esso correlate, in un luogo accessibile, conosciuto da tutti gli operatori/personale di manutenzione.
-

## Definizioni e terminologia

### Definizione di Attenzione, Avvertenza e Nota

Alcuni riferimenti importanti di questo manuale sono evidenziati e incorniciati in colori contrastanti.

#### 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 e forniscono maggiori dettagli su passaggi specifici.

### Simboli di avvertenza

Di seguito si riporta un elenco di simboli che vengono visualizzati insieme agli avvisi del TwisTorr 305 FS onboard Controller. Viene mostrato anche il pericolo che descrivono.

Un simbolo triangolare indica un'avvertenza. I significati dei simboli che possono apparire accanto alle avvertenze nella documentazione o sullo strumento stesso sono i seguenti:



Presenza di tensioni pericolose



Pericolo generico



Dichiarazione Europea di Conformità



Sito di produzione



Certificazione CSA



Waste Electrical and Electronic Equipment



Certificazione RoHS China



Marcatura UK CA

## Istruzioni per l'uso

Il seguente simbolo può essere usato sulle etichette di avvertenza attaccate allo strumento. Quando viene visualizzato questo simbolo, consultare il relativo manuale operativo o di manutenzione per la procedura corretta a cui fa riferimento tale etichetta di avvertenza.



I seguenti simboli appaiono sullo strumento per vostra informazione.

	Pericolo generico
	Presenza di tensioni pericolose
	Certificazione CE
	Certificazione CSA
	Certificazione RoHS China
	Waste Electrical and Electronic Equipment
	UK CA Marking

# Sicurezza

Questa sezione contiene le informazioni, prescritte dalla Direttiva bassa tensione 2014/35/UE, che sono essenziali per la conformità e il rispetto delle norme di sicurezza sia in generale che in relazione all'uso specifico del prodotto.

La mancata osservanza di queste istruzioni e delle altre istruzioni contenute nel presente manuale può rendere inefficaci le condizioni di sicurezza previste in fase di progettazione e causare incidenti a chi utilizza il prodotto.

Agilent Technologies declina ogni responsabilità per danni al prodotto o per la sicurezza fisica dell'operatore o di terzi, derivanti dal mancato rispetto delle norme di sicurezza indicate nella documentazione tecnica.

## Uso corretto

Questo manuale contiene importanti avvertenze e istruzioni di sicurezza da osservare affinché l'unità funzioni in sicurezza.

Il prodotto descritto in questo manuale è destinato esclusivamente all'ambito specificato nelle istruzioni. Il manuale fornisce inoltre indicazioni relative ai requisiti essenziali per l'applicazione e il funzionamento del prodotto, nonché le misure di sicurezza che possono essere adottate per garantire un funzionamento regolare. Agilent Technologies non fornisce alcuna garanzia né si assume alcuna responsabilità per applicazioni diverse da quelle descritte nel presente manuale o in cui i requisiti essenziali e le misure di sicurezza non siano rispettati. Il prodotto deve essere utilizzato solo da personale qualificato in grado di adottare le misure di sicurezza necessarie in condizioni che non causano danni o lesioni.

Gli accessori e le attrezzature utilizzati con il prodotto devono essere forniti o approvati da Agilent Technologies.

Qualsiasi operazione di regolazione o manutenzione deve essere eseguita da un tecnico professionista informato dei rischi.

Le riparazioni sul prodotto devono essere eseguite esclusivamente da personale autorizzato Agilent.

## Istruzioni per l'uso

### Uso improprio

Agilent Technologies declina ogni responsabilità derivante dall'uso improprio del TwisTorr 305 FS onboard Controller.

L'uso improprio comporterà la perdita di tutti i reclami per responsabilità e garanzie. L'uso improprio è definito come:

- installazione del dispositivo con materiale di montaggio non specificato
- funzionamento in ambiente condensante
- funzionamento in ambienti ad alta umidità fuori dal raggio specificato
- funzionamento in ambienti polverosi
- funzionamento con tensioni non specificate
- funzionamento del dispositivo in aree con radiazioni ionizzanti
- funzionamento in aree potenzialmente esplosive
- utilizzo del dispositivo in sistemi in cui stress e vibrazioni simili a impatti o forze periodiche influiscono sul dispositivo.

## Istruzioni per l'uso

### Attrezzatura di protezione

Non essendo prevista alcuna manutenzione per questo dispositivo, non sono necessari altri dispositivi di protezione.

#### **AVVERTENZA** Rischio di lesioni dovute alla caduta di oggetti



Durante il trasporto manuale dei controller, sussiste il pericolo che i carichi scivolino e cadano.

- Trasportare i controller con due mani.

### Linee guida di sicurezza per i Controller di pompe turbomolecolari

I controller della pompa turbomolecolare come descritto nel seguente manuale non devono essere aperti dall'utente per evitare il rischio di danneggiare i componenti interni.

#### **AVVERTENZA** Per evitare danni alle apparecchiature e per prevenire lesioni al personale operativo, è necessario seguire scrupolosamente le istruzioni di installazione fornite nel presente manuale!



# Trasporto & Immagazzinamento

Durante il trasporto e l'immagazzinamento del TwisTorr 305 FS onboard non devono essere superate le seguenti condizioni ambientali:

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

## Descrizione del prodotto

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 controller TwisTorr 305 FS onboard è un controller adatto ad essere utilizzato con le pompe TwisTorr 305. Ha dei connettori ausiliari tramite i quali è possibile alimentare un ventilatore aggiuntivo, comandare la valvola di vent, pilotarlo da remoto tramite un computer host collegato con linea seriale (RS232 o RS485). Nei paragrafi seguenti sono riportate tutte le informazioni necessarie a garantire la sicurezza dell'operatore durante l'utilizzo dell'apparecchiatura. Informazioni dettagliate sono fornite nel capitolo "Technical information".

## Installazione

### Preparazione per l'installazione

Il TwisTorr 305 FS onboard Controller 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 94/62 e successive modifiche.

## Set-up

**AVVERTENZA**



Il Controller (nella sua versione a tensione di rete, X3507-64131) deve essere alimentato mediante un cavo di alimentazione a tre fili (vedere tabella delle parti ordinabili) con una spina di tipo approvato a livello internazionale ai fini della sicurezza dell'utente. Utilizzare sempre questo cavo di alimentazione ed inserire la spina in una presa con un adeguato collegamento di terra onde evitare scariche elettriche e per rispettare le specifiche CE. All'interno del controller si sviluppano alte tensioni che possono recare gravi danni o la morte. Prima di eseguire qualsiasi operazione di installazione o manutenzione del controller scollegarlo dalla presa di alimentazione.

---

**AVVERTENZA**



Cavo di alimentazione: il cavo corretto per il cablaggio elettrico è un cavo a tre fili (L+N+Terra)

L = Linea

N = Neutro

La sezione del cavo deve essere almeno 1,0 mm<sup>2</sup> in caso di alimentazione 240 Vac o 2 mm<sup>2</sup> in caso di alimentazione 120 Vac per una lunghezza del cavo di riferimento pari a 3 m.

---

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 to +45 °C
- umidità relativa: 0 – 90 % (non condensante).

In ogni caso occorre che l'aria di raffreddamento possa circolare liberamente intorno all'apparato.

### ATTENZIONE

Il TwisTorr 305 FS onboard appartiene alla seconda categoria di installazione (o sovratensione) prevista dalla normativa EN 61010-1. Connettere quindi il dispositivo ad una linea di alimentazione che soddisfi tale categoria.

Il TwisTorr 305 FS onboard ha dei connettori per gli ingressi/uscite e per la comunicazione seriale che devono essere connessi ai circuiti esterni in modo che nessuna parte sotto tensione sia accessibile.

Assicurarsi che l'isolamento del dispositivo connesso al TwisTorr 305 FS onboard abbia un isolamento adeguato anche in condizione di guasto singolo come previsto dalla normativa EN 61010-1.

Per l'installazione degli accessori opzionali, vedere "Technical Information".

## Fissaggio del TwisTorr 305 FS onboard Controller

In questo paragrafo sono riportate le principali procedure operative. Prima di azionare il Controller, eseguire le connessioni elettriche e collegare il manicotto di pre vuoto facendo riferimento al capitolo "Technical Information".

### AVVERTENZA



**Per evitare danni alle persone ed all'apparato, prima di avviare la pompa, assicurarsi di avere correttamente effettuato il fissaggio della pompa al sistema da vuoto, secondo le istruzioni riportate del manuale della pompa (capitolo Installazione).**

### NOTA

Una volta collegato alla tensione di linea, per azionare il controller (e quindi per avviare la pompa) in modo semplice, collegare sul connettore J5 (Input-Output) il connettore di accoppiamento pre-cablato che viene fornito insieme al controller. Il connettore di accoppiamento pre-cablato è un modo semplice e veloce per consentire il comando del controller senza dover implementare altre connessioni esterne cablate come Serial Com o onboard I/O.

La pompa primaria e la pompa TwisTorr 305 FS possono essere attivate contemporaneamente.

# Avvio e funzionamento del Controller TwisTorr 305 FS onboard controller

Per accendere il TwisTorr 305 FS onboard è sufficiente fornire la tensione di alimentazione. Il controller riconosce automaticamente la presenza dei segnali di interlock e di START ed avvia la pompa.

Alla prima accensione, il controller avvia automaticamente la pompa in modo "soft start". Attendere il completamento dell'avvio in soft start (quindi attendere che la pompa raggiunga la massima velocità di rotazione, indicata da LED verde acceso fisso), prima di spegnere la pompa.

L'avvio della pompa in modalità soft start permette la corretta redistribuzione del lubrificante all'interno dei cuscinetti, necessaria dopo un lungo periodo di inattività della pompa).

Il periodo di fermo dopo il quale è necessario procedere ad un nuovo avvio in Soft Start è di 60 giorni e dopo il primo avvio (per il quale il controller esegue il soft start automaticamente), le successive attivazioni dell'avvio in Soft Start sono a carico dell'utente.

Per avere una partenza con "Soft Start" attivo occorre abilitare il modo suddetto tramite software o agendo sul connettore ingressi e uscite remote (onboard I/O connector).

Per la procedura di attivazione della modalità "Soft Start" fare riferimento al paragrafo "Signal Description" nel capitolo "Technical Information". Il LED verde posto sul pannello del TwisTorr 305 FS onboard indica, con la frequenza del suo lampeggio, le condizioni operative del sistema:

- acceso fisso: la pompa è in rotazione normale;
- lampeggiante lentamente (periodo di circa 400 ms): il sistema è in stato di rampa, o di Stop, o di "Waiting for interlock";
- lampeggiante velocemente (periodo di circa 200 ms): condizione di errore.

## Arresto del TwisTorr 305 FS onboard

Per arrestare il TwisTorr 305 FS onboard è sufficiente togliere la tensione di alimentazione.

## Arresto di Emergenza

Per arrestare in condizioni di emergenza il TwisTorr 305 FS onboard occorre staccare il cavo di alimentazione.

## Manutenzione

Il TwisTorr 305 FS onboard non richiede alcuna manutenzione. Qualsiasi intervento deve essere eseguito da personale autorizzato. Prima di effettuare qualsiasi intervento sul sistema scollegarlo dall'alimentazione. In caso di guasto, è possibile usufruire del servizio di riparazione Agilent o del servizio "Agilent advanced exchange service" che permette di ottenere un controller rigenerato in sostituzione di quello guasto.

**AVVERTENZA**

**Prima di effettuare qualsiasi intervento sul controller scollegare il cavo di alimentazione.**



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.



Figura 1 Logo "WEEE"

Per maggiori informazioni consultare:

<http://www.agilent.com/environment/product/index.shtml>

## Servizio Post Vendita

Nel caso in cui un cliente necessiti di un servizio di sostituzione o riparazione avanzato, si prega di contattare il distributore locale o spedire direttamente a:

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

È necessario completare il modulo "Request for Return" per restituire la pompa ad Agilent per l'assistenza (fornito alla fine di questo manuale).

## 2

# Betriebsanleitung

Informationen über diese Betriebsanleitung	28
Gültigkeit	28
Definitionen und Terminologie	29
Definition von Vorsicht, Warnung und Hinweis	29
Warnsymbole	30
Sicherheit	32
Sachgemäße Verwendung	32
Unsachgemäße Anwendung	33
Schutzausrüstung	34
Sicherheitsrichtlinien für Turbomolekularpumpen-Steuergeräte	34
Transport und Lagerung	35
Produktbeschreibung	35
Installation	36
Vorbereitung zum Einbau	36
Set-up	37
Montage von TwisTorr 305 FS onboard Controller	38
Inbetriebnahme und Betrieb des Controller TwisTorr 305 FS onboard	39
Abschalten des TwisTorr 305 FS onboard	39
Notabschaltung	39
Wartung	40
Entsorgung	41
Service	42

# Informationen über diese Betriebsanleitung

## Gültigkeit

Diese Betriebsanleitung enthält die Anweisungen für die Benutzer der TwisTorr 305 FS onboard Controller mit besonderer Bezugnahme auf die Begriffe für Sicherheit, Betrieb und Wartung auf der ersten Ebene, auf die Wartungsarbeiten beschränkt, für die der Benutzer verantwortlich ist.

Die in den spezifischen Abschnitten dargestellten Wartungsarbeiten mit spezifischen Verordnungen hinsichtlich der höheren Wartungsebene (entsprechend geschultes Personal für die Wartungsarbeiten) dürfen nicht durch den Benutzer ausgeführt werden.

Für eine korrekte Installation und Start/Stop der Absatz „Installation“ konsultieren. Für eine detailliertere technische Verwendung das Kapitel „Technical Information“ konsultieren.

### HINWEIS

- 1** Diese Betriebsanleitung enthält nützliche Informationen, damit das gesamte Personal, das die TwisTorr 305 FS onboard Controller verwendet, diese in Sicherheit verwenden und die perfekte Funktionstüchtigkeit für ihre gesamte Betriebsdauer garantieren kann.
  - 2** Diese Betriebsanleitung muss zusammen mit allen mit dieser zusammenhängenden Veröffentlichungen an einem zugänglichen Ort, der allen Bedienern/Wartungspersonal bekannt ist, aufbewahrt werden.
-

## Definitionen und Terminologie

### Definition von Vorsicht, Warnung und Hinweis

Einige wichtige Bezüge dieser Betriebsanleitung werden durch Kontrastfarben hervorgehoben und eingerahmt.

#### VORSICHT

---

Die Vorsichtshinweise werden vor Verfahren angezeigt, die Schäden am Gerät verursachen könnten, wenn sie nicht beachtet werden.

#### WARNUNG



Die Warnhinweise lenken die Aufmerksamkeit des Bedieners auf ein Verfahren oder einen spezifischen Vorgang, der schwere Verletzungen von Personen zur Folge haben könnte, wenn er nicht korrekt ausgeführt wird.

---

#### HINWEIS

---

Die Hinweise enthalten wichtige Informationen und liefern weitere Details über spezifische Arbeitsschritte.

# Warnsymbole

Das Folgende ist eine Liste von Symbolen, die in Verbindung mit den Warnungen auf dem TwisTorr 305 FS onboard Controller angezeigt werden. Die von ihnen beschriebene Gefahr wird ebenfalls angezeigt.

Ein dreieckiges Symbol weist auf eine Warnung hin. Die Bedeutung der Symbole, die neben Warnhinweisen in der Dokumentation oder auf dem Gerät selbst erscheinen können, ist wie folgt:



Allgemeine Gefahr



Allgemeine Gefahr



Europäische  
Konformitätserklärung



Herstellungsbetrieb



CSA certification



Elektro- und  
Elektronikaltgeräte



RoHS-  
Chinazertifizierung



UK CA-Kennzeichnung

## Betriebsanleitung

Das folgende Symbol kann auf am Gerät angebrachten Warnschildern verwendet werden. Wenn Sie dieses Symbol sehen, finden Sie in der jeweiligen Betriebs- oder Wartungsanleitung das richtige Verfahren, auf das sich dieses Warnschild bezieht.



Die folgenden Symbole werden zu Ihrer Information auf dem Instrument angezeigt.

	Allgemeine Gefahr
	Allgemeine Gefahr
	CE-Zertifizierung
	CSA-Zertifizierung
	RoHS-Chinazertifizierung
	Elektro- und Elektronikaltgeräte
	UK CA-Kennzeichnung

# Sicherheit

Dieser Abschnitt enthält die in der Maschinenrichtlinie 2014/35/UE vorgeschriebenen Informationen, die für die Einhaltung der Sicherheitsbestimmungen im Allgemeinen und in Bezug auf den spezifischen Einsatz der Maschine von wesentlicher Bedeutung sind.

Die Nichtbeachtung dieser Anweisungen und der anderen in diesem Handbuch enthaltenen Anweisungen kann die in der Konstruktionsphase vorgesehenen Sicherheitsbedingungen ineffizient machen und Unfälle bei den Bedienern der Maschine verursachen.

Agilent Technologies lehnt jede Verantwortung für Schäden an der Maschine oder für die physische Sicherheit des Bedieners oder Dritter ab, die sich aus der Nichtbeachtung der in den technischen Unterlagen angegebenen Sicherheitsregeln ergeben.

## Sachgemäße Verwendung

Dieses Handbuch enthält wichtige Warnungen und Sicherheitshinweise, die beachtet werden müssen, damit das Gerät sicher funktioniert.

Das in dieser Anleitung beschriebene Produkt ist ausschließlich für den in der Anleitung angegebenen Anwendungsbereich bestimmt. Das Handbuch enthält auch Angaben zu den grundlegenden Anforderungen für die Anwendung und den Betrieb des Produkts sowie zu den Sicherheitsmaßnahmen, die zur Gewährleistung eines regelmäßigen Betriebs ergriffen werden können. Agilent Technologies übernimmt keine Garantie oder Verantwortung für Anwendungen, die nicht in diesem Handbuch beschrieben sind oder bei denen die grundlegenden Anforderungen und Sicherheitsmaßnahmen nicht eingehalten werden.

