



Agilent MS120 Noise Abatement System (P/N X3701-63001)

Installation Guide



Notices

Manual Part Number

87-901-061-01
Edition A.00, July 2025

Copyright

© Agilent Technologies, Inc. 2025

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc

Via F.Ili Varian, 54
10040 Leini (TO)
ITALY

Warranty

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

Technology Licenses

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

Restricted Rights Legend

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	General Information	4
	About this guide	5
	Oil change	9
	Maintenance	9



1 General Information

About this guide	5
Oil change	9
Maintenance	9

General Information

About this guide

This guide explains the installation of the Noise Abatement System (Part Number: X3701-63001) on the MS120 pump (Part Number: G2571-64000).

The purpose of the Noise Abatement System is to reduce the noise emitted by the pump functioning.

The Noise Abatement System consists of:

2 x Noise Abatement Panel, Long

1 x Noise Abatement Panel, Front

1 x Noise Abatement Panel, Back

1 x Noise Abatement Panel Spacer

Requirements in terms of distances from the wall of the MS120 pump with the Noise Abatement System:

Table 1

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

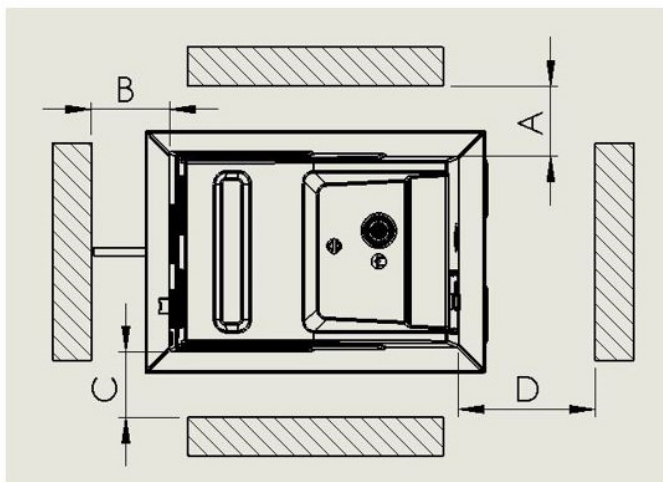


Figure 1 MS120 pump with Noise Abatement System positioning and minimum distances from the wall

General Information

WARNING



Protective gloves must be worn during installation to prevent the risk of cuts or pinching injuries.

- 1 Open the carton box and remove the Noise Abatement System.
- 2 Install the two Noise Abatement Panel, Long under the pump, as shown in figure 2:

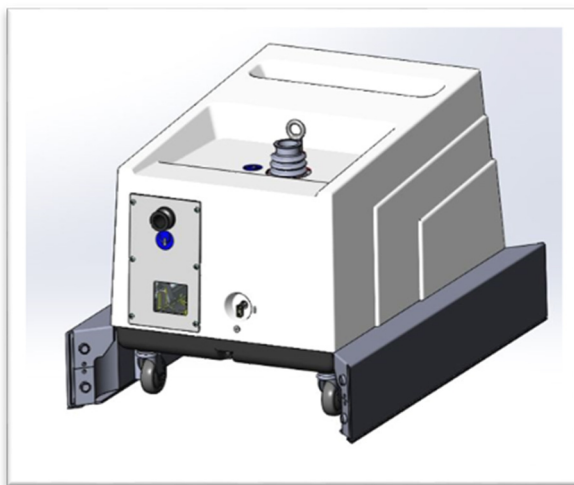


Figure 2 MS120 pump with Noise Abatement Panel, Long

General Information

- 3 Install the Noise Abatement Panel, Front (with the two grids) under the pump, opposite to the pump side with the grids, as shown in figure 3:

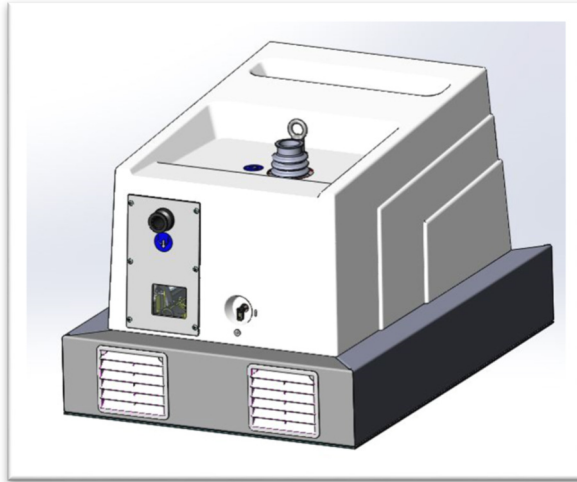


Figure 3 MS120 pump with Noise Abatement Panel, Long and Front

General Information

- 4 Screw the spacer on the Noise Abatement System, as shown in figure 4:

