

Agilent MS120 Single Stage Rotary Vane Pump (X3702-64000)

User Manual



Notices

Manual Part Number

87-900-150-01
Edition D.00, July 2024

Copyright

© Agilent Technologies, Inc. 2024

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

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

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	11
Informazioni su questo manuale	13
Validità	13
Definizioni e terminologia	14
Definizione di Attenzione, Avvertenza e Nota	14
Simboli di avvertenza	15
Sicurezza	17
Uso corretto	17
Uso improprio	19
Tabella limiti operativi	19
Attrezzatura di protezione	21
Linee guida per la sicurezza delle pompe MS120 Single Stage Rotary Vane	22
Descrizione del prodotto	24
Trasporto e immagazzinamento	25
Sollevamento	26
Installazione	27
Preparazione per l'installazione	27
Accessori forniti con la pompa	28
Assemblaggio	28
Ubicazione	29

Contents

Collegamento alla macchina utilizzatrice	30
Convogliamento aria di scarico	31
Collegamento elettrico	32
Disimballaggio	33
Utilizzo	36
Avvio e funzionamento della pompa MS120 Single Stage Rotary Vane	36
Consigli per l'utilizzo	37
Zavorratore (Gas-ballast)	37
Aspirazione vapore acqueo	38
Manutenzione	39
Sostituzione olio lubrificante	40
Sostituzione elementi disoleatori	42
Manutenzione ordinaria	44
Ricondizionamento pompa	44
Come ordinare i ricambi	44
Lubrificanti	45
Messa fuori servizio	45
Troubleshooting	46
Smaltimento	47
Servizio Post Vendita	48

Contents

2 Betriebsanleitung	49
Informationen über diese Betriebsanleitung	51
Gültigkeit	51
Definitionen und Terminologie	52
Definition von Vorsicht, Warnung und Hinweis	52
Warning Symbols	53
Sicherheit	55
Sachgemäße Verwendung	55
Unsachgemäße Verwendung	57
Tabelle mit Betriebs-Grenzwerten	57
Schutzausrüstung	59
Sicherheitsrichtlinien für die Single Stage Rotary Vane MS120 Pumpe	60
Produktbeschreibung	63
Transport und Aufbewahrung	64
Anheben	65
Einbau	66
Montagevorbereitungen	66
Mitgeliefertes Zubehör	67
Zusammenbau	67
Lage	68
Anschluss an die Maschine	69
Umleitung der Abgasluf	70
Elektrische Anschlüsse	71
Auspacken	72
Betrieb	75
Inbetriebnahme und Betrieb der Pumpe MS120 Single Stage Rotary Vane	75

Contents

Tipps für den Betrieb	76
Gas-ballast	77
Ansaugung von Wasserdampf	77
Wartung	78
Ersetzung des Schmieröls	79
Ersetzung der Ölreinigungselemente	81
Routinewartung	83
Revision der Pumpe	83
So bestellen Sie Ersatzteile	83
Schmiermittel	84
Stilllegung	84
Fehlerbehebung	85
Entsorgung	86
Service	87
3 Mode d'emploi	88
À propos de ce manuel	90
Validité	90
Définitions et terminologie	91
Signification des messages d'attention, d'avertissement et des notes	91
Symboles d'avertissement	92
Sécurité	94
Utilisation appropriée	94
Utilisation incorrecte	96
Tableau des limites opératoires	96
Équipements de protection	98
Consignes de sécurité relatives aux pompes MS120 Single Stage Rotary Vane	99

Contents

Information du produit	101
Transport et entreposage	102
Levage	103
Installation	104
Préparation pour l'installation	104
Accessoires fournis avec la pompe	105
Assemblage	105
Positionnement	106
Raccordement à la machine utilisatrice	107
Acheminement de l'air d'extraction	108
Raccordement électrique	109
Déballage	110
Utilisation	113
Démarrage et exploitation de la pompe MS120 Single Stage Rotary Vane	113
Conseils d'utilisation	114
Dispositif de lest (Gas-ballast)	115
Aspiration de vapeur d'eau	115
Entretien	116
Remplacement de l'huile lubrifiante	117
Remplacement des éléments déshuileurs	119
Entretien de routine	121
Révision pompe	121
Accessoires à commander	121
Lubrifiants	122
Mise hors service	122
Problèmes et solutions	123
Élimination	124
Service	125

Contents

4	Instructions for Use	126
	About this manual	127
	Validity	127
	Definitions and terminology	128
	Definition of Caution, Warning and Note	128
	Warning Symbols	129
	Safety	131
	Proper use	131
	Improper use	133
	Operating limits table	133
	Protective equipment	135
	MS120 Single Stage Rotary Vane Pump Safety Guidelines	136
	Product Description	139
	Transport & Storage	140
	Lifting	141
	Installation	142
	Preparation for Installation	142
	Accessories supplied with the pump	143
	Assembly	143
	Location	144
	Connection to the machine	145
	Discharge air pipe line installation	146
	Electrical connection	147
	Unpacking	148
	Use	151
	Start-up and Operation of MS120 Single Stage Rotary Vane pump	151

Contents

Tips for using	152
Gas-ballast	152
Water vapor intake	153
Maintenance	154
Oil change	155
Exhaust filters replacement	157
Ordinary maintenance	159
Pump overhaul	159
How to order spare parts	159
Lubricants	160
Decommissioning	160
Troubleshooting	161
Disposal	162
Service	163

Contents

5	Technical Information	164
	Description of the MS120 Single Stage Rotary Vane Pump	165
	Vacuum Seals	166
	Technical Specifications	167
	MS120 Single Stage Rotary Vane Pump Outline Drawing	169
	Connection to the Electric Supply	171
	Starting and Running the Pump	172
	Stopping the Pump	172
	User Interface	173
	Input/Output Communications	174
	J1 – I/O: Remote I/O signals + Serial communication	174
	Serial Communication Connection Examples	176
	How to connect the I/O port to an external cable	177
	Serial Communication Protocol	178
	Communication Format	178
	Communication Protocol	178
	Examples	181
	Window Meanings	182
	Pump Status	184
	Accessories and Spare Parts	185

1

Istruzioni per l'uso

Informazioni su questo manuale	13
Validità	13
Definizioni e terminologia	14
Definizione di Attenzione, Avvertenza e Nota	14
Simboli di avvertenza	15
Sicurezza	17
Uso corretto	17
Uso improprio	19
Tabella limiti operativi	19
Attrezzatura di protezione	21
Linee guida per la sicurezza delle pompe MS120 Single Stage Rotary Vane	22
Descrizione del prodotto	24
Trasporto e immagazzinamento	25
Sollevamento	26
Installazione	27
Preparazione per l'installazione	27
Accessori forniti con la pompa	28
Assemblaggio	28
Ubicazione	29

Istruzioni per l'uso

Collegamento alla macchina utilizzatrice	30
Convogliamento aria di scarico	31
Collegamento elettrico	32
Disimballaggio	33
Utilizzo	36
Avvio e funzionamento della pompa MS120 Single Stage Rotary Vane	36
Consigli per l'utilizzo	37
Zavorratore (Gas-ballast)	37
Aspirazione vapore acqueo	38
Manutenzione	39
Sostituzione olio lubrificante	40
Sostituzione elementi disoleatori	42
Manutenzione ordinaria	44
Ricondizionamento pompa	44
Come ordinare i ricambi	44
Lubrificanti	45
Messa fuori servizio	45
Troubleshooting	46
Smaltimento	47
Servizio Post Vendita	48

Informazioni su questo manuale

Validità

Questo manuale elenca le istruzioni per gli utenti della pompa MS120 Single Stage Rotary Vane, con particolare riferimento alle nozioni relative a sicurezza, funzionamento e manutenzione di primo livello, limitatamente alle operazioni di manutenzione di cui l'utente è responsabile.

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

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

NOTA

- 1 Questo manuale contiene informazioni utili affinché tutto il personale che utilizza la pompa MS120 Single Stage Rotary Vane possa utilizzarla in sicurezza e garantire la perfetta efficienza, per tutto il suo ciclo di vita.
 - 2 Conservare questo manuale, insieme a tutte le pubblicazioni ad esso correlate, in un luogo accessibile, conosciuto da tutti gli operatori/personale di manutenzione.
-

Definizioni e terminologia

Definizione di Attenzione, Avvertenza e Nota

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

ATTENZIONE

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

AVVERTENZA



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

NOTA

Le note contengono informazioni importanti e forniscono maggiori dettagli su passaggi specifici.

Simboli di avvertenza

Di seguito si riporta un elenco di simboli che vengono visualizzati insieme agli avvisi della pompa MS120 Single Stage Rotary Vane. Viene mostrato anche il pericolo che descrivono. Un simbolo triangolare indica un'avvertenza. I significati dei simboli che possono apparire accanto alle avvertenze nella documentazione o sullo strumento stesso sono i seguenti.

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



	Pericolo generico
	Presenza di tensioni pericolose
	Alta Temperatura
	Corrente alternata (CA)
	Messa a terra protettiva
	Pericolo di incendio
	Non disperdere nell'ambiente
	Leggere il manuale d'uso
	Attacco in aspirazione
	Attacco allo scarico

Istruzioni per l'uso

	Guanti protettivi
	Scarpe antinfortunistiche
	Attrezzatura di protezione
	Emissione di sostanze nocive
	Rischio lesioni a causa di spigoli vivi
	Caduta di oggetti
	Certificazione CSA
RoHS	Limitazione della Certificazione delle Sostanze Pericolose
CE	Certificazione CE
	Rifiuti di Apparecchiature Elettriche ed Elettroniche
UK CA	Marcatura UK CA

Sicurezza

Questa sezione contiene le informazioni prescritte dalla direttiva sulla bassa tensione 2014/35/UE, essenziali per la conformità e l'osservanza delle norme di sicurezza sia in generale sia in relazione all'uso specifico del prodotto.

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

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

AVVERTENZA



Nei casi di impiego in cui l'arresto o un guasto della pompa per vuoto possa causare danni a persone o cose, devono essere previste delle misure di sicurezza nell'impianto.

Uso corretto

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

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

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

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

AVVERTENZA



Materiali infiammabili:

Per garantire che i gas all'interno della pompa non si trovino nel rispettivo campo di infiammabilità/combustibilità, è necessario attenersi alle istruzioni e prendere nota delle precauzioni riportate qui di seguito. Quando all'interno dell'apparecchiatura sono presenti materiali potenzialmente infiammabili o combustibili, è necessario:

- Assicurarsi che il sistema sia a tenuta stagna per garantire che solo i gas previsti passino attraverso la pompa per vuoto.
- Assicurarsi che la concentrazione di tutti i gas o vapori infiammabili che entrano nella pompa (flangia di ingresso e porta di Gas Ballast) e che sono all'interno della pompa e nel condotto di scarico rimanga a meno di un quarto del limite inferiore di esplosività (LEL) ufficialmente pubblicato per i gas.
- Assicurarsi che siano presenti sistemi di sicurezza affinché la concentrazione di eventuali gas o vapori infiammabili che entrano nella pompa (flangia di ingresso e porta di Gas Ballast) e che sono all'interno della pompa e nella linea di scarico rimanga a meno di un quarto dei limiti inferiori di esplosività (LEL) ufficialmente pubblicati per i gas in qualsiasi condizione di guasto del sistema o della pompa per vuoto.

NOTA

- Prima di rimettere in funzione la pompa dopo un guasto, ispezionarla e verificare attentamente che non vi siano altri segni di danneggiamento.
- Utilizzare solo strumenti in perfetto stato di funzionamento e appositamente progettati per il lavoro da svolgere; l'utilizzo di strumenti inappropriati o inefficaci può causare gravi danni.
- Controllare sempre il lubrificante e che sia correttamente distribuito attraverso la pompa; una lubrificazione inadeguata può danneggiare seriamente la pompa.
- Dare una forma di contrassegno ai pezzi durante la rimozione per assicurarsi di rimontarli nuovamente nell'ordine corretto.
- Verificare che non siano presenti graffi o scanalature sugli alberi lavorati, nelle loro sedi all'interno della pompa o sulle superfici rettifiche della macchina. Leggeri graffi ed abrasioni possono essere eliminati con carta smerigliata molto fine o con una molatura leggera.
- Prima di assemblare un gruppo, spalmare sempre un po' di olio sulle parti interne e sulle superfici di accoppiamento. Sostituire tutte le guarnizioni con pezzi di ricambio originali prima di rimontare i componenti.

Istruzioni per l'uso

Uso improprio

Agilent Technologies declina ogni responsabilità derivante dall'uso improprio della pompa MS120 Single Stage Rotary Vane. L'uso improprio comporta la perdita di tutti i diritti a reclami per responsabilità e garanzie. Il personale responsabile del funzionamento e della manutenzione della pompa deve essere adeguatamente addestrato e deve essere a conoscenza delle norme di prevenzione degli incidenti. Le precauzioni antinfortunistiche contenute in questa sezione devono essere costantemente rispettate durante il funzionamento e la manutenzione della pompa per evitare danni agli operatori e alla pompa. Tali precauzioni sono fornite sotto forma di note di AVVERTENZA e ATTENZIONE.

AVVERTENZA È proibito aspirare attraverso la pompa:



- liquidi o sostanze solide
- gas e vapori pericolosi, esplosivi o aggressivi
- ossigeno puro o miscele d'aria arricchita di ossigeno.

AVVERTENZA



È proibito utilizzare lo scarico della pompa per creare pressioni anche limitate.

AVVERTENZA



È proibito installare la pompa in un ambiente potenzialmente esplosivo.

Tabella limiti operativi

Tab. 1

Tensione d'ingresso (V)	Stato del Controller
< 180	Guasto all'alimentazione (Errore di Sottotensione/Sovratensione – Win206 bit 5 = 1)
180 – 200	In funzione (prestazioni ridotte)
200 – 264	In funzione (prestazioni complete)
> 264	Guasto all'alimentazione (Errore di Sottotensione/Sovratensione – Win206 bit 5 = 1)

AVVERTENZA



- Il contatto con tensioni elevate può provocare la morte. Prestare sempre la massima attenzione ed osservare le norme antinfortunistiche vigenti.
- Scollegare sempre l'alimentazione elettrica alla pompa prima di effettuare interventi di manutenzione. Posizionare sopra l'interruttore generale dell'alimentazione un apposito cartello di avvertenza: **MACCHINA IN MANUTENZIONE - NON ACCENDERE**.
- La mancata dotazione della pompa con il collegamento a terra può causare gravi danni agli operatori. Assicurarsi sempre che ci sia il collegamento a terra e che sia conforme alle norme.
- Durante la pulizia della pompa e dei suoi componenti, evitare l'uso di solventi infiammabili o tossici, come benzina, benzolo, etere o alcool. Si consiglia di utilizzare una soluzione di acqua e sapone, preferibilmente in lavatrici ad ultrasuoni, avendo cura di asciugare tutte le parti pulite a temperature inferiori a 100 °C in modo da eliminare l'umidità residua.
- Sovraccarichi o guasti prolungati possono causare il surriscaldamento del motore elettrico e la fuoriuscita di fumi nocivi; togliere immediatamente l'alimentazione per precauzione e non avvicinarsi alla pompa almeno finché non si è provveduto alla ventilazione per eliminare il fumo.
- In caso di incendio non gettare acqua sulla pompa. Togliere l'alimentazione e utilizzare estintori a CO₂.
- Ispezionare attentamente le flange per assicurarsi che non siano presenti polvere, olio, sporco o difetti delle superfici di accoppiamento, prima di effettuare i collegamenti richiesti.
- Assicurarsi che tutti i giunti e gli innesti siano bloccati correttamente prima di riavviare la pompa dopo i lavori di riparazione.
- Non indossare oggetti che possano impigliarsi nei meccanismi e/o fungere da conduttori (catene, braccialetti, ecc.).
- Assicurarsi che gli attrezzi da utilizzare siano in perfetto stato di funzionamento e dotati di impugnature isolanti, ove necessario. Controllare che il materiale isolante dei cavi e che i conduttori dell'apparecchiatura di prova non presentino segni di danneggiamento.
- Non sostituire l'olio immediatamente dopo aver spento la macchina poiché l'olio potrebbe essere ancora ad alta temperatura.

NOTA

Utilizzare la molla di ritenzione per fissare il cavo di alimentazione nella presa IEC320.

Attrezzatura di protezione

L'attrezzatura di protezione degli operatori che stanno operando o eseguendo la manutenzione della pompa deve essere sempre adeguata al tipo di operazione da eseguire. Inoltre, deve soddisfare i requisiti di sicurezza della legislazione in vigore nel paese in cui viene utilizzata la macchina.

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



- A seconda della peculiarità del processo, le pompe 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 delle pompe 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.

CAUTION



Rischio di lesioni a causa di spigoli vivi

- Prima di riparare la pompa o prima di qualsiasi azione di montaggio/smontaggio della pompa del sistema, attendere l'arresto completo della pompa.
- Se necessario, indossare guanti protettivi secondo la EN 420.

Linee guida per la sicurezza delle pompe MS120 Single Stage Rotary Vane

AVVERTENZA



Nonostante le precauzioni prese in fase di progetto, esistono elementi di rischio che si presentano durante le operazioni che si eseguono in fase di uso e manutenzione.

AVVERTENZA



SUPERFICI CALDE

Durante le operazioni di manutenzione si toccano superfici che possono superare la temperatura di 80 °C. Adottare idonei mezzi di protezione in modo da evitare scottature da contatto fortuito. È buona norma, prima di effettuare qualsiasi intervento sulla pompa, attendere il suo raffreddamento.

AVVERTENZA



EMISSIONI DI SOSTANZE NOCIVE

L'aria di scarico della pompa contiene tracce di nebbie d'olio (residuo max. 2PPM/ peso equivalenti a 2,4 mg/m³). Verificare la compatibilità con l'ambiente di lavoro. Garantire un corretto ricambio d'aria oppure portare lo scarico della pompa all'esterno. Un guasto o l'usura delle tenute possono provocare perdite d'olio lubrificante. Evitare la dispersione nel terreno e l'inquinamento di altri materiali. Nel caso di aspirazione d'aria contenente sostanze pericolose (esempio agenti biologici o microbiologici), adottare dei sistemi di abbattimento prima della pompa per vuoto.

ATTENZIONE



NON DISPERDERE NELL'AMBIENTE

Gli oli esausti provenienti dalla pompa devono essere smaltiti secondo le normative vigenti nel Paese d'utilizzo della pompa.

AVVERTENZA PERICOLO GENERATO DA DEPRESSIONE



Il contatto con punti in depressione può essere causa di infortuni. Evitare il contatto con l'attacco aspirazione della pompa durante il funzionamento. Immettere aria nel circuito di aspirazione prima di ogni intervento.

AVVERTENZA PERICOLO GENERATO DA PRESSIONE



Il serbatoio della pompa è pressurizzato. Non aprire e non dimenticare aperti i tappi di carico o scarico durante il funzionamento.

AVVERTENZA Escludere sempre l'alimentazione della pompa prima di compiere operazioni di manutenzione. Apporre specifici cartelli di avvertenza: **APPARECCHIATURA IN MANUTENZIONE - NON INSERIRE L'ALIMENTAZIONE**, in corrispondenza dell'interruttore di alimentazione. Al termine ripristinare i dispositivi di sicurezza.



AVVERTENZA



SICUREZZA ELETTRICA



Nell'equipaggiamento elettrico esistono parti sottoposte a tensione che, al contatto, possono provocare gravi danni a persone e cose.

AVVERTENZA



PERICOLO DI INCENDIO



L'utilizzo della pompa per impieghi non previsti o proibiti da questo manuale, oppure la mancanza di una corretta manutenzione, possono provocare anomalie di funzionamento con rischio di surriscaldamento e incendio. In caso di incendio non usare acqua per spegnere le fiamme. Utilizzare estintori a polvere o CO₂ od altri mezzi compatibili con la presenza di equipaggiamenti elettrici ed oli lubrificanti.

Descrizione del prodotto

Questa apparecchiatura è progettata per l'uso da parte di personale addestrato. L'utente deve leggere il presente manuale di istruzioni e qualsiasi altra informazione aggiuntiva fornita da Agilent prima di mettere in funzione l'apparecchiatura.

Agilent declina ogni responsabilità per eventuali eventi che si potrebbero verificare a causa della non conformità alle presenti istruzioni, dell'uso improprio da parte di persone non addestrate, di interferenze non autorizzate con l'apparecchiatura o di qualsiasi azione contraria a quanto previsto dagli specifici standard nazionali.

I consigli rivolti al personale addetto all'installazione e alla manutenzione presuppongono che lo stesso sia esperto e preparato nell'affrontare qualsiasi problematica di manutenzione, sia meccanica che elettrica. Per qualsiasi dubbio o informazioni non riportate su questo manuale si prega di contattare il nostro servizio assistenza, comunicando sempre: modello (Model), numero di serie (Serial), anno di costruzione, riportati sulla targhetta di identificazione.

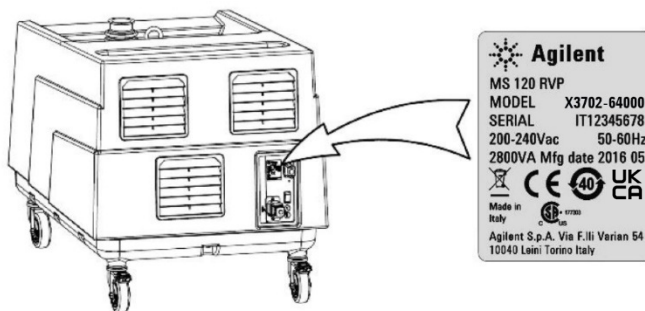


Figura 1 Targhetta di identificazione

Le Pompe MS120 Single Stage Rotary Vane sono delle pompe rotative monostadio a palette, a tenuta in bagno d'olio, azionate da motore elettrico pilotato da inverter. Queste pompe da alto vuoto sono adatte esclusivamente al pompaggio di aria e piccole quantità di vapore d'acqua. L'aspirazione di altri tipi di gas o di vapori deve essere preventivamente dichiarata ad Agilent che, se del caso, rilascerà la conformità all'impiego specifico.

Istruzioni per l'uso

Il motore elettrico flangiato è accoppiato a mezzo di giunto elastico. Il raffreddamento viene assicurato da una potente ventola centrifuga.

In aspirazione è presente un filtro a rete per proteggere la pompa da corpi solidi di diametro maggiore di 4,7 mm. Inoltre, una valvola di ritegno integrata impedisce la risalita dell'olio ed il rientro dell'aria nella camera da svuotare durante la fase d'arresto.

Nel serbatoio è inserito un sistema di separazione delle nebbie d'olio dall'aria di scarico (residuo max. 2PPM/ peso equivalenti a 2,4 mg/m³).

L'olio abbattuto viene recuperato in modo automatico dalla pompa.

Lo zavorratore impedisce la condensazione all'interno della pompa quando si aspirano piccole quantità di vapore (fare riferimento al paragrafo "Zavorratore (Gas-ballast)").

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

Trasporto e immagazzinamento

Le pompe devono essere trasportate senza olio.

Durante il trasporto e l'immagazzinamento della pompa non devono essere superate le seguenti condizioni ambientali:

- Temperatura: -20 °C ÷ 70 °C
- Umidità relativa: 0 ÷ 95% (non condensante)

Sollevamento

L'orientamento dei componenti imballati deve essere mantenuto conforme alle indicazioni fornite dai pittogrammi presenti sull'involucro esterno d'imballaggio.

Eeguire l'operazione di scarico con mezzo di sollevamento adeguato al peso della pompa. Per il sollevamento dell'imballo e della pompa fare riferimento alle figure seguenti.

Il peso dell'imballo, comprensivo della pompa, è, al massimo, di circa 120 kg, mentre il peso della sola pompa è di 100 kg.

Tab. 2 Dimensioni dell'imballo

Quota	L	B	H
Dimensione	920 mm	610 mm	790 mm

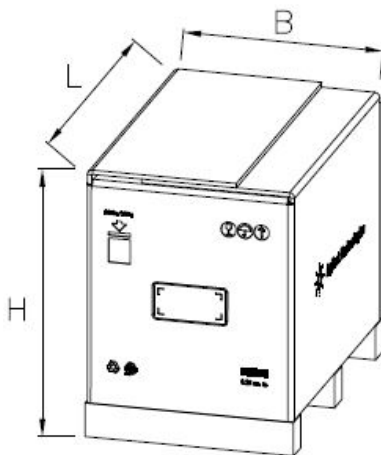


Figura 2 Pompa imballata

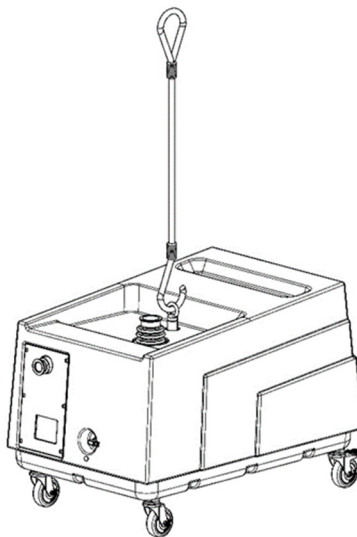


Figura 3 Pompa

Installazione

Preparazione per l'installazione

La pompa viene 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.

AVVERTENZA

Quando si disimballa la pompa, assicurarsi di non farla cadere ed evitare qualsiasi tipo di impatto improvviso o vibrazione d'urto.



Istruzioni per l'uso

Accessori forniti con la pompa

La pompa viene consegnata completa dei seguenti accessori necessari alla messa in servizio e alla manutenzione ordinaria:

- Bottiglia olio
- Beccuccio carico olio
- Tubo scarico olio
- Controconnettore

Assemblaggio

Togliere i tappi in aspirazione ed allo scarico.

La pompa viene fornita pronta all'uso, non vi sono altre parti da assemblare.

La pompa è priva di olio, vedere il paragrafo "Sostituzione olio lubrificante" per le istruzioni di carico olio.

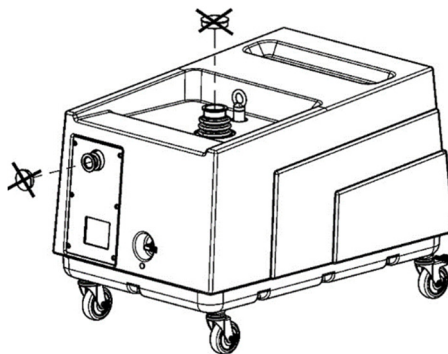


Figura 4

Istruzioni per l'uso

Ubicazione

AVVERTENZA La pompa



- deve essere bloccata agendo sui freni posti sulle ruote di appoggio, su un piano orizzontale.
- deve essere accessibile per una corretta e facile manutenzione rispettando le distanze minime da eventuali ingombri (vedere figura seguente). Dovrebbe essere inoltre accessibile con idonei mezzi di sollevamento.
- va protetta da getti o spruzzi d'acqua che potrebbero causare cortocircuiti elettrici e/o la folgorazione degli operatori vicini alla macchina.

