

Agilent IDP-4 Dry Scroll Pump

User Manual



Notices

Manual Part Number

D0129654

Edition A.00, November 2025

Copyright

© Agilent Technologies, Inc. 2025

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

Instrument Manufacturing

Agilent Technologies LDA Malaysia
Sdn Bhd Bayan Lepas
Free Industrial Zone 11900 Bayan Lepas,
Penang Malaysia

Printed in Italy

Warranty

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

Technology Licenses

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

Restricted Rights Legend

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

1 Istruzioni per l'uso	13
Informazioni su questo manuale	14
Documenti pertinenti rilevanti	14
Definizioni e terminologia	15
Definizione di Attenzione, Avvertenza e Nota	15
Simboli di avvertenza	16
Sicurezza	17
Avvertenze EMC	18
FCC	18
Utilizzo conforme	19
Uso improprio ragionevolmente prevedibile	19
Indicazioni generali sulla sicurezza per pompa per vuoto a secco Agilent	20
Personale qualificato	21
Equipaggiamento protettivo	22
Descrizione del prodotto	24
Trasporto e immagazzinaggio	24
Disimballaggio e ispezione	25
Preparazione per l'installazione	26
Procedura di avvio	30
Procedura di spegnimento	33
Smaltimento	34
Condizioni ambientali, conformità e requisiti strutturali	35

Contents

Compatibilità elettromagnetica	35
EN55011/CISPR11	35
Dichiarazione EMC di Classe A della Corea del Sud	36
ICES/NMB-001	36
Servizio Post Vendita	37
2 Betriebsanleitung	38
Informationen über diese Betriebsanleitung	39
Einschlägige unterstützende Dokumente	39
Definitionen und Terminologie	40
Definition von Vorsicht, Warnung und Hinweis	40
Warnsymbole	41
Sicherheit	42
EMV-Warnungen	43
FCC	43
Ordnungsgemäße Nutzung	44
Vernünftigerweise vorhersehbarer Missbrauch	44
Sicherheitsrichtlinien für trockene Scrollpumpe	45
Qualifiziertes Personal	46
Schutzausrüstung	47
Produktbeschreibung	49
Transport und Lagerung	49
Auspacken und Inspektion	50
Vorbereitung zum Einbau	51
Inbetriebnahme	55
Abschaltverfahren	58
Entsorgung	59

Contents

Umweltbedingungen, Konformität und Versorgungsanforderungen	60
Elektromagnetische Verträglichkeit	60
EN55011/CISPR11	60
Südkoreanische Klasse A EMV-Erklärung	61
ICES/NMB-001	61
Service	62
3 Mode d'emploi	63
À propos de ce manuel	64
Documents complémentaires pertinents	64
Définitions et terminologie	65
Signification des messages d'attention, d'avertissement et des notes	65
Symboles d'avertissement	66
Sécurité	67
Avertissement CEM	68
FCC	68
Utilisation appropriée	69
Utilisation inappropriée	69
Directive de sécurité pour la pompe à vide sèches	70
Personnel qualifié	71
Équipements de protection	72
Description du produit	74
Transport et entreposage	74
Déballage et inspection	75
Préparation pour l'installation	76
Procédure de démarrage	80

Contents

Procédure d'arrêt	83
Élimination	84
Conditions environnementales, conformité et applications pratiques	85
Compatibilité électromagnétique	85
EN55011/CISPR11	85
Déclaration CEM Classe A Corée du Sud	86
ICES/NMB-001	86
Service	87
4 Instrucciones de uso	88
Acerca del manual	89
Documentos justificativos pertinentes	89
Definiciones y terminología	90
Definición de Precaución, Advertencia y Nota	90
Símbolos de advertencia	91
Seguridad	92
Advertencias CEM	93
FCC	93
Uso adecuado	94
Uso incorrecto razonablemente previsible	94
Directrices de seguridad para bombas de vacío de desplazamiento en seco	95
Personal cualificado	96
Equipo de protección	97
Descripción del producto	99
Transporte y almacenamiento	99
Desembalaje e Inspección	100

Contents

Preparación para la instalación	101
Procedimiento de puesta en marcha	105
Procedimiento de parada	108
Eliminación	109
Condiciones medioambientales, conformidad y requisitos de los servicios públicos	110
Compatibilidad electromagnética	110
EN55011/CISPR11	110
Declaración EMC Clase A de Corea del Sur	111
CIEM/NMB-001	111
Servicio	112
5 使用说明	113
關於本手冊	114
相關證明文件	114
定義與術語	115
小心、警告和注意的定義	115
警告符號	116
安全	117
EMC 警告	118
美國聯邦通訊委員會	118
正確使用	119
合理而可預見的濫用	119
乾式渦旋幫浦安全指南	120
合資格人員	121
防護裝備	122

Contents

產品描述	124
運輸及儲存	124
開箱檢驗	125
安裝準備	126
啟動程式	130
關機程式	133
處置	134
環境條件、合規性和公用設施要求	135
電磁兼容	135
EN55011/CISPR11	135
韓國 A 類 EMC 聲明	136
ICES/NMB-001	136
服務	137
6 取扱説明書	138
この説明書について	139
関連補足資料	139
定義と用語	140
警告・注意・注の定義	140
警告の記号	141
安全性	142
EMC 警告	143
FCC	143

Contents

適切な使用	144
合理的に予見可能な誤使用	144
ドライスクロールポンプの安全ガイドライン	145
有資格者	146
保護装備	147
製品説明	149
輸送と保管	149
開梱と点検	150
設置の準備	151
接地方法	154
起動手順	155
停止手順	158
処分	159
環境条件、コンプライアンス、ユーティリティ要件	160
電磁両立性	160
EN55011/CISPR11	160
韓国クラス A EMC 宣言	161
ICES/NMB-001	161
サービス	162

Contents

7	Instructions for Use	163
	About this Manual	164
	Relevant Supporting Documents	164
	Definitions and terminology	165
	Definition of Caution, Warning and Note	165
	Warning Symbols	166
	Safety	167
	EMC Warnings	167
	FCC	167
	Proper use	168
	Reasonably Foreseeable Misuse	168
	Safety guideline for Dry Scroll Pump	169
	Qualified Personnel	170
	Protective equipment	171
	Product Description	173
	Transport & Storage	173
	Unpacking and Inspection	174
	Preparation for Installation	175
	Startup Procedure	179
	Shutdown Procedure	182
	Disposal	183
	Environmental Conditions, Compliance and Utility Requirements	184
	Electromagnetic Compatibility	184
	South Korean Class A EMC declaration	185
	ICES/NMB-001	185
	Service	186

Contents

8	Technical Information	187
	Description of the IDP-4 Dry Scroll Pump	188
	Technical Specifications	189
	IDP-4 Dry Scroll Pump Outline Drawing	191
	Performance Curves	193
	Electrical Connections	195
	Grounding Instructions	195
	Mechanical Connections	197
	Gas Ballast	199
	Accessories and Spare Parts	201
	Optional Accessories	201
	Optional Inlet Valve Kit	202
	Installation	204
	Cleaning and Preventive Maintenance	205
	External cleaning	205
	Internal cleaning as part of a Preventive Maintenance plan	206
	Verifying Performance	206
	Tip Seal Replacement	207
	Tip Seal Replacement Instructions	208
	User Maintenance and Repair	216
	Factory-Only Service	217

Contents

Service Options	218
Troubleshooting	219
Serial Connector & Pin allocation	221
Pump Serial Connector	221
Serial communication settings	222
Window Description and Protocol	223
Window Description	223
Windows Protocol	226

1 Istruzioni per l'uso

Informazioni su questo manuale	14
Documenti pertinenti rilevanti	14
Definizioni e terminologia	15
Definizione di Attenzione, Avvertenza e Nota	15
Simboli di avvertenza	16
Sicurezza	17
Avvertenze EMC	18
FCC	18
Utilizzo conforme	19
Uso improprio ragionevolmente prevedibile	19
Indicazioni generali sulla sicurezza per pompa per vuoto a secco Agilent	20
Personale qualificato	21
Equipaggiamento protettivo	22
Descrizione del prodotto	24
Trasporto e immagazzinaggio	24
Disimballaggio e ispezione	25
Preparazione per l'installazione	26
Procedura di avvio	30
Procedura di spegnimento	33
Smaltimento	34
Condizioni ambientali, conformità e requisiti strutturali	35
Compatibilità elettromagnetica	35
EN55011/CISPR11	35
Dichiarazione EMC di Classe A della Corea del Sud	36
ICES/NMB-001	36
Servizio Post Vendita	37

Informazioni su questo manuale

Il presente manuale contiene istruzioni per gli utenti e informazioni tecniche relative alla pompa per vuoto a secco IDP-4 Dry Scroll in conformità alle direttive applicabili.

Il presente documento è valido per la versione di prodotto valida al momento della pubblicazione, a condizione che il cliente o terzi non apportino modifiche al prodotto. In caso di informazioni non corrispondenti, la documentazione fornita con il prodotto deve essere considerata come documento valido.

NOTA

- 1** Il presente manuale contiene informazioni importanti che consentono a tutto il personale che opera con la pompa per vuoto a secco IDP-4 Dry Scroll di utilizzarle in modo sicuro e di garantirne le prestazioni per l'intera vita utile.
- 2** Conservare questo manuale, insieme a tutte le pubblicazioni ad esso correlate, in un luogo accessibile, conosciuto da tutti gli operatori/personale di manutenzione.

Documenti pertinenti rilevanti

X3802-90000 Dichiarazione di conformità

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 è riportato un elenco di simboli che sono visualizzati insieme alle avvertenze sulla pompa per vuoto a secco IDP-4 Dry Scroll. Viene descritto anche il pericolo correlato. Un simbolo triangolare indica un'avvertenza. I significati dei simboli che possono apparire accanto alle avvertenze nella documentazione o sul prodotto stesso sono riportati qui di seguito. Il seguente simbolo può essere usato sulle etichette di avvertenza applicate sul prodotto. Quando viene visualizzato questo simbolo, consultare il relativo manuale operativo o di servizio per la procedura corretta a cui fa riferimento l'etichetta di avvertenza.



	Certificazione CE
	Certificazione CSA
	Certificazione RoHS Cina
	Rifiuti di Apparecchiature Elettriche ed Elettroniche
	Certificazione RCM
	Sito di produzione
	Presenza di superfici surriscaldate
	Presenza di tensioni pericolose
	Pericolo generico
	Messa a terra protettiva

Istruzioni per l'uso

	Corrente alternata (CA)
	Accensione
	Spegnimento

Sicurezza

Questa sezione contiene le informazioni prescritte dalla direttiva sulla bassa tensione 2006/42/EC, che sono essenziali per la conformità e il rispetto delle norme di sicurezza in generale e in relazione all'uso specifico del prodotto.

La mancata osservanza di queste e delle altre istruzioni contenute nel presente manuale può compromettere la sicurezza 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 dall'inosservanza delle norme di sicurezza indicate nella documentazione tecnica.

Istruzioni per l'uso

Avvertenze EMC

Avvertenza EN 55011 Classe A

Questo prodotto appartiene alla Classe A. L'utilizzo di questo prodotto in ambiente domestico potrebbe causare interferenze radio; in tal caso, l'utente deve adottare misure adeguate.

FCC

Questa apparecchiatura è conforme alla Parte 15 delle disposizioni FCC. L'utilizzo è soggetto alle seguenti due condizioni: (1) l'apparecchio non deve causare interferenze nocive e (2) l'apparecchio deve accettare le interferenze ricevute, comprese le interferenze che potrebbero causare problemi di funzionamento.

NOTA

Questo apparecchio è stato testato e riscontrato idoneo ai limiti previsti per i dispositivi digitali di Classe A, conformemente alla parte 15 delle disposizioni FCC. Tali limiti sono adatti a fornire un'adeguata protezione contro le interferenze nocive se l'apparecchio viene utilizzato in ambienti commerciali. Questo apparecchio genera, utilizza e può irradiare energia a frequenza radio e pertanto, se non correttamente installato e utilizzato secondo le istruzioni del manuale, potrebbe causare interferenze nocive per le comunicazioni radio. L'utilizzo del presente apparecchio in zone residenziali potrebbe causare interferenze nocive per le comunicazioni radio e, in tal caso, l'utente deve provvedere a correggere l'interferenza a proprie spese.

Istruzioni per l'uso

Utilizzo conforme

La pompa per vuoto IDP-4 Dry Scroll è destinata esclusivamente all'uso da parte di personale specificatamente qualificato, come indicato nel presente manuale.

Tutti gli accessori e i dispositivi utilizzati con il prodotto devono essere forniti o approvati da Agilent Technologies.

Agilent Technologies non fornisce alcuna garanzia né si assume alcuna responsabilità per applicazioni diverse da quelle descritte nel presente manuale o in cui non sono rispettati i requisiti essenziali e le misure di sicurezza.

Uso improprio ragionevolmente prevedibile

Agilent Technologies declina qualsiasi responsabilità derivante dall'uso improprio e da prevedibili utilizzi non conformi della pompa per vuoto a secco IDP-4 Dry Scroll. L'uso improprio comporta la decadenza di tutte le richieste di assunzione di responsabilità e di garanzia.

L'uso improprio può riguardare, ma non limitatamente:

- collegamento della pompa alla tensione di alimentazione errata
- collegamento a massa errato o mancante
- funzionamento continuo al di sopra della pressione di ingresso consigliata
- funzionamento con sovrappressione all'ingresso della zavorra di gas
- utilizzo della zavorra di gas aperta per l'atmosfera durante il pompaggio di gas pericolosi
- funzionamento a temperatura ambiente eccessivamente elevata
- funzionamento con temperatura ambiente inferiore alle specifiche
- danni dovuti a stoccaggio e trasporto impropri della pompa
- esecuzione di manutenzione o rimozione delle coperture senza scollegare l'alimentazione di rete
- funzionamento senza elementi di flusso d'aria (copertura o ventola)
- restituzione del prodotto in imballaggi non forniti dalla fabbrica.

Indicazioni generali sulla sicurezza per pompa per vuoto a secco Agilent

Il personale responsabile del funzionamento e della manutenzione della pompa deve essere adeguatamente addestrato e deve conoscere le norme per la prevenzione degli incidenti (fare riferimento alla sezione Personale qualificato). Le precauzioni per la prevenzione degli incidenti contenute nel presente manuale devono essere rispettate costantemente durante il funzionamento e la manutenzione della pompa per evitare lesioni agli operatori e danni alla pompa o ad altre apparecchiature. Queste precauzioni sono fornite sotto forma di AVVERTENZE e di NOTE nell'intero manuale.

AVVERTENZA



Per evitare danni all'apparecchiatura e lesioni al personale operativo, attenersi scrupolosamente alle istruzioni di installazione riportate in questo manuale!

AVVERTENZA



**La pompa non deve funzionare con materiali infiammabili.
È vietato installare la pompa in un ambiente esplosivo.**

Personale qualificato

L'utilizzo della pompa per vuoto a secco IDP-4 Dry Scroll è raccomandato solo da parte di personale qualificato. Per persona qualificata si intende:

- una persona in possesso delle qualifiche appropriate, che abbia fatto esperienza o che abbia completato una formazione su come lavorare con la pompa per vuoto a secco Agilent
- una persona che adotta le misure di sicurezza necessarie e aziona la pompa in condizioni adeguate per non causare lesioni o danni
- una persona che sia ben informata e adeguatamente qualificata per seguire gli standard e le normative nazionali e locali applicabili in materia di sicurezza.

NOTA

Eventuali riparazioni su questo prodotto, oltre a quelle elencate come manutenzione standard nel presente manuale per l'utente, devono essere eseguite esclusivamente da personale autorizzato Agilent. Qualsiasi operazione di regolazione, modifica o manutenzione deve essere eseguita da un tecnico professionista informato sui rischi correlati.

Equipaggiamento protettivo

I dispositivi di protezione degli operatori che utilizzano il prodotto o eseguono la manutenzione della pompa devono essere sempre adeguati al tipo di attività da effettuare. I dispositivi di protezione utilizzati devono essere conformi ai requisiti di sicurezza secondo la normativa vigente nel paese in cui è utilizzata la macchina.

In generale l'operatore deve indossare scarpe di sicurezza durante lo spostamento, la manipolazione o l'installazione della pompa per vuoto a secco IDP-4 Dry Scroll.

AVVERTENZA



Pericolo per la salute a causa di sostanze pericolose durante la manutenzione o l'installazione.



- A seconda della peculiarità del processo, la pompa per vuoto, i componenti o i fluidi di funzionamento possono essere contaminati da sostanze tossiche, reattive o radioattive.



- Indossare attrezzatura di protezione adeguata durante la manutenzione e le riparazioni o in caso di reinstallazione.
-

AVVERTENZA

Rischio di lesioni dovute alla caduta di oggetti



Durante il trasporto manuale della pompa per vuoto, sussiste il pericolo che i carichi scivolino e cadano.



- Trasportare pompe per vuoto di piccole e medie dimensioni a due mani.
- Qualsiasi attrezzatura più pesante di 20 kg deve essere trasportata utilizzando un dispositivo di sollevamento adatto.



- Indossare scarpe antinfortunistiche con puntale in acciaio secondo la direttiva EN 347.
-

ATTENZIONE



Rischio di lesioni a causa di spigoli vivi

- Prima di riparare la pompa o prima di qualsiasi azione di montaggio/smontaggio della pompa nel sistema, attendere l'arresto completo della pompa.
 - Non operare direttamente all'interno della pompa.
 - Se necessario, indossare guanti protettivi secondo norma EN 420.
-

Descrizione del prodotto

La IDP-4 è una pompa Dry Scroll ermetica. Questa pompa è adatta al pompaggio di aria o gas inerti da un'area di bassa pressione a un'area di maggiore pressione, che non sia sostanzialmente né inferiore né superiore alla pressione atmosferica. Questa pompa non è destinata a pompare gas corrosivi, piroforici, particolati o miscele di gas infiammabili e/o combustibili e non è destinata a essere utilizzata come compressore per ottenere pressioni di scarico sostanzialmente superiori alla pressione atmosferica.

Il manuale e i documenti correlati contengono le informazioni necessarie per garantire la sicurezza dell'operatore durante l'uso dell'apparecchiatura. Informazioni dettagliate sono fornite nella sezione "Informazioni tecniche".

Trasporto e immagazzinaggio

Durante il trasporto e l'immagazzinaggio della pompa non superare i seguenti requisiti ambientali:

- Temperatura: da -20 °C a +60 °C (da -4 °F a 140 °F)
- Umidità relativa: dal 0 al 95% (senza condensa)

NOTA

È necessario considerare che durante il trasporto e lo stoccaggio della pompa per vuoto, il vapore acqueo proveniente dall'ambiente penetra all'interno della pompa indipendentemente dalla sua tenuta stagna. Per ottenere le massime prestazioni di vuoto nominale della pompa, questo vapore acqueo deve essere rimosso utilizzando le funzioni di zavorra di gas descritte nella sezione Informazioni tecniche.

Disimballaggio e ispezione

La pompa è fornita in uno speciale imballaggio protettivo. Se la confezione mostrasse segni di danni che possono essersi verificati durante il trasporto, contattare l'ufficio vendite locale. Il peso totale della pompa e dell'imballaggio è di circa 10 kg (20 libbre).

AVVERTENZA



In caso di intervento presso un centro di assistenza autorizzato Agilent, è necessario riporre la pompa nel contenitore con tutti i materiali di imballaggio. La conservazione degli imballaggi facilita qualsiasi richiesta di assistenza.

AVVERTENZA



Quando si disimballa la pompa, assicurarsi di non farla cadere ed evitare urti improvvisi o scossoni.

Preparazione per l'installazione

La pompa IDP-4 Dry Scroll è progettata e certificata per il funzionamento all'interno (categoria di installazione II) in ambienti relativamente puliti e asciutti (grado di imbrattamento 2).

Per le specifiche fare riferimento alla Tabella 1 nella sezione Informazioni tecniche. Queste pompe non sono progettate o certificate per il funzionamento all'aperto.

La pompa IDP-4 Dry Scroll non deve mai essere utilizzata in ambienti contenenti materiali esplosivi o combustibili.

La pompa IDP-4 Dry Scroll NON è intrinsecamente sicura e non deve mai essere utilizzata con gas corrosivi, piroforici, particolati o miscele di gas infiammabili e/o combustibili.

Collocare la pompa su una superficie solida e piana. Se necessario, posarla su un banco o su un tavolo, assicurandosi che è in grado di sostenere il peso della pompa e di tutte le attrezzature associate.

Collegare la pompa all'alimentazione elettrica utilizzando un cavo di alimentazione di tipo IEC-320 con una capacità di almeno 10 A.

NOTA

La causa più comune di prestazioni scarse durante l'installazione iniziale e l'avvio è il vapore acqueo assorbito sia nella pompa che nella camera da vuoto. Far funzionare la pompa con il gas ballast attivato può essere utile se la pompa è rimasta spenta per un certo periodo, permettendo al vapore acqueo di condensarsi all'interno della pompa e del sistema da vuoto. A seconda del sistema da vuoto e della quantità di vapore acqueo assorbito, possono essere necessari da circa 30 minuti fino a circa 24 ore di attivazione del gas ballast per raggiungere le prestazioni di vuoto ottimali.

ATTENZIONE

Durante il funzionamento devono essere rispettate le seguenti condizioni ambientali:

- **Temperatura:** da +5 °C a +40 °C (da 41 °F a 104 °F)
- **Umidità relativa:** dal 5 al 95% (senza condensa)

Collegare la pompa all'alimentazione elettrica utilizzando un cavo IEC-320 adatto al valore e con una tensione idonea (vedere Tabella 2).

Istruzioni per l'uso

ATTENZIONE Quando si installa la pompa per vuoto, si consiglia di posizionarla in modo tale che l'interruttore elettrico e il collegamento del cavo di alimentazione siano in una posizione facilmente accessibile per l'operatore, al fine di potere scollegare rapidamente la pompa dall'alimentazione, se necessario.

ATTENZIONE Non serrare, allentare o modificare in alcun modo i due bulloni a testa esagonale su ciascun lato della testa della pompa (vedere la figura seguente). Ciò causerà una perdita di prestazioni e/o danni alla pompa.



Figura 1

AVVERTENZA Non rimuovere o modificare alcun dispositivo di sicurezza o di isolamento dalla pompa.
In caso contrario, potrebbero insorgere seri pericoli e la garanzia potrebbe non essere più valida.



AVVERTENZA



- Questa pompa è in grado di pompare e scaricare solo aria e gas inerti. Non deve mai essere utilizzata con gas corrosivi, piroforici, particolati o miscele di gas infiammabili e/o combustibili.
 - Installare la pompa in un'area non esposta a pioggia, vapore o umidità eccessiva. Gli agenti atmosferici possono causare scosse elettriche, cortocircuiti e gravi lesioni personali.
 - Prima di procedere a ispezione, riconfigurazione o manutenzione della pompa accertarsi che l'alimentazione elettrica sia scollegata.
 - La zavorra di gas deve essere sigillata ogni volta in cui si pompano gas che non possono essere scaricati nell'atmosfera o miscelati con l'aria.
 - È disponibile un silenziatore opzionale per le applicazioni in cui è auspicabile ridurre il livello sonoro prodotto dalla pompa.
 - Per escludere l'eventualità di lesioni personali, evitare di esporre qualsiasi parte del corpo alla formazione del vuoto.
 - La pompa deve essere installata in modo che sia facile l'accesso all'interruttore di alimentazione.
 - I cavi di alimentazione approvati da Agilent per tutti i collegamenti comuni sono indicati nella sezione Collegamenti elettrici. Se fosse necessario utilizzare un cavo di alimentazione diverso, la corrente nominale del cavo di alimentazione deve essere di almeno 10A per l'impiego a 200-240V e di almeno 15A per l'impiego a 100-127V.
-

Istruzioni per l'uso

Istruzioni per la messa a terra

AVVERTENZA



Per questo prodotto deve essere effettuata la messa a terra. In caso di cortocircuito, la messa a terra riduce il rischio di scosse elettriche perché rappresenta una via di scarico della corrente elettrica. Questa pompa è dotata di un cavo di alimentazione con un connettore di messa a terra appropriato. La spina deve essere inserita in una presa correttamente installata e dotata di messa a terra in conformità a tutte le normative e le ordinanze locali.

Il connettore del cavo di alimentazione deve essere correttamente inserito nel modulo di alimentazione della pompa per evitare collegamenti difettosi o intermittenti.

ATTENZIONE

Sebbene la pompa sia in grado di pompare il particolato normalmente presente nell'atmosfera, non è progettata per processare particelle solide, prodotti chimici, polveri, solventi, condensati, liquidi o altre particelle. Questi elementi possono danneggiare l'apparecchiatura e comprometterne le prestazioni o ridurre la durata.

Si consiglia di mantenere uno spazio libero di almeno 20 cm intorno al sistema per consentire una corretta circolazione dell'aria.

ATTENZIONE

La superficie esterna della pompa deve essere pulita esclusivamente con un detergente delicato.

La pompa per vuoto a secco non contiene liquidi per eliminare polvere e detriti accumulati. Fare funzionare periodicamente la pompa con l'atmosfera per un minuto o due per detergerla.

AVVERTENZA



Verificare con un elettricista qualificato o con un addetto alla manutenzione se le istruzioni di messa a terra non fossero state completamente comprese oppure se sussistesse il dubbio che il prodotto non sia dotato di messa a terra adeguata. Non modificare la spina in dotazione. Se non si adatta alla presa, fare installare una presa adatta da un elettricista qualificato.

Collegare il prodotto solo a una presa con la stessa configurazione della spina. Non utilizzare un adattatore per questo prodotto.

Procedura di avvio

AVVERTENZA



Rimuovere i coperchi delle flange dallo scarico e dall'ingresso.



La procedura di avviamento della pompa per vuoto IDP-4 è la seguente:

1. Montare la pompa nella posizione prevista ed effettuare i collegamenti meccanici e del vuoto appropriati oppure, in alternativa, coprire l'ingresso con una flangia cieca.
2. Assicurarci che il raccordo di scarico non sia ostruito né chiuso con una flangia cieca. Se viene utilizzato un tubo di scarico o un silenziatore, le restrizioni allo scarico non devono causare una contropressione superiore a TBD 6,5 psig.
3. Assicurarci che l'interruttore ON/OFF sia nella posizione OFF '0'.
4. Collegare il cavo di alimentazione appropriato ed inserirlo nella presa di corrente. La corrente nominale del cavo di alimentazione deve essere di almeno 10A per un servizio a 200-240V e di almeno 15A per un servizio a 100-127V.
5. Premere l'interruttore principale ON/OFF sulla posizione "I".
6. Se la pompa è dotata di una valvola d'ingresso opzionale, la valvola si aprirà circa 10 secondi dopo l'accensione.