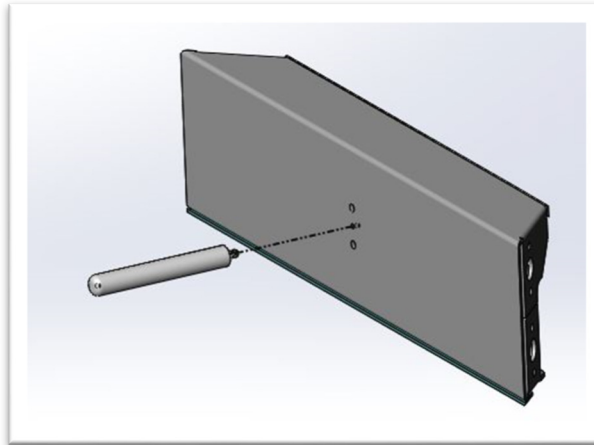


Figure 4 Installation of the spacer on the Noise Abatement Panel, Back

- 5 Install the Noise Abatement Panel, Back under the pump, as shown in figure 5:

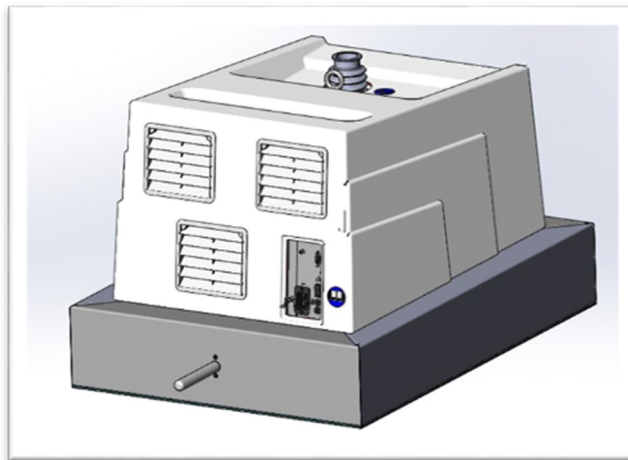


Figure 5 MS120 pump with Noise Abatement System

General Information

- 6 The installation of the system is complete.

The pump can be arranged in its work position as in figure 1: the spacer has been designed to regulate the distance B from the wall.

Oil change

WARNING



Whenever changing the oil it is recommended to use protective gloves to avoid injury caused by heat

When the MS120 Rotary Vane Pump has been operated with the noise abatement system installed it is recommended that you wait about 2 hours after turning off the MS120 before beginning the process of changing the oil to allow the internal temperature of the pump and oil to cool sufficiently.

Please refer to the section “Oil change” of the manual for the instructions on how to discharge the oil from the pump.

Maintenance

Every 100 Hours/every week, clean with soft cloth surfaces the two grids of the Noise Abatement System.

If necessary, you can use a neutral detergent.

Agilent Vacuum Products Division/Sales and Service Offices

United States

Agilent Technologies

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

Netherlands

Agilent Technologies Netherlands B.V.

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

Belgium

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

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

Brazil

Agilent Technologies Brasil

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

China

Agilent Technologies Co. Ltd (China)

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

France

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

Southeast Asia

Agilent Technologies Sales Sdn Bhd

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

India (Sales)

Agilent Technologies India Pvt. Ltd.

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

Italy

Agilent Technologies Italia S.p.A.

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

Japan

Agilent Technologies Japan, Ltd.

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

Singapore

Agilent Technologies Singapore Pte. Ltd.

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

Korea

Agilent Technologies Korea, Ltd.

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

UK and Ireland

Agilent Technologies LDA UK Limited

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

© Agilent Technologies, Inc. 2022
10/2022

India (Service)

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

Taiwan

Agilent Technologies Taiwan Limited

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

Germany and Austria

Agilent Technologies Sales & Services GmbH & Co. KG

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

Mexico

Agilent Technologies

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

Other Countries

Agilent Technologies Italia S.p.A.

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

Customer Support & Service

NORTH AMERICA:

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

EUROPE:

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

PACIFIC RIM:

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

Worldwide Web Site, Catalog and Order On- line:

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

In This Book

The manual describes the following:

- General Information

This information is subject to change without notice.

© Agilent Technologies, Inc. 2025

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



87-901-061-01

Edition 07/25
Issue A.00