Das Produkt darf nur von qualifiziertem Personal verwendet werden, das in der Lage ist, die erforderlichen Sicherheitsmaßnahmen unter Bedingungen zu treffen, die keine Schäden oder Verletzungen verursachen. Alle mit dem Produkt verwendeten Zubehörteile und Geräte müssen von Agilent Technologies geliefert oder genehmigt werden.

Einstellungs- oder Wartungsarbeiten müssen von einem Fachmann durchgeführt werden, der über die Risiken informiert ist.

Reparaturen am Produkt dürfen ausschließlich von Agilent-autorisiertem Personal durchgeführt werden.

### Unsachgemäße Anwendung

Agilent Technologies lehnt jede Verantwortung ab, die sich aus der unsachgemäßen Benutzung vom TwisTorr 305 FS onboard Controller ergibt.

Bei unsachgemäßer Verwendung verfallen alle Haftungs- und Gewährleistungsansprüche. Unsachgemäße Verwendung ist definiert als:

- Montage des Geräts unter Verwendung von nicht vorgesehenem Befestigungsmaterial
- Betrieb in Kondenswasserumgebung.
- Betrieb in einer Umgebung mit hoher Luftfeuchtigkeit über dem angegebenen Wertebereich hinaus
- Betrieb in staubiger Umgebung
- Betrieb mit anderen als den festgelegten Netzspannungen
- Betrieb des Geräts in Bereichen mit ionisierender Strahlung
- Betrieb in explosionsgefährdeten Bereichen
- Einsatz des Geräts in Anlagen, in denen stoßartige Einwirkungen und Vibrationen oder regelmäßig auftretende Kräfte auf das Gerät einwirken.

## Betriebsanleitung

### Schutzausrüstung

Da für dieses Gerät keine Wartung vorgesehen ist, ist keine weitere Schutzeinrichtung erforderlich.

#### **WARNUNG**

#### **Verletzungsgefahr durch herunterfallende Gegenstände**



Beim Tragen von Vakuum-Controllern von Hand besteht die Gefahr, dass Lasten verrutschen und herunterfallen.



- Controller mit zwei Händen transportieren.



### Sicherheitsrichtlinien für Turbomolekularpumpen-Steuergeräte

Zur Vermeidung des Risikos einer Beschädigung interner Komponenten dürfen die in dieser Bedienungsanleitung beschriebenen Turbomolekularpumpen-Steuergeräte nicht vom Benutzer geöffnet werden.

#### **WARNUNG**

Um Geräteschäden und Verletzungen des Bedienpersonals zu vermeiden, sind die in diesem Handbuch enthaltenen Installationsanweisungen unbedingt zu beachten!



# Transport und Lagerung

Beim Betrieb müssen folgende Umgebungsbedingungen eingehalten werden:

- Temperatur: von -20 °C bis +70 °C
- Relative Luftfeuchtigkeit: 0 ÷ 95% (nicht kondensierend).

# Produktbeschreibung

Diese Vorrichtung ist für eine professionelle Anwendung bestimmt. Der Benutzer muss, vor Anwendung, vorliegendes Handbuch und alle weitere von Agilent gelieferte Angaben, aufmerksam durchlesen. Agilent ist für etwaige auf teilweise oder gesamte Nichtberücksichtigung der Gebrauchsanweisungen beruhende Verantwortungen, für eine nicht geeignete Anwendung durch nicht ausgebildetes Personal, für nicht autorisierte Eingriffe oder für Anwendung unter Nichtbeachtung der nationalen Bestimmungen, nicht verantwortlich.

Der Controller TwisTorr 305 FS onboard ist für eine Anwendung mit Pumpen des Typs TwisTorr 305, geeignet. Er besitzt Hilfsverbinder, durch die ein zusätzlicher Ventilator gespeist und das Lüftungsventil betrieben werden können, eine Fernsteuerung ist durch einen, mit einer seriellen Linie verbundenen Host-Rechner, möglich (RS232 oder RS485). In den folgenden Absätzen sind alle notwendigen Informationen über die Sicherheit des Bedienungspersonals, während des Betriebs, angegeben.

Ausführliche Angaben sind im Anhang "Technical Information", enthalten.

## Installation

### Vorbereitung zum Einbau

Der TwisTorr 305 FS onboard Controller wird mit einer speziellen Schutzverpackung geliefert. Eventuelle Transportschäden müssen der zuständigen örtlichen Verkaufsstelle gemeldet werden.

Beim Auspacken vorsichtig vorgehen, damit der Controller nicht fällt oder Stößen ausgesetzt wird.

Das Verpackungsmaterial muß korrekt entsorgt werden. Es ist vollständig recyclebar und entspricht der EG-Richtlinie 94/62 für Umweltschutz.

## Set-up

**WARNUNG**



Der Controller (in seiner Netzspannungsversion, X3507-64131) muss mit einem dreidradigen Netzkabel (siehe Bestelltabelle) mit international zur Sicherheit des Benutzers zugelassenem Stecker angeschlossen werden. Es sollte immer dieses Netzkabel benutzt werden, das an eine korrekt geerdete Steckdose anzuschließen ist, um den CE Richtlinien zu entsprechen und Stromschläge zu vermeiden. Im Inneren des Controllers entstehen hohe Spannungen, die schwere Verletzungen verursachen und lebensgefährlich sein können. Vor jedem Montage- bzw. Wartungseingriff muss deshalb der Netzstecker gezogen werden.

---

**WARNUNG**



Netzkabel: Das geeignete Kabel für die Stromleitung ist ein dreidradiges (L+N+Erde) Kabel.

L = Linie

N = Neutral

Der Kabelquerschnitt muss mindestens  $1,0 \text{ mm}^2$  bei einer Netzspannung von 240 Vac oder  $2 \text{ mm}^2$  bei einer Netzspannung von 120 Vac bei einer Referenzkabellänge von 3 m betragen.

---

Beim Betrieb müssen folgende Umgebungsbedingungen eingehalten werden:

- Temperatur: von +5 °C bis +45 °C
- Relative Luftfeuchtigkeit: 0 – 90 % (nicht kondensierend).

In jedem Fall ist es erforderlich, dass die Kühlluft kann frei um das Gerät zirkulieren.

### VORSICHT

TwisTorr 305 FS onboard wird in die zweite Installationskategorie (Überspannung) der Norm EN 61010-1, eingestuft. Die Vorrichtung muss an eine Netzlinie angeschlossen werden, die dieser Kategorie entspricht.

TwisTorr 305 FS onboard verfügt außer den, für die serielle Kommunikation vorgesehenen Konnektoren, auch Ausgangs- und Eingangskonnektoren, die mit den äußeren Schaltungen verbunden werden müssen, sodass kein Teil unter Spannung zugänglich sein kann.

Sicherstellen, dass die Isolierung der mit dem TwisTorr 305 FS onboard verbundenen Vorrichtung, auch im Einzelschadenfall, so wie von Norm EN 61010-1 vorgesehen, eine angebrachte Isolierung besitzt.

Für die Installation der Options-Nebeneinrichtungen, beachten Sie bitte den Abschnitt "Technical Information"

## Montage von TwisTorr 305 FS onboard Controller

In diesem Absatz werden die wichtigsten Funktionsverfahren angegeben. Bevor Sie den Controller aktivieren, schließen Sie die Elektroanschlüsse an und schließen Sie die freie Muffe an, wie im Benutzerhandbuch beschrieben.

### WARNUNG



**Steht die Pumpe auf einem Tisch, muß auf den stabilen Stand geachtet werden, da sonst die Gefahr von Personen- und Geräteschäden besteht. Die Pumpe nie einschalten, wenn der Eingangsflansch nicht am System angeschlossen bzw. nicht mit dem Schließflansch abgedeckt ist.**

### HINWEIS

Nach Anschluss an die Netzspannung können Sie den Regler leicht bedienen (und dann um die Pumpe zu starten) indem Sie den mit dem Regler mitgelieferten vorverdrahteten Gegenstecker an den J5-Stecker (Eingang-Ausgang) anschließen.

Der vorverdrahtete Gegenstecker ist eine einfache und schnelle Möglichkeit, den Regler zu bedienen ohne dass dazu andere verdrahtete externe Verbindungen wie Serial Com oder onboard I/O benötigt werden. Die Vorpumpe und die TwisTorr 305 FS Pumpe können gleichzeitig eingeschaltet werden.

# Inbetriebnahme und Betrieb des Controller TwisTorr 305 FS onboard

Die elektrische Versorgung schaltet den TwisTorr 305 FS onboard automatisch an. Der Controller erkennt automatisch die Interlock- und Anschaltsignale und schaltet die Pumpe an.

Beim ersten Einschalten startet der Regler die Pumpe automatisch im "Softstart"-Modus. Warten Sie, bis der Softstart abgeschlossen ist (dann warten Sie, bis die Pumpe die maximale Drehzahl erreicht hat, die durch eine ständige grüne LED angezeigt wird), bevor Sie die Pumpe ausschalten.

Das Starten der Pumpe im Sanftanlauf-Modus ermöglicht die korrekte Umverteilung des Schmiermittels in den Lagern, die nach einer langen Zeit der Pumpeninaktivität erforderlich ist).

Die Stillstandszeit, nach der ein neuer Softstart erforderlich ist, beträgt 60 Tage, und nach dem ersten Start (für den die Steuerung den Softstart automatisch durchführt) ist der Benutzer verantwortlich für die nachfolgenden Softstart-Aktivierungen.

Um einen Start mit aktivem „Soft Start“ zu erzielen, ist es notwendig, diesen Modus durch die Software oder durch den Anschluss der Remote-Ein- und Ausgänge zu aktivieren) (Remote I/O connector).

Um den Modus "Soft Start" einzustellen, bitte den Absatz "Signal Description", im Kapitel "Technical Information", nachschlagen. Die grüne lichtemittierende Diode auf der Schalttafel des TwisTorr 305 FS onboard, gibt mit ihrer Leuchtfrequenz den System-Funktionszustand an:

- eingeschaltet, nicht intermittierend, die Pumpe hat eine normale Drehung;
- langsames Aufleuchten (ca. 400 ms), das System befindet sich im Ausgangs- oder Stop-Stadium oder im Stadium "Waiting for interlock";
- schnelles Aufleuchten (ca. 200 ms), gibt einen Fehler an.

## Abschalten des TwisTorr 305 FS onboard

Um den TwisTorr 305 FS onboard abzuschalten, genügt es die elektrische Versorgung abzuschalten.

## Notabschaltung

Um den TwisTorr 305 FS onboard im Notfall abzuschalten, die elektrische Verbindung unterbrechen.

### Wartung

Für den TwisTorr 305 FS onboard ist keine Wartung erforderlich. Jeder Eingriff unterliegt autorisiertem Personal. Vor jedem Eingriff, die elektrische Versorgung unterbrechen. Bei einem Störfall kann der Reparaturservice Agilent oder der Service "Agilent Advanced Exchange service" in Anspruch genommen werden, durch den man einen regenerierten Controller als Ersatz für den defekten erhält.

**WARNUNG** Trennen Sie den Controller vor jedem beliebigen Eingriff vom Netz.



Im Falle einer Verschrottung des Systems, muss diese nach den nationalen Vorschriften erfolgen.

## Entsorgung

### Bedeutung des „WEEE“ Logos auf den Schildern.

Das folgende Symbol ist in Übereinstimmung mit der EG-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 Haushalts- oder Industriegemüll 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.



Abbildung 2 Logo „WEEE“

Für weitere Informationen siehe:

<http://www.agilent.com/environment/product/index.shtml>

## Service

Sollte ein Kunde einen erweiterten Austausch- oder Reparaturservice benötigen, wenden Sie sich bitte an den örtlichen Händler oder direkt an die Email-Adresse

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

Das Ausfüllen des "Request for Return" formulars ist erforderlich, um die Pumpe zur Wartung an Agilent zurückzusenden (am Ende dieses Handbuchs angegeben).

# 3

## Mode d'emploi

À propos de ce manuel	44
Validité	44
Définitions et terminologie	45
Signification des messages d'attention, d'avertissement et des notes	45
Symboles d'avertissement	46
Sécurité	48
Utilisation appropriée	48
Utilisation inappropriée	49
Équipements de protection individuelle	50
Consignes de sécurité relatives aux contrôleurs de pompes turbomoléculaires	50
Transport et Emmagasinage	51
Description du produit	51
Installation	52
Préparation pour l'installation	52
Set-up	53
Montage du Contrôleur TwisTorr 305 FS onboard	54
Démarrage et utilisation du Contrôleur TwisTorr 305 FS onboard	55
Arrêt du TwisTorr 305 FS onboard	55
Arrêt d'urgence	55
Entretien	56
Élimination	57
Service	58

## À propos de ce manuel

### Validité

Le présent manuel contient les instructions destinées aux utilisateurs du TwisTorr 305 FS onboard Controller, notamment celles relatives à la sécurité, au fonctionnement et à la maintenance de premier niveau, dans la limite des activités de maintenance incombant à l'utilisateur.

Les opérations de maintenance décrites dans des paragraphes spécifiques qui contiennent des dispositions relevant d'un niveau de maintenance plus élevé (personnel spécialement formé aux opérations de maintenance) ne doivent pas être effectuées par l'utilisateur.

Pour une installation et une mise en marche/arrêt correctes, consultez le paragraphe « Installation ». Pour une utilisation technologique plus détaillée, consultez le chapitre « Technical Information ».

#### NOTE

- 1 Ce manuel contient des informations utiles pour que l'ensemble du personnel puisse utiliser le TwisTorr 305 FS onboard Controller en toute sécurité et pour garantir un fonctionnement parfait pendant toute sa durée de vie.
  - 2 Veuillez conserver ce manuel et tous les documents connexes dans un lieu accessible connu de tous les opérateurs et du personnel de maintenance.
-

## Définitions et terminologie

### Signification des messages d'attention, d'avertissement et des notes

Dans ce manuel, certaines informations importantes sont surlignées et encadrées avec des couleurs voyantes.

#### ATTENTION

Les messages d'attention sont affichés au début de procédures qui, si elles ne sont pas respectées, peuvent entraîner des dommages aux équipements.

---

#### AVERTISSEMENT



Les messages d'avertissement attirent l'attention de l'opérateur sur une procédure ou une pratique particulière qui, si elle n'est pas effectuée correctement, peut entraîner de graves blessures.

---

#### NOTE

Les remarques contiennent des informations importantes et fournissent des précisions sur certains passages particuliers.

---

### Symboles d'avertissement

Voici une liste de symboles qui apparaissent en conjonction avec les avertissements sur le TwisTorr 305 FS onboard Controller. Le danger qu'ils décrivent est également illustré.

Le symbole triangulaire indique un avertissement. Les significations des symboles qui peuvent apparaître à côté des avertissements dans la documentation ou sur l'appareil lui-même sont les suivantes.



Tensions dangereuses



Danger générique



Déclaration européenne de conformité



Site de fabrication



Certification CSA



Déchets d'équipements électriques et électroniques



Certification RoHS Chine










Marquage UK CA

## Mode d'emploi

Le symbole suivant peut être utilisé sur les étiquettes d'avertissement apposées sur l'appareil. Lorsque vous voyez ce symbole, reportez-vous au manuel d'utilisation ou d'entretien correspondant pour connaître la procédure correcte visée par cette étiquette d'avertissement.



Les symboles suivants apparaissent sur l'appareil pour votre information.

	Danger générique
	Tensions dangereuse
	Certification CE
	Certification CSA
	Certification RoHS Chine
	Déchets d'équipements électriques et électroniques
	Marquage UK CA

# Sécurité

Cette section contient les informations prescrites par la directive Machines 2014/35/UE, qui est essentielle à la conformité et au respect des règles de sécurité en général ainsi qu'à l'utilisation spécifique du produit.

Le non-respect de ces instructions et des autres instructions contenues dans le présent manuel peut rendre inefficaces les conditions de sécurité prévues dans la phase de conception et causer des accidents à ceux qui utilisent le produit.

Agilent Technologies décline toute responsabilité pour les dommages causés au produit ou pour la sécurité physique de l'opérateur ou des tiers résultant du non-respect des règles de sécurité indiquées dans la documentation technique.

## Utilisation appropriée

Le présent manuel contient des avertissements importants et des instructions de sécurité à respecter pour que l'appareil puisse fonctionner en toute sécurité.

Le produit décrit dans le présent manuel est destiné exclusivement au domaine d'application spécifié dans les instructions. Le manuel fournit également des indications concernant les exigences essentielles pour l'application et le fonctionnement du produit ainsi que les mesures de sécurité qui peuvent être adoptées pour garantir un fonctionnement régulier. Agilent Technologies ne fournit aucune garantie ou n'assume aucune responsabilité pour des applications autres que celles décrites dans le présent manuel ou dans lesquelles les exigences essentielles et les mesures de sécurité ne sont pas respectées.

Le produit ne doit être utilisé que par un personnel qualifié capable de prendre les mesures de sécurité nécessaires dans des conditions qui ne causent pas de dommages ou de blessures. Tous les accessoires et équipements utilisés avec le produit doivent être fournis ou approuvés par Agilent Technologies.

Toute opération de réglage ou de maintenance doit être effectuée par un technicien professionnel informé des risques.

Les réparations du produit doivent être effectuées exclusivement par le personnel autorisé d'Agilent.

## Mode d'emploi

### Utilisation inappropriée

Agilent Technologies décline toute responsabilité en raison de l'utilisation inappropriée du TwisTorr 305 FS onboard Controller.

Une utilisation inappropriée entraînera l'annulation de toutes les créances et garanties.