Potrebbero essere necessarie diverse ore affinché la pompa raggiunga la pressione finale specificata. Se la pompa non raggiunge la pressione finale prevista, ciò potrebbe essere dovuto all'accumulo di vapore acqueo all'interno della pompa. Per rimuovere il vapore acqueo, aprire il gas ballast e lasciare la pompa in funzione con il gas ballast aperto per alcune ore o durante la notte.

Istruzioni per l'uso

NOTA

La causa più comune di scarse prestazioni durante l'installazione e l'avviamento iniziali è il vapore acqueo adsorbito sia nella pompa che nella camera a vuoto. Far funzionare la pompa con il gas ballast attivato può essere utile anche se la pompa è rimasta spenta per un periodo di tempo, consentendo all'acqua di adsorbirsi sia all'interno della pompa che nel sistema a vuoto. A seconda del sistema a vuoto e della quantità di acqua adsorbita, in genere possono essere necessarie da circa 30 minuti a circa 24 ore di attivazione del gas ballast per raggiungere le massime prestazioni di vuoto.



Figura 2

NOTA

L'interruttore ON/OFF della pompa è un interruttore a bilanciere dotato di simboli conformi alla Pubblicazione IEC 417 per rappresentare le posizioni ON e OFF. La figura precedente mostra un interruttore in posizione OFF.

AVVERTENZA



Non inserire oggetti estranei attraverso le prese d'aria di raffreddamento, poiché il contatto con le pale della ventola può causare gravi lesioni personali o danni alla pompa.

AVVERTENZA



Non ostruire le prese d'aria di raffreddamento. L'ostruzione di questi fori di ventilazione può causare il surriscaldamento della pompa. Una temperatura superficiale della pompa superiore a 65 °C (150 °F) è potenzialmente dannosa. Se si osservano tali condizioni, spegnere la pompa e lasciarla raffreddare. Smontare, ispezionare la pompa per individuare eventuali danni e, se necessario, ripararla.

AVVERTENZA



Non utilizzare zavorra di gas se si utilizzano materiali potenzialmente pericolosi o che non devono essere miscelati con l'aria.

Procedura di spegnimento

Per arrestare manualmente la pompa in qualsiasi modalità di funzionamento, portare l'interruttore principale ON/OFF in posizione "0". Se la pompa è dotata di una valvola di isolamento integrata, questa si chiude rapidamente per eliminare il riflusso di gas dalla pompa ai componenti a monte.

ATTENZIONE Prima di spegnere la pompa IDP-4, assicurarsi che tutti i componenti interconnessi a monte siano opportunamente valvolati o preparati per una potenziale esposizione alla pressione atmosferica.

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 che deve essere smaltito con un apposito sistema di raccolta differenziata. Si invita pertanto l'utente finale a contattare il fornitore del dispositivo, ovvero 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 3 Logo "WEEE"

Per maggiori informazioni consultare:

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

Condizioni ambientali, conformità e requisiti strutturali

NOTA

Questa pompa richiede uno spazio libero di almeno 20 cm su tutti i lati per consentire la ventilazione, l'accesso per la manutenzione e un facile avvicinamento all'interruttore di alimentazione. Se si desidera posizionare la pompa su un banco o su un ripiano, è necessario verificare la loro taratura per sostenere il peso della pompa e di qualsiasi altra attrezzatura associata.

Compatibilità elettromagnetica

EN55011/CISPR11

Gruppo 1 apparecchi ISM: Il gruppo 1 contiene tutte le apparecchiature ISM in cui viene generata e/o utilizzata intenzionalmente energia a radiofrequenza ad accoppiamento conduttivo necessaria per il funzionamento interno dell'apparecchiatura stessa.

I **dispositivi di Classe A** sono adatti all'uso in tutte le strutture tranne quelle domestiche e quelle direttamente collegate alla rete pubblica di alimentazione a bassa tensione destinata agli edifici adibiti ad usi domestici.

Questo dispositivo è conforme ai requisiti di CISPR11, gruppo 1, Classe A come apparecchiatura radioterapica professionale. Pertanto possono sussistere potenziali difficoltà per garantire la compatibilità elettromagnetica in altri ambienti a causa di disturbi condotti e irradiati.

L'uso è soggetto alle seguenti condizioni:

- il dispositivo non deve causare interferenze nocive
- il dispositivo deve accettare le interferenze ricevute, comprese le interferenze che potrebbero causare problemi di funzionamento.

Se questa apparecchiatura causasse interferenze dannose alla ricezione radio o televisiva, che possono essere determinate spegnendo e riaccendendo l'apparecchiatura, si consiglia di adottare una o più delle seguenti misure:

- assicurarsi di utilizzare cavi appropriati per collegare il dispositivo alle periferiche
- non apportare variazioni o modifiche non espressamente approvate da Agilent Technologies, che potrebbero inficiare l'autorizzazione all'operatore per l'utilizzo dell'apparecchiatura.

Istruzioni per l'uso

Dichiarazione EMC di Classe A della Corea del Sud

Questa apparecchiatura è di Classe A, è adatta all'uso professionale ed è destinata all'uso in ambienti elettromagnetici esterni alla casa.

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자과적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

Questo dispositivo ISM è conforme alla normativa canadese ICES-001/NMB-001.

Servizio Post Vendita

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

vpt-customer@agilent.com

vpl-customer@agilent.com

È necessario completare il modulo "Richiesta di reso" per restituire il prodotto ad Agilent per l'assistenza (fornito alla fine di questo manuale).

2

Betriebsanleitung

Informationen über diese Betriebsanleitung	39
Einschlägige unterstützende Dokumente	39
Definitionen und Terminologie	40
Definition von Vorsicht, Warnung und Hinweis	40
Warnsymbole	41
Sicherheit	42
EMV-Warnungen	43
FCC	43
Ordnungsgemäße Nutzung	44
Vernünftigerweise vorhersehbarer Missbrauch	44
Sicherheitsrichtlinien für trockene Scrollpumpe	45
Qualifiziertes Personal	46
Schutzausrüstung	47
Produktbeschreibung	49
Transport und Lagerung	49
Auspacken und Inspektion	50
Vorbereitung zum Einbau	51
Inbetriebnahme	55
Abschaltverfahren	58
Entsorgung	59
Umweltbedingungen, Konformität und Versorgungsanforderungen	60
Elektromagnetische Verträglichkeit	60
EN55011/CISPR11	60
Südkoreanische Klasse A EMV-Erklärung	61
ICES/NMB-001	61
Service	62

Informationen über diese Betriebsanleitung

Dieses Handbuch enthält Anleitungen für Benutzer und technische Informationen zu den IDP-4 trockene Scrollpumpe in Übereinstimmung mit den geltenden Richtlinien.

Dieses Dokument ist für das zum Zeitpunkt der Veröffentlichung hergestellte Produkt gültig, vorausgesetzt, dass keine Änderungen am Produkt durch den Kunden oder Dritte vorgenommen werden. Im Falle widersprüchlicher Informationen ist die mit dem Produkt gelieferte Dokumentation als die gültige Dokumentation zu betrachten.

HINWEIS

- 1 Dieses Handbuch enthält wichtige Informationen, damit das Personal, das die IDP-4 trockene Scroll-Pumpe einsetzt, diese sicher betreiben und ihre Leistung über die gesamte Lebensdauer gewährleisten kann.
- 2 Bewahren Sie dieses Handbuch zusammen mit allen zugehörigen Publikationen an einem zugänglichen Ort auf, der allen Bedienern/Wartungspersonal bekannt ist.]

Einschlägige unterstützende Dokumente

X3802-90000 Konformitätserklärung

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

Im Folgenden finden Sie eine Liste von Symbolen, die in Verbindung mit Warnhinweisen auf die IDP-4 Trockene Scroll-Pumpe erscheinen. 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. 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.



	CE-Zertifizierung
	CSA-Zertifizierung
	RoHS-Chinazertifizierung
	Elektro- und Elektronikaltgeräte
	RCM-Zertifizierung
	Herstellungsbetrieb
	Heiße Oberflächen
	Hochspannung
	Allgemeine Gefahr

Betriebsanleitung

	Schutzerdungsanschluss
	Wechselstrom (AC)
	Einschalten
	Ausschalten

Sicherheit

Dieser Abschnitt enthält die von der Niederspannungsrichtlinie 2006/42/EC vorgeschriebenen Informationen, die für die Einhaltung und Befolgung der Sicherheitsvorschriften sowohl allgemein als auch in Bezug auf die spezifische Verwendung des Produkts wesentlich sind.

Die Nichtbeachtung dieser Anleitung und der anderen in diesem Handbuch enthaltenen Anleitungen kann die Sicherheit beeinträchtigen und möglicherweise zu Unfällen bei der Bedienung des Produkts führen.

Agilent Technologies lehnt jede Verantwortung für Schäden am Produkt oder für die körperliche Unversehrtheit des Bedieners oder Dritter ab, die sich aus der Nichtbeachtung der in der technischen Dokumentation angegebenen Sicherheitsvorschriften ergeben.

Betriebsanleitung

EMV-Warnungen

EN 55011 Klasse A Warnung

Dies ist ein Produkt der Klasse A. In einer häuslichen Umgebung kann dieses Produkt Funkstörungen verursachen; in diesem Fall kann der Benutzer aufgefordert werden, angemessene Maßnahmen zu ergreifen.

FCC

Dieses Gerät erfüllt die Anforderungen von Teil 15 der FCC-Vorschriften. Der Betrieb unterliegt den folgenden zwei Bedingungen: (1) Dieses Gerät darf keine schädlichen Interferenzen verursachen, und (2) dieses Gerät muss alle empfangenen Interferenzen akzeptieren, einschließlich Interferenzen, die einen unerwünschten Betrieb verursachen können.

HINWEIS

Das Gerät wurde getestet und entspricht den Grenzwerten für ein digitales Gerät der Klasse A gemäß Teil 15 der FCC-Vorschriften. Diese Grenzwerte sollen einen angemessenen Schutz gegen schädliche Störungen bieten, wenn das Gerät in einer kommerziellen Umgebung betrieben wird. Dieses Gerät erzeugt und verwendet Hochfrequenzenergie und kann diese auch ausstrahlen. Wenn es nicht in Übereinstimmung mit der Bedienungsanleitung installiert und verwendet wird, kann es schädliche Störungen des Funkverkehrs verursachen. Der Betrieb dieses Geräts in einer Wohngegend kann ebenfalls schädliche Funkstörungen verursachen; in diesem Fall ist der Benutzer verpflichtet, die Störungen auf eigene Kosten zu beheben.

Ordnungsgemäße Nutzung

Die IDP-4 Trockene Scroll-Pumpe ist ausschließlich für die Nutzung durch qualifizierte Personen als Vakuumpumpe gemäß diesem Handbuch vorgesehen.

Alle Zubehörteile und Geräte, die zusammen mit dem Produkt verwendet werden, müssen von Agilent Technologies geliefert oder genehmigt werden. Agilent Technologies übernimmt keine Garantie oder Verantwortung für Anwendungen, die von den in diesem Handbuch beschriebenen abweichen oder bei denen die grundlegenden Anforderungen und Sicherheitsmaßnahmen nicht beachtet werden.

Vernünftigerweise vorhersehbarer Missbrauch

Agilent Technologies lehnt jede Verantwortung ab, die sich aus der unsachgemäßen Nutzung und dem vorhersehbaren Missbrauch der IDP-4 Trockene Scroll-Pumpe ergibt. Unsachgemäßer Nutzung führt zum Erlöschen aller Haftungs- und Gewährleistungsansprüche.

Eine unsachgemäße Nutzung kann unter anderem sein:

- Anschluss der Pumpe an eine falsche Spannungsversorgung
- Falscher oder fehlender Erdungsanschluss
- Dauerhafter Betrieb oberhalb des empfohlenen Eingangsdrucks
- Betrieb mit Überdruck am Gasballasteingang
- Betrieb des Gasballasts mit offener Atmosphäre beim Pumpe gefährlicher Gase“
- Betrieb bei zu hoher Umgebungstemperatur
- Betrieb bei einer Umgebungstemperatur unterhalb der Spezifikation
- Beschädigung durch unsachgemäße Lagerung und Transport der Pumpe
- Durchführen von Wartungsarbeiten oder Entfernen von Verkleidungen ohne Unterbrechung der Netzstromversorgung
- Betrieb ohne Luftströmungselemente (Abdeckung oder Lüfter)
- Versand des zurückgegebenen Produkts in einer nicht werkseitigen Verpackung.

Sicherheitsrichtlinien für trockene Scrollpumpe

Das für den Betrieb und die Wartung der Pumpe zuständige Personal muss entsprechend geschult sein und die Unfallverhütungsvorschriften kennen (siehe Abschnitt Qualifiziertes Personal). Die in diesem Handbuch enthaltenen Unfallverhütungsvorschriften müssen während des Betriebs und der Wartung der Pumpe ständig beachtet werden, um Verletzungen der Bediener und Schäden an der Pumpe oder anderen Geräten zu vermeiden. Diese Vorsichtsmaßnahmen sind in Form von WARNUNG- und VORSICHT-Hinweisen im gesamten Benutzerhandbuch enthalten.

WARNUNG



Um Schäden am Gerät und Verletzungen des Bedienpersonals zu vermeiden, sind die Installationsanweisungen in diesem Handbuch genau zu befolgen.

WARNUNG



**Die Pumpe darf nicht mit brennbaren Materialien umgehen.
Es ist verboten, die Pumpe in einer explosiven Umgebung zu installieren.**

Qualifiziertes Personal

Die IDP-4 Trockene Scroll-Pumpe soll nur von qualifiziertem Personal verwendet werden. Eine qualifizierte Person ist:

- Eine Person, die über die entsprechenden Qualifikationen verfügt und Erfahrung im Umgang mit Agilent Trockenvakuumpumpe hat oder eine entsprechende Schulung absolviert hat.
- Eine Person, die die notwendigen Sicherheitsmaßnahmen befolgt und die Pumpe unter den richtigen Bedingungen betreibt, um Verletzungen oder Schäden zu vermeiden.
- Eine Person, die die geltenden nationalen und lokalen Sicherheits- und Regulierungsstandards und -vorschriften kennt und entsprechend qualifiziert ist, diese zu befolgen.

HINWEIS

Alle Reparaturen an diesem Produkt, die über die in diesem Benutzerhandbuch aufgeführten Standardwartungen hinausgehen, dürfen ausschließlich von autorisiertem Agilent-Personal durchgeführt werden. Alle Einstellungen, Änderungen oder Wartungsarbeiten müssen von einem professionellen Techniker durchgeführt werden, der über die Risiken informiert ist.

Schutzausrüstung

Die Schutzausrüstung des Bedienpersonals, das die Pumpe bedient oder wartet, muss stets der Art der ausgeführten Arbeiten angemessen sein. Die verwendete Schutzausrüstung muss den Sicherheitsanforderungen der geltenden Gesetzgebung des Landes entsprechen, in dem die Maschine eingesetzt wird.

Im Allgemeinen muss der Bediener beim Bewegen, Hantieren oder Installieren der IDP-4 Trockene Scroll-Pumpe Sicherheitsschuhe tragen.

WARNUNG



Gesundheitsgefahr durch gefährliche Stoffe bei Wartung oder Installation

- Vakuumpumpe, Komponenten oder Betriebsflüssigkeiten können je nach Prozessbesonderheit mit giftigen, reaktiven oder radioaktiven Stoffen verunreinigt sein.
- Tragen Sie bei Wartungs- und Reparaturarbeiten oder bei einer erneuten Installation angemessene Schutzausrüstung.

WARNUNG



Verletzungsgefahr durch herunterfallende Gegenstände

Beim Transport von Vakuumpumpe von Hand besteht die Gefahr, dass Lasten verrutschen und herunterfallen.

- Tragen Sie kleine und mittlere Vakuumpumpe mit zwei Händen.
- Alle Geräte, die schwerer als 20 kg sind, sollten mit einer geeigneten Hebevorrichtung transportiert werden.
- Sicherheitsschuhe mit Stahlkappe gemäß Richtlinie EN 347 tragen.

VORSICHT



Verletzungsgefahr durch scharfe Kanten

- Warten Sie vor der Wartung der Pumpe oder vor dem Ein- und Ausbau der Pumpe vom System bis zum vollständigen Stillstand der Pumpe.
 - Nicht direkt im Hochvakuumflansch betreiben.
 - Gegebenenfalls Schutzhandschuhe nach EN 420 tragen.
-

Produktbeschreibung

Die IDP-4 ist eine hermetisch abgedichtete Trocken-Scrollpumpe. Diese Pumpe eignet sich zum Fördern von Luft oder Inertgasen aus einem Bereich mit niedrigerem Druck in einen Bereich mit höherem Druck, der nicht wesentlich unter oder über dem atmosphärischen Druck liegt. Diese Pumpe ist nicht zum Fördern von korrosiven, pyrophoren, partikelbildenden Gasen oder Gasgemischen im entflammaren und/oder brennbaren Bereich vorgesehen. Sie ist auch nicht als Kompressor zum Erreichen von Förderdrücken wesentlich über dem atmosphärischen Druck vorgesehen.

Das Handbuch und die zugehörigen Dokumente enthalten die notwendigen Informationen, um die Sicherheit des Bedieners bei der Verwendung des Geräts zu gewährleisten. Ausführliche Informationen sind im Abschnitt „Technical Information“ enthalten.

Transport und Lagerung

Bei Transport und Lagerung der Pumpe sollten die folgenden Umgebungsbedingungen nicht überschritten werden:

- Temperatur: -20 °C bis +60 °C (-4 °F bis 140 °F)
- Relative Luftfeuchtigkeit: 0 bis 95% (nicht kondensierend)

HINWEIS

Es ist zu beachten, dass während des Transports und der Lagerung von Vakuumpumpe Wasserdampf aus der Umgebung in die Vakuumpumpe eindringt, unabhängig davon, wie gut die Pumpe abgedichtet ist. Um die volle Nennvakuumleistung der Pumpe zu erreichen, muss dieser Wasserdampf mit Hilfe der im Abschnitt „Technical Information“ beschriebenen Gasballastfunktionen entfernt werden.

Auspacken und Inspektion

Die Pumpe wird in einer speziellen Schutzverpackung geliefert. Wenn diese Anzeichen von Schäden aufweist, die während des Transports entstanden sein könnten, wenden Sie sich an Ihr örtliches Verkaufsbüro. Das Gesamtgewicht von Pumpe und Verpackung beträgt ca. 10 kg (20 Pfund).

WARNUNG



Der Karton und alle Verpackungsmaterialien werden im Falle einer Wartung in einem von Agilent zugelassenen Service-Center benötigt. Die Aufbewahrung dieser Verpackung erleichtert eventuelle Serviceanfragen.

WARNUNG



Achten Sie beim Auspacken der Pumpe darauf, sie nicht fallen zu lassen, und vermeiden Sie jede Art von plötzlichem Aufprall oder Stoß.

Vorbereitung zum Einbau

Die IDP-4 Pumpe ist für den Betrieb in Innenräumen (Installationskategorie II) in relativ sauberen, trockenen Umgebungen (Verschmutzungsgrad 2) ausgelegt und zertifiziert. Die technischen Daten finden Sie in Tabelle 1 im Abschnitt „Technical Information“. Diese Pumpe ist nicht für den Betrieb im Freien ausgelegt oder zertifiziert.

Die IDP-4 Pumpe darf nicht in Umgebungen eingesetzt werden, die explosive oder brennbare Materialien enthalten. Diese Pumpe ist NICHT eigensicher und dürfen niemals mit korrosiven, pyrophoren, partikelbildenden Gasen oder Gasgemischen in ihrem entflammaren und/oder brennbaren Bereich verwendet werden. Stellen Sie die Pumpe auf eine feste, ebene Fläche. Wenn sie auf einer Bank oder einem Tisch montiert werden muss, stellen Sie sicher, dass die Bank oder der Tisch für das Gewicht der Pumpe und aller zugehörigen Geräte ausgelegt ist.

Schließen Sie die Pumpe an die Stromversorgung an, indem Sie ein Netzkabel im IEC-320-Format mit einer Kapazität von mindestens 10 A verwenden.

HINWEIS

Die häufigste Ursache für eine schlechte Leistung bei der ersten Installation und Inbetriebnahme ist adsorbierter Wasserdampf sowohl in der Pumpe als auch in der Vakuumkammer. Der Betrieb der Pumpe mit aktiviertem Gasballast kann ebenfalls vorteilhaft sein, wenn die Pumpe über einen längeren Zeitraum abgeschaltet war, wodurch Wasser sowohl in der Pumpe als auch im Vakuumsystem adsorbiert werden konnte. Je nach Vakuumsystem und der Menge des adsorbierten Wassers kann es typischerweise zwischen etwa 30 Minuten und etwa 24 Stunden Gasballastbetrieb dauern, um die volle Vakuumleistung zu erreichen.

VORSICHT

Während des Betriebs müssen die folgenden Umgebungsbedingungen beachtet werden:

- **Temperatur: +5 °C bis +40 °C (41 °F bis 104°F)**
- **Relative Luftfeuchtigkeit: 5 bis 95% (nicht kondensierend)**

Schließen Sie die Pumpe mit einem Netzkabel nach IEC-320 mit der entsprechenden Strom- und Spannungsstärke an das Stromkabel an (siehe Tabelle 2).

VORSICHT

Bei der Installation der Vakuumpumpe wird empfohlen, die Pumpe so zu positionieren, dass der Netzschalter und der Anschluss des Stromkabels für den Benutzer leicht zugänglich bleiben, um die Pumpe bei Bedarf schnell vom Netz trennen zu können.

Betriebsanleitung

VORSICHT

Ziehen Sie die beiden Sechskantschrauben auf jeder Seite des Pumpekopfes nicht an, lösen Sie sie nicht und verändern Sie sie in keiner Weise (siehe folgende Abbildung). Ein solches Vorgehen führt zu Leistungseinbußen und/oder Schäden an der Pumpe.



Abbildung 4

WARNUNG



Entfernen oder verändern Sie keine Sicherheits- oder Isoliereinrichtungen an der Pumpe. Andernfalls kann dies ein ernsthaftes Sicherheitsrisiko darstellen und zum Erlöschen der Garantie führen.

WARNUNG



- Diese Pumpe ist nur zum Fördern und Absaugen von Luft und Inertgasen geeignet. Diese Pumpe ist nicht zum Fördern von korrosiven, pyrophoren, partikelbildenden Gasen oder Gasgemischen in ihrem entflammbaren und/oder brennbaren Bereich bestimmt.
 - Installieren Sie die Pumpe an einem Ort, an dem sie nicht Regen, Dampf oder übermäßiger Feuchtigkeit ausgesetzt ist. Sie können elektrische Schläge, Kurzschlüsse und schwere Körperverletzungen verursachen.
 - Vergewissern Sie sich vor der Inspektion, Neukonfiguration oder Wartung der Pumpe, dass die Stromzufuhr unterbrochen ist.
 - Der Gasballast muss versiegelt werden, wenn Gas gepumpt wird, das nicht in die Atmosphäre entweichen oder mit Luft vermischt werden kann.
 - Für Anwendungen, bei denen es wünschenswert ist, den von der Pumpe erzeugten Schallpegel zu verringern, ist ein Schalldämpfer als Option erhältlich.
 - Um Körperverletzungen zu vermeiden, darf kein Teil des menschlichen Körpers dem Vakuum ausgesetzt werden.
 - Die Pumpe muss so installiert werden, dass der Netzschalter leicht zugänglich ist.
 - Die von Agilent zugelassenen Stromkabel für alle gängigen Anschlüsse sind im Abschnitt „Elektrische Anschlüsse“ angegeben. Wenn ein anderes Stromkabel verwendet werden muss, muss die Stromstärke des Stromkabels mindestens 10A für 200-240 V und mindestens 15A für 100-127 V betragen.
-

Hinweise zur Erdung

WARNUNG



Dieses Produkt sollte geerdet werden. Im Falle eines elektrischen Kurzschlusses verringert die Erdung das Risiko eines elektrischen Schlags, indem sie einen Ableitungsweg für den elektrischen Strom schafft. Diese Pumpe ist mit einem Stromkabel ausgestattet, das mit einem Erdungsdraht und einem entsprechenden Erdungsstecker versehen ist. Der Stecker muss in eine ordnungsgemäß installierte und geerdete Steckdose eingesteckt werden, die allen örtlichen Vorschriften und Verordnungen entspricht.

Der Stecker des Stromkabels muss ordnungsgemäß in das Stromversorgungsmodul der Pumpe eingesteckt sein, um eine fehlerhafte oder unterbrochene Verbindung zu vermeiden.

VORSICHT

Obwohl die Pumpe Spuren von Partikeln fördern kann, die normalerweise in der Atmosphäre vorkommen, ist sie nicht für die Verarbeitung von Feststoffen, Chemikalien, Pulvern, Lösungsmitteln, Kondensaten, Flüssigkeiten oder anderen Partikeln ausgelegt. Diese können das Gerät beschädigen, seine Leistung beeinträchtigen oder seine Lebensdauer verkürzen.

Es wird empfohlen, rund um das System einen Freiraum von mindestens 20 cm zu lassen, um eine gute Luftzirkulation zu ermöglichen.

VORSICHT

Die Außenflächen der Pumpe dürfen nur mit milden Reinigungsmitteln gesäubert werden.

Die trockene Scroll-Pumpe enthält keine Flüssigkeit zur Reinigung von angesammeltem Staub und Ablagerungen. Lassen Sie die Pumpe regelmäßig für ein oder zwei Minuten an der Atmosphäre laufen, um sie durchzuspülen.

WARNUNG



Wenden Sie sich an einen qualifizierten Elektriker oder Servicetechniker, wenn Sie die Hinweise zur Erdung nicht vollständig verstanden haben oder wenn Sie Zweifel daran haben, ob das Gerät richtig geerdet ist. Verändern Sie den mitgelieferten Stecker nicht. Wenn er nicht in die Steckdose passt, lassen Sie die richtige Steckdose von einem qualifizierten Elektriker installieren.

Schließen Sie das Gerät nur an eine Steckdose an, die die gleiche Konfiguration wie der Stecker hat. Verwenden Sie keinen Adapter mit diesem Produkt.

Inbetriebnahme

WARNUNG



Entfernen Sie die Flanschabdeckungen von Auslass und Einlass.



Das Startverfahren für die IDP-4 Trocken-Scroll-Vakuumpumpe ist wie folgt:

1. Montieren Sie die Pumpe an ihrem vorgesehenen Standort und stellen Sie die entsprechenden mechanischen und Vakuumanlüsse her oder verschließen Sie alternativ den Einlass mit einem Blindflansch.
2. Stellen Sie sicher, dass der Auslassanschluss nicht blockiert oder verschlossen ist. Wenn ein Abluftschlauch oder Schalldämpfer verwendet wird, darf der Abgasgegendruck 6,5 psig nicht überschreiten.
3. Stellen Sie sicher, dass sich der EIN/AUS-Schalter in der AUS-Position „0“ befindet.
4. Schließen Sie das entsprechende Netzkabel an und stecken Sie das Kabel in die Netzsteckdose. Die Stromstärke des Netzkabels muss mindestens 10 A für den Betrieb bei 200–240 V und mindestens 15 A für den Betrieb bei 100–127 V betragen.
5. Drücken Sie den Hauptschalter EIN/AUS in die Position „I“.
6. Wenn die Pumpe mit einem optionalen Einlassventil ausgestattet ist, öffnet sich das Ventil etwa 10 Sekunden nach dem Einschalten.

