



## Agilent Ultra Inert Universal Press Fit

**WARNING:** Wear safety glasses to protect your eyes while handling, cutting, or installing glass or fused silica capillary columns. Use care in handling these columns to prevent puncture wounds.

This information is subject to change without notice.

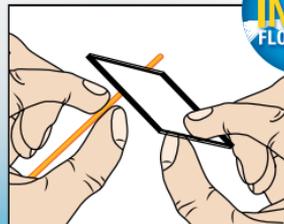
© Agilent Technologies, Inc. 2014  
Printed in USA, February 11, 2014  
5991-3944EN

## Instructions

AGILENT  
**INERT**  
FLOWPATH

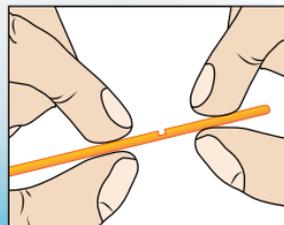
### Step 1

Lightly scribe column with wafer cutter (p/n 5181-8836).



### Step 2

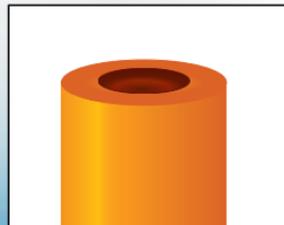
Gently break column apart.



### Step 3

Inspect column end for clean, smooth cut using magnifier (p/n 430-1020).

Wipe end of column with isopropyl alcohol and lint free cloth.



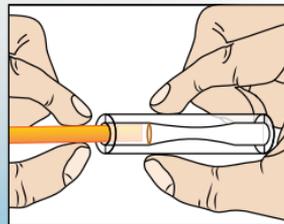
*(Continued on other side.)*



Agilent Technologies

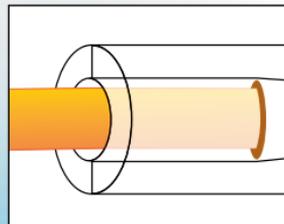
## Step 4

Grasp column near end and insert column into Press Fit.  
Push the two straight together firmly.



## Step 5

Inspect connection, using the magnifier. A brown ring completely around the end of the column indicates a leak-free seal. A tug on the column/union will assure an accurate connection. Though not required, you may also install your capillary columns using polyimide sealing resin, see [agilent.com/chem/pressfit\\_resin](https://www.agilent.com/chem/pressfit_resin)



## Step 6

Check union for leaks after your column is installed in the GC, and carrier gas is on, using an appropriate gas leak detector (p/n G3388B Agilent Electronic Leak Detector) before and after thermal cycling.