Une utilisation inappropriée est définie comme suit :

- installation de l'appareil à l'aide de matériel de montage inadapté
- utilisation dans un environnement exposé à de la condensation
- utilisation dans un environnement très humide en dehors de la plage spécifiée
- utilisation dans un environnement poussiéreux
- utilisation avec des tensions électriques non-conformes aux spécifications
- utilisation de l'appareil dans des zones exposées à des rayonnements ionisants
- utilisation dans des zones potentiellement explosives
- utilisation de l'appareil dans des systèmes soumis à des contraintes et des vibrations ou des forces périodiques qui affectent l'appareil.

## Mode d'emploi

### Équipements de protection individuelle

Comme aucune intervention d'entretien n'est prévue sur cet appareil, tout autre équipement de protection n'est pas nécessaire.

#### **AVERTISSEMENT** Risques de blessure dus à la chute d'objets



Pendant le transport manuel des contrôleur à vide, il existe un risque de glissement et de chute de la charge.

- Déplacer les contrôleurs en utilisant les deux mains.

#### Consignes de sécurité relatives aux contrôleurs de pompes turbomoléculaires

Les contrôleurs de pompes turbomoléculaires décrits dans ce manuel ne doivent pas être ouverts par l'utilisateur afin d'éviter les risques d'endommager les composants internes.

#### **AVERTISSEMENT** Pour éviter d'endommager les équipements et prévenir les risques de blessure des opérateurs, respectez rigoureusement les instructions d'installation fournies dans le présent manuel.



# Transport et Emmagasinage

Pendant le transport et le stockage du TwisTorr 305 FS onboard, veiller au respect des conditions environnementales suivantes :

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

# Description du produit

Cet appareillage a été conçu en vue d'une utilisation professionnelle. Avant toute utilisation de l'appareil, il est conseillé à l'utilisateur de lire attentivement cette notice d'instructions ainsi que toute autre indication supplémentaire fournie par Agilent qui décline par conséquent toute responsabilité en cas de non respect total ou partiel des instructions données, d'utilisation impropre par un personnel non formé, d'opérations non autorisées ou d'emploi contraire aux réglementations nationales spécifiques. Le TwisTorr 305 FS onboard est un contrôleur approprié pour être utilisé avec les pompes

TwisTorr 305. Le TwisTorr 305 FS onboard est en outre doté de connecteurs auxiliaires permettant d'alimenter un ventilateur supplémentaire, de commander la vanne de ventilation, de le piloter à distance à l'aide d'un ordinateur host connecté par ligne série. Les paragraphes suivants fournissent toutes les indications nécessaires à garantir la sécurité de l'opérateur pendant l'utilisation de l'appareillage. Des renseignements plus détaillés se trouvent dans le chapitre "Technical Information".

## Installation

### Préparation pour l'installation

Le contrôleur TwisTorr 305 FS onboard 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 les matériaux d'emballage sans autorisation. Le matériel est recyclable à 100 % et est conforme à la directive 94/62 de la CEE et aux modifications ultérieures.

## Set-up

**AVERTISSEMENT**



Le Contrôleur (dans sa version avec tension secteur, X3507-64131) doit être alimenté au moyen d'un câble d'alimentation à trois fils (voir tableau des pièces que l'on peut commander) avec une fiche du type approuvé au niveau international en vue de la sécurité de l'utilisateur. Utiliser toujours ce câble d'alimentation et introduire la fiche dans une prise pourvue d'un branchement au sol approprié à la masse, afin d'éviter toute décharge électrique et de respecter les spécifications CE. A l'intérieur du contrôleur se développent de hautes tensions qui peuvent causer de graves dommages, voire la mort. Avant d'effectuer toute opération d'installation ou d'entretien du contrôleur, le débrancher de la prise d'alimentation.

---

**AVERTISSEMENT**



Cordon d'alimentation électrique : le câble correct du raccordement électrique est un câble à trois fils (L+N+Terre).

L = Ligne  
N = Neutre

La section des fils doit être au moins de 1,0 mm<sup>2</sup> en présence d'une tension d'alimentation de 240 Vca ou de 2 mm<sup>2</sup> si la tension d'alimentation est de 120 Vca, en présence d'un câble de référence d'une longueur égale à 3 m.

---

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

- température: de +5 °C à +45 °C
- humidité relative: de 0 % à 90 % (non condensante).

Il est en tout cas nécessaire que l'air de refroidissement puisse circuler librement à l'intérieur de l'appareil.

### ATTENTION

Le TwisTorr 305 FS onboard appartient à la deuxième catégorie d'installations (ou surtension) prévue par la norme EN 61010-1. De ce fait, brancher le dispositif à une ligne d'alimentation compatible avec cette catégorie. Le TwisTorr 305 FS onboard dispose de connecteurs pour les entrées/sorties et pour la communication en série qui doivent être branchés aux circuits extérieurs de façon qu'aucune partie sous tension ne soit accessible.

S'assurer que l'isolation du dispositif branché au TwisTorr 305 FS onboard a une isolation appropriée même en condition de panne individuelle selon les termes de la norme EN 61010-1.

Pour l'installation d'accessoires en option, voir la section "Technical Information".

## Montage du Contrôleur TwisTorr 305 FS onboard

Ce paragraphe présente les principales procédures opérationnelles. Avant d'actionner le Contrôleur, procéder aux branchements électriques et brancher le manchon de prévide en se référant au Manuel d'instructions.

### AVERTISSEMENT



Pour éviter tous dommages aux personnes et à l'appareil, avant de démarrer la pompe, assurez-vous d'avoir correctement fixé la pompe au système de vide, conformément aux instructions fournies dans le manuel de la pompe (chapitre Installation).

### NOTE

Une fois connecté à la tension de ligne, pour faire fonctionner le contrôleur (et donc démarrer la pompe) de manière très simple, connectez sur le connecteur J5 (Entrée-Sortie) le connecteur homologue précâblé qui est fourni avec le contrôleur.

Le connecteur homologue précâblé est un moyen simple et rapide de permettre au contrôleur d'être utilisé sans avoir besoin de mettre en œuvre d'autres connexions externes câblées comme Serial Com ou onboard I/O.

La pompe à pré-vide et la pompe TwisTorr 305 FS peuvent être mises en marche simultanément.

# Démarrage et utilisation du Contrôleur TwisTorr 305 FS onboard

Pour allumer le TwisTorr 305 FS onboard il suffit de fournir la tension d'alimentation. Le contrôleur reconnaît automatiquement la présence de signaux d'interlock et de démarrage et il actionne la pompe.

Lors de la première mise en marche, le contrôleur démarre automatiquement la pompe en modalité « Soft Start ». Attendez que le démarrage progressif soit terminé (puis attendez que la pompe atteigne la vitesse de rotation maximale, indiquée par une DEL verte fixe), avant d'arrêter la pompe.

Le démarrage de la pompe en modalité « Soft Start » permet une redistribution correcte du lubrifiant à l'intérieur des paliers, ce qui est nécessaire après une longue période d'inactivité de la pompe).

La période d'arrêt après laquelle un nouveau démarrage Soft start est nécessaire est de 60 jours et après le premier démarrage (pour lequel le contrôleur effectue automatiquement le démarrage soft start), les activations ultérieures du démarrage soft start sont à la charge de l'utilisateur.

Pour avoir un démarrage avec « Soft Start » actif, il est nécessaire d'activer cette modalité par logiciel ou en agissant sur le connecteur des entrées et sorties à distance. (Remote I/O connector).

Pour la procédure d'activation de la modalité « Soft Start », se référer au paragraphe « SIGNAL DESCRIPTION" dans le chapitre "Technical Information)". La LED verte LD1 placée sur le panneau de la base du TwisTorr 305 FS onboard indique, par sa fréquence de clignotement, les conditions opérationnelles du système:

- allumée fixe: la pompe est en rotation normale;
- clignote lentement (période d'environ 400 ms): le système est en état de rampe, ou de stop, ou d'attente interlock (waiting for interlock).
- clignote rapidement (période d'environ 200 ms): condition d'erreur.

## Arrêt du TwisTorr 305 FS onboard

Pour arrêter le TwisTorr 305 FS onboard, il suffit de retirer la tension d'alimentation. Le contrôleur arrête immédiatement la pompe.

## Arrêt d'urgence

Pour arrêter le TwisTorr 305 FS onboard en conditions d'urgence, il faut débrancher le cordon d'alimentation.

## Entretien

Le TwisTorr 305 FS onboard n'exige aucun entretien particulier. Toute intervention doit être effectuée par un personnel agréé. En cas de panne, on peut s'adresser au service de réparation Agilent ou au service "Agilent advanced exchange service" qui permet d'obtenir un contrôleur régénéré en substitution de celui endommagé.

**AVERTISSEMENT**

**Avant d'effectuer toute opération sur le contrôleur, débrancher le câble d'alimentation.**



En cas de mise au rebut de la pompe, procéder à son élimination conformément aux réglementations nationales concernant la gestion des déchets.

## Élimination

### Signification du logo « DEEE » présent sur les étiquettes.

Le symbole représenté ci-dessous est apposé conformément à la directive CE dite « DEEE ».

Ce symbole (valable uniquement pour les pays de la Communauté européenne) indique que le produit sur lequel il est apposé NE DOIT PAS être éliminé avec des déchets ménagers ou industriels communs mais qu'il doit être confié à un centre de collecte sélective. L'utilisateur est donc invité à contacter le fournisseur du produit, qu'il s'agisse du fabricant ou d'un revendeur, pour donner lieu au processus de collecte et d'élimination, après avoir vérifié les conditions générales de vente.



Figure 3 Logo « DEEE »

Pour plus de précisions, veuillez consulter :

<http://www.agilent.com/environment/product/index.shtml>

## Mode d'emploi

### Service

Si, vous, le client avez besoin d'un service d'échange ou de réparation avancé, veuillez contacter le distributeur local ou directement par courrier à:

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

Vous devez remplir le formulaire de "Request for Return" pour retourner votre pompe à Agilent pour l'entretien (fourni à la fin du présent manuel).

## 4 Instrucciones de uso

Acerca del manual	60
Validez	60
Definiciones y terminología	61
Definición de Precaución, Advertencia y Nota	61
Símbolos de advertencia	62
Seguridad	64
Uso correcto	64
Uso incorrecto	65
Equipo de protección	66
Pautas de seguridad para Controler de bombas turbomoleculares	66
Transporte y almacenamiento	67
Descripción del producto	67
Instalación	68
Preparación para instalación	68
Puesta en marcha	69
Montaje del TwisTorr 305 FS onboard Controller	70
Puesta en marcha y funcionamiento del TwisTorr 305 FS onboard Controller	71
Parada TwisTorr 305 FS onboard	71
Parada de Emergencia	71
Mantenimiento	72
Eliminación	73
Servicio	74

## Acerca del manual

### Validez

Este manual enumera las instrucciones para los usuarios de la bomba TwisTorr 305 FS onboard Controller, con particular referencia a la información relacionada con la seguridad, la operación y el mantenimiento de primer nivel que está limitado a las operaciones de mantenimiento que son responsabilidad del usuario.

Las operaciones de mantenimiento, ilustradas en las secciones particulares, con disposiciones específicas relacionadas con el nivel más elevado de mantenimiento (personal específicamente capacitado para las operaciones de mantenimiento) no deben ser realizadas por el usuario.

Para una instalación y arranque/parada correctos, consulte el párrafo "Instalación".

Para una instalación y un encendido/apagado correctos, consulte la sección "Technical Information".

#### NOTA

- 1 Este manual contiene información útil para que todo el personal que utilice la bomba TwisTorr 305 FS onboard Controller pueda operarla de manera segura y garantizar un rendimiento perfecto, durante toda su vida útil.
- 2 Guarde este manual, junto con todas las publicaciones relacionadas, en un lugar conocido y accesible para todos los operadores/personal de mantenimiento.

## Definiciones y terminología

### Definición de Precaución, Advertencia y Nota

Algunas referencias importantes de este manual están resaltadas y enmarcadas en color contrastante.

#### PRECAUCIÓN

Los mensajes de precaución se muestran antes de los procedimientos que, si no se respetan, pueden causar daños al equipo.

---

#### ADVERTENCIA



Los mensajes de advertencia llaman la atención del operador sobre un procedimiento o práctica específica que, si no se realiza correctamente, puede causar lesiones personales graves.

---

#### NOTA

Las notas están previstas para llamar la atención sobre información importante y proporcionar más detalles en relación con pasos específicos.

---

## Símbolos de advertencia

La siguiente es una lista de los símbolos que aparecen en conjunto con las advertencias en la bomba TwisTorr 305 FS onboard Controller. También se muestra el peligro que describen.

Un símbolo triangular indica una advertencia. Los significados de los símbolos que pueden aparecer junto a las advertencias en la documentación son los siguientes:



Voltajes peligrosos



Peligro genérico



Declaración Europea de Conformidad



Lugar de fabricación



Certificación CSA



Equipo de Desecho Eléctrico y Electrónico



Certificación China RoHS



Marca UK CA

## Instrucciones de uso

El siguiente símbolo podría utilizarle en las etiquetas de advertencia fijadas al instrumento. Cuando vea este símbolo, consulte la operación relevante o el manual de servicio para el procedimiento referido por dicha etiqueta de advertencia.



Los siguientes símbolos aparecen en el instrumento para su información.

	Peligro genérico
	Voltajes peligrosos
	Certificación CE
	Certificación CSA
	Certificación China RoHS
	Equipo de Desecho Eléctrico y Electrónico
	Marca UK CA

# Seguridad

La presente sección contiene la información, prescrita por la Directiva de Maquinario 2014/35/UE, que es esencial para el cumplimiento y seguimiento de las normativas de seguridad, tanto generales como en relación con el uso específico del producto.

El incumplimiento de las presentes instrucciones y las otras instrucciones en el presente manual pueden hacer que las condiciones de seguridad previstas en la fase de diseño sean ineficaces y causen accidentes a los que operen el producto.

Agilent Technologies niega cualquier responsabilidad por daños al producto o la seguridad física del operario o terceras partes derivados del incumplimiento de las reglas de seguridad indicadas en la documentación técnica.

## Uso correcto

El presente manual contiene advertencias importantes e instrucciones de seguridad a cumplir para que la unidad funcione de manera segura.

El producto descrito en el presente manual está destinado exclusivamente al área de aplicación especificado en las instrucciones. El manual también proporciona indicaciones acerca de los requisitos esenciales para la aplicación y funcionamiento del producto, así como las medidas de seguridad que pueden ser adoptadas para garantizar el funcionamiento regular. Agilent Technologies no proporciona ninguna garantía ni asume ninguna responsabilidad por aplicaciones que no sean las descritas en el presente manual, o en las que no se respeten los requisitos y medidas de seguridad esenciales.

El producto solo puede ser utilizado por personal cualificado que pueda tomar las medidas de seguridad necesarias bajo condiciones que no causen daños o lesiones. Cualquier accesorio y equipo utilizado con el producto debe ser proporcionado o aprobado por Agilent Technologies.

Cualquier ajuste u operación de mantenimiento debe ser realizada por un técnico profesional informado sobre los riesgos.

Las reparaciones del producto deberán ser realizadas exclusivamente por personal autorizado por Agilent.

## Instrucciones de uso

### Uso incorrecto

Agilent Technologies niega cualquier responsabilidad derivada del uso incorrecto del TwisTorr 305 FS onboard Controller.

El uso incorrecto anulará todas las reclamaciones por responsabilidad y garantías.

El uso incorrecto se define como:

- instalación del dispositivo con material de montaje no especificado
- funcionamiento en un entorno de condensación
- funcionamiento en entornos de alta humedad fuera del rango especificado
- funcionamiento en entornos polvorientos
- funcionamiento con voltajes no especificados
- funcionamiento del dispositivo en áreas con radiación ionizante
- funcionamiento en zonas potencialmente explosivas
- uso del dispositivo en sistemas en los que la fuerza de impacto y las vibraciones o fuerzas periódicas afecten al dispositivo.

## Instrucciones de uso

### Equipo de protección

Como no se prevé mantenimiento para este dispositivo, no es necesario ningún otro equipo de protección.

#### **ADVERTENCIA** Riesgo de lesiones debido a la caída de objetos



Cuando se transportan los Controler de vacío manualmente, existe el peligro de que las cargas se resbalen y caigan.



- Traslade los Controler con las dos manos.



### Pautas de seguridad para Controler de bombas turbomoleculares

Los Controler de bombas turbomoleculares descritos en el siguiente manual de funcionamiento no deben ser abiertos por el usuario para evitar el riesgo de dañar los componentes internos.

#### **ADVERTENCIA** Para evitar daños al equipo y prevenir lesiones al personal operativo, se deben cumplir estrictamente las instrucciones de instalación indicadas en este manual.



# Transporte y almacenamiento

Durante el transporte y el almacenamiento del TwisTorr 305 FS onboard no pueden ser superadas las siguientes condiciones ambientales:

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

# Descripción del producto

El presente equipo está destinado a uso profesional. Antes de utilizar este sistema se aconseja al usuario que lea atentamente el manual de instrucciones así como cualquier otro tipo de información adicional que Agilent facilite. Agilent se retiene libre de eventuales responsabilidades debido a la inobservancia total o parcial de las instrucciones, a la utilización incorrecta por parte de personal no adiestrado, a intervenciones no autorizadas o a un uso contrario con las normas nacionales específicas.

El controler TwisTorr 305 FS onboard es un controler apto para ser utilizado con las bombas TwisTorr 305. Está dotado de conectores adicionales mediante los que es posible alimentar un ventilador auxiliar, dirigir la válvula de vent y pilotarlo desde control remoto mediante un ordenador host conectado con una línea serial (RS232 ó RS485). En los siguientes párrafos se han incluido todas las informaciones necesarias para garantizar la seguridad del operador durante la utilización del equipo.