Es kann mehrere Stunden dauern, bis die Pumpe den angegebenen Enddruck erreicht. Wenn die Pumpe den erwarteten Enddruck nicht erreicht, kann dies auf angesammelten Wasserdampf im Inneren der Pumpe zurückzuführen sein.

Betriebsanleitung

HINWEIS

Die häufigste Ursache für eine schlechte Leistung bei der ersten Installation und Inbetriebnahme ist adsorbierter Wasserdampf sowohl in der Pumpe als auch in der Vakuumkammer. Der Betrieb der Pumpe mit aktiviertem Gasballast kann ebenfalls vorteilhaft sein, wenn die Pumpe über einen bestimmten Zeitraum abgeschaltet war, wodurch Wasser sowohl in der Pumpe als auch im Vakuumsystem adsorbiert werden konnte. Abhängig vom Vakuumsystem und der Menge des adsorbierten Wassers kann es in der Regel zwischen etwa 30 Minuten und etwa 24 Stunden Gasballastbetrieb dauern, bis die volle Vakuumleistung erreicht ist.



Abbildung 5

HINWEIS

Der EIN/AUS-Schalter der Pumpe ist ein Wippschalter, der gemäß IEC-Publikation 417 mit Symbolen für die EIN- und AUS-Positionen versehen ist. Die folgende Abbildung zeigt den Schalter in der AUS-Position.

WARNUNG



Führen Sie keine Fremdkörper durch die Kühllüftungsöffnungen ein, da der Kontakt mit den Lüfterflügeln zu schweren Verletzungen oder zu Schäden an der Pumpe führen kann.

WARNUNG



Blockieren Sie nicht die Kühllüftungsöffnungen. Das Blockieren dieser Öffnungen kann zu einer Überhitzung der Pumpe führen. Eine Pumpeoberflächentemperatur von mehr als 65 °C (150 °F) kann potenziell schädlich sein. Wenn solche Bedingungen festgestellt werden, schalten Sie die Pumpe aus und lassen Sie sie abkühlen. Zerlegen Sie die Pumpe, prüfen Sie sie auf Schäden und führen Sie gegebenenfalls Reparaturen durch.

WARNUNG



Verwenden Sie keine Gasballastzufuhr, wenn potenziell gefährliche Materialien eingesetzt werden oder solche, die nicht mit Luft vermischt werden dürfen.

Abschaltverfahren

Um die Pumpe in jeder Betriebsart manuell abzuschalten, stellen Sie den ON/OFF-Schalter auf die Position "0". Wenn die Pumpe mit einem integrierten Absperrventil ausgestattet ist, schließt sich dieses schnell, um einen Rückfluss von Gas aus der Pumpe zu den vorgeschalteten Komponenten zu verhindern.

VORSICHT

Bevor die IDP-4 Pumpe abgeschaltet wird, stellen Sie sicher, dass alle vorgelagerten, miteinander verbundenen Komponenten ordnungsgemäß abgesperrt oder auf eine mögliche Einwirkung des Atmosphärendrucks vorbereitet sind.

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 Industiemüll 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 6 Logo „WEEE“

Für weitere Informationen siehe:

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

Umweltbedingungen, Konformität und Versorgungsanforderungen

HINWEIS

Diese Pumpe benötigt auf allen Seiten einen Freiraum von mindestens 20 cm (8 Zoll) für die Belüftung, den Zugang zur Wartung und den einfachen Zugang zum Netzschalter. Wenn die Pumpe auf einer Bank oder einem Regal aufgestellt werden soll, muss dieses so ausgelegt sein, dass es das Gewicht der Pumpe und aller anderen zugehörigen Geräte tragen kann.

Elektromagnetische Verträglichkeit

EN55011/CISPR11

Gruppe 1 ISM Geräte: Gruppe 1 umfasst alle ISM-Geräte, in denen absichtlich leitfähige Funkfrequenzenergie erzeugt und/oder verwendet wird, die für den internen Betrieb des Geräts selbst erforderlich ist.

Geräte der Klasse A sind Geräte, die zur Verwendung in allen Einrichtungen geeignet sind, die nicht zum Wohnen dienen und auch nicht direkt an ein Niederspannungsnetz angeschlossen sind, das Wohngebäude mit Strom versorgt.

Dieses Gerät erfüllt die Anforderungen von CISPR11, Gruppe 1, Klasse A als professionelles Strahlungsgerät. Daher kann es in anderen Umgebungen aufgrund von leiter- und strahlungsgebundenen Störungen zu möglichen Schwierigkeiten bei der Sicherstellung der elektromagnetischen Verträglichkeit kommen.

Der Betrieb unterliegt den folgenden zwei Bedingungen:

- Dieses Gerät darf keine schädlichen Störungen verursachen.
- Dieses Gerät muss alle empfangenen Interferenzen tolerieren, einschließlich Interferenzen, die einen unbeabsichtigten Betrieb verursachen können.

Sollte dieses Gerät den Radio- oder Fernsehempfang stören, was durch Aus- und Einschalten des Geräts festgestellt werden kann, wird dem Benutzer empfohlen, eine oder mehrere der folgenden Maßnahmen zu ergreifen:

- Sicherstellen, dass geeignete Kabel für den Anschluss des Gerätes an Zusatzgeräte verwendet werden.
- Änderungen oder Umbauten, die nicht ausdrücklich von Agilent Technologies genehmigt wurden, können dazu führen, dass der Benutzer die Berechtigung zum Betrieb des Geräts verliert.

Südkoreanische Klasse A EMV-Erklärung

Dieses Gerät entspricht der Klasse A und ist für den professionellen Einsatz und die Verwendung in elektromagnetischen Umgebungen außerhalb des Wohnbereichs geeignet.

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자파적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

Dieses ISM-Gerät entspricht den kanadischen Bestimmungen ICES-001/NMB-001.

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 care@agilent.com

vpl-customer care@agilent.com

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

3

Mode d'emploi

À propos de ce manuel	64
Documents complémentaires pertinents	64
Définitions et terminologie	65
Signification des messages d'attention, d'avertissement et des notes	65
Symboles d'avertissement	66
Sécurité	67
Avertissement CEM	68
FCC	68
Utilisation appropriée	69
Utilisation inappropriée	69
Directive de sécurité pour la pompe à vide sèches	70
Personnel qualifié	71
Équipements de protection	72
Description du produit	74
Transport et entreposage	74
Déballage et inspection	75
Préparation pour l'installation	76
Procédure de démarrage	80
Procédure d'arrêt	83
Élimination	84
Conditions environnementales, conformité et applications pratiques	85
Compatibilité électromagnétique	85
EN55011/CISPR11	85
Déclaration CEM Classe A Corée du Sud	86
ICES/NMB-001	86
Service	87

À propos de ce manuel

Le présent manuel contient les instructions et les informations techniques destinées aux utilisateurs de la pompe à vide sèches IDP-4 conformément aux directives applicables.

Ce document s'applique au produit fabriqué au moment de sa publication, sous réserve qu'aucune modification n'ait été apportée au produit par le client ou les tiers. En cas d'informations contradictoires, la documentation expédiée avec le produit doit être considérée comme celle faisant foi.

NOTE

- 1 Ce manuel contient des informations utiles pour que l'ensemble du personnel puisse utiliser la pompe à vide sèches IDP-4 en toute sécurité et pour garantir leurs performances tout au long de leur 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.

Documents complémentaires pertinents

X3802-90000

Déclaration de conformité

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

Une liste des symboles associés aux avertissements relatifs à la pompe à vide sèches IDP-4 est reportée ci-dessous. 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. 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.



	Certification CE
	Certification CSA
	Certification RoHS Chine
	Déchets d'équipements électriques et électroniques
	Certification RCM
	Site de fabrication
	Surfaces chaudes
	Tensions dangereuses
	Danger générique

Mode d'emploi

	Borne de mise à la terre de protection
	Courant Alternatif (AC)
	Mise sous tension
	Mise hors tension

Sécurité

Cette section contient les informations prescrites par la directive Basse tension 2006/42/EC, 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 consignes et des autres instructions contenues dans le présent manuel peut avoir des conséquences sur la sécurité et exposer ceux qui utilisent le produit à des accidents.

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.

Mode d'emploi

Avertissement CEM

EN 55011 Avertissement Classe A

Il s'agit d'un produit de Classe A. Dans un environnement domestique, il peut produire des interférences radio, auquel cas l'utilisateur sera tenu d'adopter les mesures appropriées.

FCC

Cet appareil est conforme à la partie 15 du règlement de la FCC. Son utilisation est soumise aux deux conditions suivantes : (1) Cet appareil ne doit pas provoquer d'interférences nocives, et (2) cet appareil doit supporter toute interférence reçue, y compris les interférences susceptibles d'entraîner un fonctionnement non désiré.

NOTE

Cet appareil a été testé et respecte les limitations d'un appareil numérique de classe A, conformément au point 15 du règlement FCC. Ces limitations sont conçues pour fournir une protection raisonnable contre les interférences nocives dans un environnement domestique. Cet équipement génère, utilise et peut émettre des fréquences radio. Il est susceptible de provoquer des interférences nocives au niveau des communications radio en cas d'installation et d'utilisation non conformes au mode d'emploi. L'utilisation de ce matériel en zone résidentielle peut également générer des interférences nuisibles au niveau des communications radio, auquel cas l'utilisateur devra corriger les interférences à ses propres frais.

Mode d'emploi

Utilisation appropriée

La pompe à spirales sèches IDP-4 est conçue pour être utilisée exclusivement par des personnes qualifiées conformément à ce manuel.

Tout accessoire et équipement utilisé avec le produit doit être fourni ou agréé par Agilent Technologies.

Agilent Technologies n'offre aucune garantie et décline toute responsabilité en cas d'utilisation dans des applications différentes de celles décrites dans ce document ou de non-respect des exigences essentielles ou des consignes de sécurité.

Utilisation inappropriée

Agilent Technologies décline toute responsabilité en cas d'utilisation inappropriée et d'usage abusif prévisible des pompes à vide sèches IDP- 3 PLUS. Une utilisation inappropriée entraînera l'annulation de toutes les créances et garanties.

Une utilisation inappropriée peut comprendre, sans s'y limiter, les situations suivantes :

- Branchement de la pompe sur une tension d'alimentation erronée
- Mise à la terre incorrecte ou manquante
- Fonctionnement continu au-dessus de la pression d'entrée préconisée
- Fonctionnement en présence de surpressions au niveau de l'entrée de lest de gaz
- Fonctionnement avec le lest de gaz ouvert vers l'atmosphère lors du pompage de gaz dangereux
- Fonctionnement à une température ambiante trop élevée
- Fonctionnement à une température ambiante inférieure aux spécifications
- Dommages dus à des modalités incorrectes de stockage et de transport de la pompe
- Exécution d'interventions d'entretien ou enlèvement des capots sans débrancher l'alimentation principale
- Fonctionnement sans éléments de circulation d'air (couverture ou ventilateur)
- Expédition du produit renvoyé dans un emballage autre que celui d'usine.

Directive de sécurité pour la pompe à vide sèches

Le personnel responsable de l'exploitation et de l'entretien des pompes doit être formé correctement et être conscient des règles de prévention des accidents (voir la section Personnel qualifié). Les précautions de prévention des accidents contenues dans ce manuel doivent être respectées en permanence pendant l'exploitation et l'entretien des pompes pour éviter les risques de blessure et d'endommagement des pompes ou d'autres équipements. Ces précautions sont données sous la forme de notes intitulées AVERTISSEMENT et ATTENTION tout au long de ce mode d'emploi.

AVERTISSEMENT



Pour éviter d'endommager les équipements et éviter de blesser le personnel préposé, les consignes d'installation reportées dans ce manuel doivent être respectées à la lettre !

AVERTISSEMENT



La pompe ne doit pas être utilisée avec des matières inflammables. Il est interdit d'installer la pompe dans un environnement explosif.

Mode d'emploi

Personnel qualifié

La pompe à spirales sèches IDP-4 ne doit être utilisée que par du personnel qualifié. Une personne qualifiée est :

- Un individu qui dispose des qualifications appropriées et de l'expérience nécessaire, ou a suivi une formation sur l'utilisation de pompes à vide sèches Agilent.
- Un individu qui respecte les consignes de sécurité nécessaires et exploite la pompe conformément aux conditions stipulées de manière à ne pas provoquer de dommages corporels ni matériels.
- Un individu qui est informé et correctement qualifié pour respecter les règlements ainsi que les normes réglementaires et consignes de sécurité en vigueur à l'échelle nationale et locale.

NOTE

Toutes les réparations effectuées sur ce produit, autres que celles considérées comme des interventions d'entretien standard dans ce mode d'emploi, doivent être réalisées exclusivement par du personnel agréé Agilent. Tout réglage, modification ou intervention d'entretien doit être effectué par un technicien professionnel ayant été informé des risques encourus.

Équipements de protection

Les équipements de protection des opérateurs qui utilisent ou entretiennent la pompe doivent toujours être appropriés au type d'opération à effectuer. Ils doivent satisfaire aux exigences de sécurité de la réglementation en vigueur dans le pays d'utilisation de l'appareil.

En règle générale, l'opérateur doit porter des chaussures de sécurité pendant la manutention, la manipulation ou l'installation des pompes à vide sèches IDP-4.

AVERTISSEMENT



Danger pour la santé en raison de la présence de substances dangereuses pendant la maintenance ou l'installation.

- En fonction des caractéristiques du processus, la pompe à vide, les composants ou les fluides de fonctionnement peuvent être contaminés par des substances toxiques, réactives ou radioactives.
- Portez des équipements de protection individuelle appropriés pendant la maintenance, les opérations de réparation ou de réinstallation.

AVERTISSEMENT

Risques de blessure dus à la chute d'objets



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

- Transportez la pompe à vide de petite et moyenne dimension avec les deux mains.
- Tout objet de plus de 20 kg doit être transporté à l'aide d'un moyen de levage adapté.
- Portez des chaussures de sécurité à pointes en acier conformes à la directive EN 347.

Mode d'emploi

ATTENTION

Risques de blessures par des arêtes vives ou tranchantes



- Avant de réparer la pompe et avant toute opération de montage/démontage de la pompe du système, attendez qu'elle soit complètement arrêtée.
 - N'intervenez pas directement à l'intérieur de la pompe.
 - Si nécessaire, portez des gants de protection conformes à la directive EN 420.
-

Description du produit

Le modèle IDP-4 est une pompe à spirales sèches hermétique. Elles sont adaptées au pompage de l'air ou de gaz inertes depuis une zone basse pression vers une zone dont la pression est supérieure qui ne se trouve pas sensiblement au-dessus ou en dessous de la pression atmosphérique. Ces pompes ne sont pas conçues pour pomper des gaz corrosifs, pyrophoriques ou formant des particules, ni des mélanges de gaz dans leur état inflammable et/ou combustible. Elles ne sont pas non plus conçues pour servir de compresseur en vue d'atteindre des pressions de refoulement sensiblement supérieures à la pression atmosphérique.

Le mode d'emploi et les documents correspondants contiennent les informations nécessaires pour garantir la sécurité de l'opérateur lorsqu'il utilise l'équipement. Des renseignements détaillés sont reportés dans la section «Technical Information».

Transport et entreposage

Lors du transport et de l'entreposage de la pompe, les conditions environnementales reportées ci-dessous ne doivent pas être dépassées :

- Température : -20 °C à +60 °C (-4 °F à 140 °F)
- Humidité relative : 0 à 95 % (sans condensation)

NOTE

Il faut savoir qu'au cours du transport et de l'entreposage de la pompe à vide, de la vapeur d'eau en provenance de l'environnement ambiant pénétrera à l'intérieur de la pompe à vide indépendamment de l'étanchéité correcte de la pompe. Pour obtenir les meilleures performances nominales à vide de la pompe, cette eau doit être éliminée à l'aide des fonctions de lest de gaz décrites dans la section «Technical Information».

Déballage et inspection

La pompe est livrée dans un emballage spécial de protection. Si celui-ci semble abîmé, suite à des dommages éventuels pendant le transport, contactez votre distributeur local. Le poids total de la pompe et de son emballage est d'environ 10 kg (20 livres).

AVERTISSEMENT



Le carton et tout le matériel d'emballage seront réutilisés si des interventions auprès d'un centre d'entretien agréé Agilent s'avèrent nécessaires. Gardez cet emballage pour faciliter les interventions d'entretien.

AVERTISSEMENT



Lors du déballage de la pompe, veillez à ne pas la laisser tomber et évitez tout impact ou choc brusque.

Préparation pour l'installation

La pompe à spirale sèches IDP-4 est conçue et certifiée pour être utilisée en intérieur (catégorie d'installation II) dans des environnements relativement propres et secs (degré de pollution 2). Voir le Tableau 1 dans la section « Technical Information » pour connaître les spécifications. Ces pompes ne sont pas conçues ni certifiées pour être utilisées à l'extérieur.

La pompe à spirale sèches IDP-4 ne doit jamais être utilisée dans des environnements ambiants qui contiennent des matières explosives ou combustibles. La pompe IDP-4 n'est PAS à sécurité intrinsèque et ne doit jamais être utilisée avec des gaz corrosifs, pyrophoriques, formant des particules ou des mélanges de gaz dans leur état inflammable et/ou combustible.

Placez la pompe sur une surface plane et robuste. Si elle doit être montée sur un banc ou une table, vérifiez que la capacité nominale du support est en mesure de soutenir le poids de la pompe et de tous les équipements associés.

Raccordez la pompe à l'alimentation électrique à l'aide d'un cordon d'alimentation de type IEC-320 d'une capacité d'au moins 10 A.

NOTE

La cause la plus fréquente de baisse de performance lors de la première installation ou de la mise en service est la présence de vapeur d'eau adsorbée à l'intérieur de la pompe et de la chambre à vide. Faire fonctionner la pompe avec le ballast de gaz activé peut s'avérer utile si la pompe est restée à l'arrêt pendant une période prolongée, permettant à l'humidité de s'y accumuler. Selon la configuration du système à vide et la quantité d'eau adsorbée, il faut généralement entre 30 minutes et 24 heures de fonctionnement avec le ballast de gaz ouvert pour atteindre les performances de vide nominales.

ATTENTION

En phase de fonctionnement, les conditions environnementales suivantes doivent être respectées :

- **Température : +5 °C à +40 °C (41 °F à 104°F)**
- **Humidité relative : 5 à 95 % (sans condensation)**

Branchez la pompe sur l'alimentation électrique à l'aide d'un câble du type IEC-320 en veillant à ce que les valeurs nominales de courant et de tension soient correctes (voir Tableau 2).

Mode d'emploi

ATTENTION Lors de l'installation de la pompe à vide, il est conseillé de positionner la pompe de manière à ce que l'utilisateur puisse accéder facilement à l'interrupteur d'alimentation et au raccordement du câble électrique pour pouvoir débrancher rapidement la pompe du réseau, si nécessaire.

ATTENTION Ne pas serrer, desserrer ni modifier en aucune façon les deux boulons hexagonaux situés de chaque côté de la tête de pompe (voir figure suivante). Cela entraînerait une perte de performance et/ou des dommages à la pompe.



Figure 7

AVERTISSEMENT



N'enlevez pas et ne modifiez pas les équipements de sécurité ou l'isolation montés sur la pompe. Dans le cas contraire, cela risque d'entraîner de graves dangers pour la sécurité et de rendre la garantie caduque.

AVERTISSEMENT



- Cette pompe est capable de pomper et d'évacuer uniquement de l'air et des gaz inertes. Cette pompe n'est pas conçue pour pomper des gaz corrosifs, pyrophoriques, ou formant des particules, ni des mélanges de gaz dans leur état inflammable et/ou combustible.
 - Installez-la dans une zone qui n'est pas exposée à la pluie, à de la vapeur ou à une humidité excessive. Cela risque de provoquer une électrocution, des courts-circuits et de graves blessures corporelles.
 - Avant d'inspecter, de reconfigurer ou d'entretenir la pompe, vérifiez que l'alimentation électrique est débranchée.
 - Le lest de gaz doit être hermétique en cas de pompage de gaz qui ne peut pas être mis à l'évent vers l'atmosphère ou mélangé à l'air.
 - Un silencieux d'échappement est disponible en option pour les applications dans lesquelles il est souhaitable de diminuer le niveau sonore émis par la pompe.
 - Pour éviter les blessures corporelles, n'exposez aucune partie du corps au vide.
 - La pompe doit être installée de manière à faciliter l'accès à l'interrupteur d'alimentation.
 - Les câbles électriques agréés par Agilent pour effectuer tous les branchements sont indiqués dans la section « Branchements électriques ». S'il est nécessaire d'utiliser un câble électrique différent, sa valeur nominale de courant doit être au moins de 10A pour un service de 200-240 V et au moins de 15A pour un service de 100-127 V.
-

Mode d'emploi

Instructions de mise à la terre

AVERTISSEMENT



Ce produit doit être mis à la terre. En cas de court-circuit électrique, la mise à la terre réduit le risque d'électrocution en fournissant un chemin de fuite du courant électrique. Cette pompe est équipée d'un câble électrique doté d'un fil de terre avec une fiche de mise à la terre appropriée. Cette fiche doit être insérée dans une prise installée correctement et mise à la terre conformément à tous les codes et règlements locaux.

Le connecteur du câble électrique doit être branché correctement sur le module d'alimentation de la pompe pour empêcher toute connexion défectueuse ou intermittente.

ATTENTION

Bien que la pompe puisse pomper les traces de particules qui se trouvent normalement dans l'atmosphère, elle n'est pas conçue pour traiter des solides, des produits chimiques, des poudres, des solvants, des condensats, des liquides ou d'autres particules. Ils peuvent endommager l'équipement, dégrader ses performances ou raccourcir sa durée de vie.

Un espace de dégagement minimum de 20 cm est conseillé tout autour du système pour permettre une circulation correcte de l'air.

ATTENTION

Les surfaces extérieures de la pompe peuvent être nettoyées avec des détergents doux uniquement.

La pompe à vide sèche ne contient pas de fluide pour nettoyer l'accumulation de poussière et de débris. Faites fonctionner régulièrement la pompe vers l'atmosphère pendant une ou deux minutes pour la nettoyer.

AVERTISSEMENT



Vérifiez avec un électricien ou un spécialiste en entretien qualifié si vous ne comprenez pas l'intégralité des instructions relatives à la mise à la terre ou si vous avez des doutes sur la mise à la terre correcte du produit. Ne modifiez pas la fiche fournie. Si elle ne s'adapte pas à la prise, faites installer une prise correcte par un électricien qualifié.

Branchez le produit uniquement sur une prise dont la configuration est identique à la fiche. N'utilisez pas d'adaptateur avec ce produit.

Procédure de démarrage

AVERTISSEMENT



Retirez les couvercles de bride du refoulement et de l'entrée.



La procédure de démarrage de la pompe à vide à spirales sèches IDP-4 est la suivante:

1. Installez la pompe à l'emplacement prévu et effectuez les connexions mécaniques et à vide nécessaires ou, à défaut, obstruez l'entrée à l'aide d'une bride aveugle.
2. Assurez-vous que la sortie d'échappement n'est pas obstruée ni fermée. Si un tuyau d'échappement ou un silencieux est utilisé, les restrictions de débit ne doivent pas générer une contre-pression supérieure à 6,5 psig.
3. Vérifiez que l'interrupteur principal ON/OFF est en position OFF (« 0 »).
4. Branchez le cordon d'alimentation approprié et connectez-le au secteur. Le cordon doit avoir une capacité minimale de 10 A pour 200 à 240 V et de 15 A pour 100 à 127 V.
5. Mettez la pompe en marche en positionnant l'interrupteur principal sur « I ».
6. Si la pompe est équipée d'une vanne d'admission optionnelle, celle-ci s'ouvrira environ 10 secondes après la mise sous tension.

Il peut s'écouler plusieurs heures avant que la pompe atteigne la pression limite spécifiée. Si cette performance n'est pas atteinte, cela peut être dû à l'humidité accumulée dans le système.

Mode d'emploi

NOTE

La cause la plus fréquente de baisse de performance lors de la première installation ou de la mise en service est la présence de vapeur d'eau adsorbée à l'intérieur de la pompe et de la chambre à vide. Faire fonctionner la pompe avec le ballast de gaz activé peut s'avérer utile si la pompe est restée à l'arrêt pendant une période prolongée, permettant à l'humidité de s'y accumuler. Selon la configuration du système à vide et la quantité d'eau adsorbée, il faut généralement entre 30 minutes et 24 heures de fonctionnement avec le ballast de gaz ouvert pour atteindre les performances de vide nominales.



Figure 8

NOTE

L'interrupteur ON/OFF de la pompe est un interrupteur à bascule, portant les symboles normalisés de la publication IEC 417 indiquant les positions Marche et Arrêt. La figure suivante illustre l'interrupteur en position OFF.

Mode d'emploi

AVERTISSEMENT



Ne jamais insérer d'objet étranger dans les ouïes de ventilation. Tout contact avec les pales du ventilateur peut entraîner des blessures graves ou endommager la pompe.

AVERTISSEMENT



Ne pas obstruer les ouvertures de ventilation. Une obstruction peut provoquer une surchauffe de la pompe. Une température de surface supérieure à 65 °C (150 °F) peut causer des dommages importants. Si une telle condition est observée, éteignez la pompe et laissez-la refroidir avant toute intervention. Procédez ensuite à un démontage et à une inspection, et effectuez les réparations nécessaires.

AVERTISSEMENT



Ne pas utiliser le ballast de gaz lors de l'emploi de substances potentiellement dangereuses ou incompatibles avec l'air.

Procédure d'arrêt

Pour arrêter manuellement la pompe quel que soit son mode de fonctionnement, mettre l'interrupteur principal ON/OFF sur "0". Si la pompe est équipée d'une vanne d'isolement intégrée, elle se fermera rapidement pour éviter le reflux de gaz depuis la pompe vers les composants en amont.

ATTENTION

Avant de procéder à l'arrêt de la pompe IDP-4, assurez-vous que tous les composants interconnectés en amont sont correctement fermés ou préparés pour une exposition potentielle à la pression atmosphérique.

