

Agilent G2630 Installing the External Flow Restrictor

Installation Guide



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2002

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.

Document Part Number

G2630-90277

Edition

First edition, April 2002

Replaces G2630-90270

Printed in USA

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

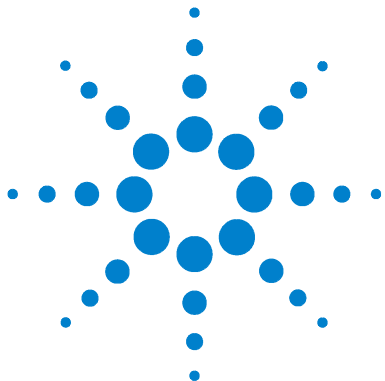
Safety Notices

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.



Installation Guide

WARNING

The external flow restrictor prevents unregulated gas flows in the event of a leak in or at the manifold. If you are using flammable gases (such as hydrogen) with your auxiliary flow module, you must install an external flow restrictor onto each flow module fitting used for a flammable gas.

Two external flow restrictors, part no. G1530-80380, are provided. Each includes a pre-installed 1/8-inch front and back ferrule set with Swagelok nut for connection to your gas line. If needed, order an additional external flow restrictor from Agilent Technologies. See [Figure 1](#).

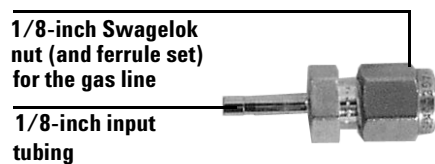


Figure 1 External flow restrictor

Tools required

7/16-inch wrenches

Table 1 Parts you may need to supply

Description	Part no.
Ferrule, 1/8-inch, stainless steel, front (10/pk)	5180-4110
Ferrule, 1/8-inch, stainless steel (10/pk)	5180-4116
Nut, 1/8-inch, stainless steel (10/pk)	5180-4104

Installation

- 1 If flammable gases are already plumbed to the auxiliary flow module, turn them off at their sources and disconnect them from the auxiliary flow module.
- 2 Loosely install a front ferrule, back ferrule, and nut on the auxiliary flow module fitting. See [Figure 2](#).

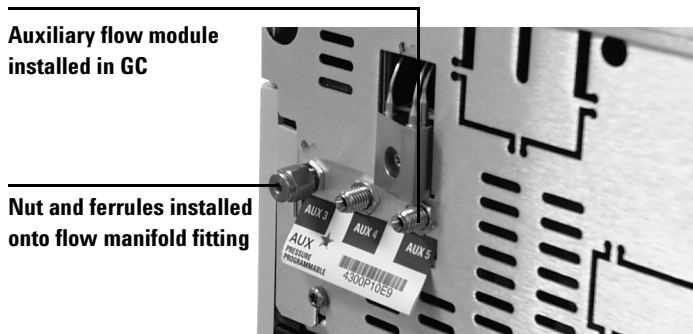


Figure 2 Loosely install nut and ferrule set

- 3 Insert the 1/8-inch tube of the external flow restrictor into the flow module fitting. Using two wrenches, tighten the nut 3/4 turn past finger-tight. See [Figure 3](#).



Figure 3 External flow restrictor installed and ready for connection to gas line

- 4 Slide the gas line into the 1/8-inch nut on the external flow restrictor.

NOTE

If a ferrule set is already installed on the gas line, first remove the 1/8-inch nut and ferrule set from the external flow restrictor.

- 5 Using two wrenches, tighten the gas line connection 3/4 turn past finger-tight.

For more information about making Swagelok connections, refer to your 6890 User Information CD-ROM.



Agilent Technologies

Agilent Technologies, Inc.

Printed in USA, April 2002



G2630-90277