Para más información consultar el capítulo "Technical information".

## Instalación

### Preparación para instalación

El TwisTorr 305 FS onboard se suministra con un embalaje especial de protección; si el embalaje muestra signos de deterioro que podrían haberse verificado durante el transporte, será necesario contactar con el ente de ventas local.

Al desempacar, tenga especial cuidado de no dejar caer el controlador ni someterlo a golpes y no deseche el embalaje en el medio ambiente.

El material es 100% reciclable y cumple con la Directiva EEC 94/62 y modificaciones posteriores.

## Puesta en marcha

**ADVERTENCIA**



El Controler (en su versión con tensión de red, X3507-64131) debe ser alimentado mediante un cable de tres hilos (ver tabla de los componentes que pueden ser ordenados) con una clavija del tipo aprobado a nivel internacional con la finalidad de respetar las normas correspondientes a la seguridad del usuario. Utilizar siempre este cable de alimentación e introducir la clavija en un enchufe con una conexión de masa adecuada para evitar descargas eléctricas. y respetar las especificaciones CE. Dentro del controler se desarrollan altas tensiones que pueden causar graves daños o la muerte. Antes de efectuar cualquier operación de instalación o mantenimiento del controler desconectarlo del enchufe de alimentación.

---

**ADVERTENCIA**



Cable de alimentación: el cable correcto para el cableado eléctrico es un cable de tres hilos (L+N+Tierra).

L = Línea

N = Neutral

La sección del cable debe ser de al menos 1,0 mm<sup>2</sup> en el caso de una fuente de alimentación de 240 Vac o 2 mm<sup>2</sup> en el caso de una fuente de alimentación de 120 Vac para una longitud de cable de referencia de 3 m.

---

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

- temperatura: de +5 °C a +45 °C
- humedad relativa: 0 – 90 % (no condensadora).

En cualquier caso, es necesario que el aire de refrigeración pueda circular libremente alrededor del aparato.

### PRECAUCIÓN

El TwisTorr 305 FS onboard pertenece a la segunda categoría de instalación (o sobretensión) prevista por la normativa EN 61010-1. Conectar, por lo tanto, el dispositivo a una línea de alimentación que cumpla con los requisitos de dicha categoría. El TwisTorr 305 FS onboard está dotado de conectores para entradas/salidas y para la comunicación que deben ser conectados a los circuitos externos de forma que ninguna parte bajo tensión sea accesible. Comprobar que el aislamiento del dispositivo conectado al TwisTorr 305 FS onboard sea el adecuado incluso ante condiciones de avería individual tal como previsto en la normativa EN 61010-1.

---

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

## Montaje del TwisTorr 305 FS onboard Controller

En este párrafo se han incluido los principales procedimientos operativos. Antes de accionar el TwisTorr 305 FS onboard, realice las conexiones eléctricas y conecte el manguito de pre-vacío siguiendo las instrucciones en la sección “Technical Information”.

### ADVERTENCIA



**Para evitar lesiones a las personas y daños al equipo, antes de poner en marcha la bomba, asegúrese de haber fijado correctamente la bomba al sistema de vacío, de acuerdo con las instrucciones proporcionadas en el manual de la bomba (capítulo Instalación).**

### NOTA

Una vez conectado a la línea de voltaje, para operar el controlador (y por lo tanto poner en marcha la bomba) de una manera muy simple, conecte en el conector J5 (Entrada-Salida) el conector de acoplamiento precableado que se suministra junto con el controlador.

El conector de acoplamiento precableado es una forma simple y rápida de permitir que el controlador sea operado sin tener la necesidad de implementar otras conexiones externas cableadas como Serial Com o onboard I/O.

La bomba pre-vacío y la bomba TwisTorr 305 FS pueden encenderse simultáneamente.

# Puesta en marcha y funcionamiento del TwisTorr 305 FS onboard Controller

Para encender el TwisTorr 305 FS onboard es suficiente suministrar la tensión de alimentación. El controler reconoce automática-mente la presencia de las señales de interlock y de puesta en marcha y activa la bomba.

Al primer encendido, el controlador arranca automáticamente la bomba en modo "soft start". Espere a que se complete el arranque en soft start (luego espere a que la bomba alcance la velocidad máxima de rotación, indicada por un LED verde fijo), antes de apagar la bomba.

El arranque de la bomba en modo soft start permite la correcta redistribución del lubricante dentro de los rodamientos (lo cual es necesario después de un largo período de inactividad de la bomba).

Después de 60 días de inactividad es necesario proceder a una nueva puesta en marcha en Soft Start y después del primer arranque (para el cual el controlador efectúa el soft start automáticamente), las posteriores activaciones de puesta en marcha en Soft Start son a cargo del usuario.

Para arrancar con "Soft Start" activo, es necesario habilitar la modalidad mencionada a través del software o actuando sobre el conector de entradas y salidas remotas (Remote I/O connector).

Para el procedimiento de activación de la modalidad "Soft Start" consultar el punto "Signal Description" del capítulo "Technical Information". El LED verde ubicado en el panel del TwisTorr 305 FS onboard indica, con la frecuencia de su centelleo, las condiciones operativas del sistema:

- luz fija: la bomba está en rotación normal;
- centelleo lento (periodo de unos 400 ms): el sistema está en estado de rampa, o de Stop, o de "Waiting for interlock";
- centelleo rápido (periodo de unos 200 ms): condición de error

## Parada TwisTorr 305 FS onboard

Para parar el TwisTorr 305 FS onboard es suficiente quitar la tensión de alimentación.

## Parada de Emergencia

Para parar el TwisTorr 305 FS onboard en condiciones de emergencia es necesario desconectar el cable de alimentación.

## Mantenimiento

El TwisTorr 305 FS onboard no necesita ningún tipo de mantenimiento. Cualquier tipo de intervención sobre el sistema deberá ser realizado por personal autorizado. Antes de efectuar cualquier intervención sobre el sistema, es necesario desconectarlo de la alimentación. En caso de avería, es posible ponerse en contacto con el servicio de reparación Agilent o el servicio "Agilent advance exchange service" que permite obtener un controlador regenerado en sustitución del averiado.

**ADVERTENCIA**

**Antes de efectuar cualquier operación en el controler desenchufar el cable de alimentación.**



En caso de que un sistema debiera ser desguazado, proceder a su eliminación respetando las normas nacionales específicas.

## Eliminación

Significado del logo "WEEE" que se encuentra en las etiquetas.

El siguiente símbolo se aplica de acuerdo con la Directiva WEEE (Residuos de Aparatos Eléctricos y Electrónicos, por sus siglas en inglés) de la CE.

Este símbolo (válido solo en los países de la Comunidad Europea) indica que el producto en el que se aplica NO debe desecharse con la basura doméstica o industrial común, sino que debe enviarse a un sistema de recolección diferenciada. Se invita al usuario final a comunicarse con el proveedor del dispositivo, ya sea directamente con la Casa Matriz o con un revendedor, para iniciar el proceso de recolección y eliminación después de verificar las condiciones contractuales de venta.



Figura 4 Logotipo "WEEE"

Para más información, consulte:

<http://www.agilent.com/environment/product/index.shtml>

## Servicio

Si un cliente necesita un servicio de intercambio avanzado o reparación, contacte con un distribuidor local o contacte por correo a:

[vpt-customer-care@agilent.com](mailto:vpt-customer-care@agilent.com)

[vpl-customer-care@agilent.com](mailto:vpl-customer-care@agilent.com)

Es obligatorio completar la "Request for Return" para devolver su bomba a Agilent para mantenimiento (proporcionada al final del presente manual).

## 5

# 使用说明

关于本手册	76
有效性	76
定义与术语	77
小心、警告和注意的定义	77
警告符号	78
安全	80
正确使用	80
使用不当	81
<b>防护装备</b>	82
涡轮分子泵控制器安全指南	82
运输及存储	83
般说明	83
安装	84
安装准备	84
设置	85
TwisTorr 305 FS 遥控器安装	86
TwisTorr 305 FS onboard Controller 的启动和运行	87
TwisTorr 305 FS <b>板载控制器的关闭</b>	87
<b>紧急停止</b>	87
维护	88
处置	89
服务	90

## 关于本手册

### 有效性

本手册列出 TwisTorr 305 FS onboard Controller 用户说明，特别注意与安全、操作和一级维护有关的注意事项，受用户负责的维护操作限制。

维护操作在特定章节中说明，对于更高等级的维护设有具体规定（接受过维护操作专门培训的人员），用户不得执行此类操作。

有关正确的安装和启动/停止，请参阅“安装”章节。

要正确安装和启动/停止，请参考“Technical Information”章节。

### 注意

- 1 本手册包含有用信息，确保所有使用 TwisTorr 305 FS onboard Controller 的人员可以安全操作，在设备整个寿命期内发挥完美效率。
- 2 将本手册和所有相关出版物放在所有操作员、维护人员方便访问的位置。

## 定义与术语

### 小心、警告和注意的定义

本手册的一些重要参考内容用背景色突出显示。

#### 小心

程序开始前，提示小心消息，如果不遵守过程指示，可能导致设备损坏。

#### 警告



警告消息提醒操作员注意特定过程或做法，如果执行错误，可能导致人员重伤。

#### 注意

注意旨在让人注意重要信息，提供具体步骤的更多详细信息。

## 警告符号

以下是在 TwisTorr 305 FS onboard Controller 上与警告一起显示的符号列表。还显示了这些符号所指的危险。

三角形符号表示警告。在文档或仪器警告旁边可能出现的符号含义如下：



表示危险电压



一般危险



欧洲符合性声明



生产现场



CSA 认证



废弃电气和电子设备



RoHS中国认证



UK CA 打標

## 使用说明

以下符号可用作附在仪器上的警告标签。当您看到此符号时，请参阅相关操作或维修手册，了解该警告标签所指的正确操作过程。



以下符号会出现在仪器上以供您参考。

	一般危险
	表示危险电压
	CE认证
	CSA 认证
	RoHS中国认证
	废弃电气和电子设备
	UK CA 打標

## 安全

本节包含机械指令 2014/35/UE 规定的信息，这些信息对于遵守常规安全法规以及机器特定用途相关安全法规至关重要。

不遵守这些说明或本手册的其他说明，将导致设计阶段设想的安全条件无效，操作员可能发生事故。

对于不遵守技术文档中的安全规定而导致的机器损坏或者操作员或第三方人身伤害，Agilent Technologies 不承担任何责任。

### 正确使用

本手册包含机器安全工作需要遵守的重要警告和安全说明。

本手册所述产品仅用于本手册指定的应用领域。手册还规定了产品应用和操作基本要求相关的说明，以及可确保正常运行的安全措施。对于将产品用于非本手册介绍的用途或者产品使用过程中不遵守基本要求和安全措施，Agilent Technologies 不承担任何责任。

必须由能够采取必要安全措施，不会导致损坏或受伤的具备资质人员使用产品。产品的配套配件和设备必须由 Agilent Technologies 提供或授权使用。

必须由了解相关风险的专业技术人员执行任何调整或维护操作。

必须由 Agilent 授权人员执行产品维修。

## 使用说明

### 使用不当

对于因不正确使用 TwisTorr 305 FS onboard Controller 而造成的任何后果，安捷伦科技公司 (Agilent Technologies) 不承担任何责任。

使用不当将导致所有责任保证和索赔保证失效。使用不当的定义为：

- 使用未指定的安装材料安装设备
- 在冷凝环境中操作
- 在规定范围外的高湿环境中操作
- 在多尘环境中操作
- 在线路电压超出规格的情况下操作
- 在电离辐射区域内操作设备
- 在潜在爆炸区域内操作
- 在类冲击应力和振动或周期性力影响设备的系统中使用设备。

## 防护装备

由于此设备无需维护，因此不需要任何其他保护设备。

### 警告

#### 坠落物体造成伤害的风险



手动运输真空控制器时，存在负载滑落和坠落的危险。



- 双手携带控制器。



## 涡轮分子泵控制器安全指南

用户不应打开以下操作手册中所述的涡轮分子泵控制器，以避免可能损坏内部组件的风险。

### 警告

为避免损坏设备并防止操作人员受伤，应严格遵守本手册中提供的安装说明！



## 运输及存储

运输和存储控制器时，应满足以下环境要求：

- 温度：-20 °C 至 +70 °C
- 相对湿度：0 ÷ 95% (无冷凝).

## 般说明

如果在使用时不遵守或不完全遵守这些说明、未经培训的人员进行不正确使用、未经授权进行设备维修或在不符合国家特定标准进行使用，安捷伦不承担任何责任。

TwisTorr 305 FS板载控制器适用于TwisTorr 305泵。它配备辅助连接器，该辅助连接器可为附加风扇供电、控制通风阀以及通过串行线路（RS232或RS485）上连接的主机对TwisTorr 305 FS板载控制器进行远程操作。以下各段包含使用设备时确保操作员安全所需的所有信息。

有关详细信息，请参阅“技术信息”附录。

## 安装

### 安装准备

控制器具有特殊的保护性包装。如果这个包装在运输过程中出现损坏的迹象，请联系当地的销售办事处。打开控制器包装时，请确保不要摔落控制器或使其受到任何形式的撞击。请勿以未经授权的方式处置包装材料。请勿以未经授权的方式处理包装材料。该材料是100%可回收的，符合EEC指令94/62号和随后颁布的修正案。

## 设置

**警告**



为了用户安全，涡轮控制器（在带有电源电压的版本中，X3507-64131）必须由 3 线电源线（请参阅可订购的零件表）和插头（国际认可）供电。始终使用此电源线并将插头插入具有接地功能的适配插座中，以避免电击并符合 EC 要求。控制器内部产生的高压会可能导致严重的人身伤害甚至死亡。在安装或维修本机之前，请务必将其与插座断开。

**警告**



电源线：正确的电线电缆是三芯 (**L+N+PE**) 电缆。

L = Line (线)

N = Neutral (中性的)

PE= Protective Earth (保护地球)

对于等于 **3 m** 的参考电缆长度，在 **240 Vac** 电源电压的情况下，导线截面必须至少为 **1.0 mm<sup>2</sup>** 或在 **120 Vac** 电源电压的情况下至少为 **2 mm<sup>2</sup>**。

在操作过程中，必须遵守以下环境条件：

- 温度：从+5°C至+45°C;
- 相对湿度：0 - 95 %（无冷凝）。

在任何情况下，冷却空气都是必要的 可以在设备周围自由循环。

小心

**TwisTorr 305 FS 板载控制器属于 EN 61010-1 标准的第二安装（或过电压）类别。因此，请将设备连接到符合上述类别的电源线。TwisTorr 305 FS 板载控制器具有输入/输出和串行通信连接器，必须将其连接到外部电路，以使电子板的带电部件无法被接触到。根据 EN 61010-1 标准，即使发生单个故障，也要确保连接到 TwisTorr 305 FS 板载控制器的设备充分绝缘。**

有关可选配件的安装，请参阅附录“Technical Information”。

## TwisTorr 305 FS 遥控器安装

本段详细介绍了主要操作程序。请参考涡轮泵说明手册，在操作涡轮控制器之前，进行所有真空歧管和电气的连接。

警告



为避免人身伤害和设备损坏，在启动泵之前，请确保已按照泵手册（安装章节）中的说明将泵正确固定到真空系统上。

注意

连接到线路电压后，要以非常简单的方式操作控制器（然后启动泵），请在 J5 连接器上连接与控制器一起提供的预接线配对连接器。

预接线的配对连接器是一种允许控制器操作的简单快捷的方法，而无需实施其他有线外部连接，如串行通信或远程输入/输出。

前置泵和 TwisTorr 305 FS 泵可以同时开启。

## TwisTorr 305 FS onboard Controller 的启动和运行

如要启动TwisTorr 305 FS板载控制器，只需提供电源即可。控制器会自动识别联锁和启动信号的存在并启动泵。

泵第一次开机时，控制器会在“慢启动”模式下自动启动泵。在关闭泵之前，请等待慢启动完成（直到LED发出稳定的绿灯光来指示泵已达到最大转速）为止。

通过“慢启动”模式启动泵，轴承内的润滑脂将以正确的方式进行重新分配（在长时间不使用泵之后必须这样做）。

如上所述，第一次启动泵时，控制器会自动执行慢启动。随后，如果泵在60天的时间内保持不活动状态，则下次启动泵时，用户必须以“慢启动”模式启动泵。

如要在激活“慢启动”的情况下启动泵，必须在软件中或通过远程I/O连接器启用慢启动模式。

有关“慢启动”模式的激活过程，请参见“技术信息”章节中的“信号说明”一节。TwisTorr 305 FS板载控制器上绿色LED灯的闪烁速度表示系统运行状况：

- 无闪烁：泵正常旋转；
- 缓慢闪烁（大约400毫秒）：系统处于突起加减速、停止或“等待互锁”的状态；
- 快速闪烁（大约200毫秒）：错误情况。

### TwisTorr 305 FS 板载控制器的关闭

如要关闭TwisTorr 305 FS板载控制器，只需断开电源即可。集成控制器会立即停止泵。

### 紧急停止

如要在紧急情况下立即停止TwisTorr 305 FS板载控制器，请拔下电源线。

## 维护

TwisTorr 305 FS板载控制器不需要任何**维护**。在系统上执行的任何工作都必须由**授权人员**执行。在系统上进行任何操作之前，必须将其与**电源断开连接**。如果发生故障，可以使用**安捷伦维修服务**。可通过**安捷伦预先更换服务**来替换控制器。

### 警告



在控制器上**进行任何操作之前，请先断开其电源**。

控制器**报废时，必须按照国家具体标准进行处理**。

## 处置

标签中的“WEEE”徽标含义。

按照 EC WEEE（废电子电气设备）指令应用以下符号。

此符号（仅在欧盟国家有效）表示所适用产品不得与家庭或工业垃圾一起处置，必须送至专门废弃物回收系统。因此，欢迎最终用户联系设备供应商（无论是母公司还是零售商），在检查销售合同条款后启动回收和处置流程。