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 9 Logo « DEEE »

Pour plus de précisions, veuillez consulter :

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

Conditions environnementales, conformité et applications pratiques

NOTE

Cette pompe requiert un espace de dégagement d'au moins 20 cm (8 pouces) de tous les côtés pour garantir la ventilation, l'accès pour l'entretien et la facilité d'accès à l'interrupteur d'alimentation. Si la pompe doit être placée sur un banc ou une étagère, ces supports doivent être conçus pour soutenir le poids de la pompe et tout autre équipement associé.

Compatibilité électromagnétique

EN55011/CISPR11

Équipements ISM du groupe 1 : le groupe 1 contient tous les équipements ISM dans lesquels est générée et/ou utilisée intentionnellement de l'énergie de fréquence radio couplée de manière conductrice indispensable au fonctionnement interne de ces équipements.

Les équipements de classe A sont adaptés à l'emploi dans tous les établissements autres que domestiques et concernent ceux qui sont connectés directement à un réseau électrique basse tension qui alimente des bâtiments à usage domestique. Ce dispositif est conforme aux exigences de CISPR11, Groupe 1, Classe A en tant qu'équipement professionnel de radiation. Par conséquent, il est possible que la compatibilité électromagnétique soit difficile à garantir dans d'autres environnements en raison des perturbations conduites et rayonnées.

L'utilisation est soumise aux deux conditions suivantes :

- Ce dispositif ne peut pas provoquer d'interférences dangereuses.
- Ce dispositif doit tolérer toutes les interférences reçues, y compris celles susceptibles de provoquer un fonctionnement indésirable.

Si cet équipement provoque des interférences au niveau de la réception d'appareils de radio ou de télévision, phénomène qui peut être déterminé en éteignant et en rallumant l'équipement, l'utilisateur est invité à prendre une ou plusieurs mesures suivantes :

- Vérifier que les câbles appropriés sont utilisés pour raccorder l'appareil à l'équipement périphérique.
- Tout changement ou modification n'ayant pas été expressément approuvé par Agilent Technologies risque d'annuler l'autorisation de l'utilisateur à exploiter l'équipement.

Mode d'emploi

Déclaration CEM Classe A Corée du Sud

Cet équipement de la Classe A est adapté à un usage professionnel et est conçu pour être utilisé dans des environnements électromagnétiques hors domiciles.

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자과적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

Ce dispositif ISM est conforme à la norme canadienne ICES-001/NMB-001.

Mode d'emploi

Service

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

vpt-customer@agilent.com

vpl-customer@agilent.com

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

4

Instrucciones de uso

Acerca del manual	89
Documentos justificativos pertinentes	89
Definiciones y terminología	90
Definición de Precaución, Advertencia y Nota	90
Símbolos de advertencia	91
Seguridad	92
Advertencias CEM	93
FCC	93
Uso adecuado	94
Uso incorrecto razonablemente previsible	94
Directrices de seguridad para bombas de vacío de desplazamiento en seco	95
Personal cualificado	96
Equipo de protección	97
Descripción del producto	99
Transporte y almacenamiento	99
Desembalaje e Inspección	100
Preparación para la instalación	101
Procedimiento de puesta en marcha	105
Procedimiento de parada	108
Eliminación	109
Condiciones medioambientales, conformidad y requisitos de los servicios públicos	110
Compatibilidad electromagnética	110
EN55011/CISPR11	110
Declaración EMC Clase A de Corea del Sur	111
CIEM/NMB-001	111
Servicio	112

Acerca del manual

Este manual contiene instrucciones para los usuarios e información técnica relacionada con la bomba IDP-4 Dry Scroll de acuerdo con las directivas aplicables.

Este documento es válido para el producto fabricado en el momento de esta publicación, siempre que no se realicen cambios en el producto por parte del cliente o de terceros. En caso de cualquier información contradictoria, la documentación enviada con el producto debe considerarse como los documentos válidos.

NOTA

- 1 Este manual contiene información importante para que todo el personal que utilice la bomba IDP-4 Dry Scroll pueda operarlas con seguridad y garantizar su rendimiento durante toda su vida útil.
- 2 Conserve este manual, junto con todas las publicaciones relacionadas, en un lugar accesible y conocido por todos los operarios/personal de mantenimiento.

Documentos justificativos pertinentes

X3802-90000 Declaración de conformidad

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 paos específicos.





Símbolos de advertencia

La siguiente es una lista de símbolos que aparecen junto con las advertencias en la bomba de espiral seca IDP-4. 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. 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.



	Certificación CE
	Certificación CSA
	Certificación China RoHS
	Equipo de Desecho Eléctrico y Electrónico
	Certificación RCM
	Lugar de fabricación
	Superficie caliente
	Voltajes peligrosos
	Peligro genérico

Instrucciones de uso

	Terminal protectora de tierra
	Corriente alterna (CA)
	Encendido
	Apagado

Seguridad

Esta sección contiene la información, prescrita por la Directiva de Baja Tensión 2006/42/EC, que es esencial para el cumplimiento y la observancia de las normas de seguridad tanto en general como en relación con el uso específico del producto.

El incumplimiento de estas instrucciones y de las demás instrucciones contenidas en este manual puede afectar a la seguridad y podría provocar accidentes a las personas que utilicen el producto.

Agilent Technologies declina toda responsabilidad por daños al producto o a la seguridad física del operador o de terceros derivados de la inobservancia de las normas de seguridad indicadas en la documentación técnica.

Advertencias CEM

EN 55011 Advertencia de clase A

Este es un producto de Clase A. En un entorno doméstico, este producto puede causar interferencias de radio, en cuyo caso el usuario puede verse obligado a tomar las medidas adecuadas.

FCC

Este dispositivo cumple la Parte 15 de las normas de la FCC. Su funcionamiento está sujeto a las dos condiciones siguientes: (1) Este dispositivo no puede causar interferencias perjudiciales, y (2) este dispositivo debe aceptar cualquier interferencia recibida, incluidas las interferencias que puedan causar un funcionamiento no deseado.

NOTA

El equipo ha sido probado y cumple los límites establecidos para los dispositivos digitales de Clase A, de conformidad con la Parte 15 de las normas de la FCC. Estos límites están diseñados para proporcionar una protección razonable contra interferencias perjudiciales cuando el equipo se utiliza en un entorno comercial. Este equipo genera, utiliza y puede irradiar energía de radiofrecuencia y, si no se instala y utiliza de acuerdo con el manual de instrucciones, puede causar interferencias perjudiciales en las comunicaciones por radio. El funcionamiento de este equipo en una zona residencial también puede causar interferencias perjudiciales en las comunicaciones de radio, en cuyo caso el usuario deberá corregir la interferencia por su cuenta.

Uso adecuado

La bomba de espiral seca IDP-4 está destinada exclusivamente al uso por parte de personas cualificadas como bomba de vacío de acuerdo con este manual.

Cualquier accesorio y equipo utilizado con el producto debe ser suministrado o aprobado por Agilent Technologies.

Agilent Technologies no ofrece ninguna garantía ni asume ninguna responsabilidad por aplicaciones distintas de las descritas en este manual o en las que no se respeten los requisitos esenciales y las medidas de seguridad.

Uso incorrecto razonablemente previsible

Agilent Technologies declina toda responsabilidad derivada del uso indebido y los usos incorrectos previsible de la bomba IDP-4 Dry Scroll. El uso indebido hará que se pierdan todas las reclamaciones de responsabilidad y garantías.

El uso indebido puede incluir, pero no se limita a:

- Conexión de la bomba a un suministro de tensión incorrecto
- Conexión a tierra incorrecta o inexistente
- Funcionamiento continuo por encima de la presión de entrada recomendada
- Funcionamiento con sobrepresión en la entrada de gas lastre
- Funcionamiento del gas lastre abierto a la atmósfera cuando se bombean gases peligrosos
- Funcionamiento con temperatura ambiente excesivamente alta
- Funcionamiento con temperatura ambiente inferior a la especificada
- Daños debidos a un almacenamiento y transporte inadecuados de la bomba
- Realización de tareas de mantenimiento o retirada de carenados sin desconectar la alimentación eléctrica
- Funcionamiento sin elementos de flujo de aire (cubierta o ventilador)
- Envío del producto devuelto en un embalaje que no es el de fábrica.

Directrices de seguridad para bombas de vacío de desplazamiento en seco

El personal responsable del funcionamiento y mantenimiento de la bomba debe estar debidamente formado y conocer las normas de prevención de accidentes (consulte la sección Personal cualificado). Las precauciones de prevención de accidentes contenidas en este manual deben respetarse continuamente durante el funcionamiento y mantenimiento de la bomba para evitar lesiones a los operarios y daños a la bomba u otros equipos. Estas precauciones se proporcionan en forma de notas de ADVERTENCIA y PRECAUCIÓN a lo largo de todo el manual del usuario.

ADVERTENCIA



Para evitar daños al equipo y lesiones al personal de servicio, deben seguirse estrictamente las instrucciones de instalación que figuran en este manual!

ADVERTENCIA



**La bomba no debe manejar materiales inflamables.
Está prohibido instalar la bomba en un entorno explosivo.**

Personal cualificado

Se recomienda que bombas IDP-4 Dry Scroll sean utilizadas únicamente por personal cualificado. Una persona cualificada es:

- Un individuo que posee las cualificaciones adecuadas y tiene experiencia, o ha completado una formación, en el trabajo con bombas de Vacío Agilent.
- Una persona que sigue las medidas de seguridad necesarias y maneja la bomba en las condiciones adecuadas para no causar lesiones o daños.
- Una persona que conozca y esté debidamente cualificada para seguir las normas y reglamentos nacionales y locales aplicables en materia de seguridad y reglamentación.

NOTA

Cualquier reparación en este producto, más allá de las indicadas como mantenimiento estándar en este manual de usuario, debe ser realizada exclusivamente por personal autorizado por Agilent. Cualquier operación de ajuste, modificación o mantenimiento debe ser realizada por un técnico profesional informado de los riesgos.

Equipo de protección

El equipo de protección de los operarios que manejan o realizan el mantenimiento de la bomba debe ser siempre el adecuado para el tipo de operación que se está ejecutando. El equipo de protección utilizado debe cumplir los requisitos de seguridad de la legislación vigente en el país en el que se utilice la máquina.

En general, el operario debe llevar calzado de seguridad mientras mueva, manipule o instale la bomba de espiral seca IDP-4..

ADVERTENCIA



Peligro para la salud debido a la presencia de sustancias peligrosas durante el mantenimiento o la instalación.

- Dependiendo de la peculiaridad del proceso, la bomba de vacío, los componentes o los fluidos de funcionamiento pueden estar contaminados con sustancias tóxicas, reactivas o radioactivas.
- Use el equipo de protección adecuado durante el mantenimiento y las reparaciones o en caso de volver a instalar el equipo.

ADVERTENCIA

Riesgo de lesiones debido a la caída de objetos



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

- Traslade la bomba de vacío de tamaño pequeño y mediano con las dos manos.
- Todo equipo de más de 20 kilos de peso debe ser transportado con un dispositivo de elevación adecuado.
- Use zapatos de seguridad con puntera de acero que cumplan con la Directiva EN 347.

PRECAUCIÓN

Riesgo de lesiones debido a la presencia de cantos vivos



- Antes de realizar el servicio técnico o antes del montaje/desmontaje de la bomba del sistema, espere a que la bomba se detenga por completo.
 - No trabaje directamente dentro de la brida de alto vacío.
 - Si es necesario, use guantes de protección que cumplan con la norma EN 420.
-

Descripción del producto

La bomba IDP-4 es una bomba de vacío de desplazamiento en seco hermética. Esta bomba es adecuada para bombear aire o gases inertes desde una zona de baja presión a una zona de mayor presión que no esté sustancialmente por debajo o por encima de la presión atmosférica. Esta bomba no está diseñada para bombear gases corrosivos, pirofóricos, formadores de partículas o mezclas de gases en su rango inflamable y/o combustible, ni está diseñada para utilizarse como compresor para alcanzar presiones de descarga sustancialmente superiores a la presión atmosférica.

El manual y los documentos relacionados contienen la información necesaria para garantizar la seguridad del operador al utilizar el equipo. Encontrará información detallada en la sección "Technical Information".

Transporte y almacenamiento

Al transportar y almacenar la bomba, no deben superarse los siguientes requisitos ambientales:

- Temperatura: -20 °C a +60 °C (-4 °F a 140 °F)
- Humedad relativa: 0 a 95% (sin condensación)

NOTA

Debe tenerse en cuenta que durante el transporte y almacenamiento de la bomba de vacío, el vapor de agua del ambiente penetrará en el interior de la bomba de vacío independientemente de lo bien sellada que esté. Para alcanzar el pleno rendimiento de vacío nominal de la bomba, este vapor de agua debe eliminarse utilizando las características de lastre de gas descritas en la sección "Technical Information".

Desembalaje e Inspección

La bomba se suministra en un embalaje protector especial. Si éste muestra signos de daños, que pueden haberse producido durante el transporte, póngase en contacto con su oficina de ventas local. El peso total de la bomba y el embalaje juntos es de aproximadamente 10 kg (20 libras).

ADVERTENCIA



La caja de cartón y todos los materiales de embalaje serán necesarios en caso de servicio en un centro de servicio autorizado por Agilent. Guardar este embalaje facilitará cualquier solicitud de servicio.

ADVERTENCIA



Al desembalar la bomba, asegúrese de no dejarla caer y evite cualquier tipo de impacto o golpe brusco.

Preparación para la instalación

La bomba de espiral seca IDP-4 está diseñada y certificada para funcionar en interiores (categoría de instalación II) en entornos relativamente limpios y secos (grado de contaminación 2).

Consulte las especificaciones en la Tabla 1 de la sección Información técnica.

Esta bomba no está diseñada ni certificadas para funcionar en exteriores.

La bomba de espiral seca IDP-4 no debe utilizarse nunca en ambientes que contengan materiales explosivos o combustibles. La bomba de espiral seca IDP-4 NO es intrínsecamente segura y nunca deben utilizarse con gases corrosivos, pirofóricos, formadores de partículas o mezclas de gases en su rango inflamable y/o combustible.

Coloque la bomba sobre una superficie sólida y nivelada. Si debe montarse sobre un banco o una mesa, asegúrese de que el banco o la mesa están preparados para soportar el peso de la bomba y de todo el equipo asociado.

Conecte la bomba a la fuente de alimentación utilizando un cable de alimentación tipo IEC-320 de al menos 10A de capacidad.

NOTA

La causa más común de un rendimiento insuficiente durante la instalación y puesta en marcha iniciales es la presencia de vapor de agua adsorbido tanto en la bomba como en la cámara de vacío. Si la bomba ha estado apagada durante un tiempo, lo que ha permitido que se acumule agua en su interior y en el sistema de vacío, puede resultar beneficioso ponerla en marcha con el lastre de gas activado. Dependiendo del sistema de vacío y de la cantidad de agua adsorbida, normalmente se necesitan entre 30 minutos y 24 horas con el lastre de gas activado para alcanzar el máximo rendimiento de vacío.

PRECAUCIÓN

Durante el funcionamiento deben respetarse las siguientes condiciones ambientales:

- **Temperatura: +5 °C a +40 °C (41 °F a 104 °F)**
- **Humedad relativa: 5 a 95% (sin condensación)**

Conecte la bomba a la fuente de alimentación utilizando un cable de alimentación de tipo IEC-320 con la intensidad y tensión nominales adecuadas (consulte la Tabla 2).

Instrucciones de uso

PRECAUCIÓN Al instalar la bomba de vacío, se recomienda colocarla de forma que el interruptor de alimentación y la conexión del cable de alimentación queden fácilmente accesibles al usuario para poder desconectarla rápidamente en caso necesario.

PRECAUCIÓN No apriete, afloje ni modifique de ninguna manera los dos tornillos de cabeza hexagonal situados a cada lado del cabezal de la bomba (véase la figura inferior). Esto provocará una pérdida de rendimiento o daños en la bomba.



Figura 10

ADVERTENCIA



No retire ni modifique ningún equipo de seguridad o aislamiento de la bomba. Si lo hace, podría crear un grave peligro para la seguridad y anular la garantía.

ADVERTENCIA



- Esta bomba sólo puede bombear y expulsar aire y gases inertes. Esta bomba no está diseñada para bombear gases corrosivos, pirofóricos, formadores de partículas o mezclas de gases en su rango inflamable y/o combustible.
 - Instálela en una zona que no esté expuesta a la lluvia, vapor o humedad excesiva. Pueden provocar descargas eléctricas, cortocircuitos y lesiones corporales graves.
 - Antes de inspeccionar, reconfigurar o dar servicio a la bomba, asegúrese de que el suministro eléctrico esté desconectado.
 - El balasto de gas debe sellarse siempre que se bombee cualquier gas que no pueda ventilarse a la atmósfera o mezclarse con aire.
 - Se dispone de un silenciador de escape opcional para aquellas aplicaciones en las que se desee disminuir el nivel sonoro producido por la bomba.
 - Para evitar lesiones corporales, evite exponer cualquier parte del cuerpo humano al vacío.
 - La bomba debe instalarse con fácil acceso al interruptor de alimentación.
 - Los cables de alimentación aprobados por Agilent para todas las conexiones comunes se enumeran en la sección de Conexiones Eléctricas. Si se requiere un cable de alimentación diferente, la corriente nominal del mismo debe ser de al menos 10 A para servicio de 200–240 V y de al menos 15 A para servicio de 100–127 V.
-

Instrucciones de uso

Instrucciones de conexión a tierra

ADVERTENCIA



Este producto debe conectarse a tierra. En caso de cortocircuito eléctrico, la conexión a tierra reduce el riesgo de descarga eléctrica al proporcionar una vía de escape para la corriente eléctrica. Esta bomba está equipada con un cable de alimentación que tiene un cable de conexión a tierra con un enchufe de conexión a tierra apropiado. El enchufe debe insertarse en una toma de corriente que esté correctamente instalada y conectada a tierra de acuerdo con todos los códigos y ordenanzas locales.

El conector del cable de alimentación debe estar correctamente enchufado en el módulo de alimentación de la bomba para evitar una conexión defectuosa o intermitente.

PRECAUCIÓN

Aunque la bomba puede bombear trazas de partículas que normalmente se encuentran en la atmósfera, no está diseñada para procesar sólidos, productos químicos, polvos, disolventes, condensados, líquidos u otras partículas. Pueden dañar el equipo, degradar su rendimiento o acortar su vida útil.

Se recomienda dejar un espacio libre mínimo de 20 cm alrededor del sistema para permitir una circulación de aire adecuada.

PRECAUCIÓN

Las superficies exteriores de la bomba sólo pueden limpiarse con detergentes suaves.

La bomba de desplazamiento en seco no contiene líquido para la limpieza del polvo y los residuos acumulados. Haga funcionar la bomba periódicamente a la atmósfera durante uno o dos minutos para enjuagarla.

ADVERTENCIA



Consulte a un electricista o técnico cualificado si no comprende completamente las instrucciones de conexión a tierra, o si tiene dudas sobre si el producto está correctamente conectado a tierra. No modifique el enchufe suministrado. Si no se adapta a la toma de corriente, solicite a un electricista cualificado que instale la toma de corriente adecuada.

Conecte el producto únicamente a una toma de corriente que tenga la misma configuración que el enchufe. No utilice adaptadores con este producto.

Procedimiento de puesta en marcha

ADVERTENCIA



Retire las tapas de las bridas del escape y de entrada.



El procedimiento de puesta en marcha de la bomba de vacío de espiral seca IDP-4 es el siguiente:

1. Instale la bomba en la posición deseada y realice las conexiones mecánicas y de vacío correspondientes o, en alternativa, cubra la entrada con una brida ciega.
2. Asegúrese que la conexión de escape no esté bloqueada ni obstruida. Si utiliza un tubo de escape o un silenciador, las restricciones de escape no deben causar una contrapresión superior a 6,5 psig.
3. Asegúrese que el interruptor de encendido/apagado esté en la posición de apagado "0".
4. Conecte el cable de alimentación correspondiente y enchúfelo a la toma de corriente. La corriente nominal del cable de alimentación debe ser de al menos 10 A para una tensión de 200-240 V y de al menos 15 A para una tensión de 100-127 V.
5. Posicione el interruptor principal de encendido/apagado a la posición "I".
6. Si la bomba está equipada con una válvula de entrada opcional, la válvula se abrirá aproximadamente 10 segundos después de encenderla.

La bomba puede tardar varias horas en alcanzar la presión máxima especificada. Si no alcanza la presión máxima esperada, puede deberse a la acumulación de vapor de agua en su interior.

Instrucciones de uso

NOTA

La causa más común de un rendimiento insuficiente durante la instalación y puesta en marcha iniciales es la presencia de vapor de agua adsorbido tanto en la bomba como en la cámara de vacío. Activar el lastre de gas puede resultar beneficioso, incluso si la bomba ha estado apagada durante un tiempo, ya que permite la adsorción del agua tanto dentro de la bomba como en el sistema de vacío. Dependiendo del sistema de vacío y de la cantidad de agua adsorbida, alcanzar el rendimiento máximo de vacío con el lastre de gas activado suele requerir entre 30 minutos y 24 horas.



Figura 11

NOTA

El interruptor de encendido/apagado de la bomba es un interruptor basculante con símbolos que, según la publicación IEC 417, representan las posiciones de encendido y apagado. La figura anterior muestra un interruptor en la posición de apagado.

ADVERTENCIA



No introduzca ningún objeto extraño a través de las rejillas de ventilación, ya que el contacto con las aspas del ventilador puede provocar lesiones personales graves o daños en la bomba.

ADVERTENCIA



No obstruya las rejillas de ventilación. Si las obstruye, la bomba podría sobrecalentarse. Una temperatura en la superficie de la bomba superior a 65 °C (150 °F) puede ser perjudicial. Si observa estas condiciones, apague la bomba y déjela enfriar. Desmonte la bomba, inspecciónela para detectar posibles daños y repárela si es necesario.

ADVERTENCIA



No utilice lastre de gas si utiliza materiales potencialmente peligrosos o que no deban mezclarse con el aire.

Procedimiento de parada

Para apagar la bomba manualmente en cualquier modo de funcionamiento, gire el interruptor principal ON/OFF a la posición «0». Si la bomba está equipada con una válvula de aislamiento integrada, se cerrará rápidamente para eliminar el reflujos de gas de la bomba a los componentes aguas arriba.

PRECAUCIÓN

Antes de apagar la bomba IDP-4, asegúrese de que todos los componentes interconectados aguas arriba estén correctamente cerrados o preparados para una posible exposición a la presión atmosférica.

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 12 Logotipo “WEEE”

Para más información, consulte:

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

Condiciones medioambientales, conformidad y requisitos de los servicios públicos

NOTA

Esta bomba requiere un espacio libre de al menos 20 cm (8 pulg.) en todos los lados para ventilación, acceso para mantenimiento y fácil acceso al interruptor de alimentación. Si se desea colocar la bomba sobre un banco o estante, éste debe estar dimensionado para soportar el peso de la bomba y de cualquier otro equipo asociado.

Compatibilidad electromagnética

EN55011/CISPR11

Equipos ISM del grupo 1: El Grupo 1 contiene todos los equipos ISM en los que se genera y/o utiliza intencionalmente energía de radiofrecuencia acoplada conductivamente que es necesaria para el funcionamiento interno del propio equipo.

Los equipos de Clase A son equipos aptos para su uso en todos los establecimientos que no sean domésticos y aquellos conectados directamente a una red de suministro de energía de bajo voltaje que alimenta edificios utilizados para fines domésticos.

Este dispositivo cumple con los requisitos de CISPR11, Grupo 1, Clase A como equipo profesional de radiación. Por lo tanto, puede haber dificultades potenciales para garantizar la compatibilidad electromagnética en otros entornos, debido tanto a las perturbaciones conducidas como a las radiadas.

La operación está sujeta a las siguientes dos condiciones:

- Este dispositivo no puede causar interferencias nocivas.
- Este dispositivo debe aceptar cualquier interferencia recibida, incluida la interferencia que pueda provocar un funcionamiento no deseado.

Si este equipo causa interferencias dañinas en la recepción de radio o televisión, lo cual se puede determinar apagando y encendiendo el equipo, se recomienda al usuario que pruebe una o más de las siguientes medidas:

- Asegúrese de utilizar los cables apropiados para conectar el dispositivo al equipo periférico.
- Los cambios o modificaciones no aprobados expresamente por Agilent Technologies podrían anular la autorización del usuario para operar el equipo.

Instrucciones de uso

Declaración EMC Clase A de Corea del Sur

Este equipo es Clase A apto para uso profesional y ha sido diseñado para su uso en entornos electromagnéticos fuera del hogar.

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자과적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

CIEM/NMB-001

Este dispositivo ISM cumple con la norma ICES-001/NMB-001 de Canadá.