Tab. 3 Minima Distanza

Quota	A	B	C	D
Dimensione	150 mm	200 mm	150 mm	500 mm

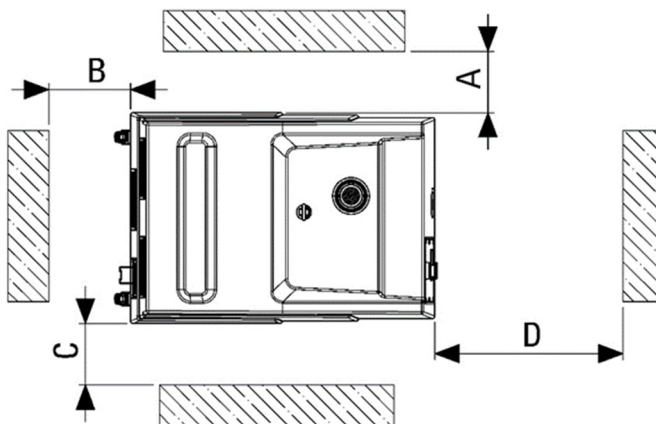


Figura 5

ATTENZIONE

La pompa non deve essere installata all'esterno e deve sempre essere protetta dagli agenti atmosferici.

Evitare che l'aria calda proveniente dallo scarico o dalle ventole di raffreddamento possa creare disagio al personale.

Non installare la pompa in una zona con polvere o altri materiali che potrebbero intasare o coprire rapidamente le superfici di raffreddamento.

Istruzioni per l'uso

Collegamento alla macchina utilizzatrice

Il collegamento della pompa alla camera da evacuare deve essere eseguito con tubazioni dello stesso diametro della bocca di aspirazione.

Il peso delle tubazioni e le eventuali dilatazioni non devono gravare sulla pompa. Effettuare il tratto finale di collegamento all'attacco aspirazione della pompa con tubo/ raccordo flessibile.

È importante che tutte le tubazioni ed i vari giunti siano a tenuta.

Tubazioni molto lunghe o di diametro piccolo diminuiscono le prestazioni della pompa.

NOTA

Questo simbolo identifica l'attacco in aspirazione.

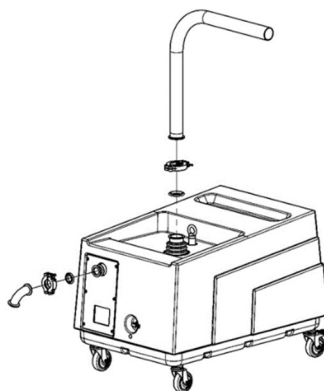


Figura 6

Istruzioni per l'uso

Convogliamento aria di scarico

Nel caso di un insufficiente ricambio d'aria nel locale pompa è possibile convogliare l'aria di scarico in altri ambienti o all'esterno.

Utilizzare tubazioni di diametro uguale alla bocca di scarico del serbatoio per una lunghezza massima di 15 m.

Per lunghezze superiori aumentare il diametro del tubo. Il peso delle tubazioni non deve gravare sulla pompa.

Utilizzare nel tratto finale raccordi o tubi flessibili.

AVVERTENZA

Questa tubazione deve essere discendente per evitare il rientro di condensa nel serbatoio della pompa.



ATTENZIONE

Non inserire rubinetti in questa tubazione. Non ostruire l'uscita.

NOTA

Questo simbolo identifica l'attacco allo scarico.



Collegamento elettrico

Verificare la tensione e la frequenza di rete con i dati riportati sulla targhetta della pompa.

Assicurarsi dell'efficienza dell'impianto di messa a terra.

AVVERTENZA

Assicurarsi che l'interruttore di accensione sia settato su off (0) onde evitare partenze indesiderate della pompa.



Eeguire l'allacciamento elettrico inserendo lo spinotto del cavo di alimentazione nella presa della pompa e bloccandolo con la molla in dotazione (vedi figura seguente).

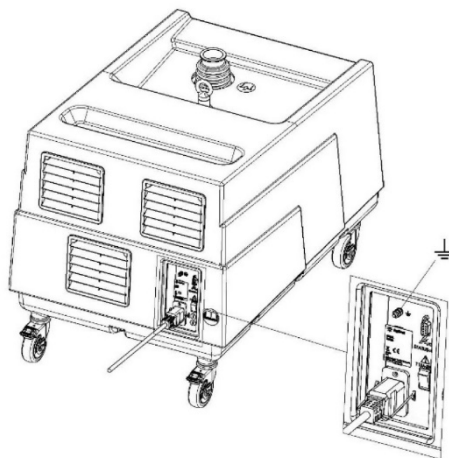


Figura 7

Istruzioni per l'uso

Disimballaggio

Durante l'operazione di disimballaggio, prestare particolare attenzione a non lasciar cadere la pompa e a non sottoporla ad urti o vibrazioni.

Non disperdere l'imballo nell'ambiente. Il materiale è completamente riciclabile e risponde alla direttiva CEE 85/399 per la tutela dell'ambiente.

NOTA

La pompa non può essere danneggiata rimanendo semplicemente esposta all'atmosfera. Si consiglia comunque di mantenerla chiusa fino al momento dell'installazione sul sistema onde evitare eventuale inquinamento da polvere.

Rimuovere le reggette che chiudono il cartone sul pallet, quindi rimuovere il cartone.

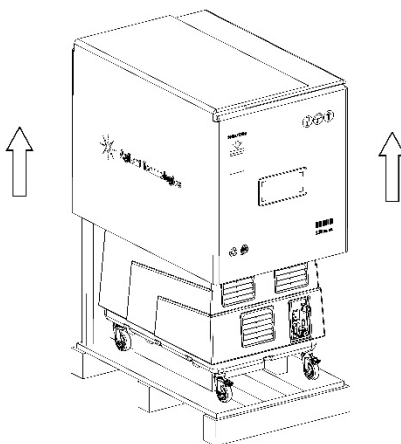


Figura 8

Istruzioni per l'uso

Rimuovere le n. 8 viti che chiudono le staffette, quindi rimuovere le n.4 staffette.

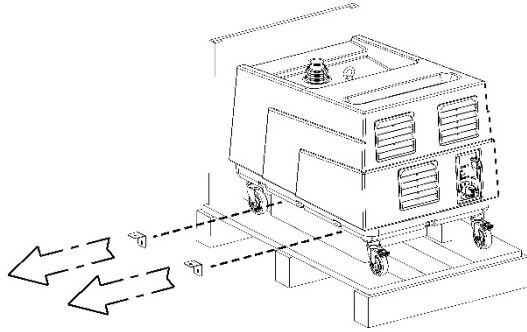


Figura 9

Sbloccare i freni di arresto delle ruote.

Fissare la rampa di discesa con n.1 vite come mostrato nella figura seguente.

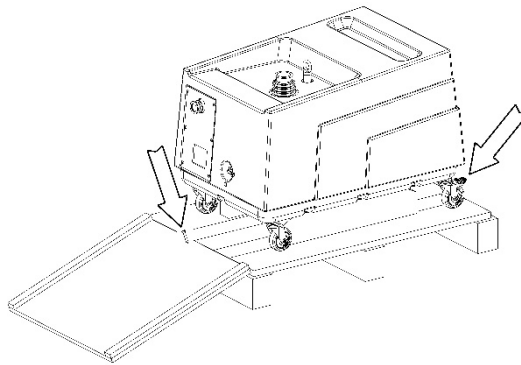


Figura 10

Istruzioni per l'uso

Spingere la pompa posteriormente per farla scivolare giù dal pallet.

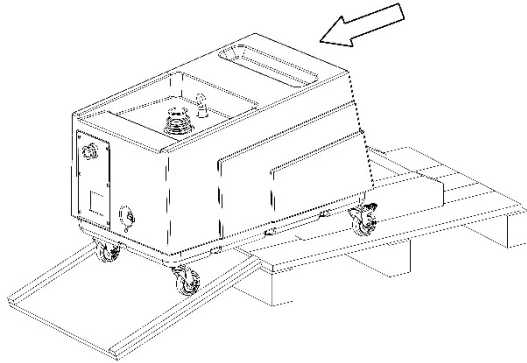


Figura 11

Utilizzo

Le pompe per vuoto descritte in questo manuale possono aspirare esclusivamente aria e piccole quantità di vapore d'acqua. L'aspirazione di altri tipi di gas o di vapori deve essere preventivamente dichiarata alla Agilent Technologies che rilascerà la conformità all'impiego specifico.

Sono adatte per l'evacuazione di sistemi chiusi o per funzionare ad un vuoto costante compreso nel seguente campo: 0,07 - 15 mbar (assoluti).

La temperatura ambiente e la temperatura di aspirazione devono essere comprese fra 12 e 35°C.

Avvio e funzionamento della pompa MS120 Single Stage Rotary Vane

Non installare e/o utilizzare la pompa in ambienti esposti ad agenti atmosferici (pioggia, gelo, neve), polveri, gas aggressivi, in ambienti esplosivi o con elevato rischio di incendio. Durante il funzionamento è necessario che siano rispettate le seguenti condizioni ambientali:

- temperatura: da +12 °C a +35 °C
- umidità relativa: 0 – 95 % (non condensante).

ATTENZIONE

Prima di avviare la pompa, occorre procedere al rifornimento di olio lubrificante, poichè la pompa viene fornita scarica.

Per maggiori informazioni relative al rifornimento di olio e per la scelta del lubrificante idoneo, fare riferimento al paragrafo "Sostituzione olio lubrificante".

NOTA

È importante che il livello dell'olio rimanga tra i valori MIN e MAX visualizzati dall'indicatore di livello posto sul lato della pompa.

Eseguire il primo riempimento attraverso il tappo (E) sino alla metà dell'indicatore di livello (F) e richiudere il tappo (E) (vedere figura seguente).

Istruzioni per l'uso

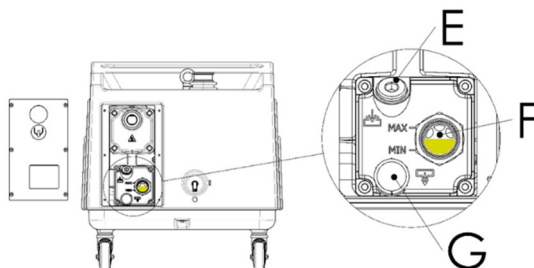


Figura 12

Accendere la pompa e portarla a vuoto massimo per almeno 2 minuti. Fermare la pompa, ricontrrollare il livello d'olio ed eseguire un eventuale rabbocco di olio ripristinando il livello corretto.

Consigli per l'utilizzo

Con temperature ambiente inferiori a 18 °C è bene riscaldare per 15 minuti la pompa facendola funzionare a pressione limite (aspirazione chiusa, senza carico).

Durante questa fase la pompa potrebbe non raggiungere i limiti di pressione dichiarati.

AVVERTENZA



La pompa non è progettata per lavorare a pressione atmosferica.

È consigliabile non superare i 5 avviamenti/ora.

Zavorratore (Gas-ballast)

La pompa MS120 è dotata di Zavorratore (Gas Ballast) e di una valvola manuale che permette il suo inserimento (I) o l'esclusione(O).

La pressione limite della pompa varia in funzione di questa scelta. Per un corretto utilizzo vedi "Aspirazione vapore acqueo.

Istruzioni per l'uso

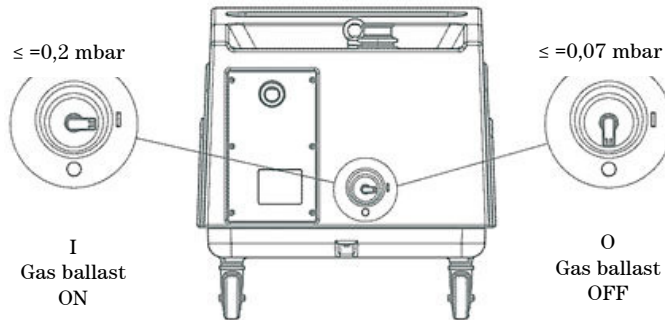


Figura 13

Aspirazione vapore acqueo

Nel caso di aspirazione di vapore acqueo è consigliato:

- Prima di aspirare vapori, portare la temperatura della pompa a regime, facendola funzionare per trenta minuti a vuoto massimo.
- Alla fine del ciclo di lavoro, nel caso di presenza di condensa nell'olio lubrificante, lasciare funzionare la pompa a vuoto massimo per almeno trenta minuti con lo zavorratore aperto.
- Effettuare questa operazione prima di fermi macchina prolungati; lo zavorratore consentirà di eliminare le condense dall'olio lubrificante.

Per la procedura d'inserimento o esclusione dello zavorratore fare riferimento all'appendice "Technical Information".

Manutenzione

Il personale responsabile del funzionamento e della manutenzione della pompa deve essere correttamente addestrato e consapevole delle norme di prevenzione infortuni.

Oltre alle prescrizioni sotto riportate, fare riferimento al paragrafo "Sicurezza".

AVVERTENZA



Prima di ogni intervento, immettere aria attraverso l'attacco in aspirazione.

AVVERTENZA



Nel caso si debba procedere ad operazioni di manutenzione della pompa al termine di un periodo di esercizio, è necessario lasciarla raffreddare, poichè la temperatura esterna può superare gli 80 °C.

AVVERTENZA



Escludere sempre l'alimentazione della pompa prima di compiere operazioni di manutenzione. Apporre specifici cartelli di avvertenza: APPARECCHIATURA IN MANUTENZIONE - NON INSERIRE L'ALIMENTAZIONE, in corrispondenza dell'interruttore di alimentazione. Al termine ripristinare i dispositivi di sicurezza.

La tabella mostra tutti gli interventi periodici necessari per mantenere in perfetta efficienza la pompa.

Manutenzioni più frequenti possono rendersi necessarie in base al tipo di utilizzo (aspirazione di vapori condensabili o sostanze inquinanti).

In questi casi solo l'esperienza diretta può suggerire i corretti intervalli di manutenzione. L'olio esausto e i pezzi di ricambio sostituiti devono essere considerati rifiuti speciali e gestiti secondo la normativa vigente nel paese d'utilizzo.

Istruzioni per l'uso

Tab. 4 Interventi di manutenzione

Intervallo di manutenzione	Descrizione intervento	Personale abilitato
24 ore / ogni giorno	Controllo livello olio prima dell'avviamento.	Operatore
100 Ore / ogni settimana	Pulire con panno morbido le superfici della pompa. Eventualmente è possibile utilizzare un detergente neutro.	Operatore
9000 Ore / ogni anno	Sostituire olio lubrificante	Tecnico qualificato
	Sostituire l'elemento disoleatore	Tecnico qualificato
14000 Ore / ogni 2 anni	Verificare i collegamenti elettrici.	Tecnico qualificato
30000 Ore / ogni 5 anni	Ricondizionamento pompa	Servizio Assistenza

Sostituzione olio lubrificante

Si raccomanda di attendere 1 ora e 30 minuti dopo lo spegnimento della pompa per effettuare l'operazione di sostituzione dell'olio lubrificante, per permettere alla temperatura della pompa e dell'olio di raffreddarsi sufficientemente.

AVVERTENZA



Utilizzare guanti protettivi per evitare scottature.

Con riferimento alla Figura 14, svitare il tappo di carico (E) e quello di scarico olio (G) solo dopo aver posizionato sotto il serbatoio un contenitore adatto (per forma e dimensioni) a raccogliere tutto l'olio della pompa (vedere Figura 15).

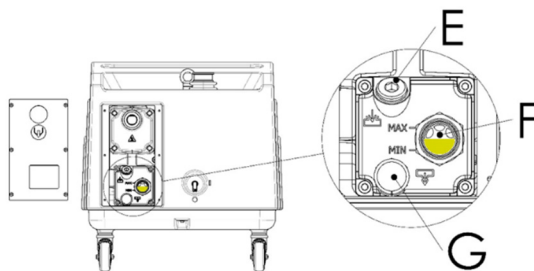


Figura 14

Istruzioni per l'uso

Avvitando il tubo scarico olio (pos. A1) sullo scarico (G) il lubrificante comincerà ad uscire. Si raccomanda di inclinare la pompa in avanti sollevando la parte posteriore della pompa di almeno 40 mm per garantire la fuoriuscita completa del lubrificante, ricordarsi di frenare le ruote per evitare che la pompa possa cadere. Quando l'olio esausto contenuto nel serbatoio sarà fuoriuscito rimontare entrambi i tappi ("E" e "G") e far girare la pompa a vuoto massimo per circa 30 secondi in modo da svuotare anche il circuito di lubrificazione dai residui d'olio esausto, quindi togliere i tappi e scaricare la rimanenza d'olio.

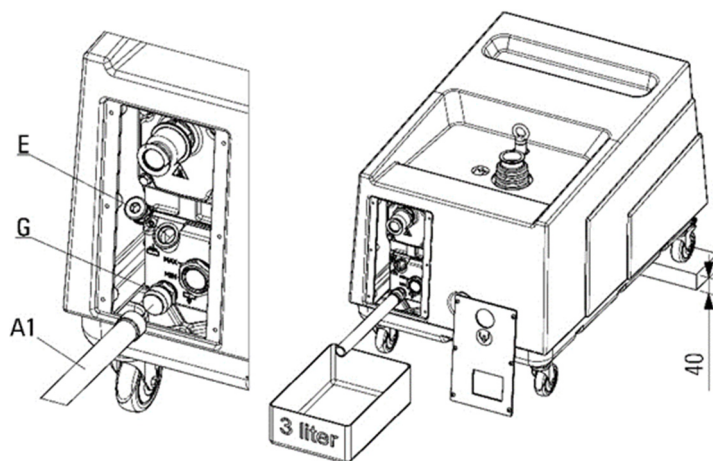


Figura 15

Se nell'olio sono presenti grosse quantità di sostanze inquinanti o si riscontra la presenza di acqua, procedere ad un lavaggio della pompa facendola funzionare a vuoto massimo per almeno 5 minuti.

Utilizzare per questa operazione una quantità di olio tale da raggiungere il livello minimo riportato sul serbatoio.

Scaricare nuovamente l'olio.

Istruzioni per l'uso

Eseguire il riempimento come segue (con riferimento alla Figura 16):

- 1 Rimuovere il tappo di carico olio (E).
- 2 Avvitare il beccuccio di carico olio (A2) sulla bottiglia di olio (A1).
- 3 Procedere al riempimento inserendo il beccuccio all'interno del foro di carico olio (Vedi "Avvio e funzionamento della pompa MS120 Single Stage Rotary Vane" e "Lubrificanti").

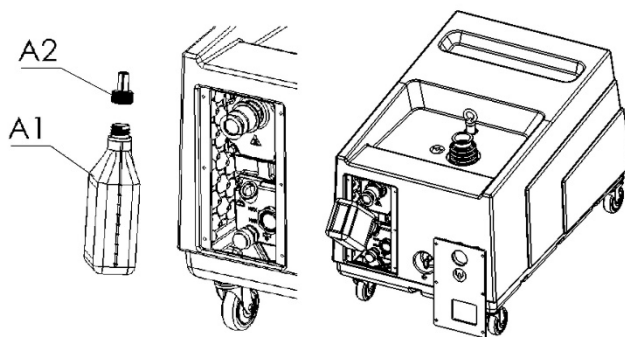


Figura 16

Sostituzione elementi disoleatori

Elementi disoleatori molto sporchi possono causare un sensibile aumento di temperatura della pompa e in casi estremi autocombustione dell'olio lubrificante.

La massima pressione ammessa nel serbatoio è di 0,6 bar misurata a portata massima (quando la pompa sta funzionando con l'aspirazione a pressione atmosferica). La pressione può essere rilevata mediante un manometro (opzionale), connesso nel foro di carico olio (Figura 16). Se è presente il manometro, verificare l'intasamento della cartuccia a pompa calda.

Istruzioni per l'uso

Per la sostituzione dell'elemento disoleatore, in riferimento a Figura 17, procedere come segue:

- 1** rimuovere il coperchio del cover (pos. C1) svitando le relative viti (C2),
- 2** togliere il coperchio serbatoio (pos. B1) svitando le relative viti (pos. B2),
- 3** estrarre e sostituire l'elemento disoleatore (pos. B4) e i relativi O.R. (pos. B5),
- 4** rimontare il coperchio di scarico (pos. B1),
- 5** se necessario sostituire la guarnizione (pos. B3),
- 6** rimontare il coperchio del cover (pos. C1).

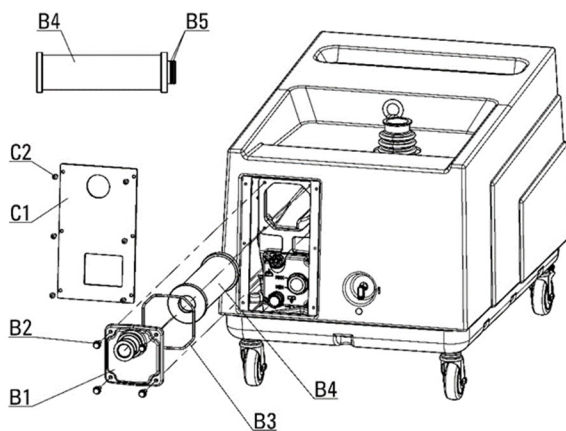


Figura 17

Istruzioni per l'uso

Manutenzione ordinaria

I ricambi necessari alla manutenzione ordinaria sono riportati nella tabella seguente. Usare solo ricambi originali. Per l'ordinazione riportare il Part Number indicato.

Tab. 5 Ricambi per la manutenzione

Part Number	Descrizione
X3760-64005	Rotary Vane Fluid AVF 60 Gold 1lt
X3702-68201	MS 120 Exhaust Filter Cartridge

Ricondizionamento pompa

Per questa operazione si consiglia di rivolgersi al Servizio Assistenza di Agilent.

Il ricondizionamento consiste nello smontaggio completo, la pulizia di tutti i particolari e la sostituzione delle parti soggette ad usura (bronzine della pompa, palette e guarnizioni). Usare solo ricambi originali. Per l'ordinazione riportare il Part Number indicato.

Tab. 6 Ricambi per ricondizionamento della pompa

Part Number	Descrizione
X3702-68202	MS 120 Major Maintenance Kit

Come ordinare i ricambi

Per ordinare i ricambi indicare sempre il modello della pompa (model), numero di matricola (serial), anno di costruzione, Part Number, descrizione e quantità richiesta.

Lubrificanti

Nella tabella seguente sono riportati gli olii consigliati per un uso generico della pompa.

Tab. 7 Lubrificanti

Temperatura ambiente	Viscosità	Olio Agilent Technologies
12 – 35 °C	ISO 53	AVF 60 Gold

Nel caso di cambio della tipologia di lubrificante, è necessario procedere ad un lavaggio della pompa facendola funzionare a vuoto massimo per almeno 5 minuti.

Utilizzare per questa operazione una quantità di olio tale da raggiungere il livello minimo riportato sul serbatoio.

Scaricare nuovamente l'olio e procedere con nuova carica olio.

Messa fuori servizio

Per la messa fuori servizio togliere l'olio dalla pompa prima della sua movimentazione. Se l'olio appare inquinato eseguire un lavaggio con olio nuovo (vedi "Sostituzione olio lubrificante").

Svuotare il serbatoio dell'olio, tappare l'aspirazione e lo scarico della pompa e immagazzinare.

NOTA

Prima di restituire la pompa al produttore per la riparazione, compilare la scheda "Request for Return" allegata a questo manuale di istruzioni e inviarla all'ufficio vendite di zona. Ricordarsi anche di inserire una copia della scheda nella scatola con cui viene restituita la pompa.

Smaltire la pompa in conformità con le normative nazionali applicabili.

Troubleshooting

NOTA

In caso di calo delle prestazioni, rumori anomali o intervento della protezione del motore, contattare il Contact Center Agilent.

Tab. 8 Troubleshooting

Inconveniente	Causa	Soluzione
Riduzione delle prestazioni	Perdite nelle tubazioni in aspirazione o sulla macchina utilizzatrice	Eliminare le perdite
	Mancanza di lubrificazione	Controllare livello e condizioni dell'olio. Ripristinare il livello o eseguire la sostituzione dell'olio
	Tensione di alimentazione troppo bassa (< 200 Vac)	Collegare la pompa ad una rete elettrica con la giusta tensione e utilizzare un cavo con la dimensione appropriata (vedi paragrafo Tabella limiti operativi)
Nebbie d'olio allo scarico	Elemento disoleatore inefficiente	Sostituire elemento disoleatore
	Elevata temperatura dovuta all'olio contaminato	Sostituire olio
	Elevata temperatura di esercizio dovuta a temperatura ambiente troppo elevata	Diminuire temperatura ambiente assicurando un migliore ricambio d'aria
	Elemento disoleatore intasato	Sostituire elemento disoleatore
La pompa non parte	Tensione di alimentazione fuori dal range previsto	Collegare la pompa a una rete con tensione idonea (vedere paragrafo Tabella limiti operativi)

Smaltimento

Significato del logo "WEEE" presente sulle etichette

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

Questo simbolo (**valido solo per i paesi della Comunità Europea**) indica che il prodotto sul quale è applicato, NON deve essere smaltito insieme ai comuni rifiuti domestici o industriali, ma deve essere avviato ad un sistema di raccolta differenziata. Si invita pertanto l'utente finale a contattare il fornitore del dispositivo, sia esso la casa madre o un rivenditore, per avviare il processo di raccolta e smaltimento, dopo opportuna verifica dei termini e condizioni contrattuali di vendita.



