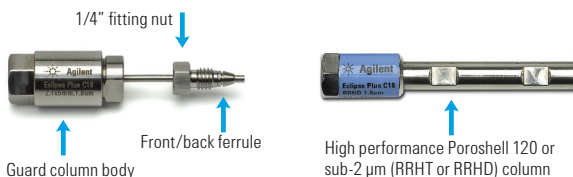


IMPORTANT INFORMATION ABOUT YOUR FAST GUARDS FOR UHPLC



Agilent **Fast Guards for UHPLC** are designed to protect high performance fast LC columns – such as RRHD, RRHT, and Poroshell 120 columns – with minimal impact on chromatographic performance. These guard columns are packed with the same particles and chemistries available in their respective Fast LC/UHPLC columns.



Installing your guard column:

Care must be taken upon installation to properly seat the guard column capillary and ferrule securely with the LC column to ensure both a leak-free and low-dispersion union. To install your guard column follow these steps:

1. Carefully remove the black rubber cap from the end of the guard column while ensuring that the fitting and ferrule (front and back) remain on the tubing.
2. Insert the guard column capillary into the inlet side of the LC column, so that the guard tubing is seated with the bottom of the column inlet.
3. Finger-tighten the 1/4" nut on the guard column into the analytical column. Once it is finger-tight, use a wrench to adjust an additional 1/4" turn to securely swage the ferrule and ensure a tight, leak-free fit. If the guard remains loose, tighten an additional 1/4" turn until the ferrule is seated and the guard is leak free.
4. Attach the tubing from the instrument and apply mobile phase flow through the guard.



Agilent Technologies

Care and use:

- Agilent Fast Guards for UHPLC are one-piece guard columns and are manufactured for high performance. After the guard column becomes clogged or the lifetime has been used, dispose of the guard column.
- Do not disassemble the guard column. Disassembly of the guard column will compromise its performance and stability.
- The guard hardware is rated to 60 °C at the maximum operating pressure. Guards are stable to their corresponding maximum column pressures:

<i>Guard Column Type</i>	<i>Pressure Limit</i>
Poroshell 120 Fast Guard	600 bar
Sub-2 µm Fast Guard	1200 bar

- Exceeding the pressure or temperature limits of the guard column will compromise the lifetime of the guard.

In order to allow for variance and flexibility in fitting choice, the ferrule on the guard column does not arrive pre-swaged.

This information is subject to change without notice.

© Agilent Technologies, Inc. 2012
Published in USA, June 25, 2012
820120-002