图 5 “WEEE” 徽标

有关更多信息，请参考：

<http://www.agilent.com/environment/product/index.shtml>

使用说明

## 服务

如果客户需要高级替换或维修服务，请联系当地经销商或直接发送邮件至

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

需要填写“Request for Return”表才能将泵退回到 Agilent 维修部门（在本手册末尾会提供）。

## 6

# 取扱説明書

この説明書について	92
有効期	92
定義と用語	93
警告・注意・注の定義	93
警告の記号	94
安全性	96
適切な使用	96
不適切な使用	97
保護装具	98
运输及存储	99
概要情報	99
取り付け	100
取り付けの準備	100
取り付け	101
TwisTorr 305 FS の取り付け	102
TwisTorr 305 FS onboard Controller コンパクトの起動と操作	103
TwisTorr 305 FS オンボードのスイッチを OFF にします	103
緊急停止	103
メンテナンス	104
処分	105
サービス	106

## この説明書について

### 有効期

本説明書には、特に安全、操作手順およびユーザーに必要なメンテナンス手順のみの簡易メンテナンスに関連する考え方を参考に、TwisTorr 305 FS onboard Controller の使用方法をユーザーに説明します。

高度なメンテナンス（メンテナンス手順の個別訓練を受けた人員）に関連した各条項とあわせて、各セクションで説明するメンテナンス手順は、ユーザーは行わないでください。

正しい取り付けと起動/停止については、「取り付け」の章をご確認ください。

正しい取付および始動/停止については、「Technical Information」セクションを参照してください。

### 注

- 1 本説明書には、TwisTorr 305 FS onboard Controller を使用する人員がすべて安全に使用でき、また機器使用寿命中に効率よく使用できるのに、役立つ情報が記載されています。
- 2 本説明書は、関連するすべての刊行物とあわせて、オペレータ/メンテナンス要員に分かりやすい、取り扱いしやすい場所に保管してください。

## 定義と用語

### 警告・注意・注の定義

このマニュアルのいくつかの重要な参照部は、対比色でハイライトされ枠で囲ってあります。

#### 注意

「注意」のメッセージは、もし監視されていない場合に装置に損害が生じる可能性を示しています。

#### 警告



「警告」のメッセージは、もし正しく操作されない場合に、重大な人身障害につながる可能性がある特定の操作や実行への操作者の注意の必要性を示しています。

#### 注

「注」は、重要な情報への注意喚起と特定の工程に関してのより詳細な情報を提供しています。

## 警告の記号

下記は、TwisTorr 305 FS onboard Controller上の警告と併せて表示される記号のリストです。危険性の説明も記載されています。

三角の記号は警告を示しています。説明書内または装置上の警告記号と併に示される記号の意味は下記の通りです：



危険な電圧



包括的な危険



CE適合宣言書



製造所



CSA 認証



廃電気・電子製品に関するEU WEEE指令



中国RoHS 指令認証











UK CAマーキング

## 取扱説明書

下記の記号は、装置に貼り付けられた警告ラベル上に使用されることがあります。もしこの記号を目視した場合、その警告ラベルに関する正しい工程のために、関連する操作またはサービスマニュアルを参照してください。



参考情報として、下記の記号は装置上に示されます。

	包括的な危険
	危険な電圧
	CE認証
	CSA 認証
	EU RoHS指令認証
	中国RoHS 指令認証
	廃電気・電子製品に関するEU WEEE指令
	UK CAマーキング

## 安全性

このセクションは、一般的、かつ機械の特定の使用に関する安全規定の遵守と監督に必須である機械指令（Machinery Directive 2014/35/UE）の情報を含みます。

これらの指示とこのマニュアル内のその他の指示に従わなかった場合、設計時に予測された安全状態が非効率に損なわれ、機械を操作する者へ事故が生じる可能性があります。

アジレント・テクノロジー株式会社は、技術的な参照書に示された安全規定の不順守により派生した、機械または操作者もしくは第三者の身体的安全への損害・損傷の責任を全て拒否します。

### 適切な使用

このマニュアルは、装置一式が安全に機能するために遵守されるべき重要な警告と安全の手引きを含みます。

このマニュアルに説明されている製品は、説明書に特記されている分野への活用に限定了目的のものです。マニュアルはまた、製品の活用と操作に必須の要件に関する指示と標準の操作を保証するために適用される安全対策を提供します。アジレント・テクノロジー株式会社は、このマニュアル内説明されているもの以外の適用、または必須の要件や安全対策に注意が払われていない場合にいかなる保証も提供せず、またはいかなる責任も負いません。

製品は、損傷や怪我を生じさせない状況下において必要な安全対策を講じることができる有資格者のみ使用することができます。製品と共に使用されたいかなる付属品や機器も、アジレント・テクノロジー株式会社により共有もしくは承認されたものでなければなりません。

いかなる調整またはメンテナンスの操作は、リスクについて承知している専門家の技術者により実行されなければいけません。

製品の修理は、専らアジレント・テクノロジー株式会社により行われなければなりません。

### 不適切な使用

アジレント・テクノロジー株式会社は、TwisTorr 305 FS onboard Controllerの不適切な使用により派生した全ての責任を拒否します。

不適切な使用は、責任と保証への全ての請求権が喪失します。不適切な使用は下記の通り定義されます：

- 指定されていない取り付け器具を使用した装置の設置
- 凝縮環境での操作
- 指定範囲外の高湿度環境での操作
- ほこりの多い環境での操作
- 仕様外の線間電圧での操作
- 電離放射線がある場所での装置の操作
- 爆発の可能性のある場所での操作
- 衝撃のような圧力や振動、または断続的な力がデバイスに影響を与えるシステム下での装置の使用。

## 保護装具

この装置にはメンテナンスは想定されていないため、他の保護装具は必要ありません。

### 警告

#### 落下物による怪我の危険性



真空コントローラーを手で運ぶ際に、荷物が滑って落下する危険性があります。



- コントローラーは両手で持ち運んでください。



#### ターボ分子ポンプ コントローラーの安全ガイドライン

以下の操作マニュアルに記載されているとおり、ターボ分子ポンプコントローラーは、内部コンポーネントが損傷するリスクを回避するために、使用者は開けないでください。

### 警告

装置の損傷と操作者の怪我を避けるために、このマニュアルに記載された取り付けの指示に厳密に従ってください！



## 运输及存储

コントローラーを輸送、保存する時は、下記の環境仕様を超過してはいけません：

- 温度範囲：-20° C から +70° C まで
- 相対湿度範囲：0 から 95 % まで（結露なし）。

## 概要情報

この装置は技術者による使用を対象としています。使用者は、この取扱説明書と Agilent により提供される他の追加情報もあわせて、装置を操作する前に全てお読みください。Agilent は、部分的であってもこれらの取扱説明に従わない場合や、訓練されていない人による不適切な使用、装置への認められていない修理、または特定の国家規格の規定に相いれないいかなる行動によって生じたいかなる事態にも責任を負いません。

TwisTorr 305 FS オンボードコントローラーは、TwisTorr 305 ポンプとの使用に適しています。追加のファンに電力を供給したり、通気弁を制御したり、シリアルライン (RS232 または RS485) に接続されたホストコンピューターによって TwisTorr 305 FS オンボードをリモート操作したりできるようにする補助コネクタと共に装備されています。

下記の章は、装置の使用時に操作者の安全を保証するために必要なすべての情報が含まれています。詳細情報は、「技術情報」の追記に記載されています。

## 取り付け

### 取り付けの準備

コントローラーは、特殊な保護梱包で提供されます。もしこの梱包に移送の間に生じた可能性のある破損が見受けられる場合、現地販売事務所に問い合わせてください。コントローラーの梱包を開ける際には、落とさない、またはいかなる形での衝撃を与えないようにしてください。不適切な方法で梱包材料を破棄しないでください。EEC指令94/62とそれに準ずる改正を遵守しています。

## 取り付け

**警告**



使用者の安全のために、ターボコントローラー（主電源電圧のバージョンで：**X3507-64131**）は。

3本のワイヤー電源コード（注文可能な部品を参照）とプラグ（国際的に認可されたもの）で電源が供給されなければいけません。電気ショックを避け、また **CE** 要件を満たすために、常にこの電源コードを使用し、適切な接地接続の電源ソケットにプラグを挿入してください。コントローラー内で発達した高電圧は、重大な損傷や死亡を引き起こす可能性があります。ユニットを取り付ける前または修理する前には常に、ソケットから接続を断ってください。

**警告**



電源コード：電気配線に適したケーブルは、3線（L+N+PE）ケーブルです。

**L = Line** (ライン)

**N = Neutral** (中性)

**PE= Protective Earth** (保護アース)

3 m と同等の長さの標準ケーブルに対して、ワイヤー部分は **240Vac** の供給電圧の場合は  $1.0\text{mm}^2$  以上、**120Vac** の供給電圧の場合は  $2\text{mm}^2$  する必要があります。

操作中、下記の環境状況は遵守されなければいけません：

- 温度：+5 °C から+45 °Cまで;
- 相対湿度：0 - 95 %（結露なし）

穴を通して自由に通気が取れるように所定の位置に設置されなければいけません。

注意

**TwisTorr 305 FS** オンボード は、**EN 61010-1** 規格の第 2 取り付け（または過電圧）カテゴリーに属します。したがって、上記のカテゴリーに準拠する主電源ラインに機器を接続してください。 **TwisTorr 305 FS** オンボードには、電子ボードの通電部分にアクセスできないように外部回路に接続する必要がある入力/出力およびシリアル通信コネクタがあります。 **EN 61010-1** 規格による単一故障の場合でも、の **TwisTorr 305 FS** オンボードに接続されている機器の分離が十分であることを確認してください。

オプションのアクセサリの取り付けについては、“Technical Information” を参照してください。

## TwisTorr 305 FS の取り付け

この章は、主な操作手順について詳しく説明します。ターボコントローラーを操作する前に、ターボポンプの取扱説明書を参照して、すべての真空マニホールドと電気接続を行ってください。

警告



人への損傷や機器への損傷を避けるために、ポンプを始動する前に、ポンプのマニュアル（「設置」の章）に記載されている指示に従って、ポンプが真空システムに正しく固定されていることを確認してください。

注

線間電圧に接続したら、非常に簡単な方法でコントローラーを操作するには、（そしてポンプを始動するために）コントローラーに付属している配線済みの相手側コネクタを J5（入出力）コネクタに接続します。

配線済みの相手側コネクタは、シリアルコムやリモート I/O などの他の配線された外部接続を行う必要なく、コントローラを操作できるようにする簡単で迅速な方法です。

フォアポンプと TwisTorr305FS ポンプを同時にオンにすることができます

## TwisTorr 305 FS onboard Controller コンパクトの起動と操作

TwisTorr 305 FS オンボードのスイッチをオンにするには、主電源を供給するだけで十分です。コントローラーは自動的にインターロックと始動信号の存在を認識し、ポンプを始動します。

ポンプが電源オンになる初回時は、コントローラーによって「ソフトスタート」モードで自動的に起動されます。ソフトスタートが完了するまで（つまり、ポンプが最大回転速度に達したことを示すLEDが一定の緑色に光るまで）待ってから、ポンプの電源を切ります。

ポンプをソフトスタートモードで起動することにより、ベアリング内の潤滑油が正しい方法で再分配されます（ポンプが長期間非稼働である後に必要です）。

前述のように、ポンプが初回に起動した時に、コントローラーはソフトスタートを自動的に始動します。その後、もしポンプが60日間非稼働のままである場合は、次回起動するときに、ユーザーはソフトスタートモードで起動しなければなりません。

「ソフトスタート」をアクティブにしてポンプを始動するには、ソフトウェアで、またはリモートI/Oコネクタでソフトスタートモードを有効にする必要があります。

「ソフトスタート」モードのアクティブ化手順については、「技術情報」の章の「信号の説明」を参照してください。

TwisTorr 305 FS オンボードのフロントパネル上の緑色のLEDの点滅速度は、システムの動作状態を示します：

- 点滅なし：ポンプは正常に回転しています。
- 低速の点滅（約400 msの周期）：システムはランプ、停止、または「インターロック待機中」の状態です。
- 高速の点滅（約200 msの周期）：エラーの状態です。

### TwisTorr 305 FS オンボードのスイッチを OFF にします

TwisTorr 305 FS オンボードのスイッチをオフにするには、主電源を切るだけで十分です。統合されたコントローラーはポンプを即座に停止します。

### 緊急停止

緊急事態にTwisTorr 305 FS オンボードを即座に停止するには、電源コードを取り外してください。

## メンテナンス

TwisTorr 305 FS オンボードはいかなるメンテナンスも必要としません。システムになされるいかなる作業も、認可された担当者により実行されなければなりません。

システムにいかなる操作を実行する前には、必ず電源を切断してください。もし故障が発生した場合は、Agilentの修理サービスを利用することもできます。交換用コントローラーは、Agilentを通じてアドバンスエクスチェンジをもとに入手できます。

**警告**



コントローラー上にいかなる作業を実行する前には、電源から接続を断ってください。

もしポンプを廃棄する場合には、特定の国の規定に基づき廃棄されなければなりません。

## 処分

ラベルに記載の“WEEE” ロゴの意味。

下記の記号は、EC WEEE（電気電子機器廃棄物）指令にしたがって適用されま  
す。

この記号（EC 諸国内のみで有効）は、この記号が適用される製品は、通常の家  
庭ゴミとあわせて処分してはいけないこと、また個別廃棄物回収システムに送付  
しなければならないことを表します。そのため、エンドユーザーは、親会社か小  
売業者かに関係なく、当該機器のサプライヤに連絡の上、契約上の販売条件の確  
認後、回収および処分プロセスを開始することが求められます。



図 6 ロゴ “WEEE”

詳細は、下記サイトを参照してください。

<http://www.agilent.com/environment/product/index.shtml>

## サービス

お客様が最新交換サービスまたは修理サービスが必要な場合は、現地代理店もしくは下記まで直接ご連絡ください：

[vpt-customer care@agilent.com](mailto:vpt-customer care@agilent.com)

[vpl-customer care@agilent.com](mailto:vpl-customer care@agilent.com)

“Request for Return”の申請の完了には、サービス提供のためにお客様のポンプをAgilentへ返送する必要があります。（このマニュアルの最後に提供されています）。

# 7

## Instructions for Use

About this manual	108
Validity	108
Definitions and terminology	109
Definition of Caution, Warning and Note	109
Warning Symbols	110
Safety	112
Proper use	112
Improper use	113
Protective equipment	114
Safety guideline for Turbomolecular Pump Controllers	114
Transport & Storage	115
Product description	115
Installation	116
Preparation for installation	116
Set-up	117
TwisTorr 305 FS onboard Controller Mounting	118
Startup and Operation of the TwisTorr 305 FS onboard Controller	119
Switching off the TwisTorr 305 FS onboard	119
Emergency Stop	119
Maintenance	120
Disposal	121
Service	122

## About this manual

### Validity

This manual lists the instructions for the users of the TwisTorr 305 FS onboard Controller, with particular reference to the notions relating to safety, operation and first level maintenance, limited to maintenance operations for which the user is responsible.

The maintenance operations, illustrated in the specific sections, with specific provisions relating to the higher level of maintenance (personnel specifically trained for maintenance operations) must not be carried out by the user.

For a correct installation and start/stop, please refer to "Installation" paragraph.

For a correct installation and start/stop, please refer to " Technical Information " section.

#### NOTE

- 1** This manual contains useful information so that all personnel using the TwisTorr 305 FS onboard Controller can operate it safely and guarantee perfect efficiency, for its entire life span.
  - 2** Keep this manual, together with all the related publications, in an accessible place known to all operators/maintenance personnel.
-

## Definitions and terminology

### Definition of Caution, Warning and Note

Some important references of this manual are highlighted and framed in contrasting color.

#### CAUTION

Caution messages are displayed before procedures which, if not observed, could cause damage to the equipment.

---

#### WARNING



Warning messages draw the operator's attention to a specific procedure or practice which, if not performed correctly, could result in serious personal injury.

---

#### NOTE

Notes are intended to call attention to important information and provide more detail regarding specific steps.

---

## Warning Symbols

The following is a list of symbols that appear in conjunction with warnings on the TwisTorr 305 FS onboard Controller. The hazard they describe is also shown.

A triangular symbol indicates a warning. The meanings of the symbols that may appear alongside warnings in the documentation are as follows:



Dangerous voltages



Generic hazard



European Declaration  
of Conformity



Manufacturing Site



CSA certification



Waste Electrical and  
Electronic Equipment



RoHS China certification






UK CA Marking

## Instructions for Use

The following symbol may be used on warning labels attached to the instrument. When you see this symbol, refer to the relevant operation or service manual for the correct procedure referred to by that warning label.



The following symbols appear on the instrument for your information.

	Generic hazard
	Dangerous voltages
	CE certification
	CSA certification
	RoHS China certification
	Waste Electrical and Electronic Equipment
	UK CA Marking

# Safety

This section contains the information, prescribed by the Low voltage Directive 2014/35/UE, which is essential for the compliance and observance of the safety regulations both generally and in relation to the specific use of the product.

Failure to comply with these instructions and the other instructions contained in this manual may render the safety conditions envisaged in the design phase inefficient and cause accidents to those operating the product.

Agilent Technologies declines all responsibility for damage to the product or for the physical safety of the operator or third parties deriving from the non-observance of the safety rules indicated in the technical documentation.

## Proper use

This manual contains important warnings and safety instructions to be observed in order for the unit to work safely.