Servicio

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

vpt-customer@agilent.com

vpl-customer@agilent.com

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

5

使用说明

關於本手冊	114
相關證明文件	114
定義與術語	115
小心、警告和注意的定義	115
警告符號	116
安全	117
EMC 警告	118
美國聯邦通訊委員會	118
正確使用	119
合理而可預見的濫用	119
乾式渦旋幫浦安全指南	120
合資格人員	121
防護裝備	122
產品描述	124
運輸及儲存	124
開箱檢驗	125
安裝準備	126
啟動程式	130
關機程式	133
處置	134
環境條件、合規性和公用設施要求	135
電磁兼容	135
EN55011/CISPR11	135
韓國 A 類 EMC 聲明	136
ICES/NMB-001	136
服務	137

關於本手冊

本手冊包含根據適用指令對 IDP-4 乾式渦旋泵的使用者說明和技術資訊。

本文件適用於發佈時生產的產品，前提是客戶或第三方未對產品進行更改。如果存在任何衝突訊息，則應將產品附帶的文件視為有效文件。

注意

- 1 本手冊包含重要訊息，以便所有使用 IDP-4 乾式渦旋泵的人員都可以操作 它們安全並保證其在整個使用壽命期間的性能。
 - 2 將本手冊以及所有相關出版物放在所有操作員/維護人員都知道的地方。
-

相關證明文件

X3802-90000 符合性聲明

定義與術語

小心、警告和注意的定義

本手冊的一些重要參考內容以背景色突出顯示。

小心

程序開始前，提示小心訊息，如果不遵守製程指示，可能導致設備損壞。

警告



警告訊息提醒操作員注意特定流程或做法，如果執行錯誤，可能導致人員重傷。

注意

注意旨在引起注意重要信息，提供具體步驟的更多詳細信息。

警告符號

以下是 IDP-4 乾燥渦旋泵上與警告一起出現的符號清單。也顯示了這些符號所指的危險。三角形符號表示警告。在文件或儀器警告旁邊可能出現的符號意義如下。以下符號可用作附在儀器上的警告標籤。當您看到此符號時，請參閱相關操作或維修手冊，以了解該警告標籤所指的正確操作流程。



	CE認證
	CSA認證
	RoHS中國認證
	廢棄電氣和電子設備
	RCM認證
	生產現場
	表面較熱
	表示危險電壓
	一般危險
	保護接地端子
	交流電 (交流)
	電源開關
	電源關閉

安全

本節包含低電壓指令 2006/42/EC 規定的信息，這些信息對於遵守一般安全法規和與產品特定用途相關的安全法規至關重要。

不遵守這些說明和本手冊中包含的其他說明可能會影響安全，並可能導致產品操作人員發生事故。

對於因不遵守技術文件中規定的安全規則而導致的產品損壞或操作員或第三方的
人身安全，Agilent Technologies 不承擔任何責任。

使用说明

EMC 警告

EN 55011 A 級警告

這是A級產品。在家庭環境中，本產品可能會造成無線電幹擾，在這種情況下，使用者可能需要採取適當的措施。

美國聯邦通訊委員會

本設備符合 FCC 規則第 15 部分的規定。操作須符合以下兩個條件：(1) 該設備不會造成有害幹擾，和(2)該設備必須接受收到的任何干擾，包括可能導致不良操作的干擾。

注意

該設備經過測試，符合 FCC 規則第 15 部分對 A 類數位設備的限制。這些限制旨在提供合理的保護，防止設備在商業環境中運作時產生有害幹擾。該設備產生、使用並可能輻射射頻能量，如果不按照說明手冊安裝和使用，可能會對無線電通訊造成有害幹擾。在住宅區操作設備也可能造成有害的無線電通訊幹擾，在這種情況下，使用者將需要自費糾正幹擾。

使用说明

正確使用

IDP-4 乾式渦旋幫浦僅供合資格人員依本手冊用作真空幫浦。

與產品一起使用的任何配件和設備必須由安捷倫科技公司提供或批准。

對於本手冊中所述的應用以外的應用或未遵守基本要求和安全措施的應用，安捷倫科技公司不提供任何保證或承擔任何責任。

合理而可預見的濫用

安捷倫科技公司不承擔因 IDP-4 乾式渦旋泵的不當使用和可預見的誤用而產生的所有責任。不當使用將導致所有責任索賠和保證被沒收。

不當使用可能包括但不限於：

- 將幫浦連接到錯誤的電源電壓
- 接地連接不當或缺失
- 在高於建議入口壓力的情況下連續運行
- 氣鎮入口超壓運行
- 泵送危險氣體時，操作氣鎮與大氣相通
- 環境溫度過高時運行
- 在低於規格的環境溫度下運行
- 泵浦儲存和運輸不當造成的損壞
- 在不斷開主電源的情況下進行維護或拆卸整流罩
- 在沒有气流元件（罩或風扇）的情況下運行。
- 使用非工廠包裝運送退回的產品。

乾式渦旋幫浦安全指南

負責泵浦操作和維護的人員必須接受適當的培訓，並且必須了解事故預防規則（請參閱合格人員部分）。在泵浦的操作和維護過程中，必須始終遵守本手冊中包含的事故預防預防措施，以避免操作人員受傷以及泵浦或其他設備損壞。這些預防措施以「警告」和「注意」註釋的形式貫穿整本使用手冊。

警告



為避免設備損壞和操作人員受傷，請嚴格遵守本手冊中的安裝說明！

警告



泵不得處理易燃材料。
禁止將幫浦安裝在爆炸性環境中。

合資格人員

IDP-4 乾式渦旋幫浦僅建議由合格人員操作。合格人員是指：

- 具有適當資格並具有使用安捷倫乾真空幫浦的經驗或已完成相關訓練的人員。
- 遵循必要的安全措施並在適當的條件下操作幫浦以免造成傷害或損壞的人員。
- 熟悉並有資格遵守任何適用的國家和地方安全和監管標準及附則的個人。

注意

除本使用手冊中列出的標準維護之外的任何對本產品的維修都必須由 Agilent 授權人員專門進行。任何調整、改造或維護作業必須由了解風險的專業技術人員進行。

使用說明

防護裝備

操作或維護幫浦的操作員的防護設備必須始終適合所執行的操作類型。使用的防護設備必須符合機器使用國現行法律的安全要求。

一般來說，操作員在移動、搬運或安裝IDP-4 幹渦旋泵時必須穿著安全鞋。

警告



維護或安裝時危險物質會導致健康危險。



- 根據製程特性，真空幫浦、組件或工作流體可能有毒、反應性或原料污染。
- 維護維修或重新安裝時，請配戴適當的防護設備。



警告



物體造成損壞的風險



手動運送真空幫浦時，有重物緩慢且不易發生的危險。



- 手工搬運中小型真空幫浦。
- 若設備重量超過 20 公斤，應使用適當的起吊設備運送。
- 穿著符合指令 EN 347，並搭配鐵鞋頭的安全鞋。

小心



鋒利邊緣導致風險受損

- 在維修幫浦之前，或在幫浦從系統執行任何安裝/操作之前，等待幫浦完全停止。
 - 不要直接在高空法蘭內操作。
 - 如有必要，請依照 EN 420 戴上防護手套。
-

產品描述

IDP-4 是一种密封式干式涡旋泵。本泵適用於將空氣或惰性氣體從低壓區輸送到高壓區，但高壓區的壓力不得顯著低於或高於大氣壓力。本泵浦不適用於輸送腐蝕性、自燃性、顆粒生成性氣體或處於易燃和/或可燃範圍內的氣體混合物，也不適用於用作壓縮機以獲得顯著高於大氣壓力的排氣壓力。

本手冊及相關文件包含保證操作人員使用設備時的安全所需的資訊。「技術資訊」部分提供了詳細資訊。

運輸及儲存

泵浦的運輸和儲存時，不得超過下列環境要求：

- 溫度： -20°C 至 +60°C (-4°F 至 140°F)
- 相對濕度： 0 至 95% (非冷凝)

注意

應該認識到，在真空幫浦的運輸和儲存過程中，無論幫浦的密封性如何，周圍環境中的水蒸氣都會滲透到真空幫浦內部。為了實現泵浦的全部額定真空性能，必須使用技術資訊部分中所述的氣鎮功能來去除水蒸氣。

開箱檢驗

該幫浦採用特殊的保護包裝提供。如果出現損壞跡象（可能是在運送過程中發生），請聯絡您當地的銷售辦事處。幫浦和包裝的總重量約為 10 公斤（20 磅）。

警告



在安捷倫認可的服務中心進行維修時，需要使用紙箱和所有包裝材料。保存此包裝將方便任何服務請求。

警告



打開幫浦包裝時，請務必不要將其跌落並避免任何突然的撞擊或震動。

安裝準備

IDP-4干式涡旋泵专为在相对清洁、干燥的室内环境中运行而设计并通过认证（安装类别III，污染等级2）。

请参阅技术信息部分的表1以获取规格说明。

这些泵未被设计或认证用于室外运行。

IDP-4干式涡旋泵绝不能在含有爆炸性或可燃物质的环境中使用。

IDP-4泵不具备本质安全性，并且绝不应与腐蚀性、自燃性、能形成颗粒的气体或处于其可燃和/或易燃范围内的气体混合物一起使用。

将泵放置在坚固、平整的表面上。

如果必须安装在工作台或桌面上，请确保该工作台或桌面能够承受泵及所有相关设备的重量。

使用额定电流至少为10安培的IEC-320型电源线将泵连接至电源。

注意

在初次安装和启动时，性能不佳的最常见原因是泵和真空腔内吸附的水蒸气。

如果泵长时间停机，导致水分在泵和真空系统内部被吸附，那么在运行时启用气镇功能也可能有助于改善性能。

根据真空系统的情况和吸附水分的多少，通常需要启用气镇约 30 分钟至约 24 小时才能达到最佳真空性能。

小心

操作期間必須遵守以下環境條件：

- 溫度：+5°C 至 +40°C（41°F 至 104°F）
- 相對濕度：5% 至 95%（非冷凝）

使用適當的 IEC-320 型電源線將幫浦連接到電源 額定電流和電壓（見表 2）。

小心

安裝真空幫浦時，建議將幫浦放置在使用者易於觸及電源開關和電源線連接的位置，以便在需要時快速斷開電源幫浦。

使用说明

小心

请勿拧紧、松开或以任何方式改动泵头两侧的两个六角螺栓（见下图）。这样做会导致性能下降和/或泵损坏。



圖 13

警告



請勿拆除或改裝幫浦上的任何安全或絕緣設備。
這樣做可能會造成嚴重的安全隱患，也可能導致保固失效。

警告



- 此幫浦只能抽吸和排出空氣和惰性氣體。此幫浦不適用於泵送腐蝕性、自燃性、形成顆粒的氣體或易燃和/或可燃範圍內的氣體混合物。
 - 安裝在不會暴露於雨水、蒸氣或過度潮濕的區域。它們可能導致觸電、短路和嚴重的人身傷害。
 - 在檢查、重新配置或維修幫浦之前，請確保電源已中斷。
 - 每當泵送任何不能排放到大氣中或與空氣混合的氣體時，必須密封氣鎮。
 - 對於需要降低幫浦產生的噪音水平的應用，可以選擇排氣消音器。
 - 為防止身體受傷，請避免將人體的任何部位暴露在真空中。
 - 幫浦的安裝位置必須能夠方便地觸及電源開關。
 - 所有常用连接所使用的 Agilent 认可电源线均在“电气连接”部分中注明。如果必须使用不同的电源线，对于 200–240V 电源，电源线的额定电流必须至少为 10A；对于 100–127V 电源，电源线的额定电流必须至少为 15A。。
-

使用说明

接地說明

警告



本產品應接地。如果發生電氣短路，接地可以為電流提供逃逸路徑，從而降低電擊風險。此幫浦配備有電源線，電源線帶有接地線和適當的接地插頭。插頭必須插入根據所有當地法規和條例正確安裝並接地的插座。

電源線連接器必須正確插入幫浦的電源模組，以防止連接故障或間歇性連接。

小心

儘管該泵可以泵送大氣中常見的微量顆粒，但它並非設計用於處理固體、化學品、粉末、溶劑、冷凝物、液體或其他顆粒。它們會損壞設備、降低其性能或縮短其使用壽命。

建議系統周圍至少留出 20 公分的自由空間，以達到適當的空氣流通。

小心

幫浦的外表面只能用溫和的清潔劑清潔。

乾式渦旋泵不含用於清除積聚灰塵和碎屑的液體。定期在大氣中運轉泵一兩分鐘以將其沖洗掉。

警告



如果您不完全理解接地說明，或者您對產品是否正確接地有疑問，請諮詢合格的電工或維修人員。請勿改裝所提供的插頭。如果與插座不匹配，請由合格的電工安裝合適的插座。

僅將產品連接至與插頭配置相同的插座。請勿將轉接器與本產品一起使用。

啟動程式

警告



移除排气口和进气口的法兰盖。



IDP-4干式涡旋真空泵的启动步骤如下:

1. 将泵安装在指定位置，并进行相应的机械和真空连接，或者用盲法兰盖封闭进气口。
2. 确保排气接头未被堵塞或封闭。若使用排气软管或消音器，排气阻力不得导致背压超过6.5 psig。
3. 确保电源开关处于关闭（“0”）位置。
4. 连接合适的电源线，并将电源线插入主电源插座。 电源线的额定电流对于200–240V电源应至少为10A，对于100–127V电源应至少为15A。
5. 将主电源开关按至“1”位置。
6. 如果泵配有可选进气阀，则在通电约10秒后阀门将自动打开。

泵可能需要数小时才能达到公布的极限真空压力规格。
如果泵未能达到预期的极限压力，原因可能是泵内积聚了水蒸气。

使用说明

注意

在初次安装和启动时，性能不佳的最常见原因是泵和真空腔内吸附的水蒸气。

如果泵长时间停机，导致水分在泵和真空系统内部被吸附，那么在运行时启用气镇功能也可能有助于改善性能。

根据真空系统的状况和吸附水分的多少，通常需要启用气镇约 30 分钟至 24 小时才能达到完整的真空性能。



圖 14

注意

泵的开/关开关为摇臂式开关，其上标有符合 IEC 第 417 号标准的符号，用于表示开启和关闭位置。下图显示的是处于关闭位置的开关。

警告



请勿将任何异物插入冷却通风口，因为与风扇叶片接触可能导致严重的人身伤害或泵损坏。

警告



请勿堵塞冷却通风口。堵塞这些通风孔可能会导致泵过热。泵表面温度超过 65 °C (150 °F) 可能会造成损坏。如果出现此类情况，请关闭泵并让其冷却。必要时应拆卸泵、检查是否损坏并进行维修。

警告



如果使用的材料具有潜在危险性或不应与空气混合，请勿使用气镇。

關機程式

若要在任何操作模式下手動關閉泵，請將主 ON/OFF 開關轉到“0”位置。如果泵浦配備了整合式隔離閥，它將快速關閉，以消除氣體從泵浦回流到上游組件。

小心

在关闭 IDP-4 泵之前，确保所有上游互连部件已正确关闭阀门，或已做好可能暴露于大气压力的准备。

處置

標籤中的「WEEE」確實意義重大

根據EC WEEE（廢電子電氣設備）指令應用以下符號。

此符號（僅在歐盟國家有效）表示所適用的產品不能與家庭或工業或一起丟棄，必須送至專門的物資回收系統。銷售合約條款後啟動回收和支付流程。



圖 15 “WEEE” 標誌

有關更多信息，請參考：

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

環境條件、合規性和公用設施要求

注意

此幫浦需要在所有側面留出至少 20 公分（8 英吋）的淨空間，以便通風、維護和方便操作電源開關。如果需要將幫浦放置在長凳或架子上，則其額定值必須能夠支撐幫浦和任何其他相關設備的重量。

電磁兼容

EN55011/CISPR11

第1組 ISM設備：第1組包含所有 ISM設備，其中包含擷取產生和/或使用設備本身內部功能所必需的連接連接RF狀態。

A類設備 適用於除住宅和直接連接到為民用建築供電的低壓供電網路的場所以外的所有場所的設備。

本設備符合 CISPR11第 1 因此，由於幹擾和輻射干擾，在其他環境中確保電磁相容性可能存在潛在障礙。

操作需滿足以下兩個條件：

- 該設備不會造成危害。
- 該設備能夠接受任何收到的干擾，其中包括可能導致意外操作的干擾。

如果該設備確實對無線電或電視接收造成乾擾（可以透過關閉和開啟設備來確定），使用者建議嘗試以下一種或多種措施：

- 確保使用合適的電纜將設備連接到外圍設備。
- 確保Agilent Technologies明確批准的變更或修改可能會使使用者操作該設備的授權無效。

使用说明

韓國 A 類 EMC 聲明

本設備適合專業用途 A類設備，適用於家庭以外的電磁環境。

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자파적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

此 ISM 設備符合加拿大 ICES-001/NMB-001 標準。

使用说明

服務

客戶如需要升級的更換或維修服務，請聯絡當地經銷商，或直接發送電子郵件至：

vpt-customer@agilent.com

vpl-customer@agilent.com

需要填寫「要求退貨」表格才能將幫浦恢復到Agilent維修部門
(在本手冊末尾會提供)。

6

取扱説明書

この説明書について	139
関連補足資料	139
定義と用語	140
警告・注意・注の定義	140
警告の記号	141
安全性	142
EMC 警告	143
FCC	143
適切な使用	144
合理的に予見可能な誤使用	144
ドライスクロールポンプの安全ガイドライン	145
有資格者	146
保護装備	147
製品説明	149
輸送と保管	149
開梱と点検	150
設置の準備	151
接地方法	154
起動手順	155
停止手順	158
処分	159
環境条件、コンプライアンス、ユーティリティ要件	160
電磁両立性	160
EN55011/CISPR11	160
韓国クラス A EMC 宣言	161
ICES/NMB-001	161
サービス	162

この説明書について

このマニュアルには、適用される指令に従った IDP-4 ドライ スクロール ポンプに関するユーザー向けの指示と技術情報が記載されています。

この資料は、この発行時点でお客様または第三者によって製品に変更が加えられていない限り、製造された製品に対して有効です。矛盾する情報がある場合は、製品と共に同封されている資料を有効な資料とみなしてください。

注

- 1 この取扱説明書には、IDP-4 ドライスクロールポンプを使用するすべての人が安全に操作し、その製品寿命全体にわたりその性能を保証するための重要な情報が含まれています。
- 2 この取扱説明書は、関連する出版物とともに、すべての操作者/メンテナンス担当者がアクセスできる場所に保管してください。

関連補足資料

X3802-90000 適合宣言

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するのより詳細な情報を提供しています。

警告の記号

以下は、IDP-4ドライスクロールポンプの警告とともに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。取扱説明書内、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。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。



	CE 認証
	CSA 認証
	中国RoHS 指令認証
	廃電気・電子製品に関するEU WEEE指令
	RCM 認証
	製造所
	高温の表面
	危険な電圧
	包括的な危険
	保護接地端子
	交流電流 (AC)
	電源オン
	電源オフ

安全性

このセクションには、低電圧指令2006/42/ECに規定されている、一般的な安全規制の遵守と製品の特定の使用に関連する安全規制の遵守に不可欠な情報が含まれています。

これらの指示、およびこの取扱説明書に記載されているその他の指示に従わない場合、安全性に影響を及ぼし、製品の操作者に事故を引き起こす可能性があります。

Agilent Technologiesは、技術文書に記載されている安全規則を遵守しなかったことに起因した製品の損傷、もしくは操作者または第三者の身体的安全について一切の責任を負いません。

EMC 警告

EN 55011 クラス A 警告

これはクラスAの製品です。この製品は、家庭環境では無線干渉を引き起こす可能性があり、その場合、ユーザーは適切な対策を講じる必要があります。

FCC

この機器は、FCC規則の第15部に準拠しています。操作は以下の2つの条件に従うものとします：(1) この機器は有害な干渉を引き起こしてはなりません。(2) この機器は、望ましくない動作を引き起こす可能性のある干渉を含む、受信したあらゆる干渉を受け入れる必要があります。

注

この機器は、FCC 規則の第 15 部に従ってテストされ、クラス A デジタル機器の制限に準拠していることが確認されています。これらの制限は、機器が商業環境で操作されるときに、有害な干渉から適切に保護するように設計されています。この機器は、無線周波数エネルギーを生成、使用、および放射する可能性があり、取扱説明書に従って設置および使用しない場合、無線通信に有害な干渉を引き起こす可能性があります。住宅地でこの機器を操作すると、有害な無線通信干渉が発生する可能性があり、その場合使用者は自己負担でそのような干渉を修正する必要があります。

適切な使用

IDP-4 ドライ スクロール ポンプは、このマニュアルに従って資格のある個人が真空ポンプとして使用することのみを目的としています。

製品に使用するアクセサリや装置は、Agilent Technologiesが提供または承認したものでなければなりません。

Agilent Technologiesは、この取扱説明書に記載されている以外の用途、または必須要件や安全対策が遵守されていない用途に対して、いかなる保証も責任も負いません。

合理的に予見可能な誤使用

Agilent Technologiesは、IDP-4 ドライ スクロール ポンプの不適切な使用および予見される誤使用に起因するすべての責任を負わないものとします。不適切な使用は、すべての責任および保証の請求権が喪失するものとします。

不適切な使用は以下が含まれますが、これらに限定されるものではありません：

- ポンプを誤った電圧電源に接続する
- 接地接続が不適切または欠落している
- 推奨される注入口圧力を超えて継続的に運転する
- ガスバラスト注入口で過圧状態で運転する
- 危険なガスをポンプで送るときにガスバラストを大気に開放して運転する
- 周囲温度が高すぎる状態で運転する
- 周囲温度が仕様より低い状態で運転する
- ポンプの不適切な保管および輸送による損傷
- 主電源を切らずにメンテナンスやカウリングの取り外しを行う
- 気流要素（カバーまたはファン）なしで運転すること
- 返品の商品を工場出荷時の梱包以外で郵送する。

ドライスクロールポンプの安全ガイドライン

ポンプの操作とメンテナンスの担当者は、適切な訓練を受け、事故防止の規則を把握する必要があります(有資格者のセクションを参照)。ポンプの操作とメンテナンスを行っている間は、操作者の負傷やポンプやその他の機器の損傷を防ぐために、この取扱説明書に記載されている事故防止の注意事項を常に遵守する必要があります。これらの注意事項は、取扱説明書の全体を通して、警告および注意の形で記載されています。

警告



機器の損傷や操作者の負傷を防ぐために、この取扱説明書に記載されている設置手順に厳密に従ってください！

警告



決してポンプで可燃性物質を扱わないでください。
爆発性のある環境にポンプを設置することは禁止されています。

有資格者

IDP-4ドライ スクロール ポンプは、資格のある担当者のみが使用することが推奨されています。

資格のある担当者とは、以下の者です：

- 適切な資格を持ち、Agilentドライ真空ポンプの使用経験があるか、トレーニングを修了した者。
- 怪我や損傷を引き起こさないように、必要な安全対策を遵守し、適切な条件下でポンプを操作する者。
- 適用される国および地方自治体の安全規制基準および条例を把握しており、それらに従い、適切な資格を有する者。

注

この取扱説明書に標準メンテナンスとして記載されている以上の本製品に対する修理は、Agilent の認定を受けた権限のある者のみが行う必要があります。いかなる調整、変更、メンテナンス作業も、危険性を熟知している専門の技術者が行う必要があります。

保護装備

ポンプを操作またはメンテナンスする操作者の保護装備は、実行する操作の種類に常に適したものでなければなりません。使用する保護装備は、機械を使用する国で施行されている法律の安全要件に準拠している必要があります。

通常、操作者は、IDP-4ドライスクロールポンプを移動、取り扱う、または設置する際には安全靴を着用する必要があります。

警告



メンテナンスまたは取り付け時に、有害物質による健康への危険。



- 過程の異常な点により、真空ポンプ、構成部分、または操作液は、有害物質、反応性物質、もしくは放射性物質に汚染される場合があります。
- メンテナンスや修理、もしくは再度取り付け時には適切で十分な保護装備を着用してください。



警告



装置の落下による損傷のリスク。



真空ポンプを手で運搬する場合、荷重が滑り落ちる危険があります。

- 両手で小型・中型の真空ポンプを運んでください。
- 20 kg 以上のいかなる装置も、適切なリフトの機器を使用して運搬されなければなりません。
- 指令 EN 347 にしたがって、爪先が鋼製の安全長靴を着用してください



注意



鋭い縁による損傷のリスク。

- ポンプを修理する前、またはシステムからポンプのいかなる取り付け/取り外しをする前に、ポンプの完全な静止をお待ちください。
 - 高真空フランジ内で直接操作しないでください。
 - 必要に応じて、指令 EN 420 に従った保護手袋を着用してください。
-

製品説明

IDP-4は、気密性のあるドライスクロールポンプです。低圧領域から大気圧より大幅に低くも高くもない高圧領域に空気または不活性ガスをポンプで送るのに適しています。このポンプは、腐食性、自然発火性、微粒子形成性のガス、または可燃性および/または燃焼性範囲のガス混合物をポンプで送るためのものではなく、大気圧より大幅に高い排出圧力を達成するためのコンプレッサーとして使用することも意図されていません。

取扱説明書および関連資料は、機器を使用する際の操作者の安全を保証するために必要な情報が含まれています。詳細情報は「技術情報」のセクションに記載されています。

輸送と保管

ポンプを輸送および保管する際は、以下の環境要件を超えないようにしてください：

- 温度： -20 °C から +60 °C (-4 °F から 140 °F)
- 相対湿度： 0 から 95% (結露なし)

注

真空ポンプの輸送および保管中は、ポンプの密閉度に関係なく、周囲の環境からの水蒸気が真空ポンプ内に浸透することを認識する必要があります。ポンプの定格真空性能を完全に発揮するには、技術情報のセクションに記載されているガスバラスト機能を使用して、この水蒸気を除去する必要があります。

開梱と点検

ポンプは特別な保護梱包で提供されます。輸送中に生じた可能性のある損傷の形跡がある場合は、最寄りの営業所にお問い合わせください。ポンプと梱包を合わせた総重量は約 10kg (20 ポンド) です。

警告



Agilent の認定サービスセンターでサービスを受ける場合、段ボールとすべての梱包材が必要になります。この梱包材を保存しておくこと、サービスの依頼がスムーズになります。

警告



ポンプを開梱する際は、落とさないように注意し、突然の衝撃やショックを避けてください。

設置の準備

IDP-4ドライスクロールポンプは、屋内での運転（設置カテゴリII）を目的として設計および認定されており、比較的清潔で乾燥した環境（汚染度2）での使用に適しています。

仕様については、「技術情報」セクションの表1を参照してください。

これらのポンプは、屋外での運転を目的として設計または認定されていません。

IDP-4ドライスクロールポンプは、爆発性または可燃性物質を含む周囲環境では絶対に使用しないでください。

IDP-4ポンプは本質安全構造ではなく、腐食性、自然発火性、粒子生成性のガス、または可燃性および／または燃焼性範囲内のガス混合物とともに使用してはなりません。

ポンプを堅固で平らな表面に設置してください。ベンチまたはテーブルに取り付ける必要がある場合は、そのベンチまたはテーブルがポンプおよび関連機器の重量を支えられるように定格されていることを確認してください。

ポンプを電源に接続する際は、少なくとも10Aの容量を持つIEC-320スタイルの電源コードを使用してください。

注

初期設置および起動時に性能が低下する最も一般的な原因は、ポンプおよび真空チャンバー内に吸着された水蒸気です。

ガスバラストを作動させた状態でポンプを運転することは、ポンプが一定期間停止しており、その間にポンプおよび真空システム内部に水分が吸着している場合にも有効です。

真空システムの構成および吸着された水分量に応じて、完全な真空性能を回復するまでにガスバラストの作動を約30分から24時間行う必要があります。

注意

操作中は、以下の環境条件を遵守する必要があります：

- 温度：+5 °C から +40 °C (41 °F から 104°F)
- 相対湿度：5 から 95% (結露なし)

適切な電流および電圧定格の IEC-320 スタイルの電源コードを使用してポンプを電源に接続してください (表 2 を参照)。

注意

真空ポンプを設置する際は、必要に応じて電源ポンプを素早く切断できるように、使用者が電源スイッチと電源コード接続に簡単にアクセスできるような位置にポンプを配置することが推奨されます。

注意

ポンプヘッドの両側にある 2 本の六角ボルトを、締め付けたり、緩めたり、いかなる方法でも変更したりしないでください（次の図を参照）。そのような操作を行うと、性能の低下やポンプの損傷を引き起こす可能性があります。