Figura 18 Logo "WEEE"

Per maggiori informazioni consultare:

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

Servizio Post Vendita

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

vpt-customer@agilent.com

vpl-customer@agilent.com

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

2

Betriebsanleitung

Informationen über diese Betriebsanleitung	51
Gültigkeit	51
Definitionen und Terminologie	52
Definition von Vorsicht, Warnung und Hinweis	52
Warning Symbols	53
Sicherheit	55
Sachgemäße Verwendung	55
Unsachgemäße Verwendung	57
Tabelle mit Betriebs-Grenzwerten	57
Schutzausrüstung	59
Sicherheitsrichtlinien für die Single Stage Rotary Vane MS120 Pumpe	60
Produktbeschreibung	63
Transport und Aufbewahrung	64
Anheben	65
Einbau	66
Montagevorbereitungen	66
Mitgeliefertes Zubehör	67
Zusammenbau	67
Lage	68
Anschluss an die Maschine	69
Umleitung der Abgasluf	70

Betriebsanleitung

Elektrische Anschlüsse	71
Auspacken	72
Betrieb	75
Inbetriebnahme und Betrieb der Pumpe MS120 Single Stage Rotary Vane	75
Tipps für den Betrieb	76
Gas-ballast	77
Ansaugung von Wasserdampf	77
Wartung	78
Ersetzung des Schmieröls	79
Ersetzung der Ölreinigungselemente	81
Routinewartung	83
Revision der Pumpe	83
So bestellen Sie Ersatzteile	83
Schmiermittel	84
Stilllegung	84
Fehlerbehebung	85
Entsorgung	86
Service	87

Informationen über diese Betriebsanleitung

Gültigkeit

Diese Betriebsanleitung enthält die Anweisungen für die Benutzer der Pumpe MS120 Single Stage Rotary Vane mit besonderer Bezugnahme auf die Begriffe für Sicherheit, Betrieb und Wartung auf der ersten Ebene, auf die Wartungsarbeiten beschränkt, für die der Benutzer verantwortlich ist.

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

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

HINWEIS

- 1** Diese Betriebsanleitung enthält nützliche Informationen, damit das gesamte Personal, das die Pumpe MS120 Single Stage Rotary Vane verwendet, diese in Sicherheit verwenden und die perfekte Funktionstüchtigkeit für ihre gesamte Betriebsdauer garantieren kann.
 - 2** Diese Betriebsanleitung muss zusammen mit allen mit dieser zusammenhängenden Veröffentlichungen an einem zugänglichen Ort, der allen Bedienern/Wartungspersonal bekannt ist, aufbewahrt werden.
-

Definitionen und Terminologie

Definition von Vorsicht, Warnung und Hinweis

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

VORSICHT

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

WARNUNG



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

HINWEIS

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

Warning Symbols

Das Folgende ist eine Liste von Symbolen, die in Verbindung mit den Warnungen auf der Pumpe MS120 Single Stage Rotary Vane angezeigt werden. Die von ihnen beschriebene Gefahr wird ebenfalls angezeigt. Ein dreieckiges Symbol weist auf eine Warnung hin. Die Bedeutung der Symbole, die neben Warnhinweisen in der Dokumentation oder auf dem Gerät selbst erscheinen können, ist wie folgt. 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.



	Allgemeine Gefahr
	Vorhandensein gefährlicher Spannungen
	Heiße Oberflächen
	Wechselstrom (AC)
	Schutzerdungsanschluss
	Brandgefahr
	Nicht in die Umwelt gelangen lassen
	Betriebshandbuch lesen
	Anschluss an die Ansaugung
	Anschluss an den Ausgang

Betriebsanleitung

	Schutzhandschuhe
	Sicherheitsschuhe
	Schutzausrüstung
	Emission von schädlichen Substanzen
	Verletzungsfahr durch scharfe Kante
	Herunterfallende Gegenstände
	CSA-Zertifizierung
RoHS	Beschränkung gefährlicher Stoffe
CE	Europäische Konformitätserklärung
	Elektro- und Elektronikaltgeräte
UK CA	UK CA-Kennzeichnung

Sicherheit

Dieser Abschnitt enthält die von der Niederspannungsrichtlinie 2014/35/EG vorgegebenen Informationen, die für die Übereinstimmung und Einhaltung der Sicherheitsvorschriften sowohl allgemein als auch in Bezug auf die spezifische Verwendung des Produkts wesentlich sind.

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

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

WARNUNG



Bei Verwendungen, bei denen ein Defekt oder ein Schaden an der Vakuumpumpe zu Personen- oder Sachschäden führen könnte, müssen an der Anlage besondere Sicherheitsvorkehrungen getroffen werden.

Sachgemäße Verwendung

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

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

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

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

WARNUNG



Entflammbare Werkstoffe:

Damit sichergestellt ist, dass die Gase in der Pumpe nicht in den entflammbaren/brennbaren Bereich gelangen, müssen die nachstehenden Anweisungen und Vorsichtsmaßnahmen befolgt werden. Wenn potenziell entflammbare oder brennbare Materialien in den Geräten vorhanden sind, müssen Sie Folgendes unternehmen:

- Sicherstellen, dass die Anlage dicht ist, um zu gewährleisten, dass nur die bestimmungsgemäßen Gase durch die Vakuumpumpe geleitet werden.
- Sicherstellen, dass die Konzentration aller brennbaren Gase oder Dämpfe, die in die Pumpe (Einlass und Gasballastanschluss) gelangen, in der Pumpe und in der Auslassleitung unter einem Viertel der für die Gase angegebenen unteren Explosionsgrenzwerte (LEL) bleibt.
- Sicherstellen, dass Sicherheitssysteme vorhanden sind, die dafür sorgen, dass die Konzentration aller brennbaren Gase oder Dämpfe, die in die Pumpe (Einlass und Gasballastanschluss) gelangen, in der Pumpe und in der Auslassleitung unter einem Viertel der für die Gase angegebenen unteren Explosionsgrenzwerte (UEG) unter jeglichen Anlagen- oder Vakuumpumpen-Störungsbedingungen beibehalten werden.

HINWEIS

- Bevor Sie die Pumpe nach einer Panne wieder in Betrieb nehmen, überprüfen Sie sie sorgfältig und prüfen Sie sie sorgfältig auf sonstige Anzeichen von Schäden.
- Benutzen Sie nur einwandfrei funktionierende und speziell für den Einsatz konzipierte Werkzeuge; Die Verwendung ungeeigneter oder unwirksamer Werkzeuge kann zu schweren Schäden führen.
- Überprüfen Sie immer den Schmierstoff und seine ordnungsgemäße Verteilung in der Pumpe; Eine unzureichende Schmierung kann die Pumpe ernsthaft beschädigen.
- Markieren Sie die Teile beim Zerlegen, um sicherzustellen, dass Sie sie in der richtigen Reihenfolge wieder zusammenbauen.
- Stellen Sie sicher, dass die bearbeiteten Wellen, ihre Sitze im Inneren der Pumpe und die maschinell geschliffenen Oberflächen keine Kratzer oder Riefen aufweisen. Leichte Kratzer und Abschürfungen können mit sehr feinem Schmirgelpapier oder durch leichtes Schleifen entfernt werden.
- Bevor Sie eine Gruppe zusammenstellen, verteilen Sie immer etwas Öl auf den Innenteilen und Passflächen. Ersetzen Sie alle Dichtungen durch Originalersatzteile, bevor Sie die Komponenten wieder zusammenbauen.

Unsachgemäße Verwendung

Agilent Technologies übernimmt keine Haftung für Schäden, die durch unsachgemäßen Gebrauch der Pumpe MS120 Single Stage Rotary Vane entstanden sind. Unsachgemäßer Gebrauch führt zum Verlust aller Haftungs- und Gewährleistungsansprüche.

Das für den Betrieb und die Wartung der Pumpe zuständige Personal muss gut ausgebildet sein und mit den Unfallverhütungsvorschriften vertraut sein. Die in diesem Abschnitt enthaltenen Unfallverhütungsvorschriften müssen bei Betrieb und Wartung der Pumpe konsequent beachtet werden, um Personen- und Pumpenschäden zu vermeiden. Diese Vorsichtsmaßnahmen werden in Form von Hinweisen wie WARNUNG und VORSICHT angezeigt.

WARNUNG



Die Ansaugung folgender Substanzen mit der Pumpe ist verboten:

- Flüssigkeiten oder Festkörper
- gefährliche, explosive oder aggressive Gase und Dämpfe
- reiner Sauerstoff oder mit Sauerstoff angereicherte Luftmischungen.

WARNUNG



Der Ausgang der Pumpe darf nicht zum Aufbau von Druck verwendet werden, auch nicht wenn es sich um geringen Druck handelt.

WARNUNG



Es ist verboten, die Pumpe in einem Raum zu installieren, in dem Explosionsgefahr herrscht.

Tabelle mit Betriebs-Grenzwerten

Tab. 9

Eingangsspannung (V)	Status des Controllers
< 180	Schaden an der Versorgung (Fehler wegen zu geringer oder zu hoher Spannung – Win206 bit 5 = 1)
180 – 200	in Betrieb (verringerte Leistung)
200 – 264	in Betrieb (volle Leistung)
> 264	Schaden an der Versorgung (Fehler wegen zu geringer oder zu hoher Spannung – Win206 bit 5 = 1)

WARNUNG



- Der Kontakt mit Hochspannung kann zum Tod führen. Seien Sie stets äußerst vorsichtig und beachten Sie die geltenden Unfallverhütungsvorschriften..
- Trennen Sie vor Wartungsarbeiten immer die Stromversorgung der Pumpe. Bringen Sie ein spezielles Warnschild über dem Leistungsschalter der Stromversorgung an: MASCHINE WIRD GEWARTET – NICHT EINSCHALTEN.
- Wenn die Pumpe nicht mit einem Erdungsanschluss ausgestattet ist, kann dies zu schweren Schäden für das Bedienpersonal führen. Stellen Sie immer sicher, dass ein Erdungsanschluss vorhanden ist und den Normen entspricht.
- Vermeiden Sie beim Reinigen der Pumpe und ihrer Komponenten die Verwendung von brennbaren oder giftigen Lösungsmitteln wie Benzin, Benzol, Ether oder Alkohol. Es wird empfohlen, eine Seifen-Wasser-Lösung zu verwenden, vorzugsweise in Ultraschallwaschmaschinen, und darauf zu achten, dass alle gereinigten Teile bei Temperaturen unter 100 °C getrocknet werden, um Restfeuchtigkeit zu entfernen.
- Längere Überlastungen oder Ausfälle können dazu führen, dass der Elektromotor überhitzt und gesundheitsschädliche Rauchgase freisetzen; Trennen Sie vorsichtshalber sofort den Strom und nähern Sie sich der Pumpe erst dann, wenn Sie für Belüftung gesorgt haben, um den Rauch auszutreiben.
- Im Brandfall kein Wasser auf die Pumpe werfen. Schalten Sie den Strom aus und verwenden Sie CO₂-Feuerlöscher.
- Überprüfen Sie die Flansche sorgfältig, um sicherzustellen, dass sich kein Staub, Öl, Schmutz oder Defekte an den Passflächen befinden, bevor Sie die erforderlichen Verbindungen herstellen.
- Stellen Sie sicher, dass alle Verbindungen und Kupplungen ordnungsgemäß verriegelt sind, bevor Sie die Pumpe nach Reparaturarbeiten wieder in Betrieb nehmen.
- Tragen Sie keine Gegenstände, die sich in den Mechanismen verfangen und/oder als Leiter dienen könnten (Ketten, Armbänder usw.).
- Stellen Sie sicher, dass die zu verwendenden Werkzeuge in einwandfreiem Zustand sind und ggf. über isolierende Griffe verfügen. Überprüfen Sie, ob das Isoliermaterial der Kabel und die Leiter des Prüfgeräts keine Beschädigungen aufweisen.
- Tauschen Sie das Öl nicht sofort nach dem Anhalten der Maschine aus, da das Öl möglicherweise noch eine hohe Temperatur hat.

HINWEIS

Verwenden Sie die Haltefeder, um das Netzkabel in der IEC320-Buchse zu sichern.

Schutzausrüstung

Die Schutzausrüstung der Bediener, die das Pumpensystem bedienen oder warten, muss immer für die Art des auszuführenden Vorgangs geeignet sein. Darüber hinaus muss es den Sicherheitsanforderungen der in dem Land, in dem die Maschine verwendet wird, geltenden Rechtsvorschriften entsprechen.

WARNUNG



Gesundheitsgefahr durch gefährliche Stoffe bei Wartung oder Installation.



- Vakuumpumpen, 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 Vakuumpumpen von Hand besteht die Gefahr, dass Lasten verrutschen und herunterfallen.



- Tragen Sie kleine und mittlere Vakuumpumpen 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 Kante



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

Sicherheitsrichtlinien für die Single Stage Rotary Vane MS120 Pumpe

WARNUNG



Trotz allen in der Planungsphase ergriffenen Vorsichtsmaßnahmen sind Betrieb und Wartung noch mit einigen Risikofaktoren behaftet.

WARNUNG



HEISSE OBERFLÄCHEN

Bei den Wartungsarbeiten müssen Oberflächen berührt werden, deren Temperatur über 80 °C betragen kann. Schützen Sie sich angemessen, um Verbrennungen durch versehentlichen Kontakt zu vermeiden. Es ist üblich, vor allen Eingriffen an der Pumpe abzuwarten, bis diese sich abgekühlt hat.

WARNUNG



EMISSIONEN SCHÄDLICHER SUBSTANZEN

Die aus der Pumpe austretende Luft enthält Spuren von Ölnebel (max. 2 PPM, entsprechend einem Gewicht von 2,4 mg/m³). Vergewissern Sie sich, dass das mit den Umständen der Arbeitsumgebung vereinbar ist. Sorgen Sie für einen geeigneten Luftaustausch oder leiten Sie die aus der Pumpe austretende Luft nach außen. Ein Defekt oder Verschleiß an den Dichtungen kann zum Austritt von Schmieröl führen. Dieses darf nicht in den Boden gelangen oder zu anderen Formen von Verschmutzung führen. Wenn Luft mit gefährlichen Substanzen angesaugt wird (beispielsweise biologische oder mikrobiologische Wirkstoffe), bauen Sie vor der Pumpe Systeme ein, die diese unschädlich machen.

VORSICHT



NICHT IN DIE UMWELT GELANGEN LASSEN

Das Altöl aus der Pumpe muss nach den in dem jeweiligen Land geltenden Vorschriften entsorgt werden.

WARNUNG



GEFAHR DURCH UNTERDRUCK

Der Kontakt mit Unterdruck kann Ursache von Unfällen sein. Vermeiden Sie den Kontakt mit der Ansaugöffnung, während die Pumpe läuft. Leiten Sie vor jedem Eingriff Luft in den Ansaugkreislauf.

WARNUNG



GEFAHR DURCH DRUCK

Der Tank der Pumpe steht unter Druck. Öffnen Sie die Deckel zum Laden und Entladen nicht während des Betriebs und lassen Sie sie nicht versehentlich unverschlossen.

WARNUNG



Klemmen Sie die Pumpe immer von der Versorgung ab, bevor Sie Wartungsarbeiten durchführen. Stellen Sie beim Stromversorgungsschalter Warnschilder auf: **GERÄT WIRD GEWARTET – STROM NICHT EINSCHALTEN!** Aktivieren Sie die Sicherheitsvorrichtungen nach dem Ende der Wartungsarbeiten wieder.

WARNUNG



ELEKTRISCHE SICHERHEIT

Einige Teile der Elektrik stehen unter Strom. Bei Kontakt kann es zu schweren Sach- und Personenschäden kommen.

WARNUNG



BRANDGEFAHR

Die Verwendung der Pumpe für andere als die in diesem Handbuch genannten oder für verbotene Zwecke sowie eine nicht ausreichende Wartung können zu Funktionsstörungen führen und bergen die Gefahr von Überhitzung und Brand. Verwenden Sie im Brandfall kein Wasser zum Löchen. Nehmen Sie Feuerlöscher mit Pulver oder CO₂ oder andere Löschmittel, die mit elektrischen Komponenten oder mit Schmieröl verträglich sind.

Produktbeschreibung

Dieses Gerät ist für den Gebrauch durch eingewiesenes Fachpersonal bestimmt. Der Benutzer muss diese Bedienungsanleitung und alle anderen von Agilent bereitgestellten zusätzlichen Informationen lesen, bevor er das Gerät in Betrieb nimmt. Agilent haftet nicht für Ereignisse, die auf die Nichteinhaltung dieser Anweisungen, die unsachgemäße Verwendung durch nicht unterwiesene Personen, nicht genehmigte Eingriffe in die Geräte oder auf Handlungen zurückzuführen sind, die im Widerspruch zu spezifischen landesweiten Normen stehen.

Die Ratschläge für das mit Installation und Wartung betraute Personal setzen voraus, dass dieses Personal kompetent und darauf vorbereitet ist, bei Problemen, wie sie bei der Wartung von Mechanik und Elektrik auftreten können, angemessen zu reagieren. Bei Unklarheiten oder wenn festgestellt wird, dass bestimmte Informationen in diesem Handbuch fehlen, wenden Sie sich bitte an unseren Kundendienst. Nennen Sie dabei immer Modell, Seriennummer und Baujahr. Sie finden diese Angaben auf dem Typenschild.

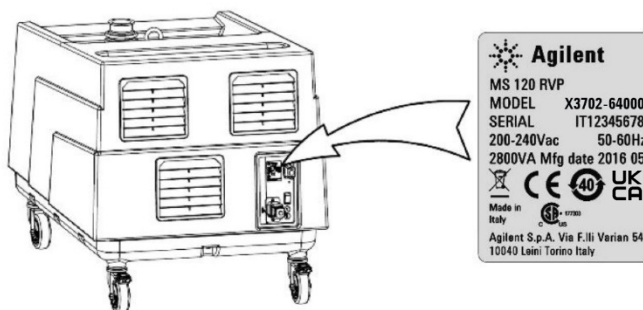


Abbildung 19 Typenschild

Die Pumpe Modell MS120 Single Stage Rotary Vane ist eine Einstadium-Rotationsflügelpumpe im Ölbad, die von einem Elektromotor inverter angetrieben. Die Pumpe Modell MS120 Single Stage Rotary Vane Pump ist eine Einstadium-Rotationsflügelpumpe im Ölbad, die von einem durch einen Inverter gesteuerten Elektromotor angetrieben wird. Die Ansaugung anderer Arten von Gas oder Dämpfen muss zuvor Agilent gemeldet werden. Agilent kann dann die entsprechende Verwendung gestatten.

Betriebsanleitung

Der geflanschte Elektromotor ist über eine elastische Verbindung angeschlossen. Die Kühlung erfolgt über ein starkes Lüfterrad.

An der Ansaugung befindet sich ein Netzfilter zum Schutz der Pumpe vor festen Fremdkörpern mit Durchmessern über 4,7 mm. Außerdem verhindert ein integriertes Rückhalteventil das Aufsteigen des Öls und das Eintreten der Luft in die beim Anhalten zu leerende Kammer.

Im Tank sitzt ein System zur Trennung der Ölnebel von der ausgestoßenen Luft (max. 2 PPM, entsprechend einem Gewicht von 2,4 mg/m³).

Das aufgefangene Öl wird von der Pumpe automatisch zurückgeführt.

Der Gas-Ballast verhindert die Kondensation in der Pumpe, wenn kleine Mengen von Wasserdampf angesaugt werden (siehe Abschnitt "Gas-ballast").

In den folgenden Abschnitten finden Sie alle Informationen, die nötig sind, um die Sicherheit während des Betriebs zu gewährleisten. Detaillierte Informationen finden Sie im Anhang "Technical Information".

Transport und Aufbewahrung

Die Pumpe muss ohne Öl transportiert werden.

Bei Transport und Einlagerung der Pumpe müssen folgende Bedingungen eingehalten werden:

- Temperatur: -20°C bis 70 °C
- relative Luftfeuchtigkeit: 0 bis 95 % (nicht kondensierend)

Anheben

Die Ausrichtung der verpackten Komponenten muss den Abbildungen auf der Außenseite der Verpackung entsprechen.

Laden Sie die Pumpe mit einer zu deren Gewicht passenden Hebevorrichtung ab. Zum Anheben der Verpackung und der Pumpe richten Sie sich nach den folgenden Abbildungen.

Das Gewicht der Verpackung mit Pumpe beträgt maximal etwa 120 kg, das Gewicht der Pumpe allein 100 kg.

Tab. 10 Abmessungen der Verpackung

Wert	L	B	H
Abmessung	920 mm	610 mm	790 mm

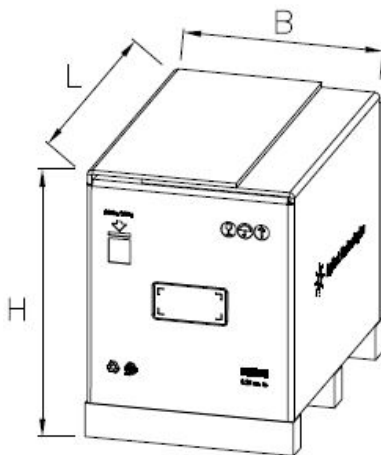


Abbildung 20 Verpackte Pumpe

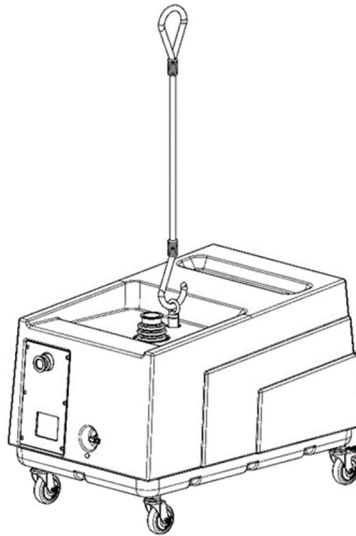


Abbildung 21 Pumpe

Einbau

Montagevorbereitungen

Die Pumpe wird in einer Spezialschutzverpackung angeliefert. Wenn diese Spuren von Beschädigungen aufweisen sollte, die während des Transports entstanden sein könnten, wenden Sie sich an Ihr örtliches Vertriebsbüro.

WARNUNG



Achten Sie beim Auspacken darauf, die Pumpe nicht fallen zu lassen und vermeiden Sie plötzliche Stöße oder Erschütterungen der Pumpe.

Betriebsanleitung

Mitgeliefertes Zubehör

Die Pumpe wird mit folgendem Zubehör geliefert, das für die Inbetriebnahme und die planmäßige Wartung gebraucht wird:

- Ölflasche
- Ölfüllschnabel
- Schlauch zum Ableiten des Öls
- I/O Mating Connector

Zusammenbau

Nehmen Sie die Deckel von der Ansaugung und vom Ausgang ab.

Die Pumpe wird betriebsbereit geliefert. Weitere Komponenten müssen nicht angebaut werden.

Die Pumpe wird ohne Öl geliefert. Sehen Sie im Absatz "Ersetzung des Schmieröls" nach, dort finden Sie die Instruktionen zum Einfüllen des Öls.

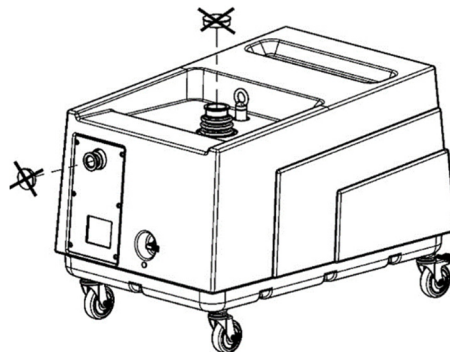


Abbildung 22

Lage

WARNUNG

Die Pumpe:



- muss blockiert werden. Stellen Sie dazu die Bremsen an den Rädern fest, wenn die Pumpe auf einer horizontalen Ebene steht.
- muss für eine korrekte und problemlose Wartung zugänglich sein, dazu müssen die korrekten Abstände zu umgebenden Hindernissen eingehalten werden (siehe folgende Abbildung). Sie muss außerdem an passenden Hebevorrichtungen befestigt werden können.
- muss vor Wasserstrahlen und Spritzern geschützt werden, die zu Kurzschlüssen und Stromschlägen führen können, durch die in der Nähe befindliches Personal verletzt werden kann.

Tab. 11 Mindestabstand

Wert	A	B	C	D
Abmessung	150 mm	200 mm	150 mm	500 mm

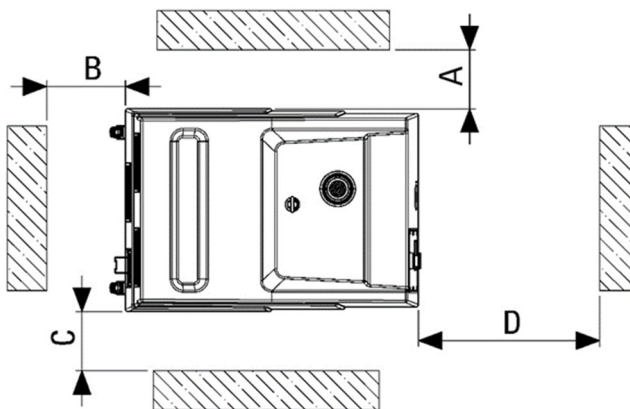


Abbildung 23

VORSICHT

Die Pumpe darf nicht außen installiert werden. Sie muss immer vor atmosphärischen Einflüssen geschützt werden.

Die heiße Luft aus dem Ausgang oder aus den Lüfterrädern kann zu Belästigungen des Personals führen.

Installieren Sie die Pumpe nicht in Bereichen, in denen sich Staub oder andere Materialien in der Luft befinden. Diese könnten die Pumpe verstopfen und die Kühlungsflächen schnell zusetzen.

Betriebsanleitung

Anschluss an die Maschine

Der Anschluss der Pumpe an den leerpumpenden Raum muss mit Leitungen erfolgen, die den gleichen Durchmesser wie die Ansaugöffnung haben.

Das Gewicht der Leitungen und der zugehörigen Erweiterung darf nicht an der Pumpe hängen. Verbinden Sie den letzten Abschnitt und die Ansaugöffnung der Pumpe mit einem Schlauch oder einem anderen flexiblen Verbindungsstück.

Alle Leitungen und Verbindungen müssen unbedingt dicht sein.

Lange Leitungen oder solche mit geringem Querschnitt verringern die Leistungsfähigkeit der Pumpe.

HINWEIS

Dieses Symbol steht für die Ansaugöffnung.

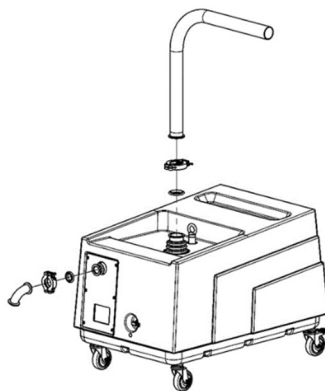


Abbildung 24

Betriebsanleitung

Umleitung der Abgasluf

Wenn der Luftaustausch im Pumpenraum unzureichend ist, kann die austretende Luft in andere Räume oder nach außen geleitet werden.

