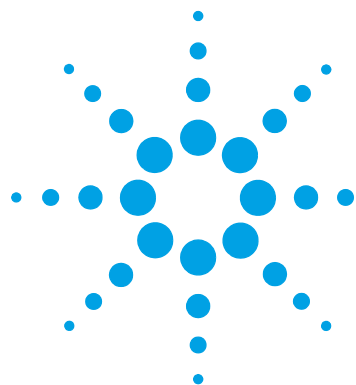




Installation of the Bio-Inert Solvent Heating Devices

Technical Note

This technical note describes the installation of the Bio-inert Solvent Heating Devices.



For bio-inert modules use bio-inert parts only!

In this note the installation of the Bio-Inert Solvent Heating Devices (9 μ L internal volume, 600 bar max pressure) (G5616-81000) into the Agilent 1290 Infinity TCC (G1316C) is described. The additional solvent heating devices can be arranged in the G1316C in various locations depending on the application needs. Some examples are shown in [Figure 1](#) on page 2.



Agilent Technologies

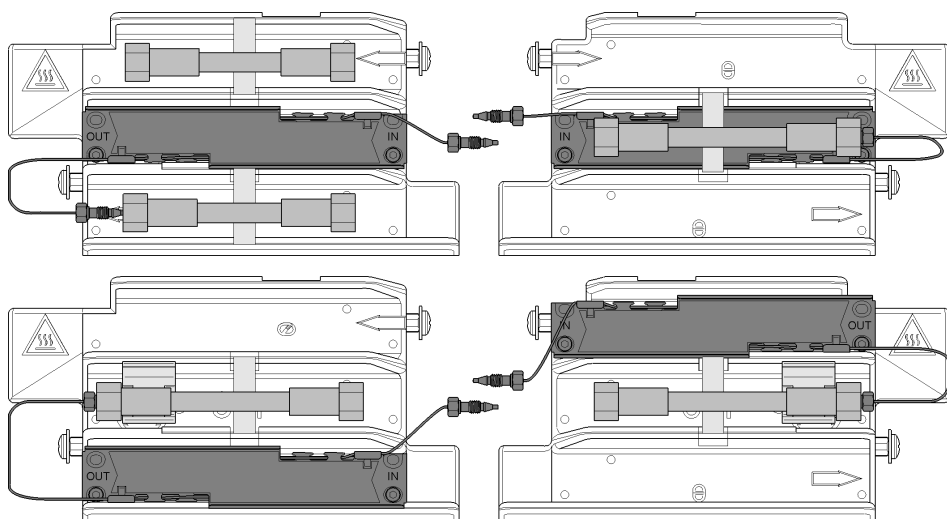


Figure 1 Arrangements of the bio-inert solvent heating devices in the 1290 Infinity TCC (G1316C)

NOTE

If the solvent heating devices are installed in the central position on the TCC's heat exchangers as shown in [Figure 1](#) on page 2, the column identification system cannot be used.

The solvent heating devices are mounted directly onto the TCC's heat exchangers.

NOTE

Don't forget to remove the protection foil from the heating transfer pad at the rear of the solvent heating device before installation.

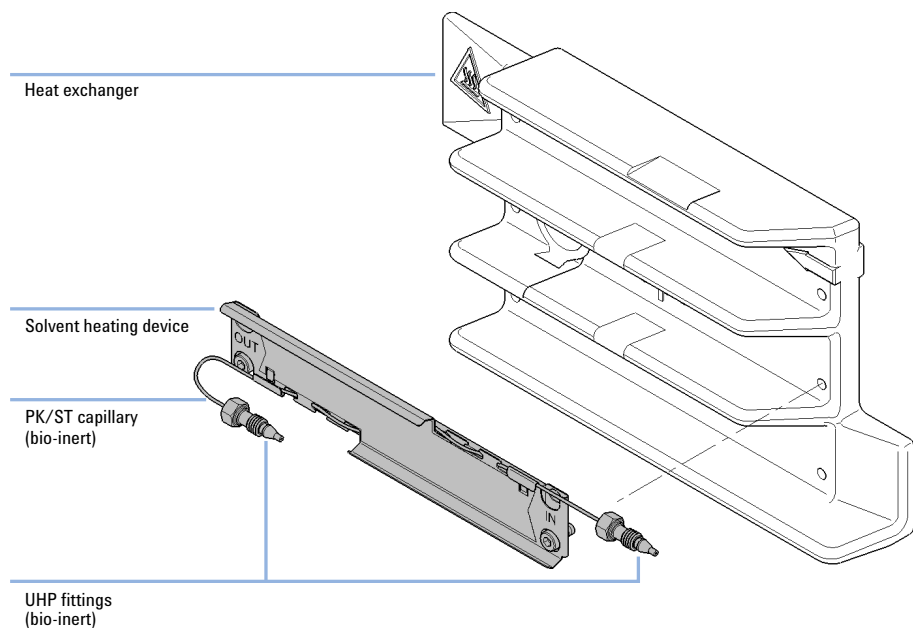


Figure 2 Installing the solvent heating device

CAUTION

Strong force/torque will damage PK/ST capillaries

Be careful when installing PK/ST capillaries.

→ Follow the instructions given in the technical note Agilent UHP-FF fittings Installation Note (G5611-90123).



G1316-90128

Part Number:
G1316-90128 Rev. B

Edition: 06/2013
Printed in Germany

© Agilent Technologies, Inc
2011-2012, 2013

Agilent Technologies, Inc
Hewlett-Packard-Strasse 8
76337 Waldbronn, Germany