

Big Universal Trap

Installation Sheet

Agilent Gas Purifiers

Part number	Fitting	Gas
RMSH-2	1/8-inch	Helium
RMSH-4	1/4-inch	Helium
RMSN-2	1/8-inch	Nitrogen
RMSN-4	1/4-inch	Nitrogen
RMSHY-2	1/8-inch	Hydrogen
RMSHY-4	1/4-inch	Hydrogen

Maximum operating pressure is 250 PSIG.

Maximum operating temperature is 100 °C.

Maximum flow rate is 8 liters per minute.

Installing the Big Universal Trap

Before installing the trap, make sure the system is free of leaks, the GC inlet and MSD vacuum chambers are clean (if using mass spectrometry detection), and the system is generally in good working order.

- 1 Shut down your GC/MSD or reduce the GC heated zone temperatures.
- 2 Set the gas source supply pressure to 10–15 psi and maintain flow in the gas source line before disconnecting it from the inlet of the old trap.
- 3 Remove the protective nut and plug from the **INLET** end of the new trap. **DO NOT** open the plug on the **OUTLET** end.
- 4 Immediately attach the new trap to the gas source tubing using one of the included ferrule sets.

CAUTION

Failure to immediately connect the trap may cause contamination of the adsorbents inside, which will reduce its adsorption capacity and contribute to elevated levels of contaminants observed by the detector. This may result in the need for additional purge time.

- 5 Insert the tubing through the nut and ferrule set until the tubing rests firmly against the shoulder in the fitting.
- 6 Finger-tighten the nut and then use a wrench to tighten the nut 3/4 turn for 1/8-inch tubing and 1-1/4 turn for 1/4-inch tubing (see [Figure 1](#)).



Figure 1 Outlet fitting closed



- 7 Wait for 3 minutes to stabilize the gas pressure inside the new trap. Increase the gas source pressure to 60 psi.
- 8 Open the outlet fitting and purge the trap for 3 minutes (see [Figure 2](#)).



Figure 2 Outlet fitting open

- 9 Adjust the gas source pressure to a normal working level. Connect the outlet fitting of the trap to the instrument tubing as described in steps 5 and 6. Wait for at least 3 minutes to purge out any air that may have entered the instrument tubing line. The trap is now ready for use (see [Figure 3](#)).

NOTE

Most instruments (such as GCs or GC-MSDs) require a waiting time for the system to stabilize and self-clean before use when the gas source or gas trap is changed. This waiting time is often 2 hours or longer. Results from analyses performed too soon after changing the trap may not be reproducible. The longer the wait time the more stable the system will be, thus protecting the instrument and the column when changing gases or the gas trap.

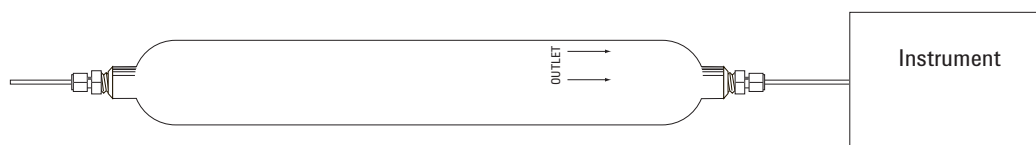


Figure 3 Completed installation

For the latest detailed MSDS information, please visit our website at www.agilent.com/chem/msds.

© Agilent Technologies, Inc. 2002

No part of this document 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.

First edition, May 2002

Printed in USA

Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610 USA



5184-3513