Verwenden Sie Leitungen, die den gleichen Durchmesser haben wie der Ausgang des Tanks bei maximaler Länge von 15 m.

Wenn die Leitungen länger sind, vergrößern Sie den Durchmesser. Das Gewicht der Leitungen darf die Pumpe nicht beschweren.

Verwenden Sie im letzten Abschnitt Schläuche oder flexible Leitungen.

WARNUNG

Diese Leitung muss ein Gefälle haben, damit kein Kondenswasser in den Tank der Pumpe gelangt.



VORSICHT

Stecken Sie diese Leitung nicht auf Hähne. Blockieren Sie den Ausgang nicht.

HINWEIS

Dieses Symbol kennzeichnet den Anschluss an den Ausgang.



Elektrische Anschlüsse

Vergewissern Sie sich, dass Spannung und Frequenz des Netzanschlusses mit den Daten auf dem Typenschild der Pumpe übereinstimmen.

Vergewissern Sie sich, dass die Erdung korrekt vorgenommen wurde.

WARNUNG

Vergewissern Sie sich, dass der Hauptschalter **wird ausgeschaltet (0) steht, um eine unerwünschte Einschaltung auszuschließen der Pumpe.**



Schließen Sie die Elektrik an, indem Sie den Stecker des Versorgungskabels an die Pumpe anschließen und mit der zugehörigen Feder blockieren (siehe Abbildung unten).

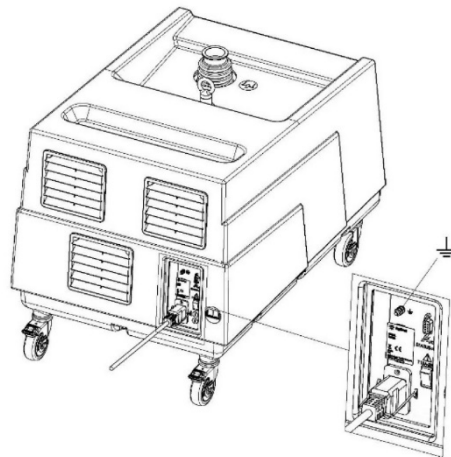


Abbildung 25

Betriebsanleitung

Auspacken

Achten Sie beim Auspacken vor allem darauf, dass die Pumpe nicht herunterfällt und keinen Stößen oder Vibrationen ausgesetzt wird.

Lassen Sie die Verpackung nicht in die Umwelt gelangen. Das Material ist vollständig recyclingfähig und entspricht den Vorgaben der Umweltschutzrichtlinie 85/399 EWG.

HINWEIS

Durch atmosphärische Einflüsse kann die Pumpe nicht beschädigt werden. Es empfiehlt sich dennoch, sie in der Verpackung zu lassen, bis sie im System installiert wird, um eine Verschmutzung durch Staub zu verhindern.

Entfernen Sie die Riemen, die den Karton auf der Palette halten. Entfernen Sie dann den Karton.

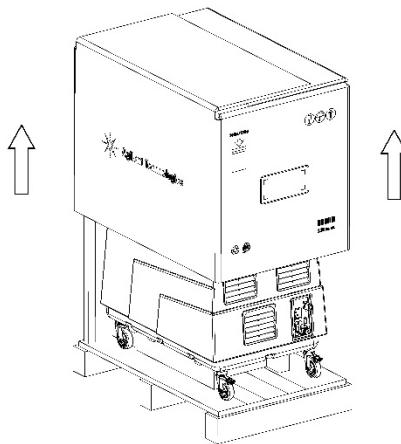


Abbildung 26

Betriebsanleitung

Entfernung der 8 Schrauben, die die Bügel halten, dann Entfernung der 4 Bügel.

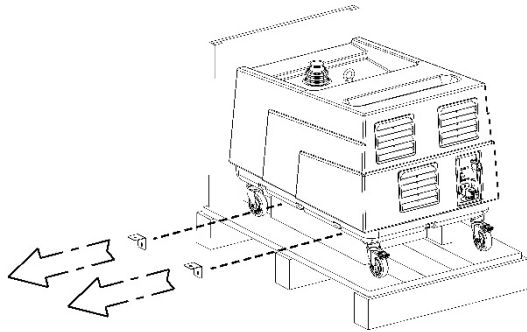


Abbildung 27

Entblocken der Radbremsen.

Befestigung der Abfahrtrampe mit einer Schraube, wie in der Abbildung gezeigt.

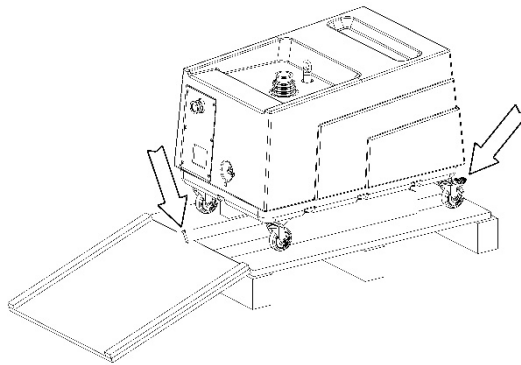


Abbildung 28

Betriebsanleitung

Pumpe rückwärts schieben, damit sie von der Palette rollt.

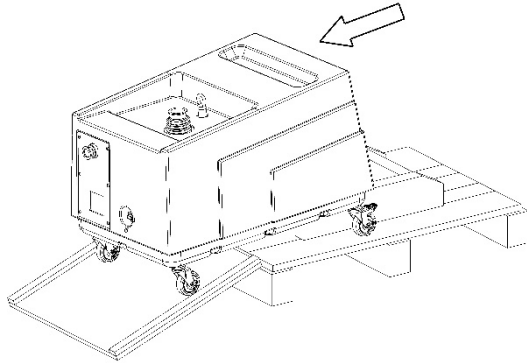


Abbildung 29

Betrieb

Die in diesem Handbuch beschriebenen Vakuumpumpen können nur Luft und kleine Mengen Wasserdampf ansaugen. Die Ansaugung anderer Arten von Gas oder Dämpfen muss zuvor der Firma Agilent Technologies gemeldet werden, die dann die entsprechende Verwendung genehmigt.

Diese Pumpen eignen sich für das Leerpumpen geschlossener Systeme oder für ein konstantes Vakuum im Bereich zwischen 0,07 und 15 mbar (absolut).

Die Umgebungstemperatur und die Temperatur der Ansaugluft müssen dabei zwischen 12 und 35°C liegen.

Inbetriebnahme und Betrieb der Pumpe MS120 Single Stage Rotary Vane

Installieren und verwenden Sie die Pumpe nicht in Räumen, die von atmosphärischen Einflüssen betroffen sind (Regen, Schnee, Frost), oder von Staub oder aggressiven Gasen. Verwenden Sie sie nicht in explosiver Umgebung oder an Orten mit Brandgefahr. Während des Betriebs müssen folgende Umgebungsbedingungen eingehalten werden:

- Temperatur: zwischen +12°C und +35 °C
- relative Luftfeuchtigkeit: 0 bis 95 % (nicht kondensierend).

VORSICHT

Bevor Sie die Pumpe starten, muss Schmieröl eingefüllt werden, da die Pumpe leer geliefert wird.

Für weitere Informationen zum Nachfüllen des Öls und zur Auswahl des geeigneten Schmiermittels schauen Sie im Abschnitt "Ersetzung des Schmieröls" nach.

HINWEIS

Der Ölpegel muss immer zwischen den Werten MIN und MAX bleiben, die an der Seite der Pumpe angezeigt sind.

Betriebsanleitung

Nehmen Sie die erste Füllung über den Stöpsel (E) vor. Füllen Sie bis zur Hälfte der Pegelanzeige (F) auf und schließen Sie den Stöpsel (E) (siehe Abbildung unten).

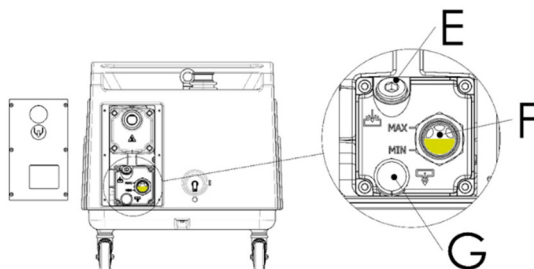


Abbildung 30

Schalten Sie die Pumpe ein und bringen Sie sie für mindestens zwei Minuten auf das maximale Vakuum. Halten Sie die Pumpe dann wieder an, prüfen Sie den Ölstand und füllen Sie dann das Öl erneut bis auf den korrekten Pegel auf.

Tipps für den Betrieb

Bei Umgebungstemperaturen unterhalb von 18 °C sollte die Pumpe 15 Minuten lang auf Temperatur gebracht werden, indem man sie mit Druck am Grenzwert (Ansaugung geschlossen, keine Last) laufen lässt.

Während dieser Phase kann es sein, dass die Pumpe nicht die angegebenen Grenzwerte für den Druck erreicht.

WARNUNG



Die Pumpe wurde nicht entwickelt, um mit atmosphärischem Druck zu arbeiten.

Es wird geraten, fünf Einschaltungen pro Stunde nicht zu überschreiten.

Gas-ballast

Die Pumpe MS120 ist mit einem Gas-Ballast und mit einem manuellen Ventil ausgestattet, das ihre Einschaltung (I) und Abschaltung (O) erlaubt.

Der Grenzwert für den Druck der Pumpe variiert je nach dieser Einstellung. Für einen korrekten Betrieb schlagen Sie unter "Ansaugung von Wasserdampf" nach.

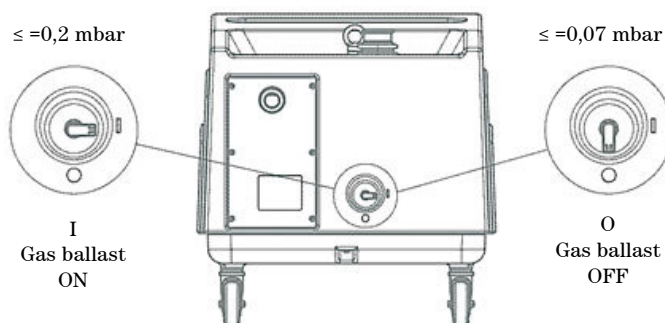


Abbildung 31

Ansaugung von Wasserdampf

Bei der Ansaugung von Dämpfen raten wir folgendes:

- Vor der Ansaugung von Dämpfen sollte die Pumpe auf Betriebstemperatur gebracht werden. Lassen Sie sie dazu eine halbe Stunde lang bei maximalem Vakuum laufen.
- Wenn nach dem Betriebszyklus Kondenswasser im Schmieröl auftritt, lassen Sie die Pumpe eine mindestens halbe Stunde lang bei maximalem Vakuum laufen. Der Gas-Ballast muss dabei offen sein.
- Führen Sie diese Schritte vor längeren Stillstandszeiten durch. Der Gas-Ballast sorgt für die Entfernung des Kondenswassers aus dem Schmieröl.

Für die Ein- und Abschaltung des Gas-Ballasts schauen Sie im Anhang "Technical Information" nach.

Wartung

Das für den Betrieb und die Wartung der Pumpe zuständige Personal muss gut ausgebildet sein und mit den Unfallverhütungsvorschriften vertraut sein.

Neben den unten wiedergegebenen Vorschriften sollten Sie auch im Abschnitt "Sicherheit" nachschlagen.

WARNUNG



Vor jedem Eingriff sollten Sie Luft in durch die Ansaugöffnung.

WARNUNG



Wenn die Pumpe nach dem Betrieb gewartet werden soll, muss sie erst abkühlen, da die Temperatur an der Außenfläche auf über 80°C steigen kann.

WARNUNG



Klemmen Sie die Pumpe immer von der Versorgung ab, bevor Sie Wartungsarbeiten durchführen. Stellen Sie beim Stromversorgungsschalter Warnschilder auf: GERÄT WIRD GEWARTET – STROM NICHT EINSCHALTEN! Aktivieren Sie die Sicherheitsvorrichtungen nach dem Ende der Wartungsarbeiten wieder.

Die Tabelle zeigt alle regelmäßig notwendigen Wartungseingriffe, damit die Pumpe in einem perfekten Zustand bleibt.

Die Notwendigkeit der häufigeren Wartungen richtet sich nach der Art der Verwendung (Ansaugung von kondensierbaren Dämpfen oder verschmutzenden Substanzen).

In diesen Fällen kann nur die Erfahrung sagen, welche Intervalle für die Wartungseingriffe richtig sind. Das Altöl und die ausgebauten Teile gelten als Sondermüll und sind nach den im jeweiligen Land geltenden Vorschriften zu behandeln.

Tab. 12 Wartungseingriffe

Wartungsintervall	Beschreibung des Eingriffs	Geeignetes Personal
24 Stunden / täglich	Kontrolle des Ölstands vor dem Start.	Bedienungspersonal
100 Stunden / wöchentlich	Reinigung der Oberfläche der Pumpe mit einem weichen Tuch. Eventuell kann ein neutrales Reinigungsmittel verwendet werden.	Bedienungspersonal
9000 Stunden / jährlich	Ersetzen des Schmieröls	qualifizierter Techniker
	Ersetzen des Ölreinigungselements	qualifizierter Techniker
14000 Stunden / alle zwei Jahre	Prüfung der elektrischen Anschlüsse	qualifizierter Techniker
30000 Stunden / alle fünf Jahre	Revision der Pumpe	Kundendienst-mitarbeiter

Ersetzung des Schmieröls

Es wird empfohlen, 1 Stunde und 30 Minuten nach dem Ausschalten der Pumpe zu warten, um den Schmierölersatzvorgang durchzuführen, damit die Temperatur der Pumpe und des Öls ausreichend abkühlen können.

WARNUNG



Verwenden Sie Handschuhe zum Schutz vor Verbrennungen.

Drehen Sie den Stöpsel (E) und den Stöpsel auf dem Ausgang (G) erst ab (Abbildung 32), nachdem Sie unter dem Tank einen nach Form und Größe passenden Behälter platziert haben, der das ganze Öl aus der Pumpe aufnehmen kann (siehe Abbildung 33).

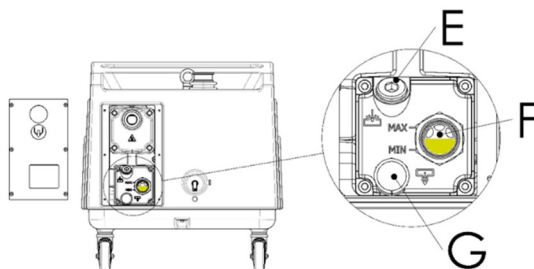


Abbildung 32

Betriebsanleitung

Nachdem Sie den Schlauch zum Ablassen des Öls (A1) auf den Ausgang (G) gesetzt haben, beginnt das Schmieröl auszulaufen. Es wird empfohlen, die Pumpe nach vorne zu kippen, indem Sie die Rückseite der Pumpe um mindestens 40 mm, damit sichergestellt ist, dass das gesamte Öl abläuft. Denken Sie auch daran, die Bremsen an den Rädern festzustellen, damit die Pumpe nicht herunterfallen kann. Wenn das alte Öl aus dem Tank abgelaufen ist, setzen Sie die Stöpsel (E und G) wieder auf und lassen Sie die Pumpe bei maximalem Vakuum etwa 30 Sekunden lang laufen, so dass auch der Schmierkreislauf sich von Ölrückständen leeren kann. Nehmen Sie dann wieder die Stöpsel ab und lassen Sie das restliche Öl ablaufen.

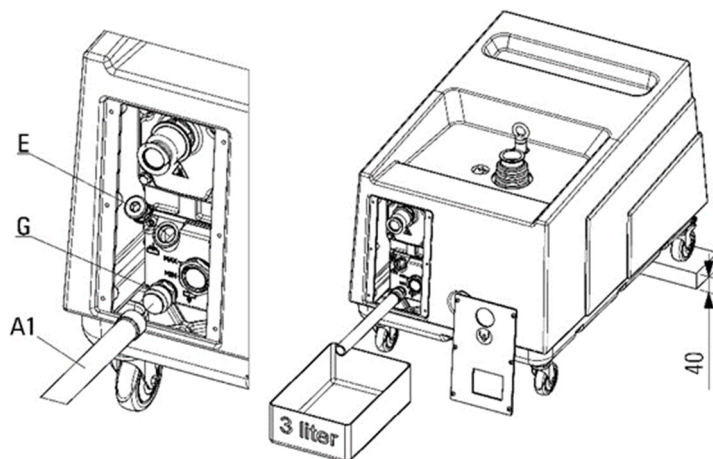


Abbildung 33

Wenn sich im Öl viele Verschmutzungen oder Wasser befinden, reinigen Sie die Pumpe, indem Sie sie bei maximalem Vakuum für etwa fünf Minuten laufen lassen.

Verwenden Sie dafür soviel Öl, dass gerade der als Minimalstand angezeigte Pegel erreicht wird.

Lassen Sie das Öl erneut ab.

Betriebsanleitung

Füllen Sie wie folgt nach (siehe Abbildung 34):

- 1 Drehen Sie den Stöpsel der Ölfüllöffnung (E) heraus.
- 2 Schrauben Sie den Öllfüllschnabel (A2) auf die Ölflasche (A1).
- 3 Füllen Sie mit Öl auf und stecken Sie dazu den Öllfüllschnabel in die Ölfüllöffnung (siehe "Inbetriebnahme und Betrieb der Pumpe MS120 Rotary Vane" und "Schmiermittel").

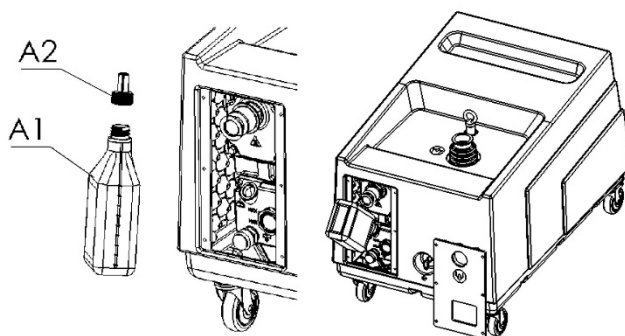


Abbildung 34

Ersetzung der Ölreinigungselemente

Stark verschmutzte Ölreinigungselemente können zu einem spürbaren Temperaturanstieg in der Pumpe und in extremen Fällen zur Selbstentzündung des Schmieröls führen.

Der maximal zulässige Druck im Tank beträgt 0,6 bar, gemessen bei maximalem Durchsatz (wenn die Pumpe mit Ansaugung bei atmosphärischem Druck arbeitet). Der Druck kann durch einen Manometer (optional) gemessen werden, der mit der Öffnung zum Einfüllen des Öls verbunden wird (Abbildung 34). Wenn der Manometer verwendet wird, prüfen Sie die Kartusche bei heißer Pumpe auf Verstopfung.

Betriebsanleitung

Zur Ersetzung des Ölreinigungselements gehen Sie wie folgt vor (Abbildung 35):

- 1** Nehmen Sie die Abdeckung (C1) ab, indem Sie die zugehörigen Schrauben (C2) herausdrehen.
- 2** Nehmen Sie den Deckel des Tanks (B1) ab, indem Sie die entsprechenden Schrauben herausdrehen (B2).
- 3** Ziehen Sie das Ölreinigungselement (B4) heraus und ersetzen Sie die O-Ringe (B5).
- 4** Setzen Sie den Deckel wieder auf den Ausgang (B1).
- 5** Ersetzen Sie wenn nötig die Dichtung (B3).
- 6** Setzen Sie die Abdeckung wieder auf (C1).

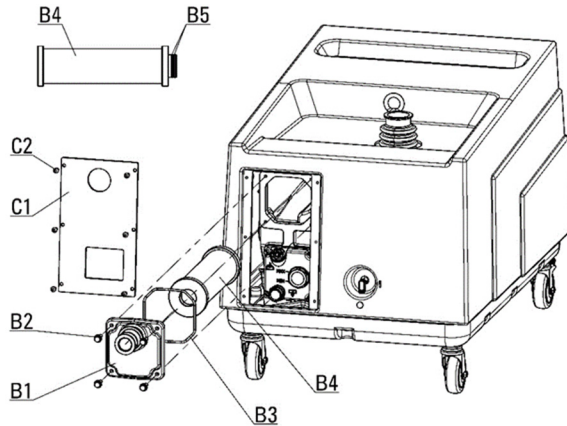


Abbildung 35

Routinewartung

Die Routinewartung finden Sie in der folgenden Tabelle. Verwenden Sie nur Original-Ersatzteile. Bei der Bestellung geben Sie bitte die Ersatzteilnummer an.

Tab. 13 Ersatzteile für die Wartung

Ersatzteilnummer	Beschreibung
X3760-64005	Rotary Vane Fluid AVF 60 Gold 1lt
X3702-68201	MS 120 Exhaust Filter Cartridge

Revision der Pumpe

Für diesen Schritt sollten Sie sich an den Kundendienst von Agilent wenden.

Die Revision besteht in einer kompletten Zerlegung, in der Reinigung aller Teile und in der Ersetzung der von Verschleiß betroffenen Teile (Lagerhülse der Pumpe, Flügel und Dichtungen). Verwenden Sie nur Original-Ersatzteile. Bei der Bestellung geben Sie bitte die Ersatzteilnummer an.

Tab. 14 Ersatzteile für die Revision der Pumpe

Ersatzteilnummer	Beschreibung
X3702-68202	MS 120 Major Maintenance Kit

So bestellen Sie Ersatzteile

Bei der Bestellung von Ersatzteilen geben Sie immer das Modell der Pumpe an, ferner die Seriennummer, das Baujahr, die Ersatzteilnummer, die Beschreibung und die gewünschte Menge.

Schmiermittel

In der folgenden Tabelle finden Sie die für den normalen Betrieb der Pumpe empfohlenen Öle.

Tab. 15 Schmiermittel

Umgebungstemperatur	Viskosität	Öl von Agilent Technologies
12 – 35 °C	ISO 53	AVF 60 Gold

Wenn auf einen anderen Typ von Schmiermittel gewechselt werden soll, muss die Pumpe gereinigt werden. Lassen Sie sie dazu für mindestens fünf Minuten bei maximalem Vakuum laufen.

Verwenden Sie dafür soviel Öl, dass gerade der als Minimalstand angezeigte Pegel erreicht wird.

Lassen Sie das alte Öl ab und füllen Sie neues nach.

Stilllegung

Für die Stilllegung müssen Sie das Öl aus dem Tank ablassen, bevor die Pumpe bewegt wird. Wenn das Öl verschmutzt erscheint, führen Sie eine Reinigung mit neuem Öl durch (siehe "Ersetzung des Schmieröls").

Leeren Sie den Tank, schließen Sie Ansaugung und Ausgang der Pumpe und lagern Sie diese ein.

HINWEIS

Bevor Sie ggf. die Pumpe zur Reparatur einschicken, muss das Formular "Request for Return" ausgefüllt und an die lokale Verkaufsstelle weitergeleitet werden. Eine Kopie dieses Formulars ist der Pumpenverpackung vor dem Versand beizulegen.

Wenn eine Pumpe entsorgt werden soll, muss sie in Übereinstimmung mit den örtlichen Gesetzen und Vorschriften entsorgt werden.

Fehlerbehebung

HINWEIS

Bei Leistungsabfall, anormale Geräuscentwicklung oder Motorschutzeingriff wenden Sie sich bitte an das Agilent Contact Center.

Tab. 16 Fehlerbehebung

Fehler	Ursache	Lösung
Leistungsmin- derung	Undichte Stellen in den Ansaugleitungen oder an der Maschine	Undichte Stellen beseitigen
	keine Schmierung	Prüfen Sie Pegel und Zustand des Öls. Füllen Sie Öl nach oder ersetzen Sie es
	Versorgungsspannung zu niedrig (< 200 V Wechselstrom)	Schließen Sie die Pumpe an eine Stromquelle mit der korrekten Spannung an und verwenden Sie ein Kabel mit passendem Querschnitt (siehe "Tabelle mit Betriebs-Grenzwerten" für den Betrieb)
Ölnebel am Ausgang	Ölreinigungselement funktionsuntüchtig	Ölreinigungselement ersetzen
	Hohe Temperatur durch verschmutztes Öl	Öl ersetzen
	Hohe Betriebstemperatur durch zu hohe Umgebungstemperatur	Umgebungstemperatur durch bessere Luftzirkulation senken
	Ölreinigungselement verstopft	Ersetzung des Ölreinigungselements
Pumpe startet nicht	Versorgungsspannung außerhalb des zulässigen Bereichs	Schließen Sie die Pumpe an ein Netz mit geeigneter Spannung an (siehe "Tabelle mit Betriebs-Grenzwerten" für den Betrieb)

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 Industriemüll 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 36 Logo "WEEE"

Für weitere Informationen siehe:

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

Service

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

vpt-customer@agilent.com

vpl-customer@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	90
Validité	90
Définitions et terminologie	91
Signification des messages d'attention, d'avertissement et des notes	91
Symboles d'avertissement	92
Sécurité	94
Utilisation appropriée	94
Utilisation incorrecte	96
Tableau des limites opératoires	96
Équipements de protection	98
Consignes de sécurité relatives aux pompes MS120 Single Stage Rotary Vane	99
Information du produit	101
Transport et entreposage	102
Levage	103
Installation	104
Préparation pour l'installation	104
Accessoires fournis avec la pompe	105
Assemblage	105
Positionnement	106
Raccordement à la machine utilisatrice	107
Acheminement de l'air d'extraction	108
Raccordement électrique	109
Déballage	110

Mode d'emploi

Utilisation	113
Démarrage et exploitation de la pompe MS120	
Single Stage Rotary Vane	113
Conseils d'utilisation	114
Dispositif de lest (Gas-ballast)	115
Aspiration de vapeur d'eau	115
Entretien	116
Remplacement de l'huile lubrifiante	117
Remplacement des éléments déshuileurs	119
Entretien de routine	121
Révision pompe	121
Accessoires à commander	121
Lubrifiants	122
Mise hors service	122
Problèmes et solutions	123
Élimination	124
Service	125

À propos de ce manuel

Validité

Le présent manuel contient les instructions destinées aux utilisateurs de la pompe MS120 Single Stage Rotary Vane, notamment celles relatives à la sécurité, au fonctionnement et à la maintenance de premier niveau, dans la limite des activités de maintenance incombant à l'utilisateur.