図 16

警告



ポンプから安全装置や絶縁装置を取り外したり改造したりしないでください。そうした場合、重大な安全上の危険が生じる可能性があり、保証が無効になる可能性もあります。

警告



- このポンプは、空気と不活性ガスのみをポンピングし、排出することができます。このポンプは、腐食性、自然発火性、微粒子形成性ガス、または可燃性および/または可燃性範囲のガス混合物をポンピングするためのものではありません。
- 雨、蒸気、過度の湿気にさらされない場所に設置してください。感電、ショート、重度の傷害を引き起こす可能性があります。
- ポンプの点検、再構成、または修理を行う前に、必ず電源を切ってください。
- ガスをポンプで汲み上げるときは、大気中に放出したり空気と混合したりできないように、ガスバラストを密閉する必要があります。
- ポンプから発生する騒音レベルを低減したい場合に、排気サイレンサーオプションを利用することができます。
- 身体の損傷を防ぐため、人体のいかなる部分も真空中にさらさないでください。
- ポンプは電源スイッチに簡単にアクセスできる場所に設置する必要があります。
- すべての一般的な接続用に認定されたアジレント製電源コードは、「電気接続」セクションに記載されています。
異なる電源コードを使用する必要がある場合は、電源コードの定格電流が **200~240V** サービスでは少なくとも **10A**、**100~127V** サービスでは少なくとも **15A** でなければなりません。

接地方法

警告



この製品は接地する必要があります。電気ショートが発生した場合、接地により電流の逃げ道が提供され、感電のリスクが軽減されます。このポンプには、適切な接地プラグ付きの接地線が付いた電源コードが装備されています。プラグは、すべての地域の法令に従って適切に設置され、接地されたコンセントに差し込む必要があります。

電源コードコネクタは、ポンプの電源モジュールに適切に差し込み、接続不良や断続的な接続を防いでください。

注意

ポンプは大気中に通常存在する微粒子をポンピングできますが、固体、化学物質、粉末、溶剤、凝縮物、液体、その他の粒子を処理するようには設計されていません。これらは機器を損傷したり、性能を低下させたり、耐用年数を短縮したりする可能性があります。

適切な空気循環を可能にするために、システムの周囲に最低 **20 cm** の空きスペースを設けることが推奨されています。

注意

ポンプの外部表面は、中性洗剤のみを使用してきれいにすることができます。

ドライスクロールポンプには、蓄積したほこりやゴミを洗浄するための液体は入っていません。定期的にポンプを大気圧で **1～2** 分間運転して、洗浄してください。

警告



接地方法が完全に理解できない場合、または製品が適切に接地されているか不明な場合は、資格のある電気技師または担当者に確認してください。付属のプラグを改造しないでください。コンセントに合わない場合は、資格のある電気技術者に適切なコンセントを取り付けてもらってください。

プラグと同じ形状のコンセントにのみ製品を接続してください。この製品にアダプターを使用しないでください。

起動手順

警告



排気口および吸気口のフランジカバーを取り外してください。



IDP-4ドライスクロール真空ポンプの起動手順は次のとおりです:

1. ポンプを所定の位置に設置し、適切な機械的および真空接続を行うか、代わりに吸入口をブランクフランジで覆ってください。
2. 排気継手が塞がれていないこと、または閉鎖されていないことを確認してください。排気ホースまたはマフラーを使用する場合、排気制限により6.5 psigを超える背圧が生じないようにする必要があります。
3. ON/OFFスイッチが「0」（OFF）位置にあることを確認してください。
4. 適切な電源コードを接続し、コードを主電源コンセントに差し込んでください。電源コードの定格電流は、200～240Vサービスの場合は少なくとも10A、100～127Vサービスの場合は少なくとも15Aである必要があります。
5. 主電源のON/OFFスイッチを「1」位置に押しってください。
6. ポンプにオプションの吸入口バルブが装備されている場合、電源投入後約10秒でバルブが開きます。

ポンプが公表されている最終圧力仕様に到達するまでに、数時間かかる場合があります。ポンプが期待される最終圧力に達しない場合は、ポンプ内部に水蒸気が蓄積している可能性があります。

NOTE

初期設置および起動時に性能が低下する最も一般的な原因は、ポンプおよび真空チャンバー内に吸着された水蒸気です。

ガスバラストを作動させた状態でポンプを運転することは、ポンプが一定期間停止しており、その間にポンプおよび真空システム内部に水分が吸着している場合にも有効です。

真空システムの構成および吸着された水分量に応じて、完全な真空性能を得るまでにガスバラストの作動を約 30 分から 24 時間行う必要があります。



図 17

注

ポンプの ON/OFF スイッチはロックスイッチ型であり、IEC 出版物 417 に準拠した記号を使用して ON および OFF の位置を示しています。

次の図は、OFF 位置にあるスイッチを示しています。

AVVERTENZA



冷却通気口に異物を挿入しないでください。ファンの羽根に接触すると、重大な人身傷害またはポンプの損傷を引き起こすおそれがあります。

警告



冷却通気口を塞がないでください。これらの通気孔を塞ぐと、ポンプが過熱するおそれがあります。

ポンプ表面温度が **65° C (150° F)** を超えると、損傷を引き起こす可能性があります。

そのような状態が見られた場合は、ポンプの電源を切り、冷却させてください。分解し、損傷がないか点検し、必要に応じて修理してください。

警告



潜在的に危険な物質、または空気と混合してはならない物質を使用する場合は、ガスバラストを使用しないでください。

停止手順

どの動作モードでも手動でポンプを停止するには、メインのON/OFFスイッチを「0」の位置に回します。ポンプに統合遮断弁が装備されている場合は、ポンプから上流のコンポーネントへのガスの逆流を防止するため、遮断弁が急速に閉じます。

注意

IDP-4 ポンプを停止する前に、上流側のすべての接続部品が適切にバルブで遮断されているか、または大気圧への曝露に備えて準備されていることを確認してください。

処分

ラベルに記載の“WEEE” ロゴの意

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 18 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>

環境条件、コンプライアンス、ユーティリティ要件

注

このポンプは、換気、メンテナンスのアクセス、および電源スイッチへの容易なアクセスのために、すべての側面に少なくとも 20 cm (8 インチ) の空きスペースが必要です。ポンプをベンチまたは棚に置く場合は、ポンプとその他の関連機器の重量を支えることが可能な定格が必要です。

電磁両立性

EN55011/CISPR11

グループ 1 ISM 装置：グループ 1 には、装置自体の内部機能に必要な伝導結合無線周波数エネルギーが意図的に生成および/または使用されるすべての ISM 装置が含まれます。

クラスAの装置は、家庭用以外のすべての施設、および家庭用の建物に供給する低電圧電源網に直接接続されている装置の使用に適している装置です。

この機器は、放射線業務用機器としてCISPR11、グループ1、クラスAの要件に適合しています。したがって、伝導および放射妨害により、他の環境での電磁両立性を確保することが困難な場合があります。

操作は以下2つの条件に従います：

- この機器は、有害な干渉を引き起こすことはありません。
- この機器は、望ましくない動作を引き起こす可能性のある干渉を含め、受信したすべての干渉を受け入れる必要があります。

この装置がラジオやテレビの受信に有害な干渉を引き起こす場合には（装置の電源を切ったり入れたりすることで判断できます）、使用者は以下の対策を1つ以上試すことを推奨します：

- 周辺機器との接続には、必ず適切なケーブルを使用してください。
- Agilent Technologies によって明示的に承認されていない変更または修正は、装置を操作する使用者の権限を無効にする可能性があります。

韓国クラス A EMC 宣言

この装置は、業務使用に適したクラス A であり、家庭外の電磁環境での使用を目的としています。

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자파적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

この ISM 機器は、カナダの ICES-001/NMB-001 に準拠しています。

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

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	164
Relevant Supporting Documents	164
Definitions and terminology	165
Definition of Caution, Warning and Note	165
Warning Symbols	166
Safety	167
EMC Warnings	167
FCC	167
Proper use	168
Reasonably Foreseeable Misuse	168
Safety guideline for Dry Scroll Pump	169
Qualified Personnel	170
Protective equipment	171
Product Description	173
Transport & Storage	173
Unpacking and Inspection	174
Preparation for Installation	175
Startup Procedure	179
Shutdown Procedure	182
Disposal	183
Environmental Conditions, Compliance and Utility Requirements	184
Electromagnetic Compatibility	184
South Korean Class A EMC declaration	185
ICES/NMB-001	185
Service	186

About this Manual

This manual contains instructions for users and technical information related to the IDP-4 Dry Scroll pump in accordance with the applicable directives.

This document is valid for the product produced at the time of this publication, provided that changes are not made to the product by the customer or third parties. In the event of any conflicting information, the documentation shipped with the product should be considered the valid documents.

NOTE

- 1 This manual contains important information so that all personnel using the IDP-4 Dry Scroll pump can operate them safely and guarantee their performance over their entire life span.
 - 2 Keep this manual, together with all the related publications, in an accessible place known to all operators/maintenance personnel.
-

Relevant Supporting Documents

X3802-90000 Declaration of Conformity

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 IDP-4 Dry Scroll pump. 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. 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.



	CE certification
	CSA certification
	RoHS China certification
	Waste Electrical and Electronic Equipment
	Regulatory Compliance Mark (RCM)
	Manufacturing site
	Hot surfaces
	Dangerous Voltages
	Generic Hazard
	Protective Ground Terminal
	Alternating current (AC)
	Power ON
	Power OFF

Safety

This section contains the information, prescribed by the Low Voltage Directive 2006/42/EC, which is essential for compliance to and observance of 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 affect safety and could potentially result in 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.

EMC Warnings

EN 55011 Class A Warning

This is a Class A product. In a domestic environment this product may cause radio interference in which case the user may be required to take adequate measures.

FCC

This device complies with Part 15 of the FCC rules. Operation is subject to the following two conditions: (1) This device may not cause harmful interference, and (2) this device must accept any interference received, including interference that may cause undesirable operation.

NOTE

The equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generated, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is also likely to cause harmful radio communications interference in which case the user will be required to correct the interference at his own expense.

Instructions for Use

Proper use

The IDP-4 Dry Scroll Pump is exclusively intended use by qualified individuals as vacuum pump as per this manual.

Any accessories and equipment used with the product must be supplied or approved by Agilent Technologies.

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.

Reasonably Foreseeable Misuse

Agilent Technologies declines all responsibility deriving from the improper use and foreseeable misuses of the IDP-4 Dry Scroll pump. Improper use will cause all claims for liability and warranties to be forfeited.

Improper use may include, but is not limited to:

- Connecting the pump to wrong voltage supply
- Improper or missing ground connection
- Operating continuously above the recommended inlet pressure
- Operating with overpressure at gas ballast inlet
- Operating gas ballast open to atmosphere when pumping dangerous gases
- Operating with excessively high ambient temperature
- Operating with ambient temperature lower than specification
- Damage due to improper storage and transportation of pump
- Performing maintenance or removing cowlings without disconnecting mains power
- Operating without airflow elements (cowling or fan).
- Shipment of returned product in non-factory packaging.

Safety guideline for Dry Scroll Pump

Personnel responsible for pump operation and maintenance must be properly trained and must be aware of the accident prevention rules (refer to Qualified Personnel section). The accident prevention precautions contained in this manual must be continuously respected during operation and maintenance of the pump to avoid injury to operators and damage to the pump or other equipment. These precautions are provided in the form of WARNING and CAUTION notes throughout the entire user manual.

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!

WARNING



**The pump must not handle flammable materials.
It is forbidden to install the pump in an explosive environment.**

Instructions for Use

Qualified Personnel

The IDP-4 Dry Scroll Pump is recommended for use only by qualified personnel. A qualified person is:

- An individual who has the proper qualifications and has experience, or has completed training on, working with Agilent Dry Vacuum Pump.
- An individual that follows the necessary safety measures and operates the pump under the proper conditions so as not to cause injury or damage.
- An individual who is knowledgeable of and properly qualified to follow any applicable national and local safety and regulatory standards and by-laws.

NOTE

Any repairs on this product, above and beyond those listed as standard maintenance in this user manual, must be carried out exclusively by Agilent authorized personnel. Any adjustment, modification or maintenance operation must be performed by a professional technician who is informed of the risks.

Protective equipment

The protective equipment of the operators who are operating or executing the maintenance of the pump must always be adequate for the type of operation being executed. Protective equipment used must comply with the safety requirements of the legislation in force in the country in which the machine is used.

In general, the operator must wear safety shoes while moving, handling, or installing the IDP-4 Dry Scroll pump.

WARNING

Danger to health by hazardous substances during maintenance or installation.



- Depending on the application, vacuum pump, components, and operating fluids can be contaminated by toxic, reactive or radioactive substances.
- Wear adequate protective equipment during maintenance and repairs or in case of reinstallation.

WARNING

Risk of injury through falling objects



When transporting vacuum pump, there is a danger of dropping the pump and of slip and falls.

- Carry small and mid-size vacuum pump with both hands.
- Any equipment heavier than 20 kg should be transported using a suitable lifting device.
- Wear safety shoes with steel toe cap according to directive EN 347.

Instructions for Use

CAUTION



Risk of injury through sharp edges and hot surfaces.

- Before servicing the pump or before any mounting or disconnecting of the pump from the system, wait for the pump to come to a complete stop and cool down to room temperature.
 - Do not place hands or fingers directly inside any parts of the pump.
 - If necessary, wear protective gloves according to EN 420.
-

Product Description

The IDP-4 is hermetic Dry Scroll pump. This pump is suitable for pumping air or inert gases from an area of low pressure to an area of higher pressure which is not substantially below or above atmospheric pressure. This pump is not intended to pump corrosive, pyrophoric, particulate-forming gases, or gas mixtures in their flammable and/or combustible range, nor is it intended to be used as a compressor to achieve discharge pressures substantially above atmospheric pressure.

The manual and related documents contain the information necessary to guarantee the safety of the operator when using the equipment. Detailed information is supplied in the "Technical Information" section.

Transport & Storage

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

- Temperature: -20 °C to +60 °C (-4 °F to 140 °F)
- Relative humidity: 0 to 95% (non-condensing)

NOTE

It should be recognized that during transportation and storage of vacuum pump, water vapor from the ambient environment will permeate inside the vacuum pump regardless of how well the pump is sealed. To achieve the full rated vacuum performance of the pump this water vapor must be removed using the gas ballast features described in the Technical Information section.

Unpacking and Inspection

The pump is supplied in a special protective packing. If this shows signs of damage which have occurred during transport, contact your local sales office. Total weight of the pump and packaging together is approximately 10 kg (20 lbs).

WARNING



The carton and all packing materials will be required in the event of service at an Agilent approved service center. Saving this packaging will facilitate any service requests.

WARNING



When unpacking the pump, be sure not to drop it and avoid any kind of sudden impact, or shock.

Preparation for Installation

The IDP-4 Dry Scroll pump is designed and certified for operation indoors (installation category II) in relatively clean, dry environments (pollution degree 2).

Refer to Table 1 in the Technical Information section for specifications. This pump is not designed or certified for operation outdoors.

The IDP-4 Dry Scroll pump should never be used in ambient environments that contain explosive or combustible materials. The IDP-4 pump is NOT intrinsically safe and should never be used with corrosive, pyrophoric, particulate-forming gases or gas mixtures in their flammable and/or combustible range.

Place the pump on a solid, level surface. If it must be mounted on a bench or table, make sure the bench or table is rated to carry the weight of the pump and all associated equipment.

Connect the pump to the power supply using an IEC-320 style power cord of at least 10 A capacity.

NOTE

The most common cause of poor performance on initial install and startup is adsorbed water vapor in both the pump and vacuum chamber. Running the pump with the gas ballast activated may also be beneficial if the pump has been shut off for a period of time, which has allowed water to adsorb inside both the pump and vacuum system. Depending on the vacuum system and amount of adsorbed water it can typically take between ~30 minutes to ~24 hours of gas ballast activation to achieve the full vacuum performance.

CAUTION

During operation the following environmental conditions must be respected:

- **Temperature:** +5 °C to +40 °C (41 °F to 104°F)
- **Relative humidity:** 5 to 95% (non-condensing)

Connect the pump to the power supply using an IEC-320 style power cord of the appropriate current and voltage rating (see Table 2).

CAUTION

When installing the vacuum pump, it is recommended that the pump be positioned in a way that the power switch and power cord connection remain easily accessible to the user in order to disconnect power pump quickly if needed.

Instructions for Use

CAUTION

Do not tighten, loosen or alter in any way the two hex head bolts on each side of the pump head (see the following figure). Doing so will cause loss of performance and/or pump damage.



Figure 19

WARNING



Do not remove or modify any safety or insulating equipment from the pump. To do so may create a serious safety hazard and could also void the warranty.

Instructions for Use

WARNING



- This pump is capable of pumping and exhausting air and inert gases only. This pump is not intended to pump corrosive, pyrophoric, particulate-forming gases or gas mixtures in their flammable and/or combustible range.
 - Install in an area that is not exposed to rain, steam, or excessive humidity. They can cause electric shock, short circuits, and severe bodily injury.
 - Before inspecting, reconfiguring, or servicing the pump, be sure the electrical supply is disconnected.
 - The gas ballast must be sealed whenever pumping any gas that cannot be vented to the atmosphere or mixed with air.
 - An exhaust silencer option is available for those applications in which it is desirable to diminish the sound level produced by the pump.
 - To prevent bodily injury, avoid exposing any part of the human body to vacuum.
 - The pump must be installed with easy access to the power switch.
 - Agilent approved power cords for all common connections are indicated at the Electrical Connections section. If it is necessary to use a different power cord, the current rating of the power cord must be at least 10A for 200-240V service and at least 15A for 100-127V service.
-

Grounding Instructions

WARNING



This product should be grounded. In the event of an electrical short circuit, grounding reduces the risk of electric shock by providing an escape path for the electric current. This pump is equipped with a power cord that has a grounding wire with an appropriate grounding plug. The plug must be inserted into an outlet that is properly installed and grounded in accordance with all local codes and ordinances.

The power cord connector must be properly plugged into the pump's power module to prevent defective or intermittent connection.

CAUTION

Although the pump can trace particulates normally found in the atmosphere, it is not designed to process solids, chemicals, powders, solvents, condensates, liquids, or other particulates. They can damage the equipment, degrade its performance, or shorten its useful life.

A minimum of 20 cm free space is recommended all around the system to allow proper air circulation.

Instructions for Use

CAUTION

The exterior surfaces of the pump may be cleaned with mild detergents only.

The dry scroll pump does not contain fluid for the cleansing of accumulated dust and debris. Run the pump periodically at atmosphere for a minute or two to flush it out.

WARNING



Check with a qualified electrician or serviceman if the grounding instructions are not completely understood, or if you are in doubt as to whether the product is properly grounded. Do not modify the plug provided. If it does not fit the outlet, have the proper outlet installed by a qualified electrician.

Connect the product only to an outlet that has the same configuration as the plug. Do not use an adapter with this product.

Startup Procedure

WARNING



Remove the flange covers from exhaust and inlet.



The startup procedure for the IDP-4 Dry Scroll Vacuum Pump is as follows:

1. Mount the pump in its intended location and make the appropriate mechanical and vacuum connections or alternatively cover the inlet with a blank off flange.
2. Ensure that the exhaust fitting is not blocked or blanked-off. If an exhaust hose or muffler is used, exhaust restrictions must not cause backpressure restriction above 6.5 psig.
3. Ensure the ON/OFF switch is in the OFF "0" position.
4. Connect the appropriate power cord and plug the cord into the mains power. The current rating of the power cord must be at least 10A for 200-240V service and at least 15A for 100-127V service.
5. Press the main ON/OFF switch to "I" position.
6. If the pump is equipped with an optional inlet valve, the valve will open approximately 10 seconds after power on.

It may take several hours for the pump to achieve published ultimate pressure specification. If the pump does not reach its expected ultimate pressure, it may be due to accumulated water vapor inside the pump.

Instructions for Use

NOTE

The most common cause of poor performance on initial install and startup is adsorbed water vapor in both the pump and vacuum chamber. Running the pump with the gas ballast activated may also be beneficial if the pump has been shutoff for a period of time, which has allowed water to adsorb inside both the pump and vacuum system. Depending on the vacuum system and amount of adsorbed water it can typically take between ~30 minutes to ~24 hours of gas ballast activation to achieve the full vacuum performance.



Figure 20

NOTE

The pump ON/OFF switch is a rocker type switch that has symbols in accordance with IEC Publication 417 to represent the ON and OFF positions. The previous figure shows a switch in the OFF position.

Instructions for Use

WARNING



Do not insert any foreign object through the cooling vents as contact with the fan blades can result in serious personal injury or damage to the pump.

WARNING



Do not block the cooling vents. Blocking these vent holes can cause pump overheating. A pump surface temperature in excess of 65 °C (150 °F) is potentially damaging. If such conditions are observed, turn the pump off and allow it to cool. Disassemble, inspect for damage, and repair, if necessary.

WARNING



Do not use gas ballast if using potentially hazardous materials or those that should not be mixed with air.

Shutdown Procedure

To shut down the pump manually in any operating mode turn the main ON/OFF switch to "0" position. If the pump is equipped with an integrated isolation valve, it will rapidly close to eliminate backflow of gas from the pump to upstream components.

CAUTION

Before shutting down the IDP-4 pump, ensure that all upstream interconnected components are properly valved off or prepared for potential exposure to atmosphere pressure.

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 21 Logo "WEEE"

For more information refer to:

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

Environmental Conditions, Compliance and Utility Requirements

NOTE

This pump requires a clear space of at least 20 cm (8 in.) on all sides for ventilation, maintenance access and easy access to the power switch. If it is desired to place the pump on a bench or shelf It must be rated to support the weight of the pump and any other associated equipment.

Electromagnetic Compatibility

EN55011/CISPR11

Group 1 ISM equipment: Group 1 contains all ISM equipment in which there is intentionally generated and/or used conductively coupled radio-frequency energy which is necessary for the internal functioning of the equipment itself.

Class A equipment is equipment suitable for use in all establishments other than domestic and those directly connected to a low voltage power supply network which supplies buildings used for domestic purposes.

This device complies with the requirements of CISPR11, Group 1, Class A as radiation professional equipment. Therefore, there may be potential difficulties in ensuring electromagnetic compatibility in other environments, due to conducted as well as radiated disturbance.

Operation is subject to the following two conditions:

- This device may not cause harmful interference.
- This device must accept any interference received, including interference that may cause undesired operation.

If this equipment does cause harmful interference to radio or television reception, which can be determined by turning the equipment off and on, the user is encouraged to try one or more of the following measures:

- Make sure that appropriate cables are used to connect the device to peripheral equipment.
- Changes or modifications not expressly approved by Agilent Technologies could void the user's authority to operate the equipment.

Instructions for Use

South Korean Class A EMC declaration

This equipment is Class A suitable for professional use and is for use in electromagnetic environments outside of the home.

A 급 기기
(업무용 방송통신기자재)

이 기기는 업무용 (A 급) 전자과적합기기로서 판매자 또는 사용자는 이 점을 주의하시기 바라며, 가정외의 지역에서 사용하는 것을 목적으로 합니다.

ICES/NMB-001

This ISM device complies with Canadian ICES-001/NMB-001.

Service

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

vpt-customer@agilent.com

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 IDP-4 Dry Scroll Pump	188
Technical Specifications	189
IDP-4 Dry Scroll Pump Outline Drawing	191
Performance Curves	193
Electrical Connections	195
Grounding Instructions	195
Mechanical Connections	197
Gas Ballast	199
Accessories and Spare Parts	201
Optional Accessories	201
Optional Inlet Valve Kit	202
Installation	204
Cleaning and Preventive Maintenance	205
External cleaning	205
Internal cleaning as part of a Preventive Maintenance plan	206
Verifying Performance	206
Tip Seal Replacement	207
Tip Seal Replacement Instructions	208
User Maintenance and Repair	216
Factory-Only Service	217
Service Options	218
Troubleshooting	219
Serial Connector & Pin allocation	221
Pump Serial Connector	221
Serial communication settings	222
Window Description and Protocol	223
Window Description	223
Windows Protocol	226

Description of the IDP-4 Dry Scroll Pump

The Agilent IDP-4 Dry Scroll pump is an oil-free, hermetic compact, quiet, vacuum pump with speed control. The IDP-4 utilizes digital motor control technology, providing uniform vacuum performance at all global mains frequencies (50 Hz and 60 Hz) and input voltages (100-127 VAC and 200-240 VAC).

