

Agilent External Sample Filter for Micro GC

Installation Instructions

Using the External Filter Unit

The external sample filter, p/n CP736729, is recommended as a prefilter for the Agilent Micro GC. Whenever possible, remove moisture from samples introduced to the Micro GC. More information about sample gas requirements can be found in the *Agilent 490 Micro Gas Chromatograph User Manual*, part number G3581-90001.

The filter connection depends on the instrument's sample inlet type, [Figure 1](#), and [Figure 2](#) on page 2 show the filter components and how it is installed. [Figure 1](#) highlights the unheated front sample inlet. Connection for the (un)heated rear inlet is given in [Figure 2](#) on page 2.

Agilent advises replacement of the external filter unit at regular intervals. A package of 5 replacement filters is p/n CP736467. To install the filter element, the male part of the filter (CP736737) must be hand-tightened into the female part (CP736736), followed by a 1/8 turn with a 7/16-inch wrench. Orient the arrow on the female part of the filter towards the sample inlet.

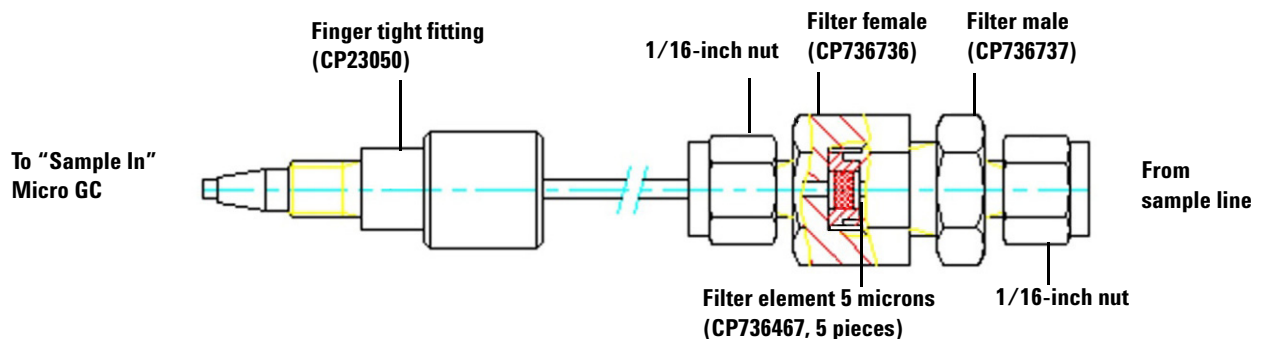


Figure 1 Unheated sample line (front) connection



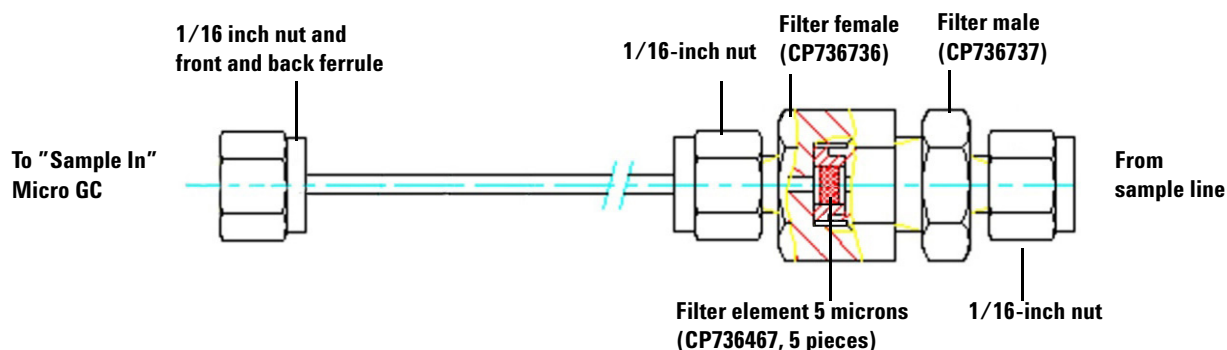


Figure 2 (Un)heated sample line (rear) connection

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.

© Agilent Technologies, Inc. 2014
 Printed in USA
 First edition, October 2014
 Agilent Technologies, Inc.
 5301 Stevens Creek Boulevard
 Santa Clara, CA 95051 USA



G3581-90011