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

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

NOTE

- 1 Ce manuel contient des informations utiles pour que l'ensemble du personnel puisse utiliser la pompe MS120 Single Stage Rotary Vane, en toute sécurité et pour garantir un fonctionnement parfait pendant toute sa durée de vie.
 - 2 Veuillez conserver ce manuel et tous les documents connexes dans un lieu accessible connu de tous les opérateurs et du personnel de maintenance.
-

Définitions et terminologie

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

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

ATTENTION

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

AVERTISSEMENT



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

NOTE

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

Symboles d'avertissement

Voici une liste de symboles qui apparaissent en conjonction avec les avertissements sur la pompe MS120 Single Stage Rotary Vane.


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.



	Danger générique
	Tensions dangereuses
	Surfaces chaudes
	Courant Alternatif (AC)
	Borne de mise à la terre de protection
	Danger d'incendie
	Ne pas déverser dans l'environnement
	Lire le manuel d'utilisation
	Raccord d'aspiration
	Raccord d'évacuation

Mode d'emploi

	Gants de protection
	Chaussures de sécurité
	Equipements de protection
	Emission de substances dangereuses
	Risques de blessures par des arêtes vives
	Chute d'objets
	Certification CSA
RoHS	Restriction de la certification des substances dangereuses
CE	Déclaration européenne de conformité
	Déchets d'équipements électriques et électroniques
UK CA	Marquage UK CA

Sécurité

Cette section contient les informations prescrites par la Directive basse tension 2014/35/EU, qui est essentielle à la conformité et au respect des règles de sécurité en général ainsi qu'à l'utilisation spécifique de la machine.

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

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

AVERTISSEMENT



Prévoir des mesures de sécurité sur l'installation lorsque l'arrêt ou une défaillance de la pompe à vide risquent de causer des dommages corporels et matériels.

Utilisation appropriée

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

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

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

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

AVERTISSEMENT



Matières inflammables :

Pour éviter que les gaz à l'intérieur de la pompe n'atteignent leur plage d'inflammabilité/de combustion, vous devez respecter les consignes et prendre note des précautions reportées ci-dessous. Lorsque des matières potentiellement inflammables ou combustibles sont présentes dans l'équipement, vous devez :

- Vérifier que le système est étanche pour vous assurer que seuls les gaz prévus passent par la pompe à vide.
- Vérifier que la concentration de n'importe quels gaz ou vapeurs inflammables qui entrent dans la pompe (bride d'entrée et orifice "Gas Ballast"), à l'intérieur de la pompe et dans le tuyau d'évacuation reste inférieure à un quart de la Limite inférieure d'explosivité (LIE) des gaz.
- Vérifier que les systèmes de sécurité sont en place pour maintenir la concentration de n'importe quels gaz ou vapeurs inflammables qui entrent dans la pompe (bride d'entrée et orifice "Gas Ballast"), à l'intérieur de la pompe et dans le tuyau d'évacuation à une valeur inférieure à un quart de la Limite inférieure d'explosivité (LIE) des gaz quel que soit le circuit ou dans toutes les conditions de défaillance de la pompe.

NOTE

- Avant de remettre la pompe en service après une panne, inspectez-la et vérifiez soigneusement tout autre signe de dommage.
- Utiliser uniquement des outils en parfait état de fonctionnement et spécialement conçus pour le travail ; l'utilisation d'outils inappropriés ou inefficaces peut causer de graves dommages.
- Vérifiez toujours le lubrifiant et sa bonne répartition dans la pompe ; une lubrification inadéquate peut endommager sérieusement la pompe.
- Donnez aux pièces une forme de marquage lorsque vous les démontez pour vous assurer de les remonter dans le bon ordre.
- Vérifier qu'il n'y a pas de rayures ou de rainures sur les arbres usinés, dans leurs sièges à l'intérieur de la pompe ou sur les surfaces rectifiées par la machine. De légères rayures et abrasions peuvent être éliminées avec du papier émeri très fin ou par un léger meulage.
- Avant de constituer un groupe, étalez toujours un peu d'huile sur les pièces intérieures et les surfaces de contact. Remplacez tous les joints par des pièces de rechange d'origine avant de remonter les composants.

Utilisation incorrecte

Agilent Technologies décline toute responsabilité liée à une utilisation incorrecte de la pompe MS120 Single Stage Rotary Vane.

Une utilisation incorrecte provoquera une annulation de la garantie et des droits de recours en responsabilité. Le personnel responsable de l'exploitation et de l'entretien de la pompe doit avoir été formé correctement et est tenu de connaître les règles de prévention des accidents. Les consignes de prévention des accidents contenues dans cette section doivent être respectées tout au long de l'exploitation et des interventions d'entretien de la pompe afin d'éviter de blesser les opérateurs et d'endommager la pompe. Ces précautions sont indiquées par les notes AVERTISSEMENT et ATTENTION.

AVERTISSEMENT



Il est interdit d'aspirer avec la pompe:

- des liquides ou des substances solides
- des gaz et des vapeurs dangereux, explosifs ou agressifs
- de l'oxygène pur ou des mélanges d'air enrichi d'oxygène.

AVERTISSEMENT



Il est interdit d'utiliser l'évacuation de la pompe pour créer des pressions même limitées.

AVERTISSEMENT



Il est interdit d'utiliser la pompe dans des milieux potentiellement explosifs.

Tableau des limites opératoires

Tab. 17

Tension d'entrée (V)	État du Controller
< 180	Défaut d'alimentation (Erreur de Sous-tension/Surtension – Win206 bit 5 = 1)
180 – 200	En fonction (performances réduites)
200 – 264	En fonction (performances complètes)
> 264	Défaut d'alimentation (Erreur de Sous-tension/Surtension – Win206 bit 5 = 1)

AVERTISSEMENT



- Le contact avec des tensions élevées peut entraîner la mort. Soyez toujours extrêmement prudent et respectez les règles de prévention des accidents en vigueur.
- Débranchez toujours l'alimentation électrique de la pompe avant les travaux de maintenance. Placez un panneau d'avertissement spécial sur le disjoncteur d'alimentation électrique : MACHIN EN ENTRETIEN - NE PAS ALLUMER.
- Le fait de ne pas fournir à la pompe une connexion à la terre peut causer de graves dommages aux opérateurs. Assurez-vous toujours qu'il y a une connexion à la terre et qu'elle est conforme aux normes.
- Lors du nettoyage de la pompe et de ses composants, évitez d'utiliser des solvants inflammables ou toxiques, tels que l'essence, le benzol, l'éther ou l'alcool. Il est recommandé d'utiliser une solution d'eau et de savon, de préférence dans des machines à laver à ultrasons, en prenant soin de sécher toutes les pièces nettoyées à des températures inférieures à 100 °C afin d'éliminer l'humidité résiduelle.
- Des surcharges ou des pannes prolongées peuvent provoquer une surchauffe du moteur électrique et un dégagement de fumées nocives ; coupez immédiatement l'alimentation électrique par mesure de précaution et ne vous approchez pas de la pompe au moins jusqu'à ce que vous ayez prévu une ventilation pour chasser la fumée.
- En cas d'incendie, ne jetez pas d'eau sur la pompe. Coupez l'alimentation électrique et utilisez des extincteurs à CO₂.
- Inspectez soigneusement les brides pour vous assurer qu'il n'y a pas de poussière, d'huile, de saleté ou de défauts des surfaces de contact, avant d'effectuer les connexions requises.
- Assurez-vous que tous les joints et accouplements sont correctement verrouillés avant de redémarrer la pompe après des travaux de réparation.
- Ne portez aucun objet susceptible de s'emmêler dans les mécanismes et/ou de servir de conducteur (chaînes, bracelets...).
- S'assurer que les outils à utiliser sont en parfait état de fonctionnement et disposent de poignées isolantes, si nécessaire. Vérifier que le matériau isolant des câbles et que les conducteurs de l'équipement de test ne présentent aucun signe de dommage.
- Ne remplacez pas l'huile immédiatement après l'arrêt de la machine car l'huile peut encore être à haute température.

NOTE

Utilisez le ressort de rétention pour fixer le câble secteur dans la prise IEC320.

Équipements de protection

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

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, les pompes 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 les pompes 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.

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.
- Si nécessaire, portez des gants de protection conformes à la directive EN 420.

Consignes de sécurité relatives aux pompes MS120 Single Stage Rotary Vane

AVERTISSEMENT



Malgré les précautions prises en phase de projet, il existe des éléments de risque qui se présentent durant les opérations effectuées durant l'utilisation et l'entretien.

AVERTISSEMENT



SURFACES CHAUDES

Durant les opérations d'entretien, on touche des surfaces dont la température peut dépasser 80°C. Adopter des moyens de protection appropriés pour éviter les brûlures dues à un contact fortuit. On recommande d'attendre le refroidissement de la pompe avant toute intervention sur cette dernière.

AVERTISSEMENT



ÉMISSIONS DE SUBSTANCES NOCIVES

L'air d'extraction de la pompe contient des traces de brouillards d'huile (résidu max. 2PPM/ poids équivalant à 2,4 mg/m³). Vérifier la compatibilité avec le milieu de travail. Garantir un renouvellement correct de l'air ou amener l'évacuation de la pompe à l'extérieur. Une défaillance ou l'usure des garnitures peut provoquer des fuites de huile lubrifiante. Éviter le déversement sur le terrain et la pollution d'autres matériaux. En cas d'aspiration d'air contenant des substances dangereuses (comme des agents biologiques ou microbiologiques), prévoir des systèmes de captage avant la pompe à vide.

AVERTISSEMENT



NE PAS DÉVERSER DANS L'ENVIRONNEMENT

Les huiles usées provenant de la pompe doivent être éliminées conformément aux réglementations en vigueur dans le pays d'utilisation.

AVERTISSEMENT DANGER GÉNÉRÉ PAR LA DÉPRESSION



Le contact avec des points en dépression peut provoquer des accidents. Éviter le contact avec le raccord d'aspiration de la pompe durant le fonctionnement. Injecter de l'air dans le circuit d'aspiration avant toute intervention.

AVERTISSEMENT DANGER GÉNÉRÉ PAR LA PRESSION



Le réservoir de la pompe est pressurisé. Ne pas ouvrir et ne pas oublier de fermer les bouchons de remplissage et de vidange durant le fonctionnement.

AVERTISSEMENT Toujours exclure l'alimentation de la pompe avant d'effectuer des opérations d'entretien. Appliquer des pancartes spécifiques d'avertissement : ENTRETIEN EN COURS SUR L'APPAREIL - NE PAS METTRE SOUS TENSION, au niveau de l'interrupteur d'alimentation. Rétablir les dispositifs de sécurité au terme de l'opération.



AVERTISSEMENT



SÉCURITÉ ÉLECTRIQUE

L'équipement électrique comprend des pièces sous tension qui, en cas de contact, peuvent provoquer des dommages corporels et matériels graves.

AVERTISSEMENT



DANGER D'INCENDIE

L'utilisation de la pompe pour des usages non prévus ou interdits dans ce manuel ou un entretien incorrect peuvent provoquer des anomalies de fonctionnement et générer un risque de surchauffe et d'incendie. Ne pas éteindre les flammes avec de l'eau en cas d'incendie. Utiliser des extincteurs à la poudre ou au CO₂ ou d'autres moyens compatibles avec la présence d'équipements électriques et d'huiles lubrifiantes.

Information du produit

Cet équipement est destiné à être utilisé par un personnel formé. L'utilisateur doit lire ce manuel d'instructions et toute autre information supplémentaire fournie par Agilent avant d'utiliser l'équipement. Agilent ne sera pas tenu responsable en cas d'événement se produisant en raison de la non-conformité à ces instructions, de l'utilisation inappropriée par des personnes non formées, de l'interférence non autorisée avec l'équipement ou de toute action contraire à ce qui est énoncé par les normes nationales spécifiques.

Les conseils s'adressent à un personnel préposé à l'installation et à l'entretien expert et préparé pour affronter tout problème d'entretien, mécanique et électrique. En cas de doute ou d'informations ne figurant pas dans ce manuel, veuillez contacter notre service d'assistance en communiquant toujours le modèle (Model), le numéro de série (Serial), l'année de construction indiqués sur la plaque d'identification.

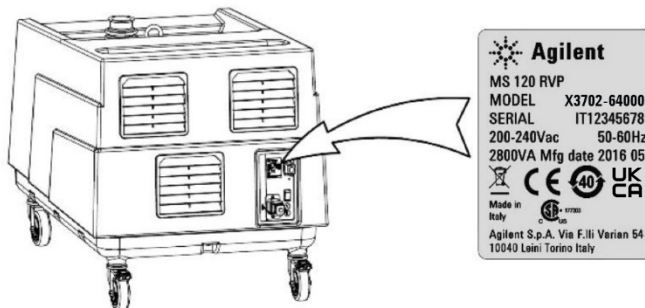


Figure 37 Targhetta di identificazione

Les MS120 Single Stage Rotary Vane Pumps sont des pompes rotatives à simple étage à palettes, étanches à bain d'huile, actionnées par un moteur électrique commandé par inverter. Ces pompes à vide poussé sont exclusivement indiquées pour le pompage de l'air et de petites quantités de vapeur d'eau. L'aspiration d'autres types de gaz ou de vapeur doit être préalablement déclarée à la société Agilent qui délivrera au besoin la conformité à l'usage spécifique.

Mode d'emploi

Le moteur électrique bridé est accouplé par l'intermédiaire d'un joint élastique. Le refroidissement est assuré par un ventilateur centrifuge puissant.

À l'aspiration, un filtre à crépine protège la pompe contre les corps solides d'un diamètre supérieur à 4,7 mm. En outre, un clapet de retenue intégré empêche la remontée de l'huile et le retour de l'air dans la chambre à vide au cours de la phase d'arrêt.

Dans le réservoir, un système de séparation des brouillards d'huile de l'air d'extraction (résidu max. 2PPM/ poids équivalant à 2,4 mg/m³).

L'huile éliminée est automatiquement récupérée par la pompe.

Le dispositif de lest empêche la condensation à l'intérieur de la pompe lorsqu'on aspire de petites quantités de vapeur (voir le paragraphe "Dispositif de lest (Gas-ballast)").

Les paragraphes suivants fournissent toutes les informations nécessaires pour garantir la sécurité de l'opérateur durant l'utilisation de l'appareil. On fournit des informations détaillées dans l'appendice "Technical Information".

Transport et entreposage

Les pompes doivent être transportées sans huile.

Respecter les conditions ambiantes suivantes durant le transport et le stockage:

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

Mode d'emploi

Levage

L'orientation des composants emballés doit rester conforme aux indications fournies par les pictogrammes présents à l'extérieur de l'emballage.

Effectuer l'opération de déchargement avec un moyen de levage compatible avec le poids de la pompe. Pour le levage de l'emballage et de la pompe voir les figures suivantes.

Le poids de l'emballage, comprenant la pompe, atteint un maximum de 120 kg, tandis que le poids de la pompe seule est de 100 kg.

Tab. 18 Dimensions de l'emballage

Cote	L	B	H
Dimension	920 mm	610 mm	790 mm

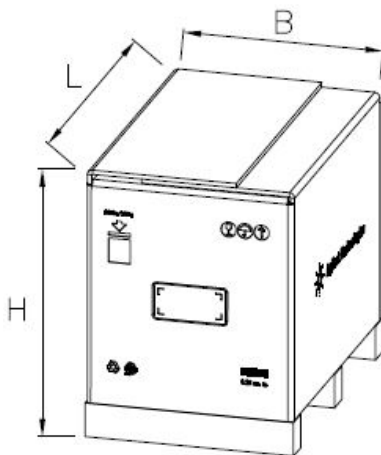


Figure 38 Pompe emballée

Mode d'emploi

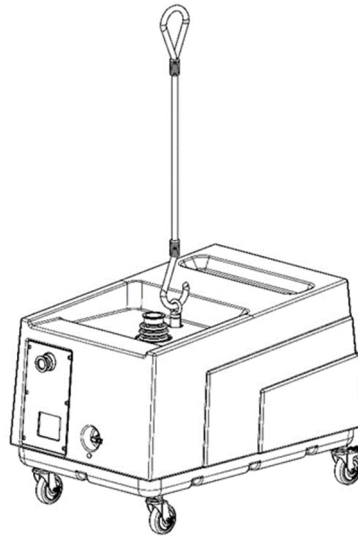


Figure 39 Pompe

Installation

Préparation pour l'installation

La pompe est livrée dans un emballage de protection spécial. En cas de détection de signes de dommages, susceptibles de s'être produits pendant le transport, contactez votre service commercial local.

AVERTISSEMENT



Lors du déballage de la pompe, veillez à ce qu'elle ne tombe pas et évitez de l'exposer à des chocs brusques ou des vibrations liées à un impact.

Mode d'emploi

Accessoires fournis avec la pompe

La pompe est livrée avec les accessoires suivants nécessaires pour la mise en service et l'entretien ordinaire:

- Bouteille d'huile
- Embout de remplissage d'huile
- Type de vidange de l'huile
- I/O Mating Connector.

Assemblage

Enlever les bouchons sur l'aspiration et la vidange.

La pompe est fournie prête à l'emploi, sans aucune partie à assembler.

La pompe est sans huile, voir le paragraphe "Remplacement de l'huile lubrifiante" pour les instructions du remplissage en huile.

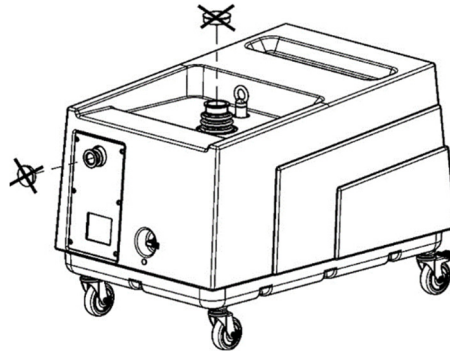


Figure 40

Positionnement

AVERTISSEMENT La pompe:



- doit être bloquée au moyen des freins situés sur les roues d'appui, sur un plan horizontal.
- doit être accessible pour un entretien correct et aisé en respectant les distances minimales par rapport aux encombrements éventuels (voir la figure suivante). Elle devrait également être accessible avec des moyens de levage appropriés.
- doit être protégée contre les jets ou les éclaboussures d'eau qui risquent de provoquer des courts-circuits électriques et / ou l'électrocution des opérateurs à proximité de la machine.

Tab. 19 Distance minimale

Cote	A	B	C	D
Dimension	150 mm	200 mm	150 mm	500 mm

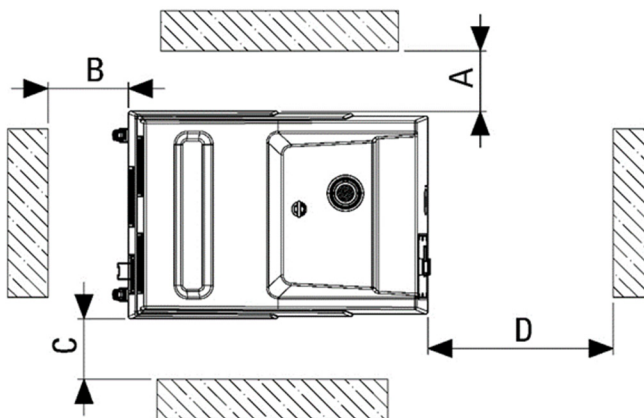


Figure 41

La pompe ne pas être installée à l'extérieur et elle doit toujours être protégée contre les agents atmosphériques.

ATTENTION

Éviter que l'air chaud provenant de l'évacuation ou des ventilateurs de refroidissement ne gênent le personnel.

Ne pas installer la pompe dans une zone où la poussière ou d'autres matériaux risquent d'engorger ou de couvrir rapidement les surfaces de refroidissement.

Mode d'emploi

Raccordement à la machine utilisatrice

Le raccordement de la pompe à la chambre à évacuer doit être réalisé avec des conduites du même diamètre que la bouche d'aspiration.

Le poids des conduites et les éventuelles dilatations ne doivent pas peser sur la pompe. Réaliser le segment final de connexion au raccord d'aspiration de la pompe avec un tube / raccord flexible.

Il est important que toutes les conduites et les différents joints soient étanches.

Des conduites très longues ou de petit diamètre réduisent les performances de la pompe.

NOTE

Ce symbole identifie le raccord d'aspiration.

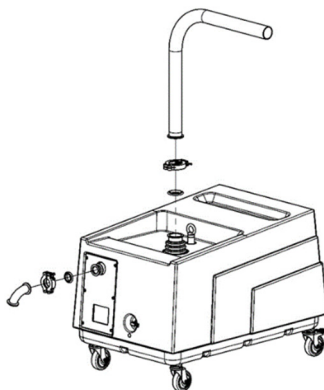


Figure 42

Mode d'emploi

Acheminement de l'air d'extraction

En cas de renouvellement insuffisant de l'air dans le local de la pompe, on peut acheminer l'air d'extraction vers d'autres locaux ou à l'extérieur.

Utiliser des conduites du diamètre de la bouche d'évacuation du réservoir et d'une longueur maximale de 15 m.

Pour des longueurs supérieures, augmenter le diamètre du tube. Le poids des conduites ne doit pas peser sur la pompe.

Utiliser des raccords ou des tuyaux flexibles pour réaliser le segment final.

AVERTISSEMENT

Cette conduite doit être inclinée pour éviter le retour de la condensation dans le réservoir de la pompe.



ATTENTION

Ne pas installer de robinets sur cette conduite. Ne pas obstruer la sortie.

NOTE

Ce symbole identifie le raccord d'évacuation.



Raccordement électrique

Vérifier la tension et la fréquence de réseau avec les données indiquées sur la plaque de la pompe.

S'assurer que le circuit de mise à la terre fonctionne.

AVERTISSEMENT

S'assurer que l'interrupteur d'allumage est mis sur le hors tension (0) pour éviter les démarrages indésirables de la pompe.



Réaliser le raccordement électrique en introduisant la broche du câble d'alimentation dans la prise de la pompe et en la bloquant avec le ressort fourni (voir la figure suivante).

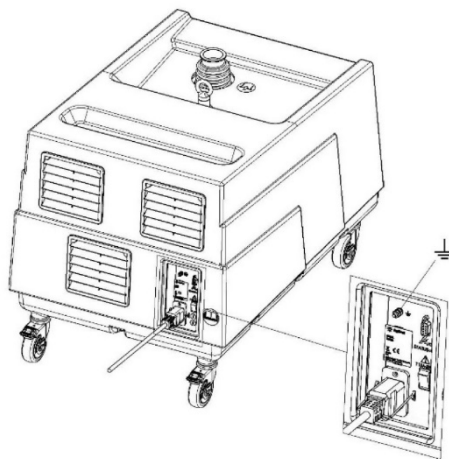


Figure 43

Mode d'emploi

Déballage

Durant l'opération de déballage, faire particulièrement attention à ne pas laisser tomber la pompe et à ne pas la soumettre à des chocs ou à des vibrations.

Ne pas abandonner l'emballage dans l'environnement. Le matériel est complètement recyclable et il est conforme à la directive CEE 85/399 pour la protection de l'environnement.

NOTE

La pompe ne se détériore passivement lorsqu'elle est exposée à l'atmosphère. On conseille toutefois de ne pas la débiller jusqu'au moment de l'installation sur le système pour éviter toute contamination éventuelle due à de la poussière.

Enlever les feuilards qui ferment le carton sur la palette puis enlever le carton.

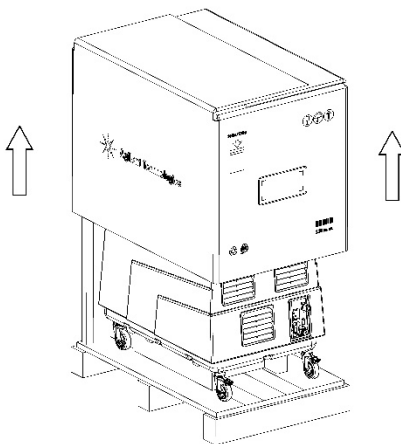


Figure 44

Mode d'emploi

Extraire les 8 vis qui ferment les étriers, puis enlever ces derniers.

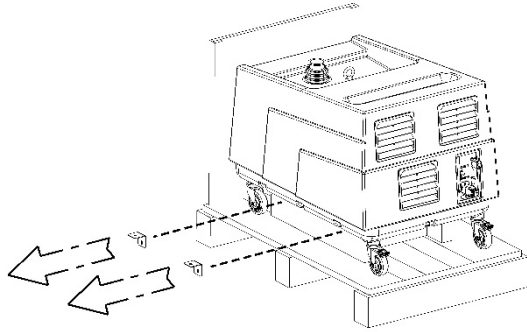


Figure 45

Débloquer les freins d'arrêt des roues.

Fixer la rampe de descente avec 1 vis d'après la figure.

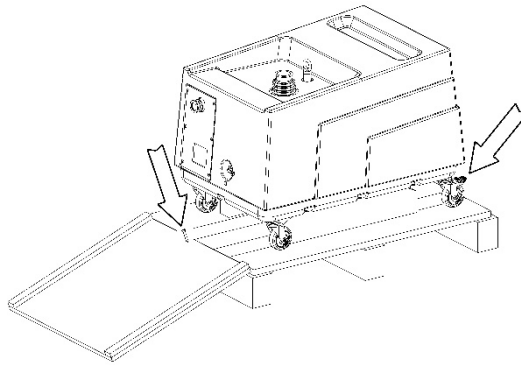


Figure 46

Mode d'emploi

Pousser la pompe par l'arrière pour la faire glisser et la descendre de la palette.

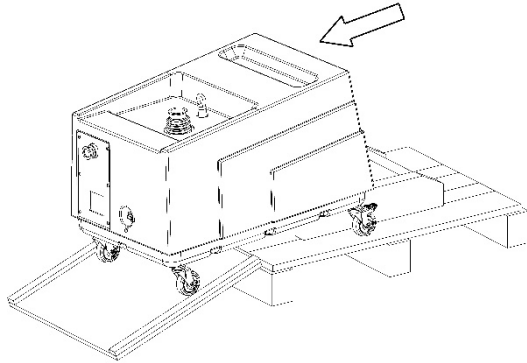


Figure 47

Utilisation

Les pompes à vide décrites dans ce manuel peuvent exclusivement aspirer de l'air et de petites quantités de vapeur d'eau. L'aspiration d'autres types de gaz ou de vapeurs doit être préalablement déclarée à la société Agilent Technologies qui délivrera au besoin la conformité à l'usage spécifique.