The product described in this manual is intended exclusively for the area of application specified in the instructions. The manual also provides indications regarding the essential requirements for the application and operation of the product as well as the safety measures that can be adopted to guarantee regular operation. Agilent Technologies does not provide any guarantee or assume any responsibility for applications other than those described in this manual or in which the essential requirements and safety measures are not respected.

The product must only be used by qualified personnel who are able to take the necessary safety measures under conditions that do not cause damage or injury. Any accessories and equipment used with the product must be supplied or approved by Agilent Technologies.

Any adjustment or maintenance operation must be performed by a professional technician informed about the risks.

Repairs on the product must be carried out exclusively by Agilent authorized personnel.

## Instructions for Use

### Improper use

Agilent Technologies declines all responsibility, deriving from the improper use of the TwisTorr 305 FS onboard Controller.

Improper use will cause all claims for liability and warranties to be forfeited. Improper use is defined as:

- installation of the device with unspecified mounting material
- operation in condensing environment
- operation in high humidity environment out of the specified range
- operation in dusty environment
- operation with line voltages out of specifications
- operation of the device in areas with ionizing radiation
- operation in potentially explosive areas
- use of the device in systems in which impact-like stress and vibrations or periodic forces affect the device.

## Instructions for Use

### Protective equipment

As no maintenance is foreseen for this device, any other protective equipment is not needed.

#### **WARNING**

#### **Risk of injury through falling objects**



When transporting vacuum controllers by hand, there is a danger through loads slipping and falling down.



- Carry controllers two-handed.



### Safety guideline for Turbomolecular Pump Controllers

Turbomolecular pump controllers as described in the following operating manual should not be opened by the user to avoid the risk of damaging internal components.

#### **WARNING**

To avoid damage to equipment and to prevent injuries to operating personnel the installation instructions as given in this manual should be strictly followed!



### Transport & Storage

When transporting and storing the TwisTorr 305 FS onboard, the following environmental conditions should not be exceeded:

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

### Product description

This equipment is destined for professional use. Before operating the equipment, the user should read this instruction manual and any other additional information supplied by Agilent. Agilent declines any responsibility for non-compliance, even partial, with these instructions, improper use by untrained persons, unauthorized repairs of the equipment or use not complying with specific national standards.

The TwisTorr 305 FS onboard controller is suitable for use with TwisTorr 305 pumps. It is equipped with auxiliary connectors that make it possible to power an additional fan, to control the vent valve and for remote operation of the TwisTorr 305 FS onboard by a host computer connected on a serial line (RS232 or RS485).

The following paragraphs contain all the information necessary to guarantee operator safety when using the equipment.

For detailed information, refer to the "Technical Information" section.

## Installation

### Preparation for installation

The TwisTorr 305 FS onboard Controller is supplied in a special protective packing. If this shows signs of damage which may have occurred during transport, contact your local sales office.

When unpacking the controller 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 94/62 and subsequent amendments.

## Set-up

**WARNING**



For user safety, the Controller (in grid voltage version, X3507-64131) must be powered by a 3-wire power cord (see orderable parts table) and plug (internationally approved). Always use this power cord and insert the plug in a socket with a suitable ground connection to avoid electrical shock and to comply with EC requirements. The high voltage developed inside the controller can cause severe injury or death. Before installing or servicing the unit, always disconnect it from the socket.

---

**WARNING**



Power supply cord: the correct cable for electrical wiring is a three wires (L+N+PE) cable.

L = Line

N = Neutral

PE = Protective Earth

The wire section has to be at least 1.0 mm<sup>2</sup> in case of 240 Vac supply voltage or 2 mm<sup>2</sup> in case of 120 Vac supply voltage for a reference cable length equal to 3 m.

---

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 complied with:

- temperature: from +5 °C to +45 °C
- relative humidity: 0 – 90 % (non-condensing).

In any case it is necessary that the cooling air can circulate freely around the apparatus.

### CAUTION

The TwisTorr 305 FS onboard belongs to the second installation (or overvoltage) category of the EN 61010-1 standard. Therefore, connect the device to a mains line that complies with the above category. The TwisTorr 305 FS onboard has Input/Output and serial communication connectors that must be connected to external circuits in such a way that no live parts of electronic board are accessible. Make sure that the insulation of the device connected to the TwisTorr 305 FS onboard is adequate even in the case of a single fault as per the EN 61010-1 standard.

---

For installation of optional accessories, see "Technical Information" section.

## TwisTorr 305 FS onboard Controller Mounting

This paragraph details the main operating procedures. Prior to operating the Controller, make all vacuum manifold and electrical connections, referring to the Turbo pump instruction manual.

### WARNING



**To avoid damage to people and damage to the equipment, before starting the pump, make sure that you have correctly fastened the pump to the vacuum system, according to the instructions given in the pump manual (chapter Installation).**

### NOTE

Once connected to the line voltage, to operate the controller (and then to start the pump) in a very simple way, connect on the J5 connector (Input-Output) the pre-wired mating connector that is supplied together with the controller.

The pre-wired mating connector is a simple and quick way to allow the controller to be operated without having the need to implement other wired external connections like Serial Com or onboard I/O.

The forepump and TwisTorr 305 FS pump can be switched on at the same time.

# Startup and Operation of the TwisTorr 305 FS onboard Controller

To switch on the TwisTorr 305 FS onboard, it is sufficient to provide the mains supply. The controller automatically recognizes the presence of the interlock and start signals and starts the pump.

The first time the pump is powered on, it is started automatically in “Soft Start” mode by the controller. Wait until the Soft Start completes (that is, until the LED emits a fixed green light indicating that the pump has reached its maximum rotation speed), before switching off the pump.

By starting the pump in Soft Start mode, the grease inside the bearings is redistributed in the correct manner (necessary after the pump has been inactive for a long period).

As stated, the first time the pump is started, the controller executes Soft Start automatically. Subsequently, if the pump remains inactive for a period of time of 60 days, the next time it is started the user must start it in Soft Start mode.

To start the pump with “Soft Start” activated, Soft Start mode must be enabled in the software or through the Remote I/O connector.

For the “Soft Start” mode activation procedure, see the “Signal Description” paragraph in the “Technical Information” chapter. The flashing rate of the green LED on the TwisTorr 305 FS onboard front panel indicates system operating conditions:

- with no flashing: the pump is rotating normally;
- slow flashing (period of about 400 ms): the system is in ramp, or in Stop, or in “Waiting for interlock” status;
- fast flashing (period of about 200 ms): error condition.

## Switching off the TwisTorr 305 FS onboard

To switch off the TwisTorr 305 FS onboard, it is sufficient to remove the mains supply. The integrated controller immediately stops the pump.

## Emergency Stop

To stop the TwisTorr 305 FS onboard immediately in an emergency condition, remove the power cord.

## Maintenance

The TwisTorr 305 FS onboard does not require any maintenance. Any work performed on the system must be carried out by authorized personnel. Before carrying out any operation on the system, it must be disconnected from the power supply. If a fault occurs, it is possible to use the Agilent repair service. Replacement controllers are available on an advance exchange basis through Agilent.

**WARNING**



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

If a controller is to be scrapped, it must be disposed off 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.



**Figure 7** Logo "WEEE"

For more information refer to:

<http://www.agilent.com/environment/product/index.shtml>

## Service

Should a customer need an advanced exchange or repair service, please contact local distributor or directly mail to

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

Completion of the "Request for Return" form is required to return your pump to Agilent for service (provided at the end of this manual).

## 8 Technical Information

Description of the TwisTorr 305 FS onboard Controller	124
Technical specification	125
Controller Outline	127
TwisTorr 305 onboard Controller Installation	129
Bottom Mounting	129
Side Mounting	130
Interconnections	131
P3 – Vent	131
P5 – External Fan	132
J5 – In-Out	132
Signal Description	134
How to Connect the Open Collector Inputs of the Controller	137
How to Connect the Outputs of the Controller	141
J6 – Serial	141
Input power connection on the model X3507-64130	143
RS 232/RS 485 communication description	143
Communication Format	143
Communication Protocol	144
Examples	146
Window Meanings	148
Air Cooling Kit Installation	152
Use in the Presence of Magnetic Fields	155
Accessories and Spare Parts	155

## Description of the TwisTorr 305 FS onboard Controller

Two different versions of the controller are available:

X3507-64130      TwisTorr 305 FS onboard controller 24V

X3507-64131      TwisTorr 305 FS onboard controller 110/230V

The controller is a solid-state frequency converter with the following features:

- Drives the pumps of the TwisTorr 305 FS family.
- Powers the pump cooling fan.
- Drives the vent valve.
- Remote I/O compatible with the previous version.
- Serial communication compatible with the previous RS 232 and 485 versions; due to some new features which are present in the software, some minor changes to the serial protocol have been made.

### NOTE

With reference to the previous 304 and 301 pump families, the serial “window protocol” differences implemented for the 305 family are the following:

- Added WIN001, WIN143, WIN144, WIN167, WIN157
- Removed WIN106, WIN107

Refer to the “Technical Information – window meanings” chapter to get the relevant details.

---

The dedicated controller is a solid-state frequency converter which is driven by a single chip microcomputer and consists of two PCBs which include power supply and 3-phase output, analog and input/output section, microprocessor and digital section. The power supply, together with the 3-phase output, converts the single phase AC mains supply or 24Vdc supply into a 3-phase, low voltage, medium frequency output which is required to power the pump.

The controller can be operated by a remote host computer via the serial connection. A Windows-based software is available (optional).

## Technical specification

The following table lists the main technical data of the TwisTorr 305 FS onboard Controller.

**Table 1 TwisTorr 305 FS onboard Controller Technical Specification**

TwisTorr 305 FS onboard Controller Technical data	
Pwr supply (mains):	
Input voltage:	100 - 240 Vac
Input freq.:	50 - 60 Hz
Max input pwr:	300 VA
Stand-by pwr:	10 W
Max oper. power:	150 W with water or air cooling
Protection fuse	4 A
Pwr supply (24 Vdc):	
Input voltage:	24 Vdc
Max input pwr:	200 W
Stand-by pwr:	10 W
Max oper. power:	150 W with water or air cooling
Protection fuse	8 A
Compliance with:	EN61010-1 (2010)/A1 (2019) EN 61326-1 (2013)
Power cable	With European or NEMA plug 3 meters long (optional)
Serial communication (TwisTorr 305 kit)	RS232 cable with a 9-pin D type male connector and a 9-pin D type female connector, and TwisTorr software (optional)
Storage temperature	-20 ° C to +70 ° C
Input	75 Vac, three phase, 1010 Hz
Internal use only	Max. altitude 2000 m
Installation category	II

## Technical Information

TwisTorr 305 FS onboard Controller Technical data	
Pollution degree	2
Operating ambient temperature	5 ° C to 40 ° C
Weight	1 kg (2,2 lbs)

### NOTE

When the TwisTorr 305 FS onboard has been stored at a temperature less than 5°C, wait until the system has reached the above mentioned temperature.

## Controller Outline

The outline dimensions for the TwisTorr 305 FS onboard Controllers are shown in the following figure.

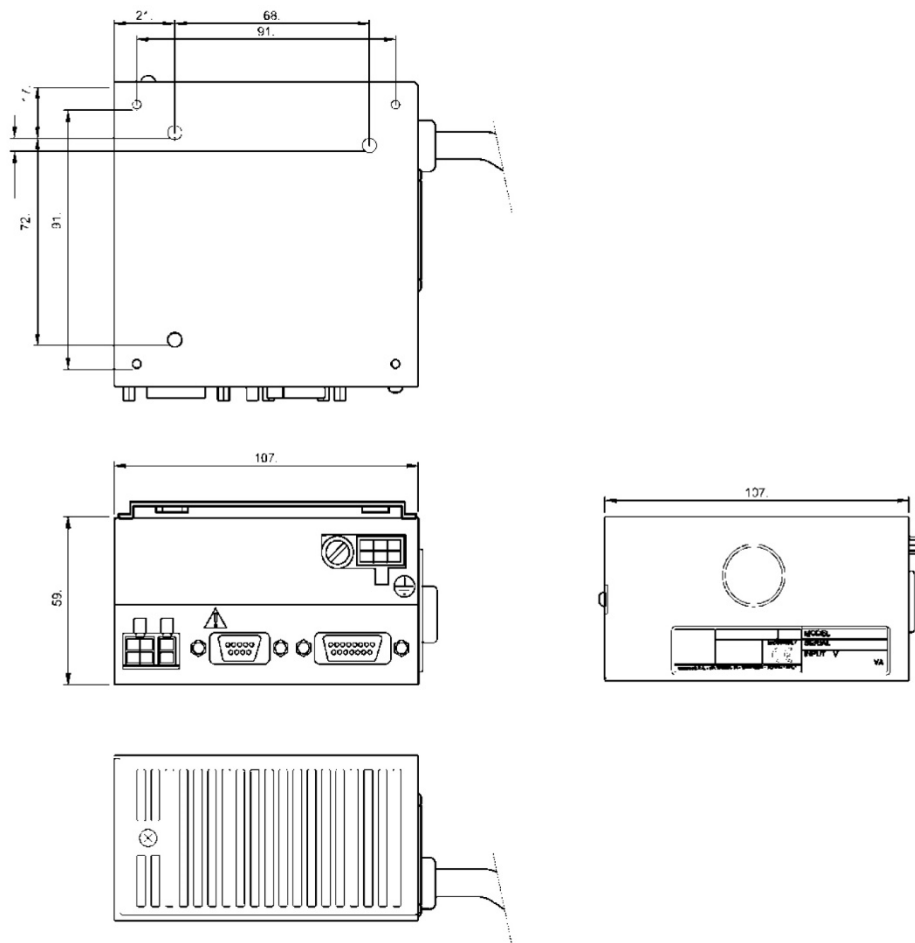


Figure 8 Outline 24 V – X3507-64130

## Technical Information

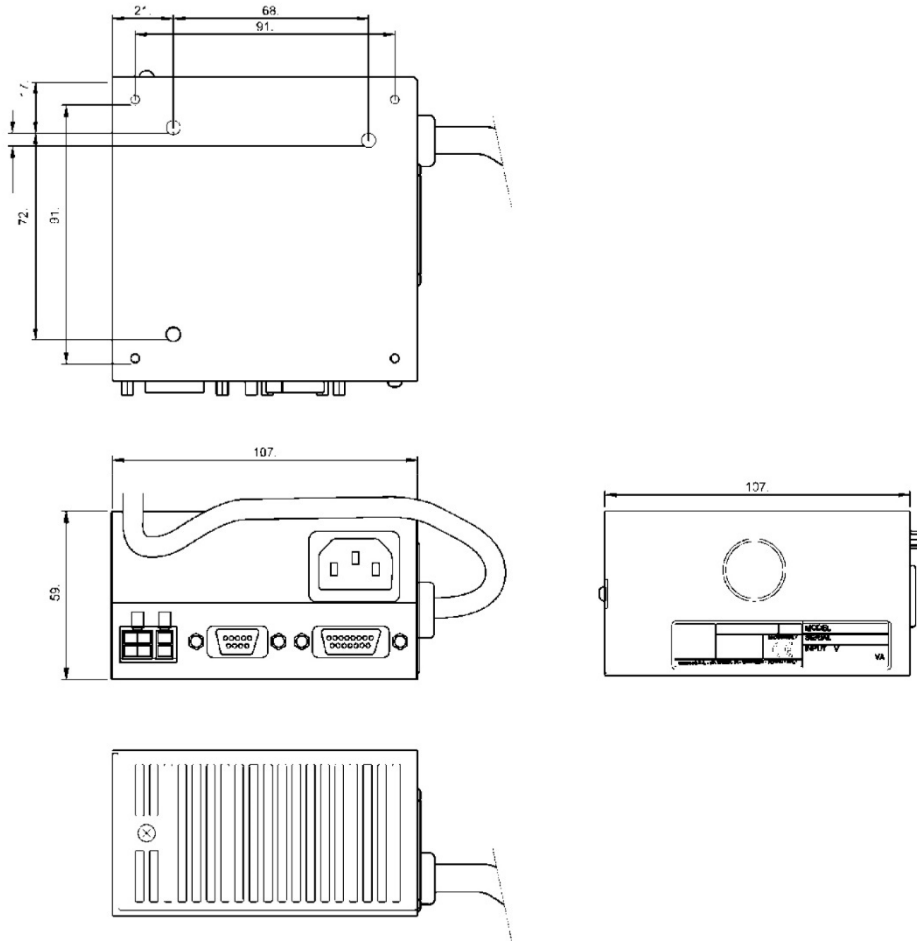


Figure 9 Outline X3507-64131

## TwisTorr 305 onboard Controller Installation

The controller can be mounted in two positions:

- bottom mounting (as per the complete system)
- side mounting

To install the controller execute the following procedures.

### Bottom Mounting

See the following figure.

- 1 Turn the pump upside-down;
- 2 Screw the 4 fixing studs (provided with the accessories bag) in the holes on the bottom of pump;
- 3 Place the controller on the studs aligning the 4 rubber holes on the top of the controller case;
- 4 Press the controller towards the pump;
- 5 Turn the pump again;
- 6 Plug the line cord and connect the pump cable to the pump;
- 7 Connect the mating connector with the jumper on the interlock signal to start the pump.

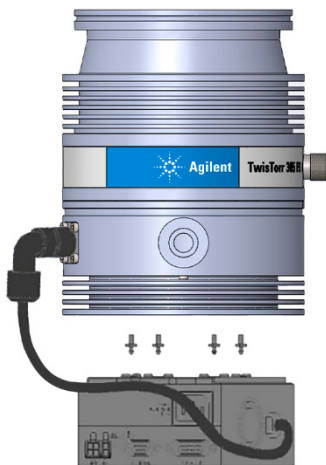


Figure 10 Mounting 1

## Technical Information

### Side Mounting

#### NOTE

The L-shaped bracket (P/N X3500-68012) is available as an option.