The IDP-4 Dry Scroll Pump models include:

- X3802-64000 IDP-4 Dry Scroll Pump, Hermetic
- X3802-64010 IDP-4 Dry Scroll Pump, Hermetic, Inlet Valve
- X3802-64400 IDP-4 Dry Scroll Pump, Hermetic, Cryo

Technical Specifications

Table 1 IDP-4 Technical Specifications

Model	IDP-4 Dry Scroll Pump
Interface dimensions	H=181 mm x W=139 mm x L= 288 mm (7.1X5.5X11.3") standard pump H=237 mm x W=139 mm x L= 288 mm (VPI Valve) H=237 mm x W=139 mm x L= 338 mm (with all accessories)
Peak pumping speed (ISO 21360-1: 2020) (ISO 21360-2: 2020)	65 L/min, (4 m3/h, 2.35 cfm) at full speed setting Includes average peak pumping speed for both N2 and He without inlet valve configuration
Media	No corrosive, explosive or particulate forming gases
Ultimate Pressure* (ISO 21360-1: 2020) (ISO 21360-2: 2020)	1.8 +/- 0.2 x 10 ⁻¹ Torr (2.4 +/- 0.27 x 10 ⁻¹ mbar) at 3300 rpm (default setting)
Leak Rate	<1 x 10 ⁻⁶ sccs helium
Maximum recommended continuous inlet pressure	1.0 atmosphere (0 psig)
Maximum inlet pressure at nominal voltage	1.0 atmosphere (0 psig)
Maximum outlet pressure***	0.45 bar (6.5 psig)
Inlet connection	NW16
Exhaust connection	NW16
Gas ballast	Female 1/8" National Pipe Thread (adapter provided)
Ambient operating temperature	5 °C to 40 °C (41 °F to 104 °F)
Ambient Relative Humidity	5-95%; non-condensing
Storage temperature	-20 °C to 60 °C (-4 °F to 140 °F)
Motor rating: EN 60034-1-2:2007 EN 60034-30 (GB 18613-2020)	0.16 HP (0.12Kw)
Operating voltages**	Single phase 100-127VAC 50/60Hz 200-240 50/60Hz
Motor protection	Thermal protection/over current protection/fuse

Technical Information

Table 1 IDP-4 Technical Specifications (continued)

Model	IDP-4 Dry Scroll Pump
Operating speed	Variable Speed (2200-3300rpm), default setting 3300rpm +/-10%
Cooling system	Air-cooled
Weight	Pump only: 7.7kg (17 lbs)
Shipping weight:	Approximately 9.8 kg (22 lbs) with accessories
Sound Pressure Levels (EN ISO 2151:2008) (ISO 4871:1996) (ISO 11201:2010)	50 to 55 +/- 2.5 dba (min to max RPM setting) with default configuration (gas ballast closed)
Fuse type	5A slow blown 250V 5X20 mm
Installation	Indoor use, Equipment Class 1, Overvoltage Category II, Pollution Degree 2
Altitude	2000 m (6562 ft)
Emissions and Immunity	Electromagnetic Compatibility 2014/30/EU:2014 EN 61326-1: (2013) as a Class A device EN 61326-1: (2013) as a Class A device EN 61000-3-2:2019 Class A EN 61000-3-3:2014/A2:2022 Class A EN 61000-6-2:2020
Safety	Machinery Directive 2006/42/EC: 2006 EN 1012-2:1996+A1: 2010 EN ISO 12100:2012 Low Voltage Directive 2006/95/EC: 2007 EC 61010-1 Ed. 3.1 b:2018 CSA certified for US and Canada CE, ROHS

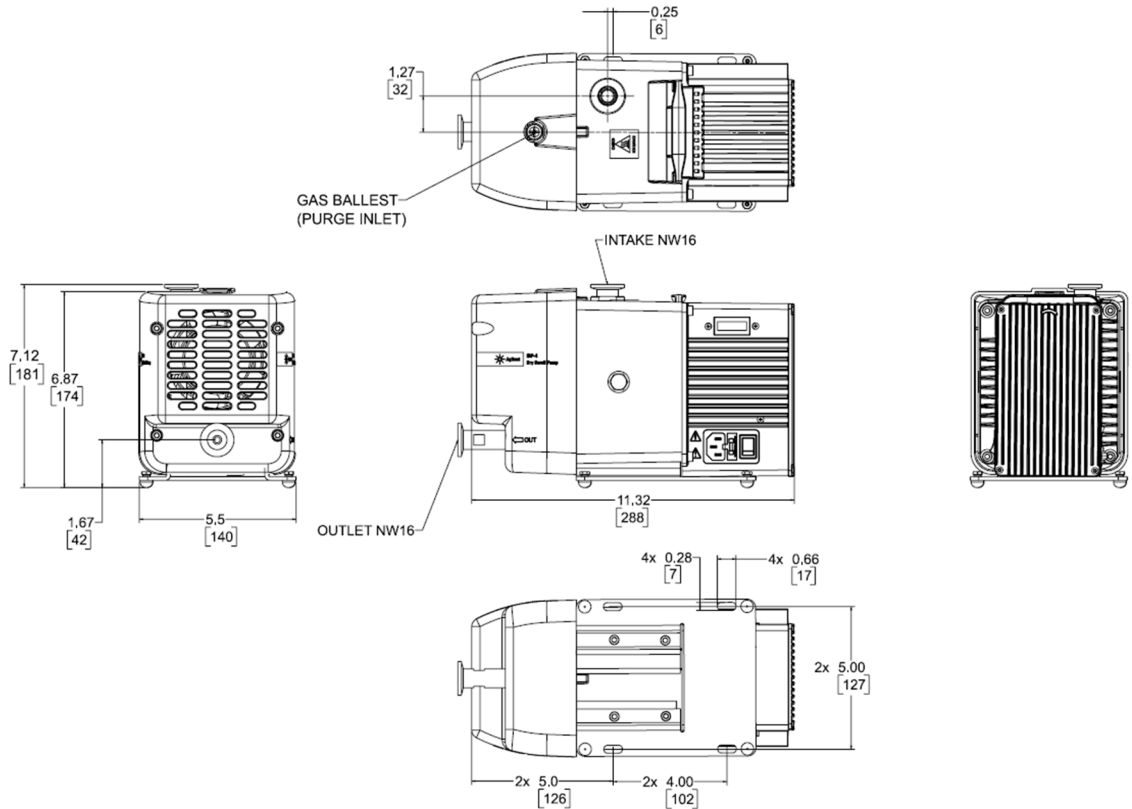
* If the pump is new or has been exposed to atmosphere for an extended period of time, it is possible that water can be present which can affect the pump's ability to achieve ultimate pressure. To remove water from the pump and system, activate the gas ballast using the procedure in the gas ballast paragraph on page 199. Ultimate pressure will vary for different configuration and RPM setting.

** Mains supply voltage fluctuations shall not exceed $\pm 10\%$ of the nominal voltage. For operation in Japan, the pump will function at 100VAC, but may offer reduced peak pumping speed at maximum load conditions. For operation in Japan at high inlet pressures and high pumping loads, use of 200VAC is recommended.

*** For customers with outlet pressure might experience different sound quality and slower pumping speed.

Technical Information

IDP-4 Dry Scroll Pump Outline Drawing



Note:

The IDP-4 mounting holes are backward compatible with the IDP-3.

Figure 22 IDP-4 Outline Drawing

Technical Information

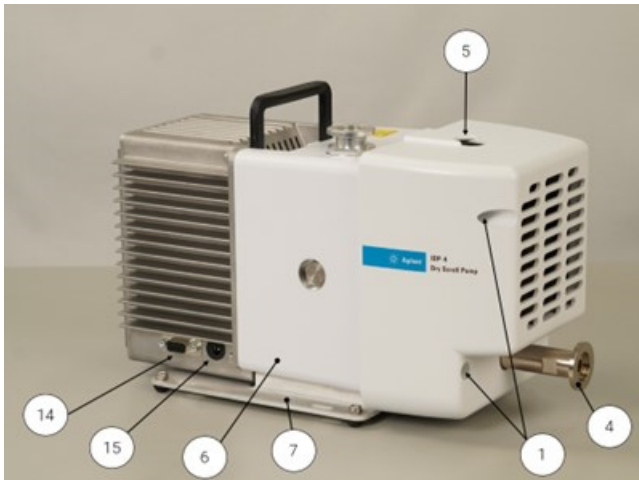
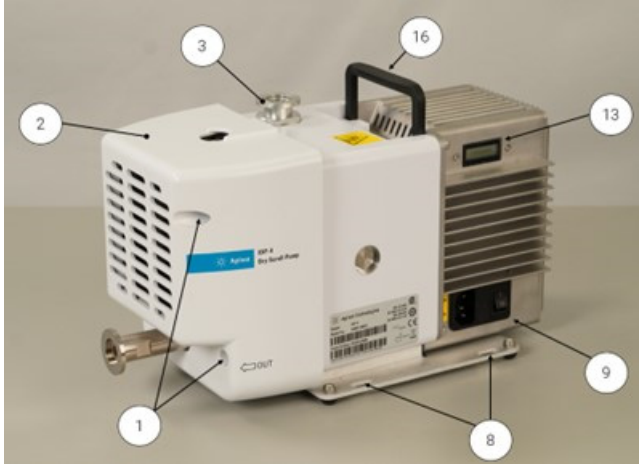


Figure 23 Principal Items

1. Front Cowling Bolts; M5 (4)
2. Front Cowling
3. Inlet Flange & Screen (NW16)
4. Exhaust Flange (NW16)
5. Gas Ballast Port
6. Frame
7. Base Plate
8. Mounting Holes; (4)
9. Power Entry Module
10. Power Cable Connection
11. Fuse Holder
12. Power Switch
13. Hour Meter
14. Serial communication
15. Inlet Valve connector
16. Handle

Performance Curves

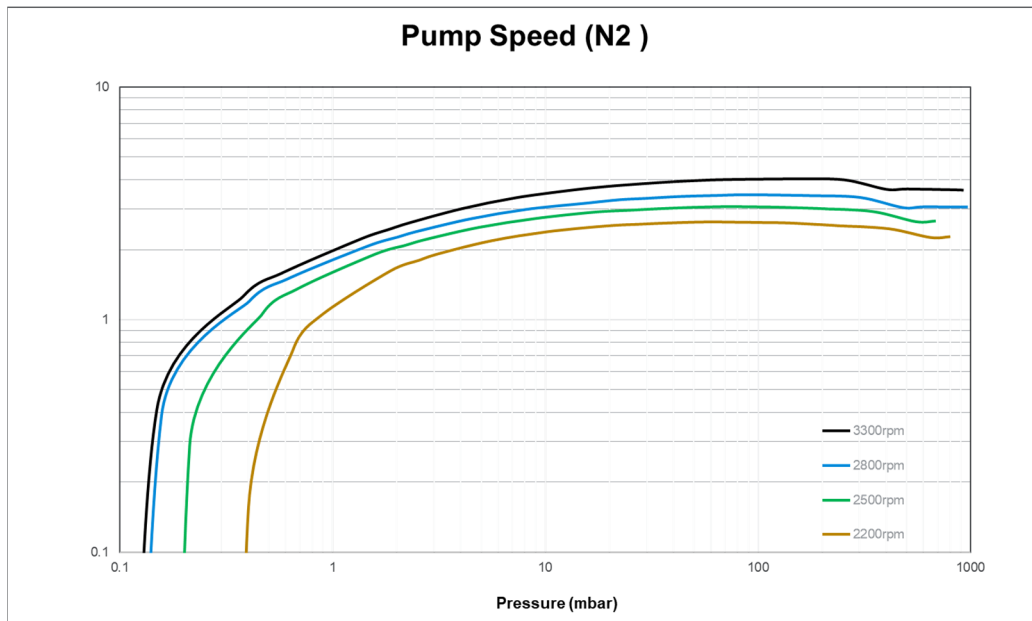


Figure 24 IDP-4 Pumping Speed (Average) – without inlet valve

Technical Information

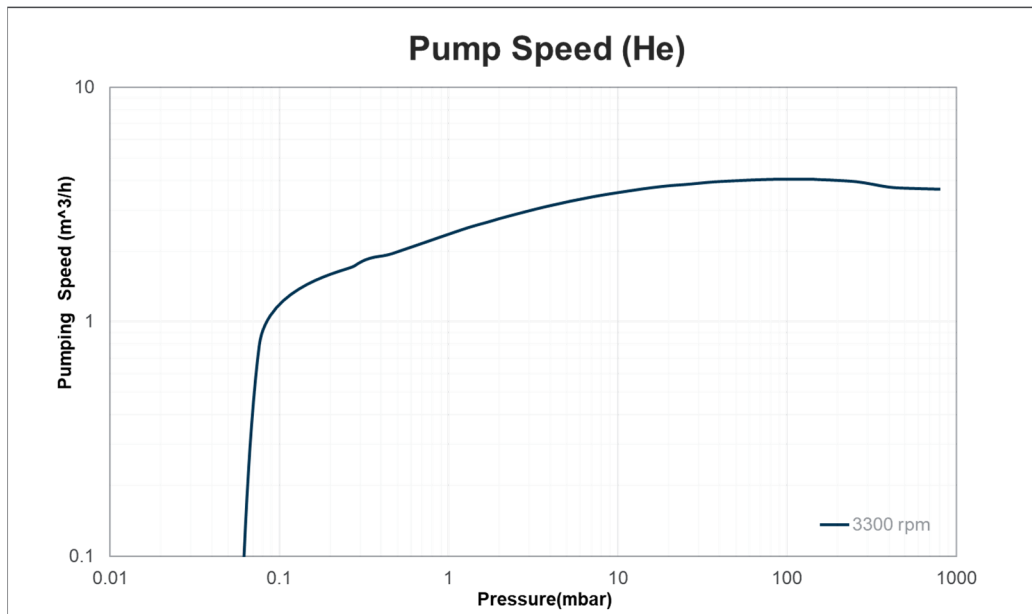


Figure 25 IDP-4 Power Draw Curves

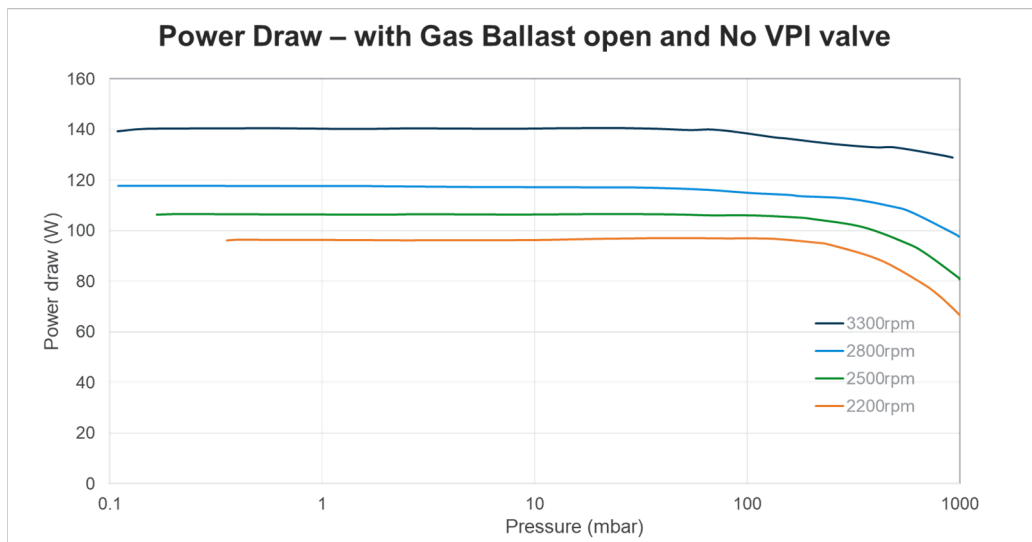


Figure 26 IDP-4 Power draw without valve (average)

Technical Information

Electrical Connections

Power Cord

Several power cord options equipped with an IEC-320 connector are available from your Agilent dealer. Descriptions of the available power cords and their ordering numbers are given in the following table.

Table 2 Standard IDP-4 Power Cord Options

Country	Power Cord Specification	Part Number
Europe	10 A / 220-230 VAC, 2.5 m	656494220
Denmark	10 A / 220-230 VAC, 2.5 m	656494225
Switzerland	10 A / 230 VAC, 2.5 m	656494235
UK/Ireland	13 A / 230 VAC, 2.5 m	656494250
India	10 A / 220-250 VAC, 2.5 m	656494245
Israel	10 A / 230 VAC, 2.5 m	656494230
Japan	12 A / 100 VAC, 2.3 m	656494240
North America	10 A / 230 VAC, with C13 2.5 m (NEMA 6-15)	656494255
North America	13 A / 125 VAC, 2.0 m	656458203

Grounding Instructions

WARNING



This product should be grounded. In the event of an electrical short circuit, grounding reduces the risk of electric shock by providing an escape path for the electric current. This pump is equipped with a power cord that has a grounding wire with an appropriate grounding plug.

WARNING



If the product must be reconnected for use on a different type of electric circuit, the connector should be replaced by qualified service personnel.

Technical Information

Extension Cords

In general, extension cords are not recommended for use with this product.

If you must use an extension cord with this product:

- Agilent recommends using only extension cords with a minimum of 16-gage wire and a maximum length of 25 feet (7.6 m).
- Use only a 3-wire extension cord that will accept the plug.
- Make sure your extension cord is in good condition.
- Be sure the extension cord is rated high enough to carry the current your product will draw. An undersized cord will cause a drop in line voltage resulting in loss of power and overheating.

The I/O and Serial Ports should only be connected to external control circuitry which is properly isolated from line voltage.

Proper installation will reduce the influence of external noise and help to meet the EMC requirements. Cables for remote I/O or serial communication must not exceed 30 m length. Contact Agilent technical support if assistance is needed to make special communication cables.

NOTE

Take care to have good, secure contact between the metallic case and the external shield of the wire. This operation must be done at least on the controller side.

Inrush Currents

The IDP-4 Dry Scroll pump limits inrush currents by utilizing digital motor control. This pump takes in AC power and delivers DC power to a brushless DC motor. At pump start, the drive frequency is ramped up over several seconds. This provides for smooth starting and limits the inrush current to the maximum steady state running current.

Technical Information

Mechanical Connections

Pump Inlet Connection

For optimized vacuum performance use NW16, or larger, clean vacuum hardware with as short a length as practical between the pump and vacuum chamber to reduce conductance losses. Inserting a bellows type hose between the pump and vacuum chamber is recommended as it will provide both vibration isolation and strain relief.

CAUTION

Do not connect heavy items directly to the inlet or exhaust fittings of the pump as it could result in fatigue failures in the pump's inlet or exhaust fittings. It is recommended instead to use a flexible bellows type of hose with a separate mechanical support for heavy items.

Pump Exhaust Connection

The pump is fitted with a NW16 exhaust fitting. To avoid overheating the pump, do not restrict the exhaust flow with long lengths of small diameter tubing. Use as short as practical lengths of 10 mm ID, or larger, diameter hardware. If an exhaust filter is used, make sure that it is cleaned or replaced on a regular basis.

WARNING



This product should only be used to pump air and inert gases only. This pump is not intended for use with corrosive, pyrophoric, particulate-forming gases or gas mixtures in their flammable and/or combustible range.

WARNING



Do not use gas ballast if using potentially hazardous materials.

Technical Information

CAUTION Operate the pump in ambient temperatures only between 5 °C to 40 °C (41 °F to 104 °F), otherwise damage to the pump or shortened operating life may result.

CAUTION Do not block the fan ducts. Blocking these ducts can cause pump overheating. Note that the pump will automatically shut off and could potentially impact the service life of the pump.

CAUTION Note that the pump will automatically shut off in the event of a discharge pressure in excess of the maximum recommended 6.5 psig back pressure or in the event of an elevated pump head temperature caused by an excessive exhaust back pressure.

CAUTION Provision, such as an Integrated Isolation Valve, should be incorporated in the vacuum system upstream to the rouging pump to protect damage to components (i.e., a turbomolecular pump) from a sudden venting to atmosphere, such as in the event of an over temperature or excess exhaust pressure which would cause the pump to shut off.

Gas Ballast

The standard configuration of this pump is shipped with the gas ballast port sealed off by a 1/4-20 thread screw (item 5 on Principal Items Figure 28). This configuration is suitable for relatively dry applications, however, in presence of water vapor, solvents, etc., are being pumped, remove the seal screw and replace it with Part Number IDP3GBKIT, which includes an adapter and 1/8" NPT sintered filter plug. Do not exceed 25 in-lb of torque when installing the adapter into the manifold.

For applications where the ingress of atmospheric air is undesirable, dry nitrogen at a flow rate of approximately 5 l/min can be provided to the gas ballast.



Figure 27

NOTE

Water handling capability is 4 g/hr under typical conditions, with up to 10 g/hr in higher pressure ranges.

CAUTION

Even very low amounts of water or condensable gases could require the use of gas ballast, as liquid will slowly, but continuously accumulate in the pump eventually resulting in loss of vacuum performance.

WARNING



The gas ballast must be sealed whenever pumping any gas not intended to be vented to atmosphere.

Technical Information

CAUTION

Pumping high water vapor loads can cause a temporary increase in ultimate pressure. Pumping water vapor loads higher than the water vapor handling capability of the gas ballast will temporarily compromise the vacuum performance and can reduce tip seal life.

CAUTION

Operating the pump continuously without periodically activating the gas ballast on a regular basis will in most applications result in a buildup of water in the pump resulting in a severe but recoverable loss of vacuum performance and in addition a reduced tip seal life.

The most common cause of poor performance on initial install and startup is adsorbed water vapor in both the pump and vacuum chamber. Running the pump with the gas ballast activated may also be beneficial if the pump has been shut off for a period of time, which has allowed water to adsorb inside both the pump and vacuum system. Depending on the vacuum system and amount of adsorbed water it can typically take between ~30 minutes to ~24 hours of gas ballast activation to achieve the full vacuum performance.

CAUTION

To achieve the published vacuum performance the pump must be run with the gas ballast activated for a period of time after the initial installation and startup or if the pump has been in storage in order to remove any water from both the pump and vacuum chamber. Not activating the gas ballast is the most common reason for unnecessary downtime.

Accessories and Spare Parts

Optional Accessories

The accessories listed in the following table are available for use with the IDP-4 Dry Scroll Pump.

Table 3 Accessories

Part Number	Description
IDP3GBKIT	Gas Ballast kit
VPI16IDP24DC	VPI Valve-24VDC, IDP-4
EXSLRIDP3	Exhaust Silencer Kit
IDP3GBKIT	Breather Vent and Adapter
X3802-68001	Serial communication cable kit

Technical Information

Optional Inlet Valve Kit

The IDP-4 Scroll Pump is equipped with an exhaust valve to prevent rapid leak-up to atmosphere when turned off. However, any vacuum pump generates a small increase in inlet line pressure when turned off. A small amount of gas or pump debris can be carried back into a vacuum line or chamber when the pump is turned off. To prevent this in applications where the process is sensitive to pump debris, install a fast acting, automatic, normally closed valve.

For your convenience, Agilent offers an optional inlet valve and operator kit for the IDP-4 (see following figure).



Figure 28 IDP-4 with optional inlet valve

- The opening of this valve occurs between 8 - 10 seconds after pump startup.
- Valve closing occurs ≤ 25 msec after pump shut off.

Technical Information

The inlet valve kit can be obtained separately for existing standard IDP-4 pump. Refer "Installation" section for detailed procedures.

The inlet valve kit can be obtained separately for existing standard IDP-4 Dry Scroll pump. Refer to kit part number VPI16IDP24DC for upgrade or purchase model X3802-64010 (IDP-4 Std Valve).

The IDP-4 Inlet valve technical specifications are shown in the following table.

Table 4

Leak Rate:	
Body	< 1X10 ⁻⁹ atm cc/sec He
Seal	<1X10 ⁻⁹ atm cc/sec He
Closing time (after solenoid valve power loss)	25 ms
Closing burst	0.01 torr-liter
Materials:	
Body	Aluminum (6061-T6)
Seal	Viton
Temperature Range:	
Valve Body	0~100 °C
Solenoid	0~50 °C

Installation

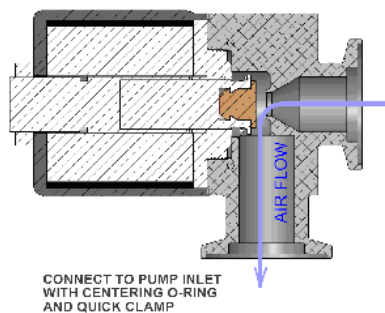


Figure 29 Mount Inlet Valve with IDP-4 Pump

To install:

1. Connect the solenoid valve to the pump inlet with a centering o-ring and quick clamp as shown in Figure 26. Plug the connector at the end of the solenoid valve to the 24 V DC output as shown in Figure 27.

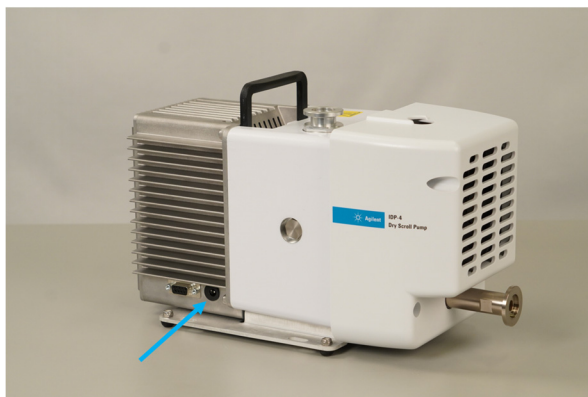


Figure 30

WARNING



The Integrated Isolation Valve should never be used for long term isolation of the vacuum chamber, and in no circumstances should vulnerable equipment such as turbomolecular pump or diffusion pump be operated when the backing pump is off.

Cleaning and Preventive Maintenance

Personnel responsible for pump operation and maintenance must be properly trained and aware of accident prevention rules.

WARNING



- **Death may result from contact with high voltages. Always take extreme care and observe accident prevention regulations in force.**
 - **When pump is powered up, be careful of moving parts and high voltages.**
 - **If you must perform maintenance on the pump, allow it to cool to ambient temperature as the temperature may be in excess of 60 °C.**
 - **Always disconnect the power supply to the pump and the power cable unplugged from pump before beginning any maintenance work.**
 - **Pump has auto-restart features. There is a risk of finger damage from the cooling fan after a thermal trip and auto restart. Always disconnect the pump from mains power before doing any service operation.**
-

External cleaning

The exterior surfaces of the IDP-4 pump may be cleaned with a cloth slightly moistened with water. If necessary, isopropyl alcohol or a mild detergent may be used instead. Do not pour or spray any liquids onto the pump.

Do not use solvents or cleaning products that could damage the plastic cowlings.

Internal cleaning as part of a Preventive Maintenance plan

Unlike conventional oil filled vacuum pump, Agilent Dry Scroll pumps do not contain oil which requires periodic replacement. During the typical operation of the pump, the pump will accumulate a minor buildup of tip seal dust and condensed water within the pumping mechanism – especially if being used to pump atmospheric air. Proactive gas ballasting and venting of tip seal dust described below are important aspects of a preventive maintenance plan as the combination of tip seal dust and trapped water has been known to result in premature tip seal wear.

To remove condensed water vapor buildup, periodically activate the gas ballast. It is recommended that the pump's gas ballast be activated at the initial installation of the pump, after maintenance, and as often as needed to remove water buildup which can result in increased base pressure. Until experience is gained with the specific application, the schedule found in Table 5 can be used as a starting point. Once the rate of buildup of water vapor in the pump is characterized for the application, this process may be required more or less frequently.

For the cleansing of accumulated tip seal dust and debris inside the pump head, run the pump periodically with the inlet cycled to atmospheric pressure multiple times for one or two minutes each. Until experience is gained with the specific application, vent the pump regularly as above, then adjust the venting schedule according to the conditions observed.

Verifying Performance

Prior to initiating any service requests on IDP-4 pump, the first step is to ensure that any loss of vacuum performance is not caused by a buildup of water in the pump. Initiating the gas ballast procedures described in the Technical Information section for ~12hrs (or typically less time) will in many cases restore the pump to its full rated vacuum performance and eliminate an unnecessary repair and associated downtime.

Agilent provides standard operating procedures for measuring base pressure and determining leak rate in the IDP-4 Preventative Maintenance Manual which should be followed whenever evaluating the vacuum performance of the pump.

NOTE

Warranty repair requests found to be the result of water build up in the pump or incorrect base pressure or leak rate measurements will not be covered.

Tip Seal Replacement

The life of a tip seal is highly variable depending on the application as well as inlet pressure(s) and accumulated water, but Agilent recommends that until experience is gained on the application, the schedule found in the table below can be used as a starting point. Replacing the tip seal on a regular basis will avoid unplanned downtime. In general, it is recommended to change the tip seals before they need to be changed.

Table 5 Suggested maintenance based on duty level

	Light-Duty	Medium-Duty	Heavy-Duty
Description	Intermittent use Pump is shut down between experiments.	Continuous low-pressure use. Ex. turbo backing Low gas load	Continuous high inlet pressure use. High gas load
Vent to atmosphere	Upon start up and preparation for the experiment, vent the inlet to atmosphere for 5 seconds once the pump has reached full speed. Ensure the exhaust is not restricted.	1x per month if the application permits. Properly isolate any sensitive equipment prior to venting. Run the pump at full speed with the inlet fully open for 5 seconds. Ensure the exhaust is not restricted.	1x per year Properly isolate any sensitive equipment prior to venting. Run the pump at full speed with the inlet fully open for 5 seconds. Ensure the exhaust is not restricted.
Gas Ballast Use	Upon start up and preparation for the experiment, run the pump with the inlet capped and the gas ballast open for 4 hours to dry out any water accumulated during a preceding idle period.	Continuous gas ballast use is recommended if the application permits.	Periodic use of the gas ballast is likely required. Continuous use of the gas ballast is recommended if pumping a steady load of water vapor.
Tip Seal Replacement	Perform a minor maintenance on the pump every 10,000 hours, or when the pump's ultimate pressure cannot be reached with overnight gas ballast use.	Perform minor maintenance on the pump every 18 - 24 months. Ensure that performance degradation is not due to water vapor prior to completing maintenance.	Perform minor maintenance once per year.