Elles sont indiquées pour l'évacuation de systèmes fermés ou pour fonctionner avec un vide constant compris dans le champ suivant : 0,07 - 15 mbar (absolus).

La température ambiante et la température d'aspiration doivent être comprises entre 12 et 35°C.

Démarrage et exploitation de la pompe MS120 Single Stage Rotary Vane

Ne pas installer et/ou utiliser la pompe dans des milieux exposés aux agents atmosphériques (pluie, gel, neige), aux poussières, aux gaz agressifs, dans des milieux explosifs ou présentant un risque d'incendie important. Respecter les conditions environnementales suivantes durant le fonctionnement :

- température : de +12 °C à +35 °C
- humidité relative : 0 – 95 % (non condensante).

ATTENTION

La pompe est fournie sans huile lubrifiante ; ainsi, avant de la démarrer la pompe, faire le plein en huile.

Pour de plus amples informations relatives à l'approvisionnement en huile et au choix du lubrifiant approprié, voir le paragraphe "Remplacement de l'huile lubrifiante".

NOTE

Il est important que le niveau d'huile se stabilise entre les valeurs MIN et MAX affichées par l'indicateur de niveau situé sur le côté de la pompe.

Procéder au premier remplissage à travers le bouchon (E) jusqu'à la moitié de l'indicateur de niveau (F) et refermer le bouchon (E) (voir la figure suivante).

Mode d'emploi

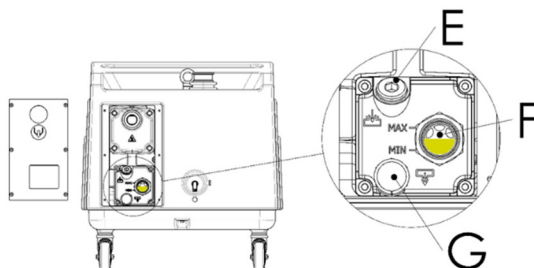


Figure 48

Allumer la pompe et l'amener au vide maximum pendant au moins 2 minutes. Arrêter la pompe, recontrôler le niveau d'huile et éventuellement faire l'appoint en rétablissant le niveau correct.

Conseils d'utilisation

À des températures inférieures à 18 °C, on conseille de chauffer la pompe pendant 15 minutes en la faisant tourner à la pression limite (aspiration fermée, sans charge).

Au cours de cette phase, la pompe pourrait ne pas atteindre les limites de pression déclarées.

AVERTISSEMENT



La pompe n'a pas été conçue pour fonctionner à la pression atmosphérique.
On conseille de ne pas dépasser 5 démarrages/heure.

Mode d'emploi

Dispositif de lest (Gas-ballast)

La pompe MS120 est équipée d'un dispositif de lest (Gas Ballast) et d'une vanne manuelle qui permet son activation (I) ou son exclusion (O).

La pression limite de la pompe varie en fonction de ce choix. Pour une utilisation correcte, voir "Aspiration de vapeur d'eau".

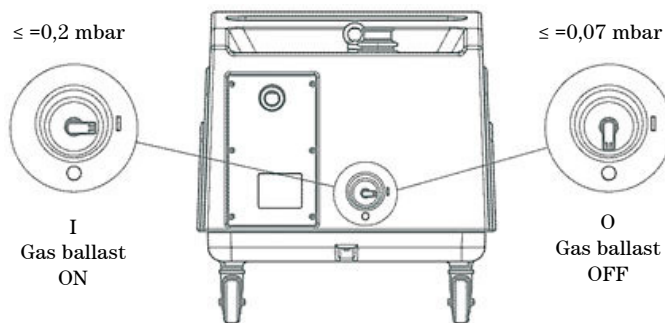


Figure 49

Aspiration de vapeur d'eau

Si l'on aspire de la vapeur d'eau, on conseil :

- d'aspirer d'abord les vapeurs, d'amener la température de la pompa à plein régime, en la faisant fonctionner pendant 30 minutes au vide maximum.
- au terme du cycle de fonctionnement, en présence de condensation dans l'huile lubrifiante, laisser fonctionner la pompe au vide maximum pendant au moins 30 minutes avec le dispositif de lest ouvert.
- effectuer cette opération avant les arrêts prolongés de la machine; le dispositif de lest permettra d'éliminer les condensations de l'huile lubrifiante.

Pour la procédure d'activation ou d'exclusion du dispositif de l'Est, consulter l'appendice "Technical Information".

Entretien

Le personnel préposée à l'utilisation et à l'entretien de la pompe doit être bien formé et posséder une connaissance approfondie des normes de prévention des accidents.

Outre les prescriptions indiquées ci-après, consulter le paragraphe "Sécurité".

AVERTISSEMENT Avant toute intervention, introduire de l'air par le raccord d'aspiration.



AVERTISSEMENT S'il faut procéder à des opérations d'entretien de la pompe au terme d'une période de fonctionnement, il faut la laisser refroidir car la température externe peut dépasser 80 °C.



AVERTISSEMENT Toujours exclure l'alimentation de la pompe avant d'effectuer des opérations d'entretien. Appliquer des pancartes spécifiques d'avertissement : **ENTRETIEN EN COURS SUR L'APPAREIL - NE PAS METTRE SOUS TENSION**, au niveau de l'interrupteur d'alimentation. Rétablir les dispositifs de sécurité au terme de l'opération.



Le tableau indique toutes les interventions périodiques nécessaires pour garder la pompe en parfait état de marche.

Des interventions d'entretien plus fréquentes peuvent être nécessaires selon le type d'utilisation (aspiration de vapeurs condensables ou de substances polluantes).

Dans ces cas, seule l'expérience directe peut suggérer les intervalles d'entretien corrects. L'huile usée et les pièces de rechange remplacées doivent être considérées comme des déchets spéciaux et gérées conformément à la réglementation en vigueur dans le pays d'utilisation.

Mode d'emploi

Tab. 20 Interventions d'entretien

Intervalle d'entretien	Description intervention	Personnel habilité
24 heures / par jour	Contrôle du niveau d'huile avant le démarrage.	Opérateur
100 heures / par semaine	Nettoyer les surfaces de la pompe avec un chiffon doux. On peut éventuellement utiliser un détergent neutre.	Opérateur
9000 heures / par an	Remplacer l'huile lubrifiante	Technicien qualifié
	Remplacer l'élément déshuileur	Technicien qualifié
14000 heures / tous les 2 ans	Vérifier les raccordements électriques.	Technicien qualifié
30000 heures / tous les 5 ans	Révision pompe	Service d'Assistance

Remplacement de l'huile lubrifiante

L'est recommandé d'attendre 1 heure et 30 minutes après l'arrêt de la pompe pour effectuer l'opération de remplacement de l'huile de lubrifiant pour permettre à la température de la pompe et de l'huile de refroidir suffisamment.

AVERTISSEMENT



Utiliser des gants de protection pour éviter de se brûler.

D'après la Figure 50, dévisser le bouchon de remplissage (E) et le bouchon de vidange d'huile (G) uniquement après avoir positionné sous le réservoir un récipient indiqué (au niveau de la forme et des dimensions) afin de récupérer toute l'huile de la pompe (voir Figure 51).

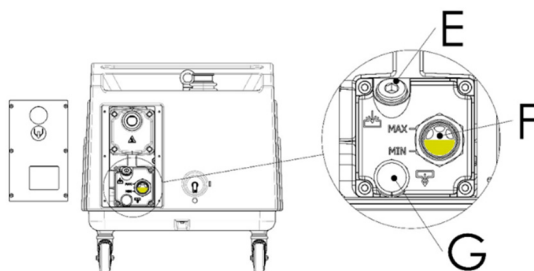


Figure 50

Mode d'emploi

Lorsqu'on visse le tube de vidange d'huile (rep. A1) sur l'orifice d'évacuation (G), le lubrifiant commence à s'écouler. L'est recommandé d'incliner la pompe vers l'avant en soulevant l'arrière de la pompe d'au moins 40 mm pour garantir l'écoulement complet du lubrifiant; ne pas oublier de freiner les roues pour éviter une chute de la pompe. Quand l'huile usée contenue dans le réservoir se sera complètement écoulee, remonter les deux bouchons ("E" et "G") et faire tourner la pompe au vide maximum pendant environ 30 secondes de manière à éliminer également du circuit de lubrification les résidus d'huile usée, puis enlever les bouchons et vidanger l'huile restante.

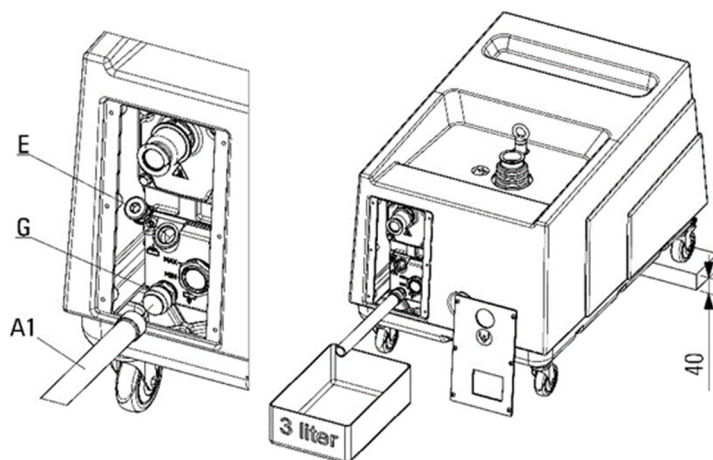


Figure 51

Si l'huile contient de grandes quantités de substances polluantes ou si l'on relève la présence d'eau, procéder à un lavage de la pompe en la faisant fonctionner au vide maximum pendant au moins 5 minutes.

Pour cette opération, utiliser une quantité d'huile suffisante pour atteindre le niveau minimum indiqué sur le réservoir.

Vidanger de nouveau l'huile.

Mode d'emploi

Procéder au remplissage comme suit (voir la Figure 52) :

- 1 Enlever le bouchon de remplissage d'huile (E).
- 2 Visser l'embout de remplissage d'huile (A2) sur la bouteille d'huile (A1).
- 3 Procéder au remplissage en insérant l'embout à l'intérieur du trou de ravitaillement en huile (voir "Démarrage et exploitation de la pompe MS120 Single Stage Rotary Vane" et "Lubrifiants").

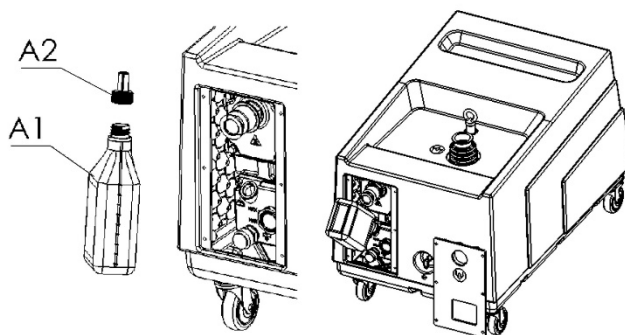


Figure 52

Remplacement des éléments déshuileurs

Des éléments déshuileurs très encrassés peuvent provoquer une augmentation sensible de la température de la pompe et dans des cas extrêmes, une autocombustion de l'huile lubrifiante.

La pression maximale admise dans le réservoir est de 0,6 bars mesurée au débit maximum (quand l'aspiration fonctionne à la pression atmosphérique). On peut relever la pression avec un manomètre (en option), raccordé à l'orifice de remplissage en huile (Figure 52). En présence du manomètre, vérifier l'engorgement de la cartouche lorsque la pompe est chaude.

Mode d'emploi

Pour le remplacement de l'élément déshuileur, voir Figure 53, et procéder comme suit :

- 1** enlever le couvercle du cover (rep. C1) en desserrant les vis correspondantes (C2),
- 2** enlever le couvercle du réservoir (rep. B1) en desserrant les vis correspondantes (rep. B2),
- 3** extraire et remplacer l'élément déshuileur (rep. B4) et les joints toriques correspondants (rep. B5),
- 4** remonter le couvercle de vidange (rep. B1),
- 5** au besoin, remplacer le joint (rep. B3),
- 6** remonter le couvercle du cover (rep. C1).

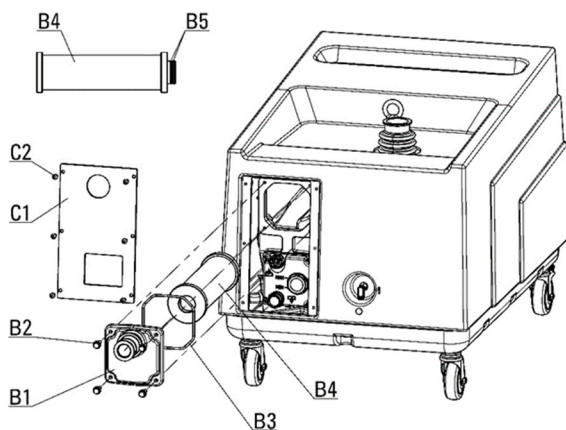


Figure 53

Mode d'emploi

Entretien de routine

Les pièces de rechange nécessaires pour l'entretien de routine sont indiquées dans le tableau suivant. Utiliser exclusivement des pièces de rechange originales. Au moment de la commande, indiquer le Part Number fourni.

Tab. 21 Pièces de rechange pour l'entretien

Part Number	Description
X3760-64005	Rotary Vane Fluid AVF 60 Gold 1lt
X3702-68201	MS 120 Exhaust Filter Cartridge

Révision pompe

Pour cette opération, on conseille de s'adresser au Service d'Assistance de la société Agilent.

La révision consiste à démonter complètement, à nettoyer toutes les pièces et à remplacer les pièces d'usure (coussinets de la pompe, palettes et joints). Utiliser exclusivement des pièces de rechange originales. Au moment de la commande, indiquer le Part Number fourni.

Tab. 22 Pièces de rechange pour la révision de la pompe

Part Number	Description
X3702-68202	MS 120 Major Maintenance Kit

Accessoires à commander

Pour commander les pièces de rechange, toujours indiquer le modèle de la pompe (model), le numéro de série (serial), l'année de construction, le Part Number, la description et la quantité nécessaire.

Lubrifiants

Le tableau suivant indique les huiles préconisées pour un usage générique de la pompe.

Tab. 23 Lubrifiants

Température ambiante	Viscosité	Huile Agilent Technologies
12 – 35 °C	ISO 53	AVF 60 Gold

Lorsqu'on change de type de lubrifiant, il faut procéder à un lavage de la pompe en la faisant fonctionner au vide maximum pendant au moins 5 minutes.

Pour cette opération, utiliser une quantité d'huile suffisante pour atteindre le niveau minimum indiqué sur le réservoir.

Vidanger de nouveau l'huile et procéder à un nouveau remplissage.

Mise hors service

Pour la mise hors service, vidanger l'huile de la pompe avant sa manutention. Si l'huile est contaminée, procéder à un lavage avec de l'huile neuve (voir "Remplacement de l'huile lubrifiante").

Vider le réservoir d'huile, boucher l'aspiration et vider la pompe avant le stockage.

NOTE

Avant de renvoyer la pompe à l'usine pour réparation, vous devez remplir le formulaire « Request for Return » joint à ce manuel d'instruction et l'envoyer au bureau commercial. Vous devez également joindre une copie de ce document au colis contenant la pompe avant l'expédition.

Mettez la pompe au rebut conformément aux lois et réglementations locales.

Problèmes et solutions

NOTE

En cas de détérioration des performances, de niveau de bruit anormal ou de intervention de la protection du moteur, veuillez prendre contact avec Agilent Contact Center.

Tab. 24 Problèmes et solutions

Problèmes	Cause	Solution
Réduction des performances	Fuite au niveau des conduites d'aspiration ou sur la machine utilisatrice	Éliminer les fuites.
	Manque de lubrification	Contrôler le niveau et l'état de l'huile Rétablir le niveau ou remplacer l'huile.
	Tension d'alimentation trop faible (< 200 Vca)	Raccorder la pompe à un réseau électrique suffisamment puissant et utiliser un câble d'une dimension appropriée (voir paragraphe "Tableau des limites opératoires").
Brouillards d'huile à l'échappement	Éléments déshuileur inefficace	Remplacer l'élément déshuileur
	Haute température due à une huile contaminée	Remplacer l'huile
	Haute température d'exercice due à une température ambiante trop élevée	Réduire la température ambiante en assurant un meilleur renouvellement de l'air
	Élément déshuileur engorgé	Remplacer l'élément déshuileur
La pompe ne démarre pas	Tension d'alimentation hors de la gamme prévue	Raccorder la pompe à un réseau présentant une tension appropriée (voir paragraphe "Tableau des limites opératoires").

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 54 Logo « DEEE »

Pour plus de précisions, veuillez consulter :

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

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-care@agilent.com

vpl-customer-care@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 Instructions for Use

About this manual	127
Validity	127
Definitions and terminology	128
Definition of Caution, Warning and Note	128
Warning Symbols	129
Safety	131
Proper use	131
Improper use	133
Operating limits table	133
Protective equipment	135
MS120 Single Stage Rotary Vane Pump Safety Guidelines	136
Product Description	139
Transport & Storage	140
Lifting	141
Installation	142
Preparation for Installation	142
Accessories supplied with the pump	143
Assembly	143
Location	144
Connection to the machine	145
Discharge air pipe line installation	146
Electrical connection	147
Unpacking	148
Use	151
Start-up and Operation of MS120 Single Stage Rotary Vane Pump	151
Tips for using	152
Gas-ballast	152
Water vapor intake	153
Maintenance	154
Oil change	155
Exhaust filters replacement	157
Ordinary maintenance	159
Pump overhaul	159
How to order spare parts	159
Lubricants	160
Decommissioning	160
Troubleshooting	161
Disposal	162
Service	163

About this manual

Validity

This manual lists the instructions for the users of the MS120 Single Stage Rotary Vane Pump, with particular reference to the notions relating to safety, operation and first level maintenance, limited to maintenance operations for which the user is responsible.

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

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

NOTE

- 1 This manual contains useful information so that all personnel using the MS120 Single Stage Rotary Vane Pump can operate it safely and guarantee perfect efficiency, for its entire life span.
 - 2 Keep this manual, together with all the related publications, in an accessible place known to all operators/maintenance personnel.
-

Definitions and terminology

Definition of Caution, Warning and Note

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

CAUTION

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

WARNING

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

NOTE











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

Warning Symbols

The following is a list of symbols that appear in conjunction with warnings on the MS120 Single Stage Rotary Vane 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.



	Generic hazard
	Dangerous voltages
	Hot surfaces
	Alternate current (AC)
	Protective Ground Terminal
	Fire hazard
	Do not dispose into the environment
	Read the operating instructions
	Inlet port
	Exhaust port

Instructions for Use

	Protective gloves
	Safety shoes
	Protective equipment
	Hazardous substances
	Risk of injury through sharp edges
	Falling objects
	CSA certification
RoHS	Restriction of Hazardous Substances certification
CE	CE certification
	Waste Electrical and Electronic Equipment
UK CA	UK CA Marking

Safety

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

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

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

WARNING



In case of applications where the pump stop or failure can cause damages to people or things, safety measures for the system must be adopted.

Proper use

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

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

The product must only be used by qualified personnel who are able to take the necessary safety measures under conditions that do not cause damage or injury. Any accessories and equipment used with the product must be supplied or approved by Agilent Technologies. Any adjustment or maintenance operation must be performed by a professional technician informed about the risks.

WARNING



Flammable Materials:

To ensure that gases inside the pump do not enter their flammable/combustible range you must obey the instructions and take note of the precautions given below. When potentially flammable or combustible materials are present within the equipment you must:

- Ensure the system is leak tight to safeguard that only the intended gases are passing through the vacuum pump.
 - Ensure that the concentration of any flammable gases or vapors entering the pump (inlet and gas ballast port), inside the pump, and in the exhaust pipeline remain at less than one quarter of the gases' published lower explosive limit (LEL).
 - Ensure that safety systems are in place to maintain the concentration of any flammable gases or vapors entering the pump (inlet and gas ballast port), inside the pump, and in the exhaust pipeline remain at less than one quarter of the gases' published lower explosive limits (LEL) in any system or vacuum pump fault conditions.
-

NOTE

- Before putting the pump back into operation after a breakdown, inspect it and check carefully for any other signs of damage.
 - Use only tools that are in perfect working order and specially designed for the job; use of inappropriate or ineffective tools may cause serious damage.
 - Always check the lubricant and that it is properly distributed through the pump; inadequate lubrication may damage the pump seriously.
 - Give the parts some form of marking as you strip them down to ensure that you reassemble them again in the proper order.
 - Check that there are no scratches or grooves on the machined shafts, in their seats inside the pump or on machine-ground surfaces. Slight scratches and abrasions may be eliminated with very fine emery paper or by a little light grinding.
 - Before putting a group together, always spread a little oil over inner parts and mating surfaces. Replace all seals with original spare parts before reassembling components.
-

Improper use

Agilent Technologies declines all responsibility, deriving from the improper use of the MS120 Single Stage Rotary Vane Pump. Improper use will cause all claims for liability and warranties to be forfeited. Personnel responsible for pump operation and maintenance must be well-trained and must be aware of the accident prevention rules. The accident prevention precautions contained in this section must be continuously respected during operation and maintenance of the pump to avoid damage to operators and to the pump. These precautions are provided in the form of WARNING and CAUTION notes.

WARNING

The pump must not handle:



- liquids or solid substances
- dangerous, explosive or aggressive gases and vapors
- pure oxygen or air mixtures enriched with oxygen.

WARNING



It is prohibited to use the discharge of the pump to create even limited pressures.

WARNING



It is forbidden to install the pump in a potentially explosive environment.

Operating limits table

Tab. 25

Input Voltage (V)	Controller Status
< 180	Power fail (Undervoltage/Overvoltage Error – Win206 bit 5 = 1)
180 – 200	Operative (reduced performance)
200 – 264	Operative (full performance)
> 264	Power fail (Undervoltage/Overvoltage Error – Win206 bit 5 = 1)

Instructions for Use

WARNING



- Death may result from contact with high voltages. Always take extreme care and observe the accident prevention regulations in force.
- Always disconnect the power supply to the pump before maintenance work. Place a special warning sign over the power supply breaker switch: MACHINE UNDERGOING MAINTENANCE - DO NOT POWER ON.
- Failure to provide the pump with an earth connection may cause serious damage to operators. Always ensure that there is an earth connection and that it complies with the standards.
- When cleaning the pump and its component parts, avoid the use of flammable or toxic solvents, such as petrol, benzol, ether or alcohol. The recommendation is to use a soap and water solution, preferably in ultrasound washing machines, taking care to dry all the cleaned parts at temperatures under 100 °C in order to eliminate residual moisture.
- Prolonged overloads or breakdowns may cause the electric motor to overheat, and to release noxious smoke; remove the power immediately as a precaution and do not approach the pump at least until you have provided ventilation to drive out the smoke.
- In case of fire, do not throw water on the pump. Switch the power off and use CO₂ extinguishers.
- Carefully inspect the flanges to ensure that there is no dust, oil, dirt or defects of the mating surfaces, before making the required connections.
- Ensure that all joints and couplings are locked correctly before starting the pump again after repair work.
- Do not wear any objects that may become entangled in the mechanisms and/or act as conductors (chains, bracelets, etc.).
- Ensure that the tools to be used are in perfect working condition and have insulating grips, where necessary. Check that the insulating material of the cables and that the conductors of the test equipment do not show any signs of damage.
- Do not replace the oil immediately after stopping the machine as the oil may still be at high temperature.

NOTE

Use the Retention Spring to secure the mains cable into the IEC320 socket.

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 performed. Furthermore, it must comply with the safety requirements of the legislation in force in the country in which the machine is used.

WARNING

Danger to health by hazardous substances during maintenance or installation.



- Depending on the process peculiarity, vacuum pumps, components or 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 pumps by hand, there is a danger from loads slipping and falling.

- Carry small and mid-size vacuum pumps two-handed.
- 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

- Before servicing the pump or before any mounting/ dismounting action of the pump from the system, wait for the complete standstill of the pump.
 - If necessary, wear protective gloves according to EN 420.
-

MS120 Single Stage Rotary Vane Pump Safety Guidelines

WARNING



Despite of all the precautions adopted when designing the equipment, there are some risk elements that arise during operation and servicing.

WARNING



HOT SURFACES

The pump surfaces may exceed the temperature of 80°C.
Adopt the necessary safeguards to avoid burns due to chance contact.
Before carrying out any maintenance on the pump, be sure the pump is cool.

Instructions for Use

WARNING



HARMFUL SUBSTANCES EMISSIONS

The discharged air contains part of traces of oil mist (maximum residual 2PPM/weight corresponding to 2.4 mg/m³). Check the compatibility with the environment. Make sure a correct air change is allowed otherwise convey the pump discharge outside. A failure or the seals wear can cause an oil leakage. Avoid the dispersion to the ground and the pollution of other materials. In case that any air containing dangerous substances must be pumped down (for example, biological or microbiological agents), make sure to adopt filtering systems before introducing air in the work environment.

CAUTION



DO NOT DISPOSE INTO THE ENVIRONMENT

Used discharged oil from the pump must be disposed in accordance with the regulations in force in the Country of use.

WARNING



HAZARD CAUSED BY VACUUM

Any contact with parts under vacuum can cause injuries. Avoid any contact with the pump inlet port during the pump operation. Introduce air in the inlet circuit before every operation cycle.

WARNING



HAZARD CAUSED BY PRESSURE

The pump tank is pressurized. Do not open the oil filling and discharge plugs during operation.

Instructions for Use

WARNING



Always disconnect the power supply to the pump before starting maintenance work. Place a special warning signs over the power supply breaker switch: **MACHINE UNDERGOING MAINTENANCE - DO NOT POWER ON**. When finished, remove the safety warning.

WARNING



ELECTRIC SAFETY

Some components of the electric equipment are electrically charged during operation. Any contact may cause serious injuries to persons or objects.

WARNING



FIRE HAZARD

The use of the pump in situations unforeseen or not recommended by this manual, as well as lack of correct maintenance, may create high risks for overheating or fire. In case of a fire do not use water to extinguish but use a powder CO₂ extinguisher or other means compatible with the electrical equipment and lubricating oil.

Product Description

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

The suggestions given to the staff engaged in the installation and servicing assumes that the personnel is expert and prepared in facing any problem of servicing, both mechanical and electrical. For any questions or information not included in this manual, please contact our Service Department, always providing: model (Model), serial number (Serial), year of manufacture, stated on the pump name plate.