See the following figure.

- 1 Screw the 4 fixing studs (provided with the accessories bag) in the holes on the L-shaped bracket;
- 2 Place the controller on the studs aligning the 4 rubber holes on the top of the controller;
- 3 Press the controller towards the bracket;
- 4 Turn the pump upside-down;
- 5 Place the bracket on the pump bottom;
- 6 Screw a little bit the three screws M4 in the holes on the pump bottom.

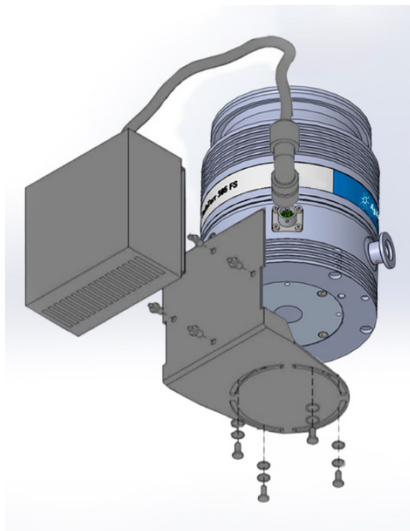


Figure 11 Side Mounting

## Interconnections

The following figure shows the TwisTorr 305 FS onboard interconnections.

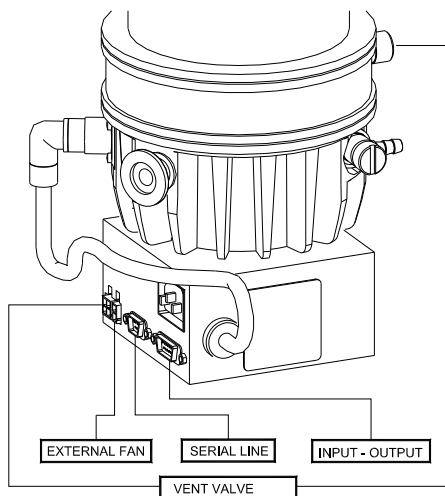


Figure 12 TwisTorr 305 FS onboard interconnection

### P3 – Vent

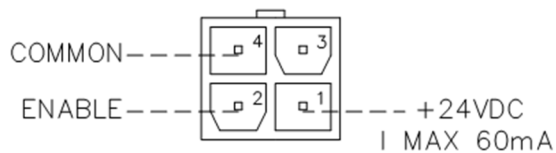
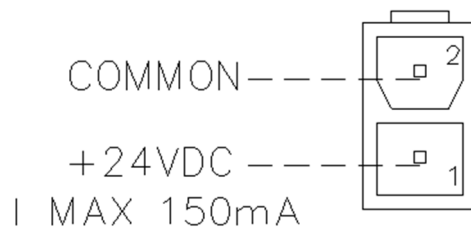


Figure 13 P3 Vent

This is a dedicated 24 Vdc connector to control the optional vent valve.

## Technical Information

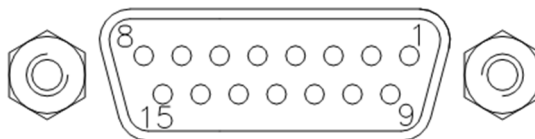
### P5 – External Fan



**Figure 14** External Fan

This is a dedicated 24 Vdc connector to supply the optional external fan.

### J5 – In-Out



**Figure 15** In-Out

This connector carries all the input and output signals to remote control the TwisTorr 305 FS onboard.

It is a 15-pins D type connector; the available signals are detailed in the table, the following paragraphs describe the signal characteristics and use.

## Technical Information

**Table 2** Signal characteristics and use

Pin N.	Signal Name	IN-/OUTPUT
1	Start/Stop (+)	In
2	Start/Stop (-)	In
3	Interlock (+)	In
4	Interlock (-)	In
5	Speed setting (+)	In
6	Speed setting (-)	In
7	Soft start (+)	In
8	Soft start (-)	In
9	+ 24 vdc	Out
10	Spare	
11	Programmable set point (+)	Out
12	Contact: normally open	
13	Fault output	Out
14	Programmable analog signal (+)	Out
15	<ul style="list-style-type: none"><li>• Ground</li><li>• Programmable analog signal (-)</li></ul>	Out

## Signal Description

**Start/Stop:** input signal to start or stop the pump. With the prewired mating connector, supplied with the unit, the interlock (+) signal is connected to the +24Vdc pin and the START/STOP (-) signal to the GROUND pin: in this condition the pump automatically starts as soon as the controller recognises the input supply ("Plug & Pump").

**Interlock:** input signal to control the pump rotation. With the supplied cover connector the interlock (+) signal is connected to the +24 Vdc pin and the interlock (-) signal to the GROUND pin.

**Soft start:** this input is used to provide a "soft start" to the pump; in this condition the ramp-up time could be up to 45 min.

**Speed setting:** PWM input signal to set the pump speed. The PWM signal characteristics must be the following:

- frequency: 100 Hz +/-20%
- amplitude: from 5 to 24 V
- duty cycle range: from 25% to 75% (toff/T) corresponding to a rotational frequency from 800 Hz to 1010 Hz linearly. With duty cycle <25% rotational frequency = 800 Hz, duty cycle >75% rotational frequency = 1010 Hz.

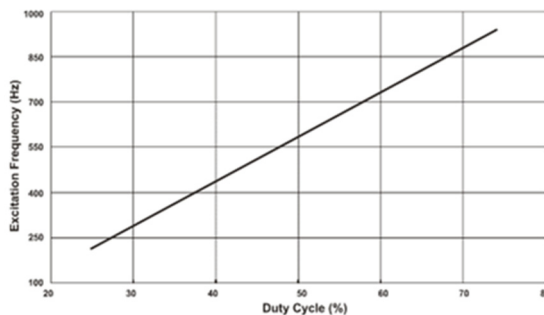


Figure 16

### NOTE

The duty cycle percentage is referred to the low level portion of the PWM signal.

**Programmable analog signal:** this output signal is a voltage (from 0 to 10 Vdc) proportional to a reference quantity (frequency or power) set by the user. The default setting is the frequency (see the following example diagram).

## Technical Information

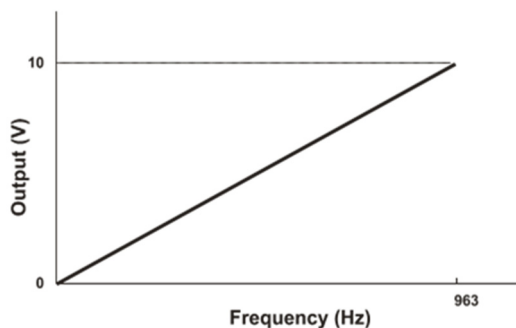


Figure 17

**Fault:** this open collector output signal is ON when a system fault condition is detected.

**Programmable set point:** this open collector output signal is enabled when the reference quantity chosen (frequency, current or time) is higher than the set threshold. The signal can be "high level active" (that is the output is normally at 0 Vdc and becomes 24 Vdc when activated), or "low level active" (that is the output is normally at 24 Vdc and becomes 0 Vdc when activated). Moreover, if the reference quantity is the frequency or the current drawn, it is possible to set the hysteresis (in % of the threshold value) to avoid bouncing.

For example:

- reference quantity: frequency
- threshold: 500 Hz
- hysteresis: 1%
- activation type: "high level"

## Technical Information

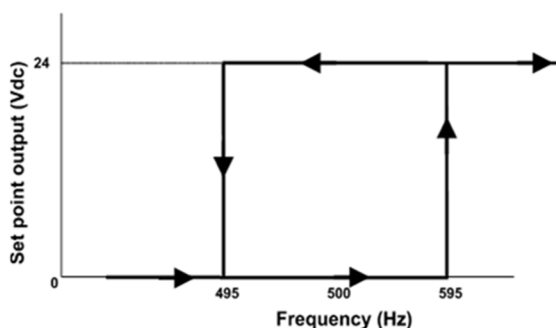


Figure 18 Set point output

The set point output stays at 0 Vdc until the frequency becomes higher than 505 Hz (that is 500 Hz + 1% of 500 Hz), then the output goes at 24 Vdc and stays at 24 Vdc until the frequency becomes lower than 495 Hz (that is 500 Hz – 1% of 500 Hz).

It is possible to delay the set point checking for a programmable delay time.

The PROGRAMMABLE SET POINT signal has the following default settings:

- reference quantity: frequency
- threshold: 909 Hz
- hysteresis: 2 %
- activation type: high level
- delay time: 0 second

### NOTE

The communication S/W between the controller and the pump (A-Plus) allows the operator to set all the programmable feature.

When no external input-output device is available this connector must be closed with the supplied mating connector that short-circuits the START and INTERLOCK inputs with the GROUND input (see the following figure).

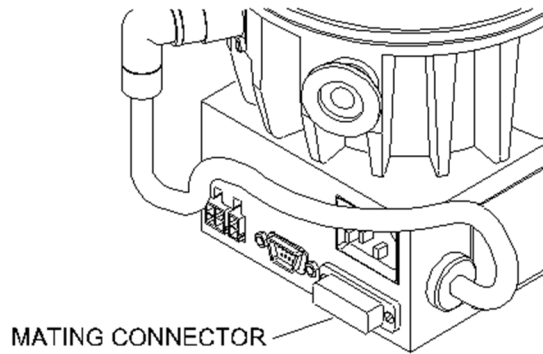


Figure 19

## How to Connect the Open Collector Inputs of the Controller

Here below there are the typical connections of the open collector input of the integrated controller to an external system. Two cases are considered:

- 1 the customer supplies the 24 Vdc
- 2 the customer does not supply the 24 Vdc

Please note that on the connector a 24 Vdc, 60 mA voltage, a GROUND signal and the open collector pin are available.

## Technical Information

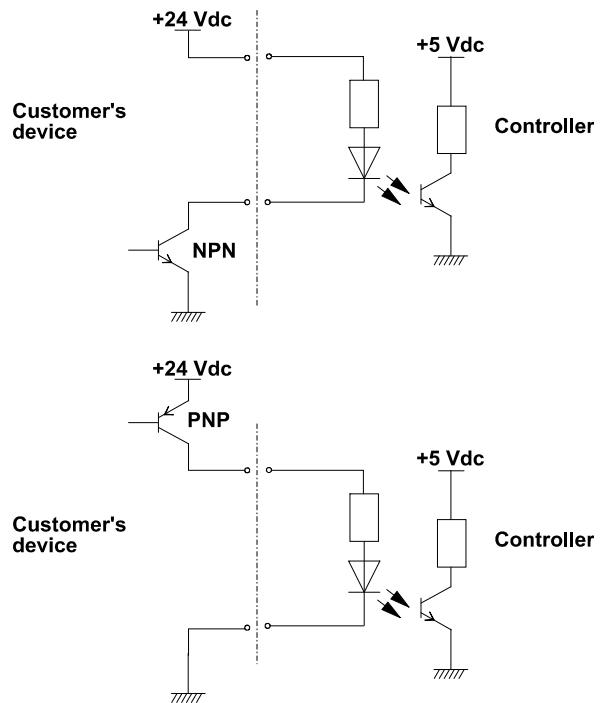


Figure 20 Case 1

## Technical Information

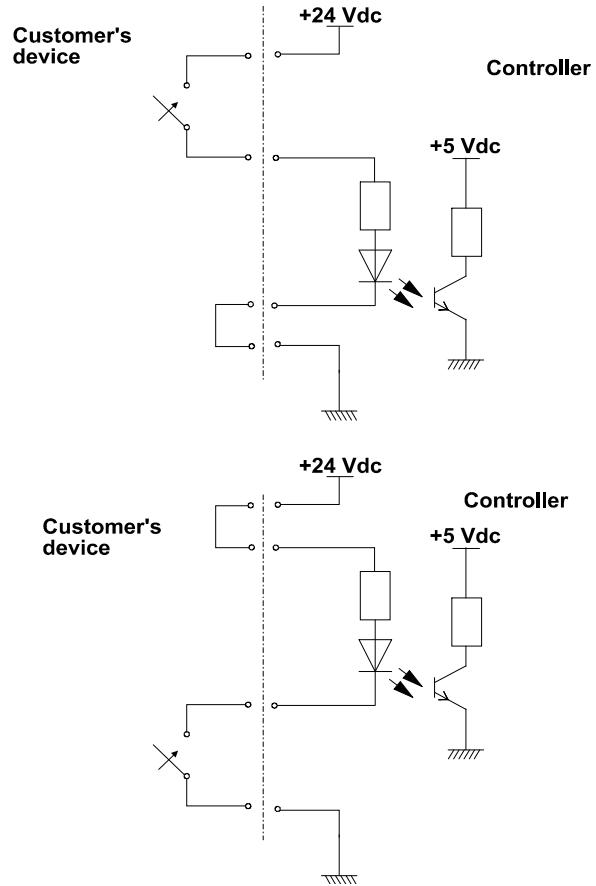


Figure 21 Case 2 with relay utilisation

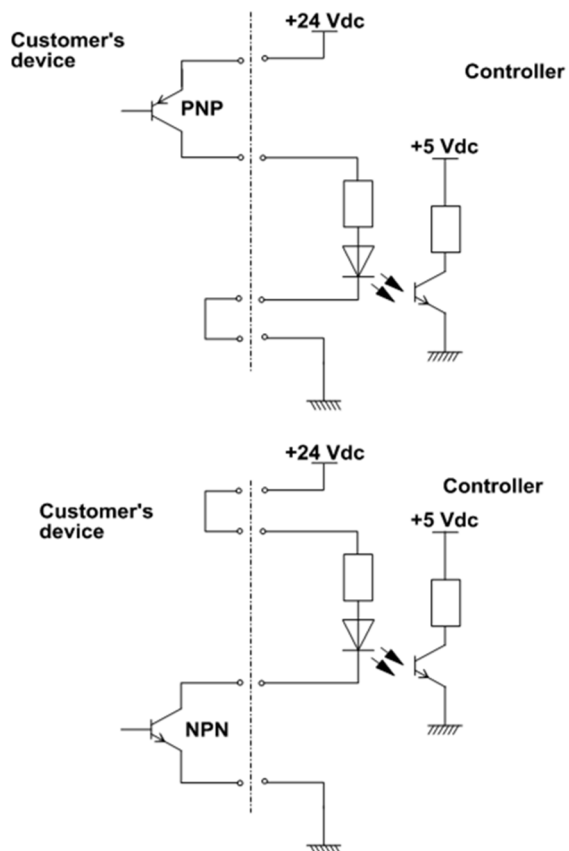


Figure 22 Case 2 with transistor utilisation

## How to Connect the Outputs of the Controller

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. The figure example refers to the programmable set point signal on pins 11 and 9.

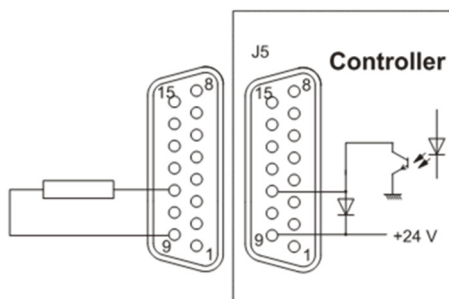


Figure 23

### J6 – Serial

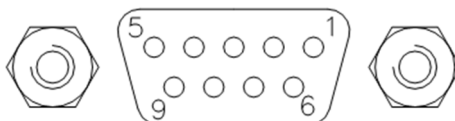


Figure 24 Pin connector

This is a 9 pin D-type serial input/output connector to control via an RS 232 or RS 485 connection the TwisTorr 305 FS onboard.

## Technical Information

**Table 3** Signal characteristics and use

Pin N.	Signal Name
1	+5 V (OUT)
2	TX (RS232)
3	RX (RS232)
4	A+ (RS422 Optional)
5	GND
6	A + (RS485)
7	B - (RS422 Optional)
8	B - (RS485)
9	RESERVED

### NOTE

Note that the vent valve can also be controlled by means of the serial connection.

A serial communication kit with a serial cable and the A-Plus software is available (optional).

## Input power connection on the model X3507-64130

The following figure shows the input power connection relevant to the model n. X3507-64131.

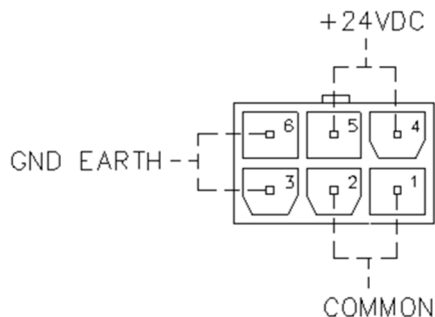


Figure 25

## RS 232/RS 485 communication description

Both the RS 232 and the RS 485 interfaces are available on the connector J6.

The communication protocol is the same (see the structure below), but only the RS 485 manages the address field. Therefore to enable the RS 485 is necessary to select the type of communication as well as the device address by means of the TwisTorr software.

### Communication Format

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

## Technical Information

### 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:

Table 4

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

## Technical Information

- <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>.

