



MS120 Noise Abatement System

P/N X3701-63001

Installation Guide

Notices

Manual Part Number

87-901-061-01

Edition 1, October 2019

Copyright

© Agilent Technologies, Inc. 2018

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.lli Varian, 54

10040 Ieini (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.

1 General Information

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:

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

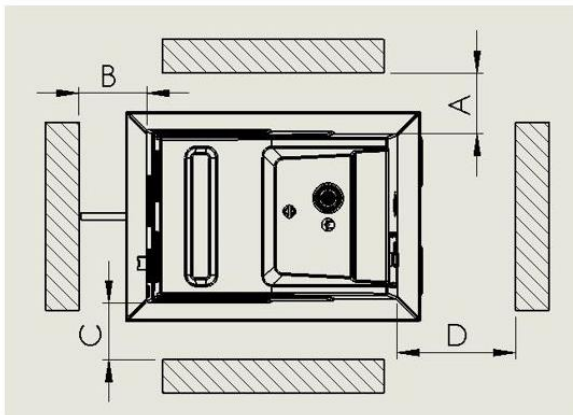


Figure 1. MS120 pump with Noise Abatement System positioning and minimum distances from the wall

- 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:

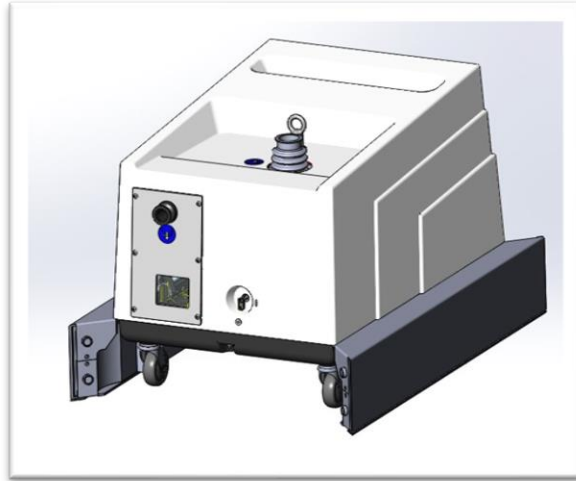


Figure 2. MS120 pump with Noise Abatement Panel, Long

- 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:

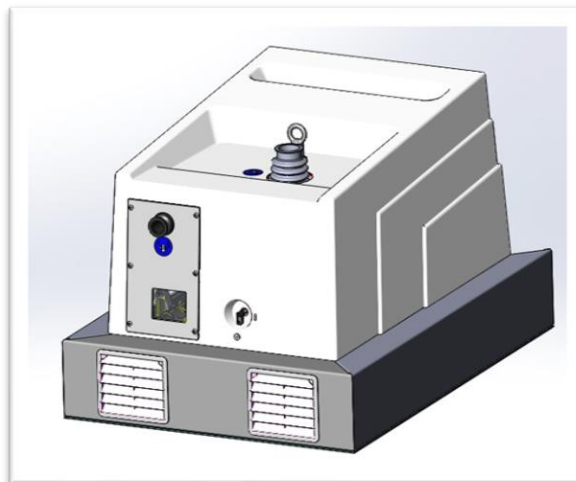


Figure 3. MS120 pump with Noise Abatement Panel, Long and Front

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

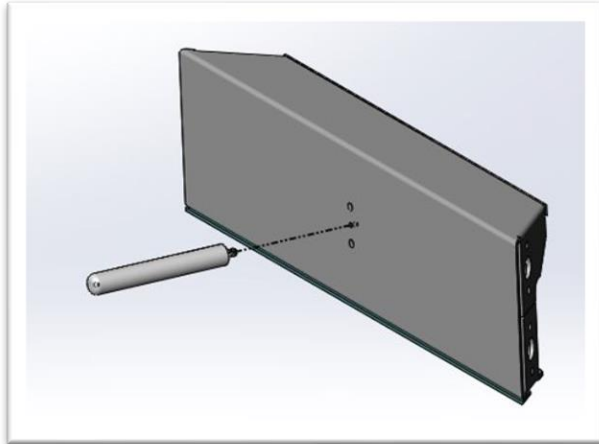


Figure 4. Installation of the spacer on the Noise Abatement Panel, Back

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

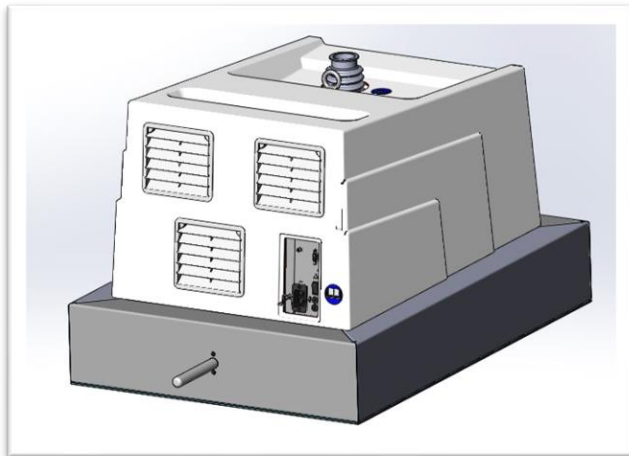


Figure 5. MS120 pump with Noise Abatement System

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.

www.agilent.com

© Agilent Technologies, Inc. 2018

Edition 1/2019

87-901-061-01