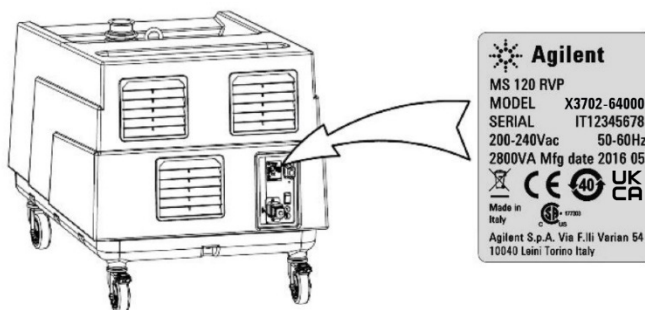


Figure 55 Plate

The MS120 Single Stage Rotary Vane Pumps are single-stage, rotary vane pumps oil sealed, driven by an electric motor controlled by an inverter. These high vacuum pumps are suitable for pumping air and small quantity of water vapor only. Pumping of other types of gas or vapors must be declared in advance to Agilent Technologies that will give the conformity to the specific use.

Instructions for Use

The flanged electric motor is coupled by means of an elastic coupling. The cooling is ensured by a powerful centrifugal fan.

At the pump inlet there is a mesh filter in order to protect it from solid parts having diameter larger than 4.7 mm. Furthermore, an integrated non-return valve prevents the oil coming back and the return of air in the chamber to be pumped down during the stop phase.

In the tank there is a system of oil vapor separation from the discharged air (maximum residual 2PPM/weight corresponding to 2.4 mg/m³).

The separated oil is recovered automatically by the pump.

The gas ballast valves prevent condensation inside the pump when pumping down small quantity of vapor (see paragraph "Gas ballast valves").

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

Transport & Storage

The pumps must be stored or transported without oil.

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

- Temperature: from -20 °C to 70 °C
- Relative humidity: 0 to 95% (non-condensing)

Lifting

The orientation of the packed components must correspond to the instructions given by the pictograms on the external covering of the packaging.

For unloading use a lifting equipment suitable for the pump weight. For lifting the packaging and the pump, refer to the following figures.

Total weight of the pack, including the pump, is max. 120 kg. Weight of the pump only is 100 kg.

Table 26 Packing size

Quote	L	B	H
Dimension	920 mm	610 mm	790 mm

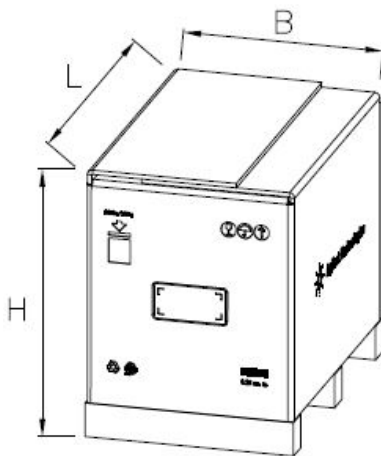


Figure 56 Pump packaged

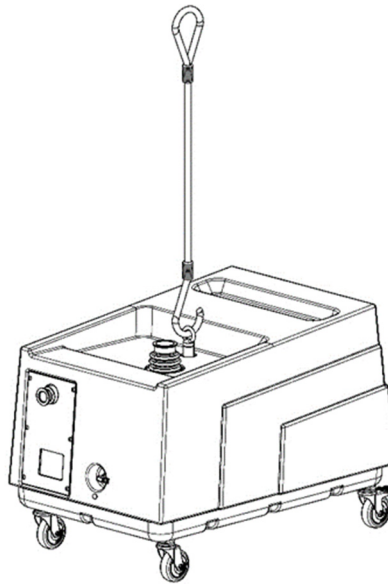


Figure 57 Pump only

Installation

Preparation for Installation

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

WARNING



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

Instructions for Use

Accessories supplied with the pump

The pump is delivered with the following accessories required for commissioning and maintenance:

- Oil Tank
- Nozzle cap load oil
- Oil drain pipe
- I/O Mating Connector
- Spare fuses.

Assembly

Remove inlet and exhaust plastic caps.

The pump is supplied ready for use, there are not other parts to be assembled.

Pump is without oil, see paragraph "Oil change" for the instructions on how to fill in the oil.

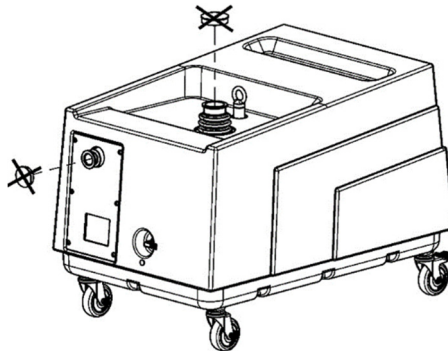


Figure 58

Instructions for Use

Location

WARNING



The pump

- may be blocked from movement by engaging the brakes on the wheels, on a horizontal plane.
- must be accessible for correct and easy maintenance, by respecting the minimum distances from possible obstructions (see following figure). It should be accessible to suitable lifting equipment.
- it must be protected from jets or sprays of water which could cause electrical shorts and / or to the electrocution of operators near to the machine.

Table 27 Minimum distance

Quote	A	B	C	D
Dimension	150 mm	200 mm	150 mm	500 mm

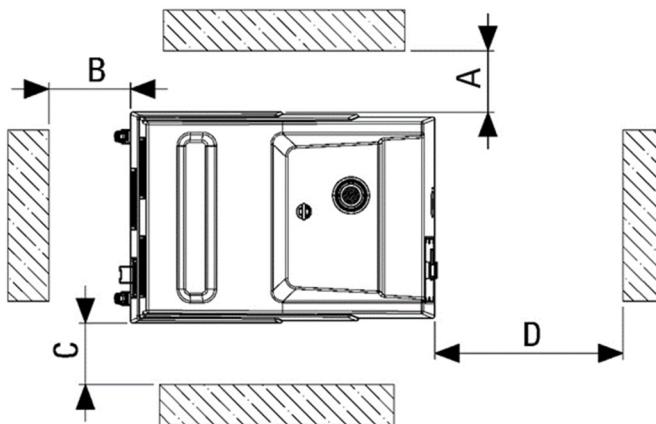


Figure 59

CAUTION

The pump must not be installed outdoors and must always be protected from the weather.

Avoid warm air coming from the exhaust or the cooling fans causing discomfort to the personnel.

Do not install the pump in a dusty area or where other materials may block or cover the cooling surfaces quickly.

Instructions for Use

Connection to the machine

The connection to the chamber to be pumped down must be carried out by means of pipes of the same diameter as the inlet port.

Pipe weights and expansions, if any, must not rest on the pump. It is advisable to make the final connection to the pump inlet port with flexible pipes or fittings.

It is important that all the pipes and the different fittings are tight.

Very long or small diameter pipes decrease the pump performances.

NOTE

This symbol identifies the inlet port.

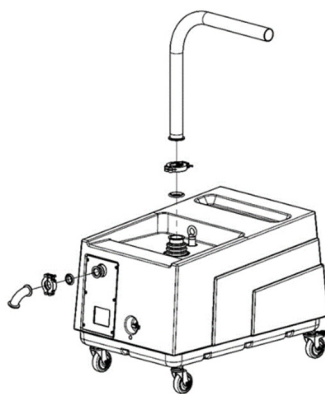


Figure 60

Instructions for Use

Discharge air pipe line installation

If the pump has been installed in a room with poor air exchange, it is possible to pipe the pump discharge air to other rooms or outside.

Use pipes with the same diameter as the tank discharge port with a maximum length of 15 m.

For longer pipes increase pipe diameter. Pipe weights must not rest on the pump. In the final length use flexible pipes or pipe fittings.

WARNING



This pipe must be descending, to avoid the condensate going back to the tank

CAUTION

Do not connect ball valves to this pipeline. Do not clog the outlet.

NOTE

This symbol identifies the exhaust port.



Electrical connection

Check the main voltage and frequency in use to correspond to the data stamped on the pump name plate.

Make sure the grounding is correctly done.

WARNING

Make sure the power switch is set to off (0) in order to avoid unintended initialization of the pump.



Perform the electrical connection by inserting the plug of the power cord into the outlet of the pump and block it with the supplied spring (see following figure).

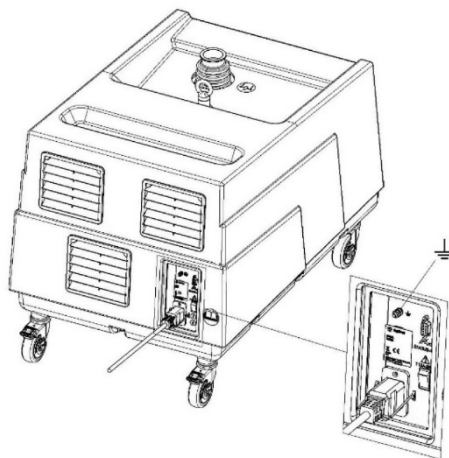


Figure 61

Instructions for Use

Unpacking

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

Do not dispose of the packing materials in an unauthorized manner. The material is 100% recyclable and complies with EEC Directive 85/399.

NOTE

Normal exposure to the environment cannot damage the pump. Nevertheless, it is advisable to keep it closed until it is installed in the system, thus preventing any form of pollution by dust.

Remove the strips that close the carton box on the pallet. Remove the carton box.

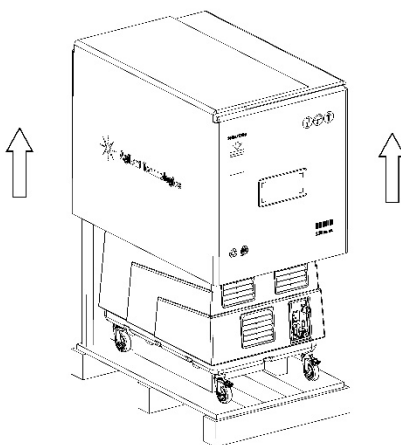


Figure 62

Instructions for Use

Remove 8 screws which close the clamps and then remove 4 clamps.

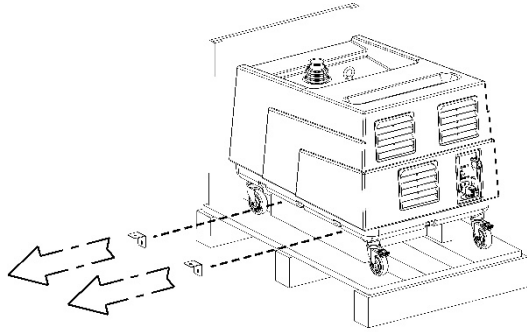


Figure 63

Unlock the brakes of the wheels.

Fasten the drop ramp with one screw as shown in the following figure.

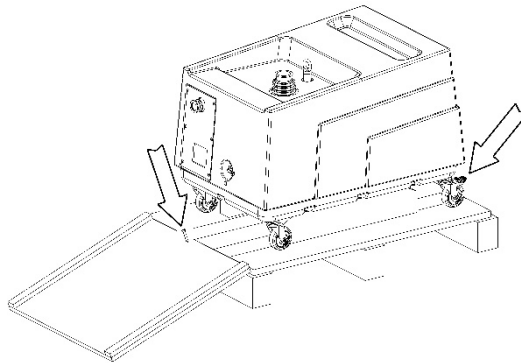


Figure 64

Instructions for Use

Push the pump behind to let it slide down from the pallet.

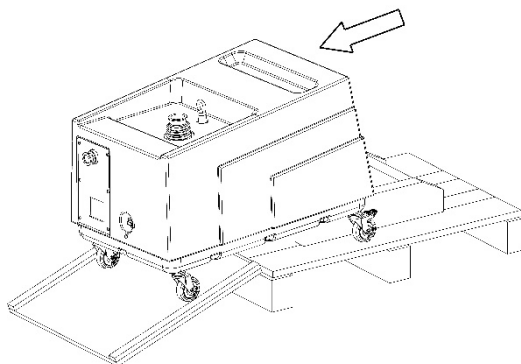


Figure 65

Use

These vacuum pumps have been designed to handle air and small quantity of water vapor only. Handling of other types of gas or vapors must be declared in advance to Agilent Technologies that will give the conformity to the specific use.

They are suitable to evacuate closed systems or to operate at a constant vacuum within the following vacuum range: 0.07 - 15 mbar (abs.).

The ambient temperature and the inlet temperatures must be included between 12° and 35 °C.

Start-up and Operation of MS120 Single Stage Rotary Vane Pump

Do not install or use the pump in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk.

During operation, the following environmental conditions must be respected:

- temperature: from +12 °C to +35 °C
- relative humidity: 0 – 95% (non-condensing).

CAUTION

Before starting the pump, fill up with lubricating oil as the pump is delivered empty.

Further information for the oil refurbishment and the choice of the suitable oil, refer to paragraph "Oil change".

NOTE

It is important that the oil level stays within the range MIN and MAX that is indicated by the level gauge on the side of the pump.

Carry out the first filling up through the plug (E) up to the half of the sight glass (F) and close the plug (E) (see following figure).

Instructions for Use

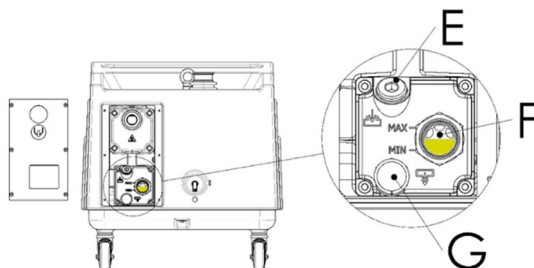


Figure 66

Start the pump and take it to the maximum vacuum level for at least 2 minutes. Stop the pump, check again the oil level and add the lacking oil, if necessary, in order to get the correct oil level.

Tips for using

When the room temperature is lower than 18°C, it is a good practice to let the pump operate at the ultimate pressure (Inlet port closed, without load) for about 15 minutes.

During this period the pump may not reach the stated pressure limits.

WARNING



Avoid operating the pump with inlet port vented to atmospheric pressure.

It is recommended not to exceed 5 pump starts per hour.

Gas-ballast

The pumps of the MS series are equipped with gas ballast valve and with a manual valve which permits its use or its exclusion.

The limit pressure of the pump depends on this selection. For a correct use, please refer to paragraph "Water vapor suction".

Instructions for Use

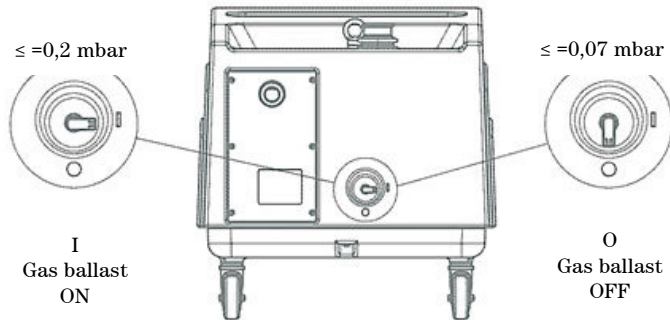


Figure 67

Water vapor intake

In order to pump down any water vapor, the following operations are recommended:

- Before introducing water vapor, let the pump run for thirty minutes at the maximum vacuum in order to bring the pump temperature to its operating value.
- At the end of the working cycle, in case of further presence of condensate in the oil, let the pump run for at least thirty minutes at the maximum vacuum with Gas-ballast open.
- It is advisable to carry out this operation before stopping the pump for a long time. The gas ballast valve will allow the elimination of water condensate from the lubricating oil.

See "Technical Information" for the insertion procedure or exclusion of the gas ballast.

Maintenance

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

In addition to the requirements listed below, please refer to the paragraph "Safety".

WARNING



Introduce air through the inlet port before every maintenance operation.

WARNING



If you have to perform maintenance on the pump after a considerable time in operation, leave it to cool as temperature of the outer surface may be in excess of 80 °C.

WARNING



Always disconnect the power supply to the pump before starting maintenance work. Place a special warning signs over the power supply breaker switch: MACHINE UNDERGOING MAINTENANCE - DO NOT POWER ON. When finished, remove the safety warning.

In order to keep the pump operating at a high efficiency level, it is mandatory to follow all periodical service points listed in the table below.

However, more frequent service operations may be necessary depending on what the pump is used for (suction of condensable vapors, suction polluting substances).

For such cases, only direct experience can indicate the correct service frequency needed. The exhausted oil and the replaced spare parts must be considered as special waste products and handled according to the local regulations in the Country of use.

Instructions for Use

Table 28 Service operations

Servicing frequency	Description of the operation	Authorized personnel
24 Hours/every day	Check oil level before starting.	Operator
100 Hours / every week	Clean with soft cloth surfaces of the pump. If necessary, you can use a neutral detergent.	Operator
	Change the lubricating oil.	Skilled Worker
9000 Hours / every year	Change the lubricating oil.	Skilled Worker
	Replace the exhaust filter.	Skilled Worker
14000 Hours / every 2 years	Check the electrical connections.	Skilled Worker
30000 Hours / every 5 years	Pump overhaul.	Customer Service

Oil change

It is recommended to wait a hour and a half after turning off the pump before beginning the process of changing the lubricant oil to allow the internal temperature of the pump and oil to cool sufficiently.

WARNING



Use protective gloves to avoid injury caused by heat.

Please refer to Figure 68 Unscrew the oil filling plug (E) and the discharge plug (G) only after having placed underneath the pump tank a suitable container suitable container (see Figure 69 proper size and shape) for collecting the total quantity of oil.

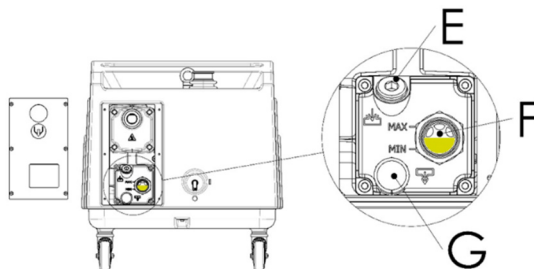


Figure 68

Instructions for Use

Screwing the oil drain pipe on the discharge (G), the lubricant will start to drain. It is recommended to tilt the pump forwards by raising the back 40 mm to ensure the complete draining of the lubricant, remember to brake the wheels to prevent the pump from falling. Once the oil in tank has completely been discharged from the tank, reassemble both plugs ("E" and "G") and let the pump run under vacuum for about one minute, so that the lubricating line gets emptied and any oil residual keeps inside the pump. Then remove the plugs and discharge the rest of the oil.

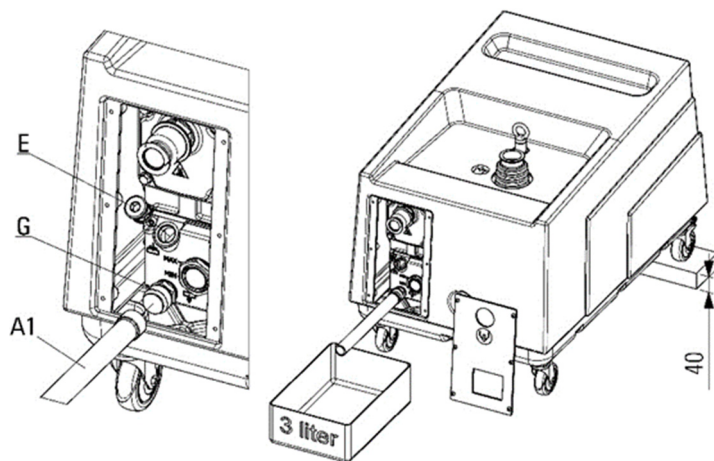


Figure 69

If the oil is polluted or if some water is in the oil, clean the pump by letting it run with a suitable quantity of oil (up to the minimum level shown on the tank) at maximum vacuum level for at least 5 minutes.

Drain the oil again.

Instructions for Use

Fill in the oil following these steps (see Figure 70):

- 1 Unscrew the oil charge plug (E).
- 2 Screw the oil nozzle cap (A2) on the oil bottle (A1).
- 3 Proceed with the new oil filling inserting the oil nozzle cap in the oil charge hole (see "Start-up and Operation of MS120 Single Stage Rotary Vane Pump" and "Lubricants").

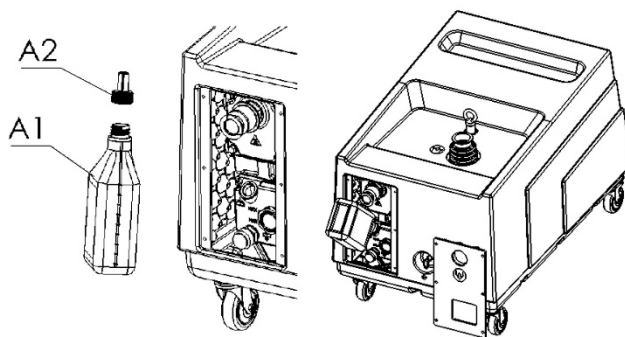


Figure 70

Exhaust filters replacement

Very dirty exhaust filters may cause a considerable pump temperature increase and in extreme cases oil lubricant spontaneous ignition.

Maximum allowed pressure in the tank is 0.6 bar measured at the maximum capacity (when the pump is working with the inlet open to atmospheric pressure). If a pressure gauge (optional) has been fitted to the oil filling hole (Figure 70) check the exhaust filter blockage with the pump warm.

Instructions for Use

For to replace the exhaust filter (see Figure 71):

- 1** Remove the cover (pos. C1) by unscrewing its screws (C2).
- 2** Remove the tank cover (pos. B1) unscrewing its screws (pos. B2).
- 3** Extract and replace the exhaust filter (pos. B4) and its O Ring (pos. B5).
- 4** Reassemble the fixing cartridge disk (pos. B1).
- 5** If necessary, replace the gasket (pos. B3).
- 6** Reassemble the discharge cover (pos. C1).

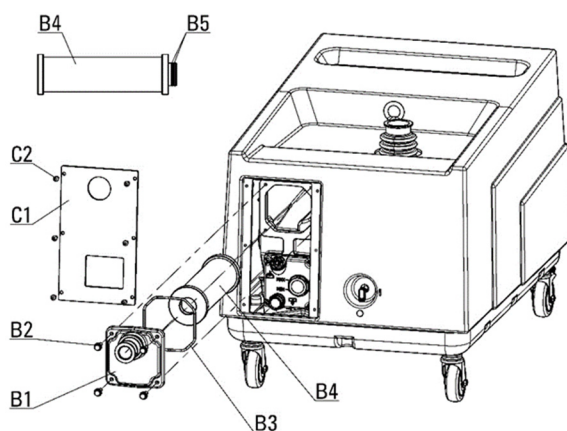


Figure 71

Instructions for Use

Ordinary maintenance

The spares necessary for the ordinary maintenance are shown in the table below. Use only genuine spare parts. When placing the order, please specify the P/N.

Table 29 Normal servicing Spares

Part Number	Description
X3760-64005	Rotary Vane Fluid AVF 60 Gold 1lt
X3702-68201	MS 120 Exhaust Filter Cartridge

Pump overhaul

This operation can be performed only by the Agilent Contact Center.

The overhaul consists of a complete disassembly, cleaning of all components as well as replacement of parts that are subject to wear (pump bushings, vanes and gaskets). Use only genuine spare parts.

When placing the order, please specify the P/N.

Table 30 Pump overhaul spares

Part Number	Description
X3702-68202	MS 120 Major Maintenance Kit

How to order spare parts

When ordering spare parts, always state the pump model, serial number, year of production, Part number, description and quantity needed.

Lubricants

In the following table are reported the oil recommended for generic use.

Table 31 Lubricants

Ambient temperature	Viscosity	Agilent Technologies oil
12 – 35 °C	ISO 53	AVF 60 Gold

In case of change of the type of lubricant, it is necessary to proceed with a pump by letting it run at maximum vacuum for at least 5 minutes.

For this operation use a quantity of oil such as to reach the minimum level indicated on the tank.

Discharge again the oil and proceed with new oil charge.

Decommissioning

Drain the oil from the pump prior to the removal. If the oil is polluted, flush the pump with fresh oil (see "Oil change").

Drain the oil from the tank, plug the inlet and the discharge ports and store the pump without oil.

NOTE

Before returning the pump to the factory for repair, the "Request for Return" sheet attached to this instruction manual must be completed and sent to the local sales office. A copy of the sheet must be inserted in the pump package before shipping.

If a pump is to be discarded, it must be disposed of in compliance with local laws and regulations.

Troubleshooting

NOTE

In case of drop in performance, anomalous noise or motor protection tripping, please get in contact with Agilent Contact Center.

Table 32 Troubleshooting

Trouble	Cause	Remedy
Reduction in performances	Inlet pipes or machine are leaking	Identify leaking point and seal it
	No lubrication	Check oil level and oil conditions. Fill with oil to the right oil level or change the oil
	Supply Voltage too low (< 200 Vac)	Connect the pump to a mains with the correct voltage and use a cable with the appropriate size (see Operating limits table)
Discharge oil mist	Inefficient exhaust filters	Replace exhaust filters
	High temperature due to polluted oil	Oil change
	High operating temperature due to high ambient temperature	Decrease room temperature by allowing a better air exchange
	Blocked exhaust filters	Replace exhaust filters
Pump doesn't start	Supply Voltage out of the range	Connect the pump to a mains with the correct voltage (see Operating limits table)

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 72 Logo "WEEE"

For more information refer to:

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

Service

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

vpt-customer@agilent.com

vpl-customer@agilent.com

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

5 Technical Information

Description of the MS120 Single Stage Rotary Vane Pump	165
Vacuum Seals	166
Technical Specifications	167
MS120 Single Stage Rotary Vane Pump Outline Drawing	169
Connection to the Electric Supply	171
Starting and Running the Pump	172
Stopping the Pump	172
User Interface	173
Input/Output Communications	174
J1 – I/O: Remote I/O signals + Serial communication	174
Serial Communication Connection Examples	176
How to connect the I/O port to an external cable	177
Serial Communication Protocol	178
Communication Format	178
Communication Protocol	178
Examples	181
Window Meanings	182
Pump Status	184
Accessories and Spare Parts	185

Description of the MS120 Single Stage Rotary Vane Pump

The MS120 Single Stage Rotary Vane Pump model is: **X3702-64000**.

The MS120 Single Stage Rotary Vane Pumps is a rotary vane lubricated vacuum pump, with oil recirculation.