The controller can answer with the following response types:

**Table 5**

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 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	

## Technical Information

### Command: SOFT-START (ON)

Source: PC

Destination: Controller

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

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

### Command: SOFT-START (OFF)

Source: PC

Destination: Controller

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

Source: Controller

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

## Window Meanings

Table 6

WIN#	Read/ Write	Type	Meaning	Admitted Values
000	R/W	L	Start/Stop (in remote mode the window is a read only)	Start = 1; Stop = 0 (default = 0)
001	R/W	L	Low Speed Activation command	Active = 1; Non Active = 0 (default = 0)
008	R/W	L	Remote or Serial configuration	Remote = 1; Serial = 0 (default = 1)
100	R/W	L	Soft Start activation command (write only in Stop condition)	Active = 1; Non Active = 0 (default = 0)
101	R/W	N	Set Point type	0 = Frequency; 1 = Current; 2 = Time (default = 0)
102	R/W	N	Set Point threshold (expressed in Hz, mA or s)	(default = 909)
103	R/W	N	Set Point delay: time between the pump start and the set point check (seconds)	0 to 65535 (default = 0)
104	R/W	L	Set Point signal activation type: the signal can be "high level active" or "low level active"	0 = high level active; 1 = low level active (default = 0)
105	R/W	N	Set point hysteresis (in % of threshold)	Set point hysteresis, in % of threshold (default = 2)
108	R/W	N	Baud Rate	0 = 600 bit/s 1 = 1200 bit/s 2 = 2400 bit/s 3 = 4800 bit/s 4 = 9600 bit/s (default = 4)
109	W	L	Pump life/Cycle time/Cycle number reset	To reset the counters, write '1'
110	R/W	L	Interlock type. The interlock can be checked at start or continuously	Impulse = 0 Continuous = 1 (default = 1)
111	R/W	N	Analog output type: output voltage signal proportional to frequency or power	0 = frequency; 1 = power (default = 0)
120	R/W	N	Rotational frequency setting (Hz)	700 to 1010 (default = 1010)
121	R/W	N	Maximum rotational frequency in Hz (active only in Stop condition)	700 to 1010 (default = 1010)

## Technical Information

WIN#	Read/ Write	Type	Meaning	Admitted Values
122	R/W	L	Set vent valve on/off (on = closed)	700 to 1010 (default = 1010)
125	R/W	N	Set the vent valve operation	Automatic = 0 (ref. to note 1) On command = 1 (ref. to note 2) (default = 0)
126	R/W	N	Vent valve opening delay (expressed in 0.2 sec)	0 to 65535 (corresponding to 0 to 13107 sec) (default = 15)
140	R/W	L	Vent Valve selection (N.C./N.O.)	0 = N.C. 1 = N.O. (default = 1)
143	R/W	N	External fan configuration (always on or driven by serial command)	0 = Always On 1 = Serial (default = 0)
144	R/W	L	External Fan Setting (ON/OFF serial command)	On = 1; Off = 0 (default = 0) Not active if W143=0
157	R/W	N	Gas Type Selection	0 = Ar 1 = N2 2 = He (default = 0) Ref. to item 4 of the note
167	R	L	Stop Speed Reading (it reads the pump speed while in stop mode)	On = 1; Off = 0 (default = 0)
200	R	N	Pump current in mA dc	
201	R	N	Pump voltage in Vdc	
202	R	N	Electrical power delivered to the pump in W (pump current x pump voltage duty cycle)	
203	R	N	Driving frequency in Hz	
204	R	N	Pump body temperature n °C	
205	R	N	Pump status	0 = Stop; 1 = Interlock; 2 = Ramp; 3 = Autotuning; 4 = Braking; 5 = Normal; 6 = Fail
206	R	N	Error code	Bit description: refer to the picture which is located on the bottom of this table
300	R	N	Cycle time in minutes (set to zero by a reset command)	0 to 999999
301	R	N	Cycle number (set to zero by a reset command)	0 to 9999
302	R	N	Pump life in hours (set to zero by a reset command)	0 to 999999

## Technical Information

WIN#	Read/ Write	Type	Meaning	Admitted Values
400	R	A	CRC EPROM (QE)	QE8XXXX (where "XXXX" are variable chars)
402	R	A	CRC Param. (PA)	PA8XXXX (where "XXXX" are variable chars)
404	R	A	CRC Parameter structure	"XXXX"
503	R/W	N	RS485 Address	0 to 31 (default = 0)
504	R/W	N	Serial type select	0 = RS 232; 1 = RS 485 (default = 0)
724	R/W	N	Run Up Time (in minutes)	0 to 30 (ref. to note 3) (default = 15)
725	R/W	L	Run Up Time activation command	On = 1; Off = 0 (default = 1)

### NOTE

All the serial windows which are not mentioned in the relevant table are reserved for future use or reserved to Agilent Service.

These windows are not intended to be used by the end user.

## Technical Information

### NOTES

- 1 Automatic means that when the controller stops, the vent valve is opened with a delay defined by window n. 126; when the controller starts, the vent valve is immediately closed.
- 2 On command means that the vent valve is opened or closed by means of window n. 122.
- 3 Run Up Time is the maximum amount of time for which the pump is allowed to spin at a speed value less than the normal operating speed. The pump is not intended to spin for an indefinite time at a speed value lower than 700Hz. If the pump, due to the gas flow applied to it, spins at a speed value less than 700Hz, the Run Up timer starts to count; when the counter value exceeds the Run Up Time set value, the pump will stop and the error condition "Run Up Time" will be issued by the control unit.
- 4 Select the gas type which is appropriate for the customer process in order to achieve the full performance of the pump.

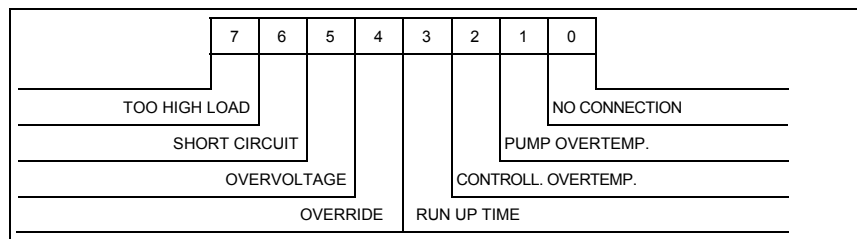
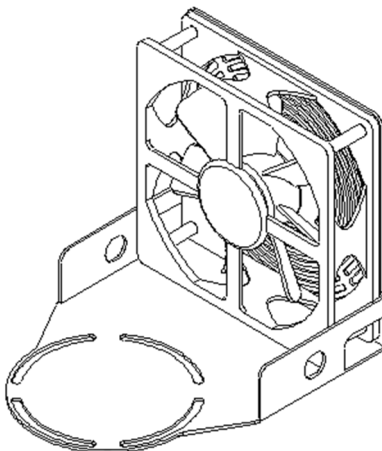


Figure 26 Window N. 206 Bit Description

## Air Cooling Kit Installation

An air cooling kit (mod. X3500-68010) is available to improve the TV 301 cooling during heavy operational conditions (optional).



**Figure 27**

Fan specifications:

- air flow: 147 m<sup>3</sup>/h
- input voltage: 24 Vdc
- dimensions: 119 x 119 x 32 mm
- power: 2.6 W

The fan bracket is shaped so that it can be mounted close to the pump.

In case the pump is located far from the control unit and the control unit is connected to the pump via an extension cable, the cooling kit is already provided with a specific extension cable (which is included in its package) in order to connect it to the controller.

To fix the fan to the TwissTorr 305 FS onboard case execute the following procedure (see the following figure).

## Technical Information

- 1 Fix the fan to the suitable bracket by means of the furnished screws;
- 2 Fix the bracket to the pump body between the pump and the controller;
- 3 Connect the fan supply to the P4 connector of the controller.

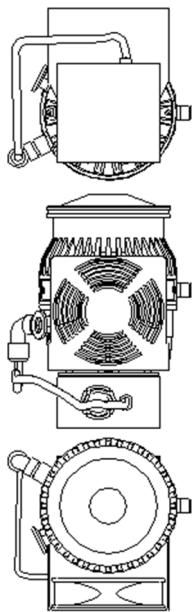


Figure 28

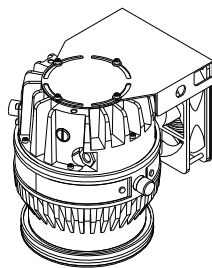


Figure 29

## Technical Information

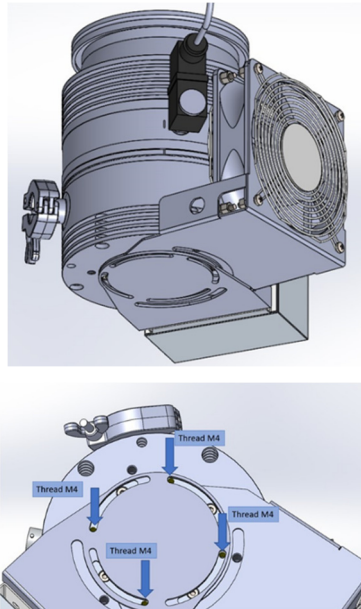


Figure 30

## Use in the Presence of Magnetic Fields

Magnetic fields induce eddy currents in the rotor of a turbomolecular pump that tend to prevent rotation. This causes increased electrical power consumption by the motor, most of which is dissipated in the rotor. In general, therefore, an increase in pump current can be expected.

## Accessories and Spare Parts

**Table 7** Accessories and spare parts

Description	Part Number
Mains cable NEMA Plug, 3m long	9699958
Mains cable European Plug, 3m long	9699957
Serial cable	9699883
RS232 Serial to B/T Adapter	X3514-68003
Air cooling kit for TwisTorr 305 FS Controller	X3500-68010
Vent valve for TwisTorr 305 FS Controller, N.O. orifice 1.2 mm	9699834
Vent valve for TwisTorr 305 FS Controller, N.O. orifice 0.5 mm	9699834M006
Vent valve for TwisTorr 305 FS Controller, N.C. orifice 1.2 mm	9699834M010
Vent valve for TwisTorr 305 FS Controller, N.C. orifice 0.5 mm	9699834M011
L-shaped bracket	X3500-68012

**NOTE**

The communication S/W between the controller and the pump (A-Plus) is available on the Agilent website:

<https://www.agilent.com/en/products/vacuum-technologies/vacuum-leak-detection-software/a-plus-software>

The B/T adapter (X3514-68003) can be used to connect the Vacuum Link app to the controller.

<https://www.agilent.com/en/product/vacuum-technologies/vacuum-leak-detection-software/vacuum-link-app>



## ***Vacuum Products Division***

*Dear Customer,*

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

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

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

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

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

*Sincerely,*

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

**Giampaolo LEVI**

*Vice President and General Manager  
Agilent Vacuum Products Division*

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

## 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.Ili Varian, 54 – 10040 Leini (TO) – Italy

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

NAME	COMPANY	FUNCTION
<p>ADDRESS:</p> <p>TEL. N° : <span style="margin-left: 200px;">FAX N° :</span></p> <p>E-MAIL:</p>		
<p>PROBLEM / SUGGESTION :</p>          <p>REFERENCE INFORMATION (model n°, serial n°, ordering information, time to failure after installation, etc.):</p>          <p style="text-align: right;">DATE</p>		
CORRECTIVE ACTION PLAN / ACTUATION (by AGILENT VPD)		LOG N°

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



**Vacuum Products Division  
Instructions for returning products**

Dear Customer,

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

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

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

**Product preparation**

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

**Shipping**

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

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

**EUROPE:**

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

[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

**NORTH AMERICA:**

Fax: 1 781 860 9252  
Toll Free: 800 882 7426

[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)

**PACIFIC RIM:**

please visit our website for individual  
office information

<http://www.agilent.com>



**TERMS AND CONDITIONS**

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

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



**Vacuum Products Division  
Request for Return Form**

Customer information		
Company :		Contact Name:
Address:		Tel: <input type="text"/> Fax: <input type="text"/>
		Email: <input type="text"/>

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

Type of return
<input type="checkbox"/> Non Billable <input type="checkbox"/> Billable <input checked="" type="checkbox"/> New PO # (hard copy must be submitted with this form): _____ <input type="checkbox"/> Exchange <input type="checkbox"/> Repair <input type="checkbox"/> Upgrade <input type="checkbox"/> Consignment/Demo <input type="checkbox"/> Calibration <input type="checkbox"/> Evaluation <input type="checkbox"/> Return for Credit

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

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

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

**NOTE:** If a product is received at Agilent which is contaminated with a toxic or hazardous material that was not disclosed, **the customer will be held responsible** for all costs incurred to ensure the safe handling of the product, and is liable for any harm or injury to Agilent employees as well as to any third party occurring as a result of exposure to toxic or hazardous materials present in the product.

## Agilent Vacuum Products Division/Sales and Service Offices

### United States

#### **Agilent Technologies**

121 Hartwell Avenue  
Lexington, MA 02421 - USA  
Ph. +1 781 861 7200  
Fax: +1 781 860 5437  
Toll-Free: +1 800 882 7426  
[vpl-customer@agilent.com](mailto:vpl-customer@agilent.com)

### Netherlands

#### **Agilent Technologies Netherlands B.V.**

Customer Contact Center  
Laan van Langerhuize 1, toren A-8  
1186 DS Amstelveen  
Tel. +31 020 547 2600  
Fax +31 020 654 5748  
[customer@agilent.com](mailto:customer@agilent.com)

### Belgium

#### **Agilent Technologies Belgium S.A./N.V.**

Customer Contact Center  
Pegasus Park  
De Kleetlaan 12A bus 12  
B-1831 Diegem  
Tel. +32 2 404 92 22  
Fax +32 2 626 46 30  
[customer@agilent.com](mailto:customer@agilent.com)

### Brazil

#### **Agilent Technologies Brasil**

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

### China

#### **Agilent Technologies (China) Co. Ltd**

No.3, Wang Jing Bei Lu, Chao Yang District  
Beijing, 100102, China  
Tel: +86 (0)10 64397888  
Fax: +86 (0)10 64392765  
Toll free: 400 8206778 (mobile)  
Toll free: 800 8206778 (landline)  
[vacuum.cnmarketing@agilent.com](mailto:vacuum.cnmarketing@agilent.com)  
[vpc-customerservice@agilent.com](mailto:vpc-customerservice@agilent.com)

### France

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

### Southeast Asia

#### **Agilent Technologies Sales Sdn Bhd**

Unit 201, Level 2 uptown 2,  
2 Jalan SS21/37, Damansara Uptown  
47400 Petaling Jaya, Selangor, Malaysia  
Ph. +603 7712 6181  
Fax: +603 7727 1239  
Toll free: 1 800 880 805  
[vps-customerservice@agilent.com](mailto:vps-customerservice@agilent.com)

### India (Sales)

#### **Agilent Technologies India Pvt. Ltd.**

Unit Nos 110- 116, & Part of 101 & 109  
First Floor, Splendor Forum,  
Plot No.-3, District Centre, Jasola  
New Delhi-110025  
Ph. +91 11 4623 7100  
Fax: +91 4623 7105  
Toll Free: 1 800 180 1517

### Italy

#### **Agilent Technologies Italia S.p.A.**

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

### Japan

#### **Agilent Technologies Japan, Ltd.**

9-1 Takakura-cho Hachioji-city,  
Tokyo, Japan  
Tel.: +81- 3-5232-1253  
Fax: +81-120-565-154  
Toll-Free: +81-120-477-111  
[jp-vvt-sales.pdl-ext@agilent.com](mailto:jp-vvt-sales.pdl-ext@agilent.com)

### Singapore

#### **Agilent Technologies Singapore Pte. Ltd.**

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

### Korea

#### **Agilent Technologies Korea, Ltd.**

Ilshin Building 4F  
Yongsan-gu Hannam-daero  
Seoul Korea 04418  
Tel: +82 (0)2 2194 9449  
Fax: +82 (0)2 2194 9853  
Toll free: 080 222 2452  
[vpc-customerservice@agilent.com](mailto:vpc-customerservice@agilent.com)

### UK and Ireland

#### **Agilent Technologies LDA UK Limited**

Lakeside Cheadle Royal Business Park  
Cheadle, Cheshire SK8 3GR,  
United Kingdom  
Ph. +44 01865291570  
Fax +44 01865291571  
Toll free: 00 800 234 234 00  
Toll free fax: 00 800 345 345 00  
[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

© Agilent Technologies, Inc. 2021  
02/2021

### India (Service)

Agilent Technologies India Pvt. Ltd.  
C-Block, RMZ Centennial Plot Number- 8A, 8B, 8C,  
8D,  
Doddanakundi Industrial Area, ITPL Road,  
Mahadevapura Post, Bangalore- 560048  
Ph. +91 80 40614000  
Fax: +91 80 40148991

### Taiwan

#### **Agilent Technologies Taiwan Limited**

No. 20 Gao-shuang Road, Ping-zhen Dist  
Tao-Yuan City  
32450 Taiwan, R.O.C.  
Tel: +886 3 4959004  
Toll free: 0800 018 768  
[vpw-customerservice@agilent.com](mailto:vpw-customerservice@agilent.com)

### Germany and Austria

#### **Agilent Technologies Sales & Services GmbH & Co. KG**

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

### Mexico

#### **Agilent Technologies**

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

### Other Countries

#### **Agilent Technologies Italia S.p.A.**

Via F.lli Varian, 54  
10040 Leini, (Torino) - Italy  
Tel.: +39 011 997 9111  
Fax: +39 011 997 9350  
Toll-Free: 00 800 234 234 00  
[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

### Customer Support & Service

#### **NORTH AMERICA:**

Toll Free: 800 882 7426  
[vpl-ra@agilent.com](mailto:vpl-ra@agilent.com)  
[Lexington-service@agilent.com](mailto:Lexington-service@agilent.com)

#### **EUROPE:**

Toll Free: 00 800 234 234 00  
[vpt-customer@agilent.com](mailto:vpt-customer@agilent.com)

#### **PACIFIC RIM:**

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

#### **Worldwide Web Site, Catalog and Order On- line:**

[www.agilent.com/chem/vacuum](http://www.agilent.com/chem/vacuum)  
Representatives in most countries

## In This Book

The manual describes the following:

- Istruzioni per l'uso
- Bedienungshandbuch
- Notice de mode d'emploi
- Manual de instrucciones
- 用户手册
- ユーザーマニュアル
- Instruction for Use
- Technical information

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies, Inc. 2022

Edition A.00, 06/22



87-901-060-01