Technical Information

The tip seal replacement can be performed using the kit below. It contains the necessary tip seals, the main O-ring.

Tip Seal Replacement Instructions

Table 6 Parts & Tools Required

Qty	Part Number	Description
1	IDP3TS	Replacement Tip Seal Set
1	---	4mm Allen Wrench
1	---	Razor blade or side-cutting pliers

WARNING



If dangerous gases were being pumped, flush the pump with air or inert gas for at least 10 minutes prior to disassembly.

Technical Information

The following figure shows the various components involved in a tip seal replacement procedure. Refer to the figure 28 as you follow the procedure (Video is available at Agilent.com).

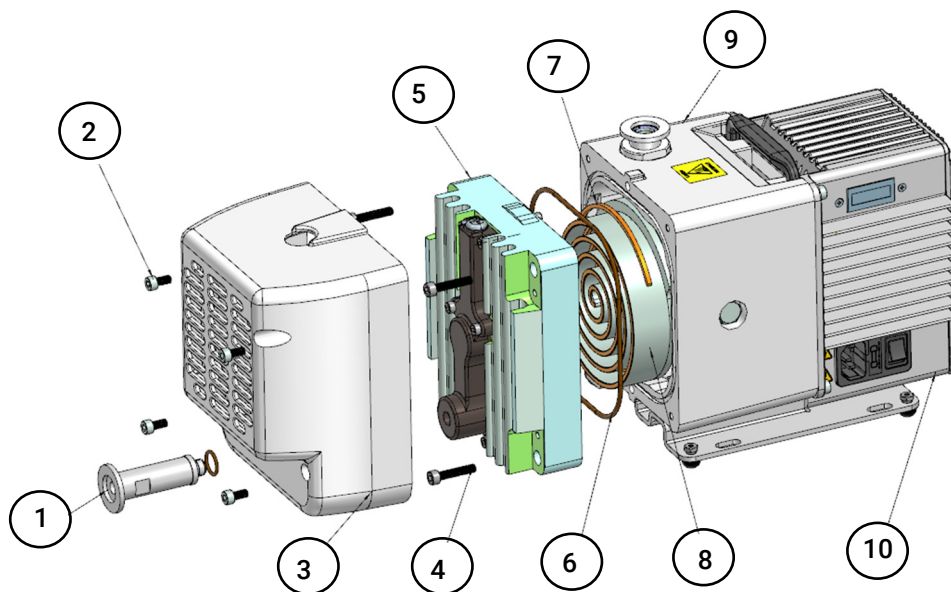


Figure 31 Exploded Diagram

1. Exhaust Fitting
2. Front Cowling Bolts M5(4)
3. Front Cowling
4. Frame Bolt, M5(4)
5. Outboard Hosing
6. Main O-ring
7. Tip Seals
8. Orbiting Plate
9. Frame
10. Motor and controller

Technical Information

Disassembling the pump and removing the worn tip seals

1. Power off the pump and disconnect it from electrical power.



2. Remove the exhaust fitting by unscrewing it from the front of the pump.



3. Remove the (4) 4mm cap head bolts fastening the cowling to the pump.



4. Remove the front cowling and disconnect the fan connector – set the cowling aside.

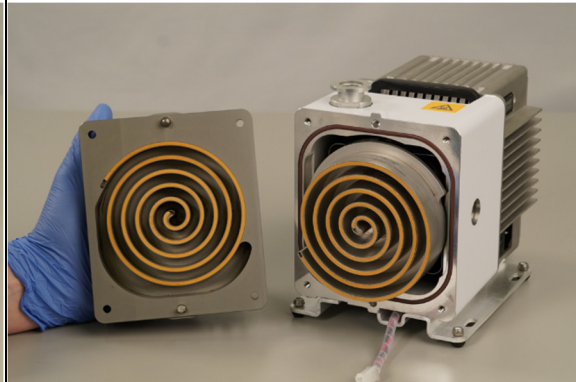


Technical Information

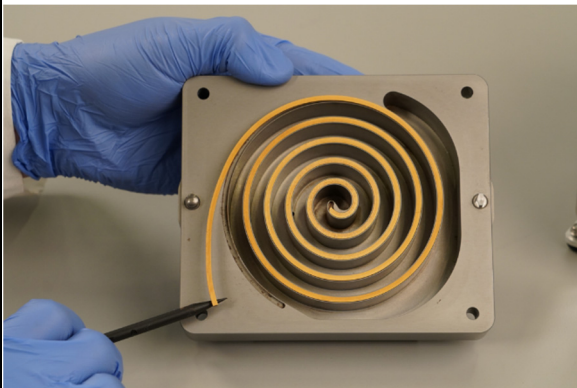
5. Remove the (4) 4mm cap-head bolts fastening the Outboard housing to the pump frame.



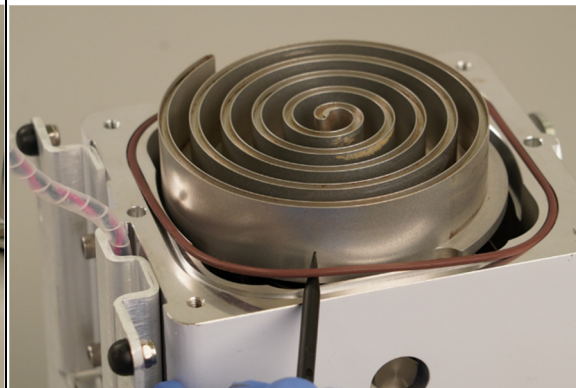
6. Remove the outboard housing axially off the frame.



7. Using a soft tool (wood or plastic pick) remove the tip seals from the Outboard housing and the outer stage of the Orbiting plate. Discard the used tip seals.



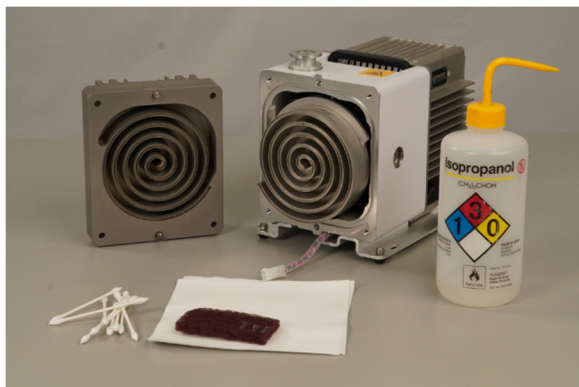
8. Remove and discard the main O-ring.



Technical Information

Cleaning the scroll surfaces and sealing surfaces:

1. Clean both working surfaces of the scrolls and remove any build up or dust from the pumping mechanism. This is best done with a soft abrasive pad, isopropyl alcohol, and lint free cloths.



NOTE Ensure that the main o-ring groove on the frame and the outboard housing sealing face are clean and free of debris. Use a lint free cloth and isopropyl alcohol to clean them if necessary.

Allow the surfaces of the scrolls to air dry before proceeding.

WARNING Do not use compressed air to blow tip seal dust off parts – this will lead to a large amount of airborne dust.



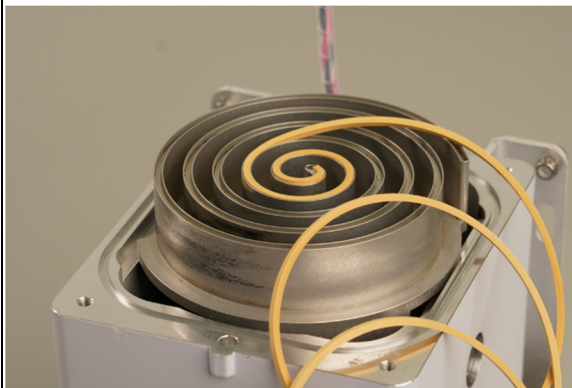
Technical Information

Replacing the tip seals and reassembling the pump:

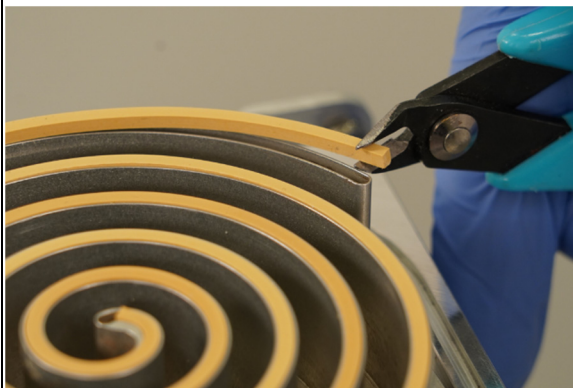
NOTE

Two tip seals are provided in the kit; one fits the orbiting scroll and the other fits the outboard housing scroll.

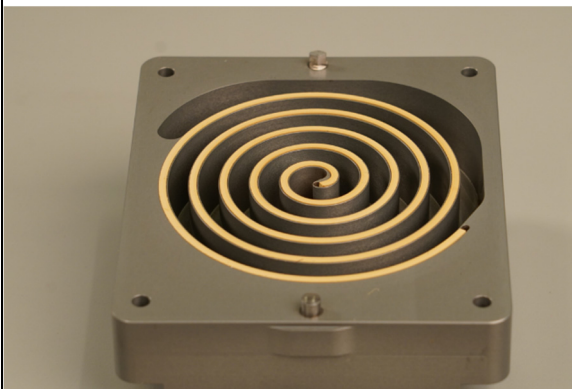
1. Starting at the center of the orbiting plate, install the tip seal into the groove, ensuring that the curl memory of the tip seal matches the curvature of the tip seal groove.



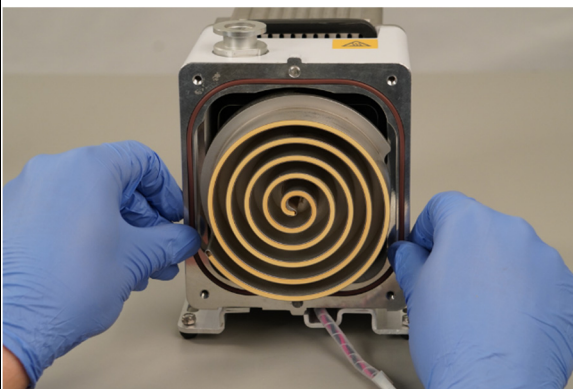
2. Once the tip seal groove has been filled, cut the tip seal so that there is a 3mm space between the end of the tip seal and the outer end of the tip seal groove.



3. Install the tip seal on the outboard housing - ensuring that the curl of the tip seal matches the curvature of the scroll. Once the tip seal groove has been filled, trim the tip seal in the same manner as step 2.

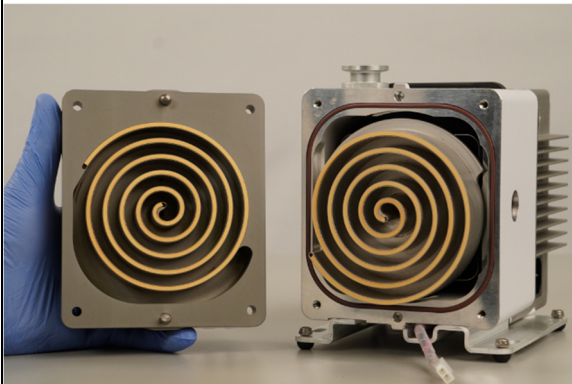


4. Place the new main O-ring into the groove in the frame.

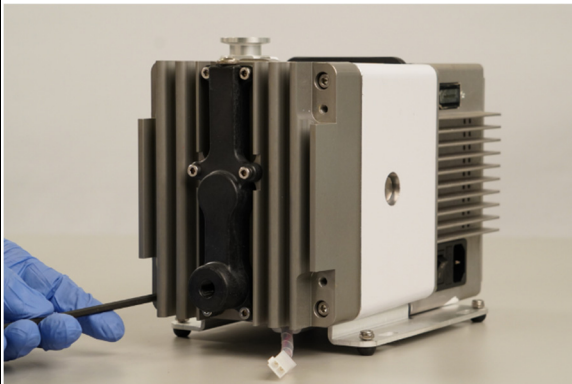


Technical Information

5. Carefully replace the outboard housing by lining up the locating pins. Ensure the tip seal has not fallen out of its groove.



6. Reinstall the (4) M5 bolts and torque to 3.4 N-m (30 in-lb).



7. Reconnect the fan/cowling to the electrical connector on the pump.



8. Reinstall the front cowling and tighten the (4) M5 bolts.



9. Reinstall the exhaust fitting.



10. Reconnect the pump to mains power and power on.



NOTE

Newly installed Tip Seals may require several hours of run time to seat properly and enable the pump to meet speed and base pressure specifications.

Testing the Pump:

Power on the pump and run it for about 5 seconds. Verify that the front fan is running.

If you hear loud noises or observe labored operation, this indicates that the Tip Seal or main O-ring are possibly out of place.

Disassemble and reassemble the pump following the relevant steps in the tip seal replacement instructions to ensure all parts are assembled correctly.

User Maintenance and Repair

The following table provides the list of part numbers available for purchase to maintain and repair your IDP-4 Dry Scroll Pump. Additional video references related to user maintainable and replaceable parts can be found on the product page on Agilent.com.

Table 7 Maintenance Intervals and Part Numbers

Service Operation	Interval	Parts Needed	
		Part Number	Description
Regular Maintenance			
Gas Ballast Activation	See Table 5 for suggested intervals based on application	None	
Vent pump to atmospheric pressure		None	
Minor Maintenance: Tip Seal & Main O-ring replacement		IDP3TS	IDP-3, 4 Tip Seal Kit
Replace Exhaust filter element	Yearly, if exhaust silencer is installed.	REPLSLRFILTER2	FILTER ELEMENT, IDP3,4 EXHAUST
If-Required Component Replacement			
<i>Front Cowling and cooling fan</i>	If damaged or fan not functioning	X3802-67300	IDP-4 Fan & Cowling Kit
<i>Feet</i>	If broken or worn	IDP3VIBISOKIT	IDP-3, 4 Vibration Isol. Feet Kit
<i>Plastic exhaust Manifold</i>	If cracked, broken, or leaking.	X3803-67005	IDP-3, 4 Std Exhaust mfl'd Kit
<i>VPI Valve</i>	If leaking or damaged	VPI16IDP24DC	Sub-Assy, VPI Valve-24VDC, IDP3, 4
<i>Standard Exhaust Flange</i>	If damaged	X3802-67120	IDP-4 Std Exhaust Flange Kit

Technical Information

Factory-Only Service

Some maintenance tasks and repairs can only be performed by Agilent service personnel. See the next section for service options.

Table 8 Factory Only Service

Service Operation	Interval
Major Maintenance: Includes a tip seal replacement plus the replacement of all bearings, O-rings, and poppet valves.	Every 5 years of continuous operation, or on audible bearing noise. May be required more frequently for aggressive applications.
Motor replacement	Every 10 years of continuous operation or if required.
Drive Train Components Repair (shaft, scrolls, bearings).	If damaged or noisy.

Service Options

In the event of performance degradation or loss of function of the IDP-4 pump there are several options to facilitate the repair or exchange of the vacuum pump.

Factory Repair

If more extensive repairs on the motor or drive train are required (bearings, crankshaft, scrolls), Agilent offers factory repair services. Factory repair entails sending the pump to Agilent for a complete rebuild and resolution of performance issues. Once complete, the pump will be tested and return shipped to the customer. Contact your local Agilent representative for more information.

Advance Exchange

Agilent also may offer Advance Exchange service for some models. If available, fully rebuilt and tested pumps are kept in stock worldwide ready to ship. Once the rebuilt unit is received by the customer, they must package the defective module or full pump in the advance exchange packaging provided and return it to Agilent. Contact your local Agilent representative for more information and for active exchange part numbers.

NOTE

For further service options, please contact your local service center listed in the back of this manual and ask for your local representative.

Should a customer need service, please contact local distributor or directly e-mail to:

vpt-customer@agilent.com

vpl-customer@agilent.com

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

Troubleshooting

Use the Troubleshooting chart (Table 9) to assist in defining a problem, determining a possible cause, and defining action steps to remedy the situation.

Table 9 Troubleshooting Chart

Problem	Possible Causes	Corrective Actions
Will not start.	Motor or controller thermal protector open	Allow the pump to cool. Identify the cause of overload.
	Excessive voltage drop in the power cord.	Check the size and length of the cable.
	Defective motor	Inspect. Contact Agilent.
	Pump jammed	Inspect. Contact Agilent.
Poor ultimate pressure	System leak	Locate and repair the leak.
	Water vapor in the pump or system	Flush pump with air or dry nitrogen.
	Water in the pump	Activate the Gas Ballast.
	Gas ballast plugged	Replace breather vent. Contact Agilent.
	Solvent in the pump	Activate the Gas Ballast. Install trap or filter.
	Worn tip seal	Replace tip seals.
	Poor conductance to pump	Replumb with shorter and/or larger diameter tubing.
Hammering noise	Low pumping speed.	Check pump RPM setting. Set RPM to expected value.
	Pump overheated	Check the ventilation to pump. Check ambient temperature.
	Check the ambient temperature	Check the intake screen. Flush pump. Disassemble the pump head and inspect.
Debris in pump	Debris in pump	Check intake screen. Flush pump. Disassemble pump head and inspect.
	Pump runs intermittently	Motor or controller thermal protector is cycling open and closed. *

Technical Information

Table 9 Troubleshooting Chart (continued)

Problem	Possible Causes	Corrective Actions
Motor speed intermittent slowing	Power limit to protect from high power draw due to high torque on the motor,	Identify the cause of overload.
Pump stops	Overcurrent protection is triggered*	Identify the cause of overload.
Pump does not pump for first 30 seconds.	The Optional Isolation Valve is closed until the pump ramps to a minimum rpm to achieve vacuum.	Check the connection of the isolation valve.
Hour meter display is blank.	Internal pump controller failure or no AC power to the pump.	Check branch circuit breaker plug AC power cord into unit and/or wall outlet. Contact Agilent.
Grinding noise which gets slowly louder over time.	Possible Bearing Failure	Pump Module rebuild or exchange

- * The IDP-4 pump is equipped with an auto-reset thermal motor protector. This protector will automatically shut down the pump when it detects an overload condition and will automatically restart the pump when the motor has cooled to within an acceptable temperature range.

Serial Connector & Pin allocation

Pump Serial Connector

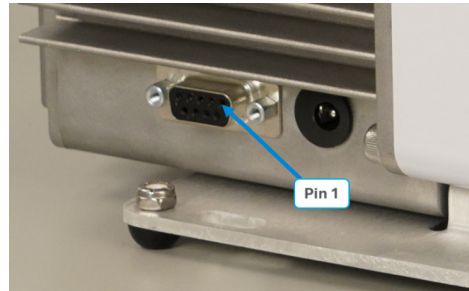
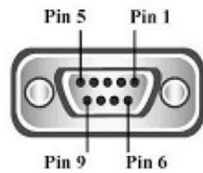


Figure 32 Female DSUB9 Pin Connections

Table 10 DSUB9 Serial Pin Connections

Pin #	Function
1	+5V 50mA
2	Not Connected
3	Not Connected
4	For Factory Use Only
5	0V
6	RS485 (A+)
7	For Factory Use Only
8	RS485 (B-)
9	Not Connected

CAUTION

Pin1 & Pin0: Contains an auxiliary output 5VDC 50mA power supply. Make sure not to exceed the 50mA limit.

Serial communication settings

The serial communication uses the RS485 protocol.

To read/change RS485 address read/write WIN503.

See WINDOW protocol description for further details.

- Baud Rate: 4800, 9600 (default), 19200, 38400 bps
- Data Bits: 8
- Parity: None
- Stop Bits: 1
- Flow Control: None

Window Description and Protocol

Window Description

Table 11 Window meanings

Win #	Read	Write	Type	Name	Meaning	Default (Factory Set)	Unit
0	Always	Always	L	Start/Stop	Start/Stop command	-	-
108	Always	Always	N	Baud Rate	It defines the serial communication baud rate: 3 = 4800 bit/s 4 = 9600 bit/s 5 = 19200 bit/s 6 = 38400 bit/s	4 = 9600 bit/s	Index
127	Always	Always	N	Rotation Frequency Setting	It defines the target rotation frequency in Hz	3300	Rpm
200	Always	Off	N	Bus current	It reads the actual value of the bus current supplied by controller to electric motor	-	0.1 Ampere
201	Always	Off	N	3phase voltage	It reads the voltage applied to the motor windings	-	Volt
202	Always	Off	N	Output power	It reads the actual value of power supplied by controller to electric motor	-	W
205	Always	Off	N	Pump status	Real time pump status: 0 = Stop 2 = Start 5 = Normal 6 = Fail	-	Index
206	Always	Off	N	Error code	This location has a bit significant: 31-13 – RESERVED 12 – Locked Rotor 11 – RESERVED 10 - RESERVED 9 - RESERVED 8 - RESERVED 7 - RESERVED 6 - Bus Overvoltage 5 - Hall Error 4 - Motor Overtemperature 3 - Heatsink Overtemperature 2 - RESERVED 1 - RESERVED 0 - System Fault	-	bitfield
216	Always	Off	N	Board Temperature	It reads the Heatsink Assembly temperature	-	°C

Technical Information

Table 11 Window meanings (continued)

Win #	Read	Write	Type	Name	Meaning	Default (Factory Set)	Unit
225	Always	Always	L	Inlet valve control	Control inlet valve open or close command This function is disabled during the 10-second startup delay that occurs when power is supplied to the pump or the power switch is turned ON.	-	-
227	Always	Always	L	Gas ballast valve control	Control gas ballast open or close command	-	-
232	Always	Off	N	Rotational Frequency	Rotational Reading Hz	-	0.1Hz
233	Always	Off	N	Rotational Frequency	Rotational Reading RPM	-	RPM
234	Always	Off	N	Bus Voltage	It reads the Bus Voltage	-	Volt
236	Always	Off	N	Torque	Estimate Mechanical Torque	-	N-mm
300	Always	Off	N	Cycle time	It indicates the time of the last complete cycle power on to power off	-	Minutes
301	Always	Off	N	Cycle number	It indicates the number of cycles start/stop during all the pump life.	-	-
302	Always	Off	N	Pump life	It indicates the total hours of pump life. The motor is running.	-	Hours
303	Always	Off	N	Total Error Counts	Total errors counts for life.	-	-
319	Always	Always	A	Controller Part Number	It is the Part Number of the controller	-	-
320	Always	Always	A	Pump Part Number	It is the Part Number of the pump	-	-
321	Always	Always	A	Pump Special Model	It indicates the Part Number of special pumps	-	-
322	Always	Always	A	Pump Serial Number	It indicates the Serial Number of the pump	-	-
323	Always	Always	A	Controller Serial Number	It indicates the Serial Number of the controller	-	-
324	Always	Always	A	Pump Modification Level	It indicates the upgrading level of the pump	-	-
325	Always	Always	A	Controller Modification Level	It indicates the upgrading level of the controller	-	-
358	Always	Always	N	Pump life since last service	It indicates the days after the last resetting	-	Days
399	Always	Off	N	Controller life	It indicates the total hours of Controller life. The total time the controller is powered on.	-	Hours

Technical Information

Table 11 Window meanings (continued)

Win #	Read	Write	Type	Name	Meaning	Default (Factory Set)	Unit
400	Always	Off	A	Firmware CRC	It defines the CRC of the actual embedded Firmware	-	-
401	Always	Off	A	Bootloader CRC	It defines the CRC of the embedded bootloader program	-	-
406	Always	Off	A	Firmware Revision	It indicates the Firmware Revision	-	-
503	Always	Always	N	RS485 Address	It defines the serial address in RS485 communication (accessible with every serial address)	-	-
710	Always	Always	N	Fail history win -ROW number	This is a LOG function. Indicates the row containing the error. Refer to Note below.	-	-
711	Always	Always	N	Fail history win - COLUMN number	This is a LOG function. 1 is time, 2 is error code. Refer to Note below.	-	mins
712	Always	Off	N	Fail history win - CELL VALUE	This is a LOG function. This is the error code or timestamp. Refer to Note below.	-	Mins/bitfield

NOTE

The IDP-4 retains a log of past error codes (from WIN#206) in its memory even after loss of power. Past error codes can be viewed by navigating the Error Log Table with WIN#710 and WIN#711 and displaying the values with WIN#712. These windows can be used to navigate the Error Log Table as described below.

- WIN#710 – Row Selector (range 0 – 22)
- When the table is full, new entries overwrite rows starting from 0.
- WIN#711 – Column Selector (1 = Timestamp, 2 = Error Code)
- WIN#712 – Value Display

Examples:

1. After setting WIN#710 to 5 and WIN#711 to 2, WIN#712 will show the error code in row 5.
2. After setting WIN#710 to 5 and WIN#711 to 1, WIN#712 will show the time stamp of the error in row 5.

Windows Protocol

Refer to Agilent Technical Support for details on programming.

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
Toll Free: 00 800 234 234 00

vpt-customer@agilent.com

NORTH AMERICA:

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

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:	Fax:
	Email:	

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

Type of return	
<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"/> 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"/>	

**I declare that the above information is true and complete to the best of my knowledge and belief.
I understand and agree to the terms and conditions on page 2 of this document.**

Name:	Authorized Signature:
Position:	
Date:	

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

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies

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

Brazil

Agilent Technologies Brasil

Avenida Marcos Penteadro 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 Co. Ltd (China)

No.3, Wang Jing Bei Lu, Chao Yang District
Beijing, 100102, China
Tel: +86 (0)10 64397888
Fax: +86 (0)10 64397666
Toll free: 400 8203278 (mobile)
Toll free: 800 8203278 (landline)
vacuum_cnmarketing@agilent.com
vpc-customerservice@agilent.com

France

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

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

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

Japan

Agilent Technologies Japan, Ltd.

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

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

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

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-care@agilent.com

© Agilent Technologies, Inc. 2022
10/2022

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
Tel: +886 3 4959004
Toll free: 0800 018 768
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.Ili 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-care@agilent.com

Customer Support & Service

NORTH AMERICA:

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

EUROPE:

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

PACIFIC RIM:

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

Worldwide Web Site, Catalog and Order On- line:

www.agilent.com/chem/vacuum
Representatives in most countries

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

This information is subject to change without notice.

© Agilent Technologies, Inc. 2025

Agilent Technologies LDA Malaysia
Sdn Bhd Bayan Lepas
Free Industrial Zone 11900 Bayan Lepas,
Penang Malaysia



D0129654

Edition A.00
November 2025