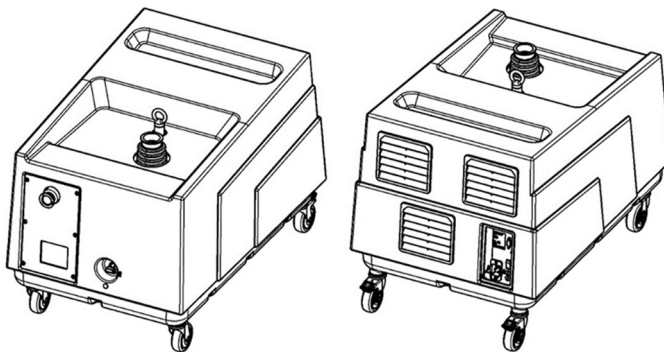


Figure 73

This vacuum pump is suitable for pumping only air and small quantity of water vapor. Pumping of other types of gas or vapors must be declared in advance to Agilent Technologies that will give the conformity to the specific use.

The main features are:

- all parts in direct contact with the fluid pumped are free of copper alloys;
- all materials are carefully selected to provide extended life;
- due to its design features and low number of gaskets, the pump requires little maintenance, disassembly and reassemble are easy and require minimal time.

Technical Information

The oil guarantees perfect sealing of the discharge valves, enters the pump to ensure lubrication and sealing of the parts inside, facilitates heat dissipation and reduces pump noise.

The pump is equipped with a special anti-suckback device which automatically isolates the vacuum system when the pump stops. This avoids rises in pressure or oil flow in the vacuum system while air is allowed back into the stator chambers.

The air entering the pump after the anti-suckback device has closed prevents the oil in the casing from filling the stator chambers.

Vacuum Seals

A special feature of this pump is the low number of gaskets that are employed. The seals in the circuit are obtained by means of FKM gaskets. Sealing of the rotor shaft is guaranteed by a rotating gasket with dust-guard lip. The suction flange and duct are sealed by mean of OR gaskets.

Technical Specifications

Table 33 Technical Specifications

	Units	Value
Nominal speed* (at 1500 rpm)	m ³ /h	156
	cfm	92
Effective capacity at 5 mbar abs. (3.75 Torr)	m ³ /h	≥ 120
	cfm	≥ 71
Ultimate total pressure with gas ballast valve open*	mbar	≤ 0,2
	Torr	≤ 0.150
Ultimate total pressure with gas ballast valve closed*	mbar	≤ 0,07
	Torr	≤ 0.0525
Nominal Input Voltage and frequency (~1)	V	200 – 240
	Hz	50 - 60
Max input power	VA	3000
Factory Set Frequency	Hz	50 (factory setting)
Touch current	mA	3.0 maximum
Maximum Ambient temperature	°C	35
Sound pressure level (without gas ballast valve) **	dB(A)	60 ±1
Water vapor tolerance	mbar	15
	Pa	1500
Water vapor pumping capacity	kg/h	1
	Lbs/h	2.2
Operating temperature range	°C	+12 to + 35
	°F	+54 to +95
Ground stud max current	Amp	15

Technical Information

Table 33 Technical Specifications (continued)

	Units	Value
Inlet flange		ISO KF DN40
		ISO KF NW40
Outlet flange		ISO KF DN25
		ISO KF NW25
Oil capacity	l	1,8
	Gal US	0.47
Type lubricating fluid		AVF 60 Gold
Total weight	kg	100
	lbs	220
IP Value		20
Installation category		II
Pollution degree		2
Max altitude	m	2000

* According to PNEURO standard 6602

** At ultimate total pressure, 20 °C (68 °F) room temperature

CE marks

EN61326-1 2013 Electrical equipment for measurement control and laboratory use EMC Requirements

Emission: EN55011 Class B; EN61000-3-2; EN61000-3-3

Immunity Industrial Levels:

EN61000-4-2 EN61000-4-8

EN61000-4-3 EN61000-4-11

EN61000-4-4 EN61000-4-12

EN61000-4-5 EN61010-1

EN61000-4-6

EN61010-1, 2010 (Safety requirements for electrical equipment for measurement control laboratory use)

MS120 Single Stage Rotary Vane Pump Outline Drawing

The following figures show the pumps layout and dimensions:

Outline Drawings

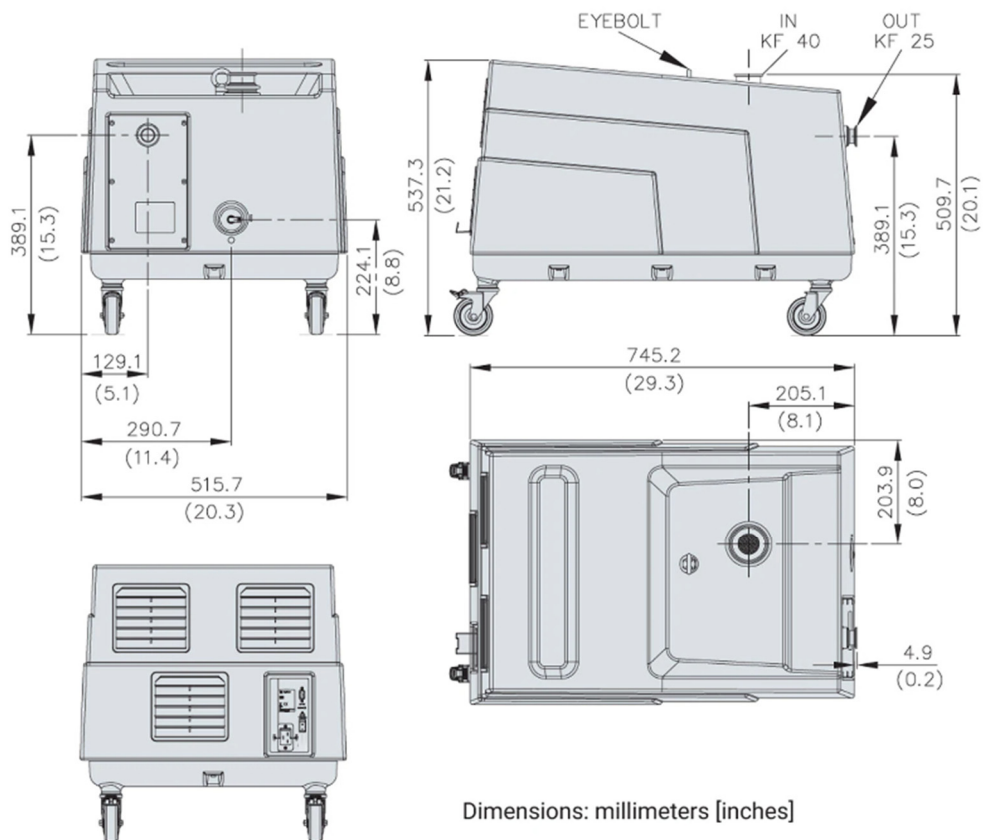


Figure 74 MS120 Single Stage Rotary Vane Pumps dimensions

Technical Information

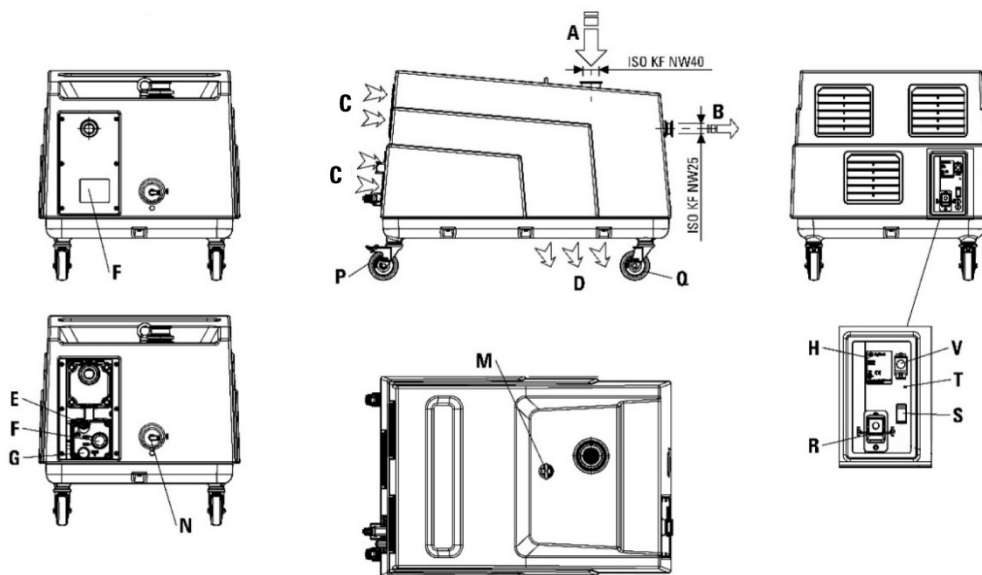


Figure 75 MS120 Single Stage Rotary Vane Pumps layout

Table 34 Identification of components

Reference (previous figure)	Component
A	Inlet
B	Exhaust
C	Cooling air inlet
D	Cooling air outlet
E	Oil filling plug
F	Oil level sight glass
G	Oil drain (with extension)
H	Pump name plate
M	Lifting eyebolt
N	Gas ballast valve (ON-OFF)
P	Swivel caster with brake
Q	Swivel caster
R	Mains
S	ON-OFF Switch
T	Status Led
V	Serial port

Pumping speed curve at 1500rpm

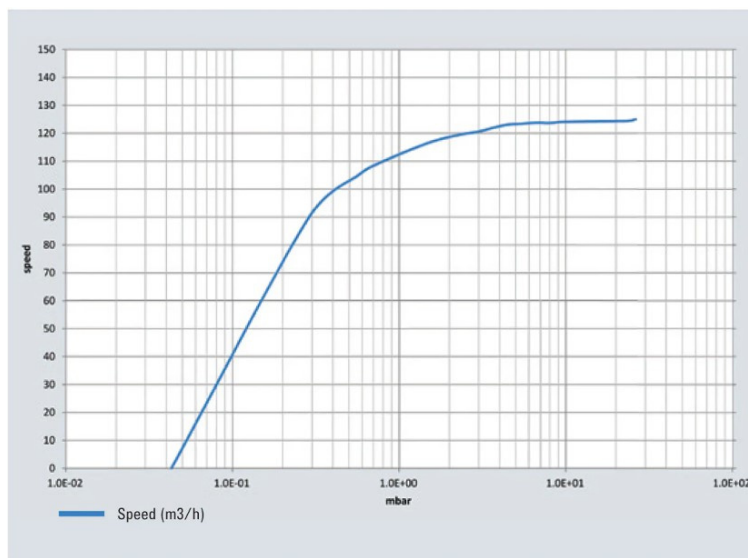


Figure 76 MS120 Single Stage Rotary Vane Pumps Pumping Speed

Connection to the Electric Supply

CAUTION

It is recommended to connect the pump to the power supply through a dedicated safety switch on the main electrical panel of the installation, or in proximity of the power supply connection point.

WARNING



The pump must be installed in a way that allows an easy interruption of the line voltage.

Power supply cord: The correct cable for electrical wiring is a three wires (Ph+N+Earth) cable. The wire section has to be at least 2 mm² (AWG14).

Starting and Running the Pump

The pump can be operated in REMOTE mode through the Remote I/O signals or in SERIAL mode through the RS232 or RS485 Serial communication.

The default configuration is REMOTE mode.

NOTE

To allow the pump starting you have to wire properly the START pins by connecting the mating connector provided with the pump or by connecting pin 7 to pin 9 on J1 – I/O connector.

Stopping the Pump

There are no special procedures for switching the pump off; it needs only to be disconnected from the electric power by means of the bipolar switch. When the pump is stopped, the anti-suckback device makes it possible to maintain vacuum in the vessel connected on the inlet flange of the pump if the tank is hermetically sealed.

User Interface

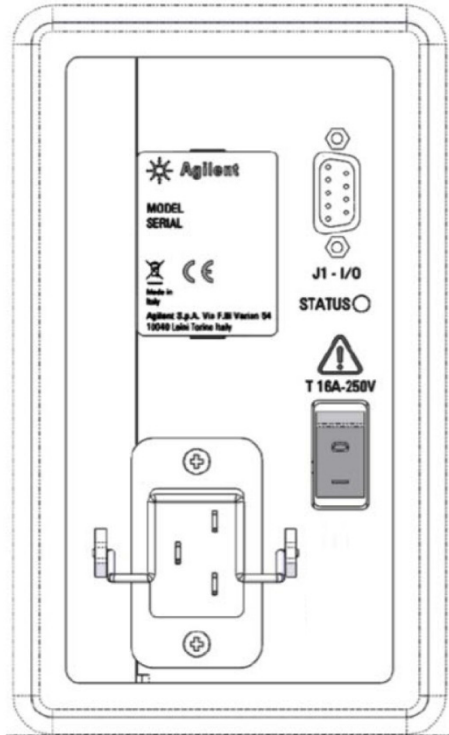


Figure 77

NOTE

Use the Retention Spring to secure the mains cable into the IEC320 socket.

Input/Output Communications

J1 – I/O: Remote I/O signals + Serial communication

The connector J1 (D-SUB 9 pins Female connector type) hosts both Remote I/O signals and RS232/485 communication ports.

The RS232 communication port has standard pin out. It can be used with a standard of the shelf RS232 cable.

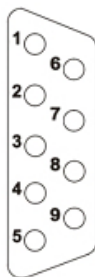


Figure 78

Pin assignment

Table 35 Pin assignment

Pin N.	Signal Name	IN / OUT
1	Status (N.O. relay contact)	out
2	TX (RS232)	
3	RX (RS232)	
4	Status (N.O. relay contact)	out
5	Ground	
6	A+ (RS485)	
7	START/STOP (Interlock in Serial Mode)	in
8	B- (RS485)	
9	+24V (50 mA max)	out

Technical Information

- Interlock/Status: N.O. Solid State Relay contact – It is closed as soon as the rotational frequency exceed the threshold defined with Win102 (plus histeresys defined by Win105 as % of Win102).
- START/STOP: If the inverter is managed by remote signals, it commands the pump starting. This is a 0 / 24V digital input active high. To activate the input (and starting the pump) it should be tied to 24Vdc (external 24Vdc or pin 9 on the I/O connector), a non active state could be obtained either leaving the input not connected or connecting it to GND.

More in detail:

- Active state: $V_{input} \geq 12V$ or $I_{sink} \geq 1.6\text{ mA}$

- Non Active state: $V_{input} \leq 4V$ or $I_{sink} \leq 0.1\text{ mA}$

Here below there is a typical connection of the START/STOP input to an external system output.

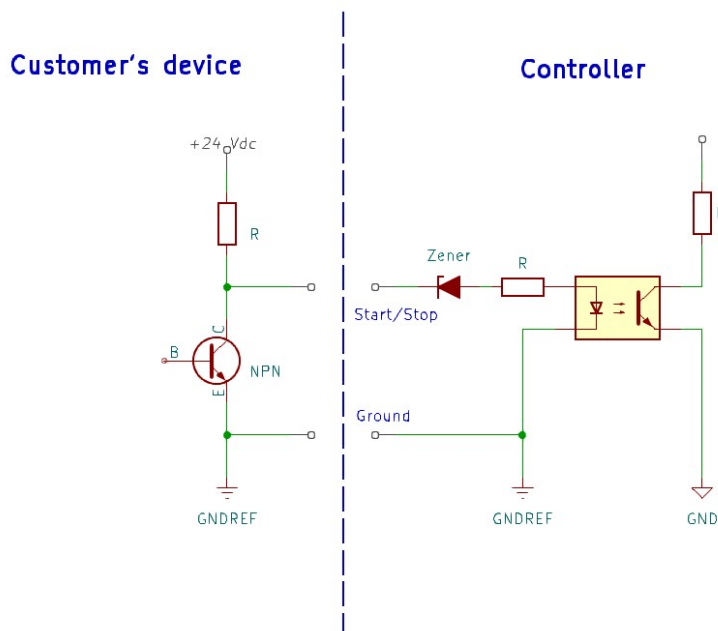


Figure 79

NOTE

The pump operation is I/O signals dependent so you must always plug-in the provided mating connector to start the pump.

Technical Information

Serial Communication Connection Examples

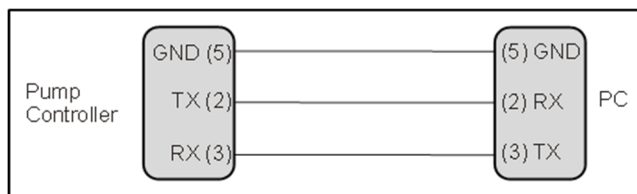


Figure 80

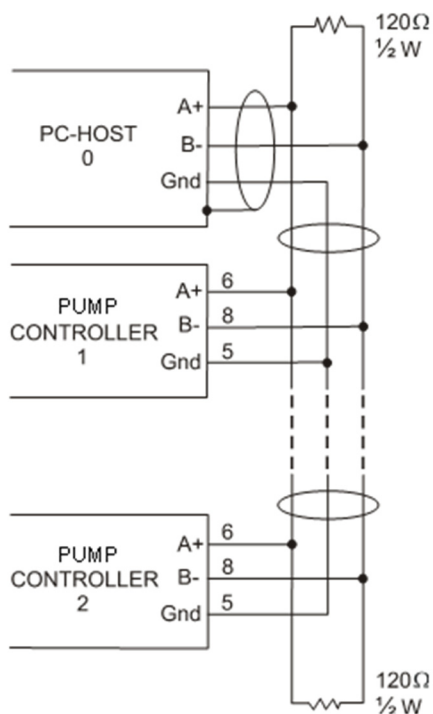


Figure 81

The communication port mating connector is supplied with the unit. See the host computer's instruction manual for further details on serial port connections.

Technical Information

How to connect the I/O port to an external cable

The following picture shows the right procedure to connect a cable to the I/O port connector.

A shielded cable of 30 m maximum length has to be utilized for both serial and I/O port connection.

NOTE

Take care to have a good contact (soldered) between the metallic connector case and the external shield of the cable. Moreover, this connection has to be assured at least on the controller side.

In this way, you will be sure to reduce the influence of the external noise and to accomplish the EMC requests.

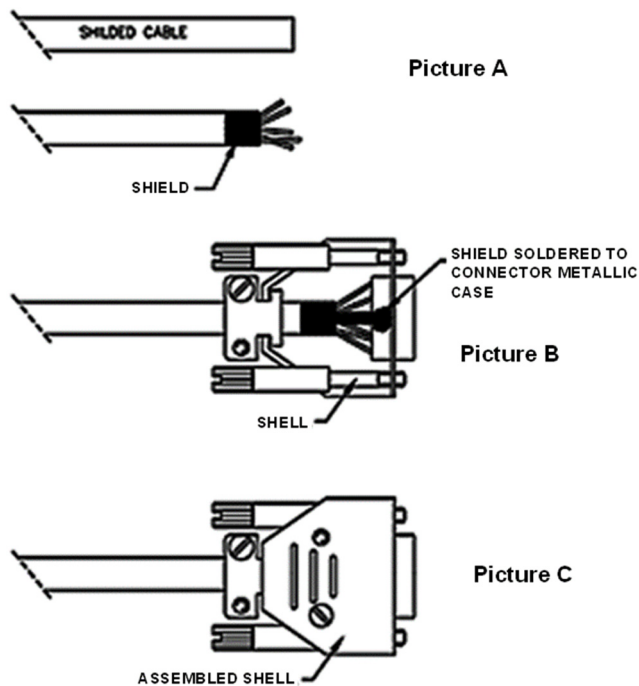


Figure 82

Technical Information

Serial Communication Protocol

Both RS-232 and RS-485 interfaces are available on the J1 I/O connector.

Communication Format

- 8 data bit
- no parity
- 1 stop bit
- baud rate: 600/1200/2400/4800/9600/19200/38400 programmable

Communication Protocol

The communication protocol is a MASTER/SLAVE type where:

- Host = MASTER
- Controller = SLAVE

The communication is performed in the following way:

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

The MESSAGE is a string with the following format:

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

Where:

NOTE

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

Technical Information

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

Table 36

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

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

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

NOTE

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

Technical Information

The controller can answer with the following response types:

Table 37

Type	Length	Value	Description
Logic	1 byte	-	after a read instruction of a logic window.
Numeric	6 bytes	-	after a read instruction of a numeric window.
Alphanumeric	Variable 1÷14 bytes	-	after a read instruction of an alphanumeric window.
ACK	1 byte	(0x6)	the command execution has been successfully completed.
NACK	1 byte	(0x15)	the command execution has been failed.
Unknown Window	1 byte	(0x32)	the specified window in the command is not a valid window.
Data Type Error	1 byte	(0x33)	the data type specified in the command (Logic, Numeric or Alphanumeric) is not accorded with the specified window.
Out of Range	1 byte	(0x34)	the value expressed during a write command is out of the range value of the specified window.
Win Disabled	1 byte	(0x35)	the specified window is Read Only or temporarily disabled (for example you can't write the Soft Start when the Pump is running).

Examples

Command: START

Source: PC

Destination: Pump

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

Source: Pump

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

Command: STOP

Source: PC

Destination: Pump

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

Source: Pump

Destination: PC

02	80	06	03	38	35
STX	ADDR	ACK	ETX	CRC	

Command: READ PUMP STATUS

Source: PC

Destination: Pump (with address = 3)

02	83	3	3	3	30	03	3	37
ST	ADD	WINDOW			R	ET	CRC	
X	R				D	X		

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

Destination: PC

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

Window Meanings

The following table describes the windows available to the user. The columns R/W indicates if the window is read-only (R) or if it can also be modified (RW).

The windows default values are shown in bold.

Table 38 Window Meaning

WIN	TYPE	R/W	DESCRIPTION
000	L	RW	START/STOP (1= START; 0= STOP)
008	L	RW	REMOTE/SERIAL CONFIGURATION (1= Remote; 0= Serial)
102	N	RW	REMOTE I/O SET POINT VALUE [Hz] (0 – 127) 40
105	N	RW	REMOTE I/O SET POINT HYSTERESIS [%] (0 – 100) 2
108	N	RW	SERIAL COMMUNICATION BAUD RATE (0-6) [0=600, 1=1200, 2=2400, 3=4200, 4=9600, 5=19200 bit/s]
120	N	RW	ROTATIONAL FREQUENCY SETTING [Hz] (40 - 60) 58
201	N	R	3PHASE VOLTAGE [V] 3phase Voltage supplied to the motor
202	N	R	POWER [W] Electrical power supplied to the motor
203	N	R	DRIVING FREQUENCY [Hz] Reads the actual electrical driving frequency
205	N	R	STATUS (0=Stop; 1=Wait interlock; 2=Start; 3=Autotuning; 5=Normal; 6=Fail)
206	N	R	ERROR CODE: [bit field] Bit 7: Rotor locked Bit 6: Short Circuit Bit 5: Vline Under/Over Voltage Bit 4: Motor Overtemp Bit 3: Line Fault Bit 2: Controller Overtemp Bit 1: Vdc BUS Under Voltage Bit 0: Over Current
211	N	R	CONTROLLLER POWER OUTPUT HEATSINK TEMPERATURE [°C]
216	N	R	CONTROLLER TEMPERATURE [°C] Controller inlet air temperature
222	N	R	CONTROLLLER POWER SUPPLY HEATSINK TEMPERATURE [°C]
234	N	R	BUS VOLTAGE [V] Controller internal DC BUS Voltage
235	N	R	LINE VOLTAGE [Vpk] Supply line peak Voltage

Technical Information

Table 38 Window Meaning (Continued)

WIN	TYPE	R/W	DESCRIPTION
300	N	R	CYCLE TIME [min] It indicates the time of the ongoing cycle (from the last Start) or the time of the last cycle if the pump is in Stop status
301	N	R	CYCLE NUMBER Indicates the total number of Start/Stop cycles carried by the pump
302	N	R	PUMP LIFE [h] Indicates the total hours of pump life in running condition
307	N	R	PUMP STAND-BY [h] Indicates the total hours of pump life in stand-by (pump switched-on but not in running condition)
320	A	R	PUMP MODEL NUMBER
321	A	R	PUMP SPECIAL MODEL NUMBER
322	A	R	PUMP SERIAL NUMBER
406	A	R	SW PROGRAM LISTING CODE & REVISION
407	A	R	SW PARAMETER LISTING CODE & REVISION
503	N	RW	RS485 Address [0 - 31] It defines the RS485 serial communication address
504	L	RW	Serial communication standard [0 = RS232; 1 = RS485]

Pump Status

The Pump Status indicates the operating status of the pump.

The possible Status are the following:

- Stop = the pump is waiting for a START command
- Ramp = the pump has been started and is increasing the rotational frequency in order to reach the target set point
- Normal Operation = the pump is running at the target frequency
- Fail = the pump is not running because has been detected an error condition. The error condition can be discovered reading the W206 value through serial communication interface
- Autotuning = the pump rotational speed is not able to reach (or maintain) the target value due to a too high gas load.

The Pump Status can be detected by looking at the STATUS LED or reading the W205 value through serial communication interface.

Table 39

Led Status	Controller Status
Green blinking slowly (once every 2 seconds)	Stop
Green blinking slowly (twice every 1 second)	Ramp – Autotuning
Green fix	Normal Operation
Red fix	Fail

NOTE

When you switch the controller on an electronic self-test is performed, during such phase you'll see the led Orange. During this phase the pump is not able to be started.

Accessories and Spare Parts

Table 40 Accessories and Spare Parts

Part Number	Description
X3701-63001	MS 120 Noise Abatement System
X3760-64005	Rotary Vane Fluid AVF 60 Gold 1lt
X3702-64001	Power Cord USA Plug-IEC320 15A
X3702-64006	Power Cord UE Plug-IEC320 15A
X3702-64003	Power cord UK PLUG-IEC320 13A
9699883	Serial cable and A-PLUS
X3702-68001	MS 120 Remote I/O Adapter Cable
X3702-68201	MS 120 Exhaust Filter kit
X3702-68202	MS 120 major maintenance kit

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:
Address:
Contact Name:
Tel:
Fax:
Email:

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

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

Health and safety
The product has been exposed to the following substances:
Substances (please refer to MSDS forms)
Trade name
Chemical name
Chemical Symbol
CAS Number
Toxic
Harmful
Corrosive
Reactive
Flammable
Explosive (*)
Radioactive (*)
Biological (*)
Oxidizing
Sensitizer
Other dangerous substances

Goods preparation
If you have replied YES to one of the above questions. Has the product been purged?
If yes, which cleaning agent/method:
Has the product been drained from oil?
I confirm to place this declaration on the outside of the shipping box.

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

Name:
Position:
Date:
Authorized Signature:

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 018 786
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
- Instruction for Use
- Technical information

This information is subject to change without notice.

© Agilent Technologies, Inc. 2024

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



87-900-150-01

Edition 07/24
Issue D.00