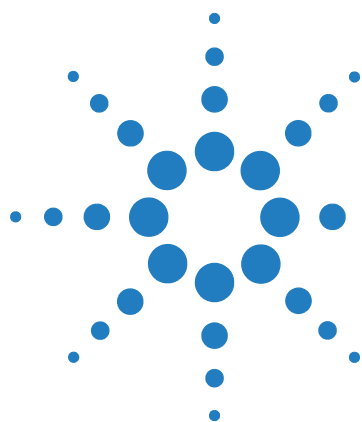




Installation of the Low Dispersion Heat-Exchanger Double Assemblies in the 1290 Series Thermostatted Column Compartment (G1316C)

In this note the installation of the additional Low Dispersion Heat Exchanger Double Assembly into the Agilent 1290 Series thermostatted column compartment (G1316C TCC) is described.

The device is typically mounted into the center location of either the left or the right heater element where it can support two columns.

The additional heater can be arranged in the G1316C in various locations depending on the application needs. Some examples are shown in [Figure 1](#).

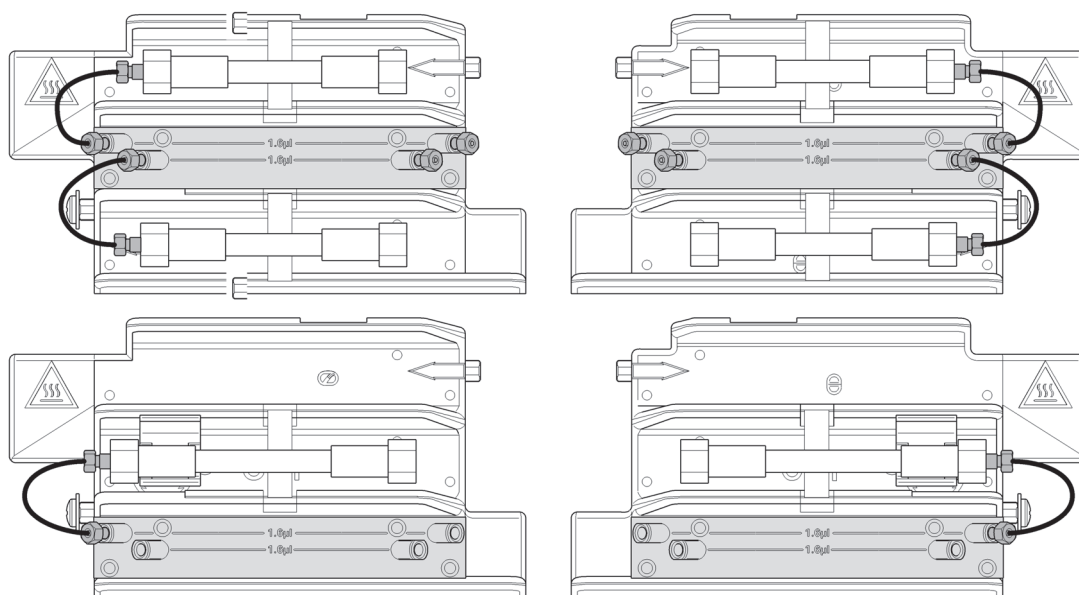


Figure 1 Arrangements of Heater and Cooling Devices (G1316C)

NOTE

If the additional heater and cooling devices are used as shown in [Figure 1](#) (top), the column identification system cannot be used. If the column identification system is required, fix the heater and cooling devices in the upper or lower locations or fix them right/left of the current location.

NOTE

The maximum flow rate to be used with the Low Dispersion Heatexchangers is 2.5 ml/min at 100 °C and 20 °C ambient.



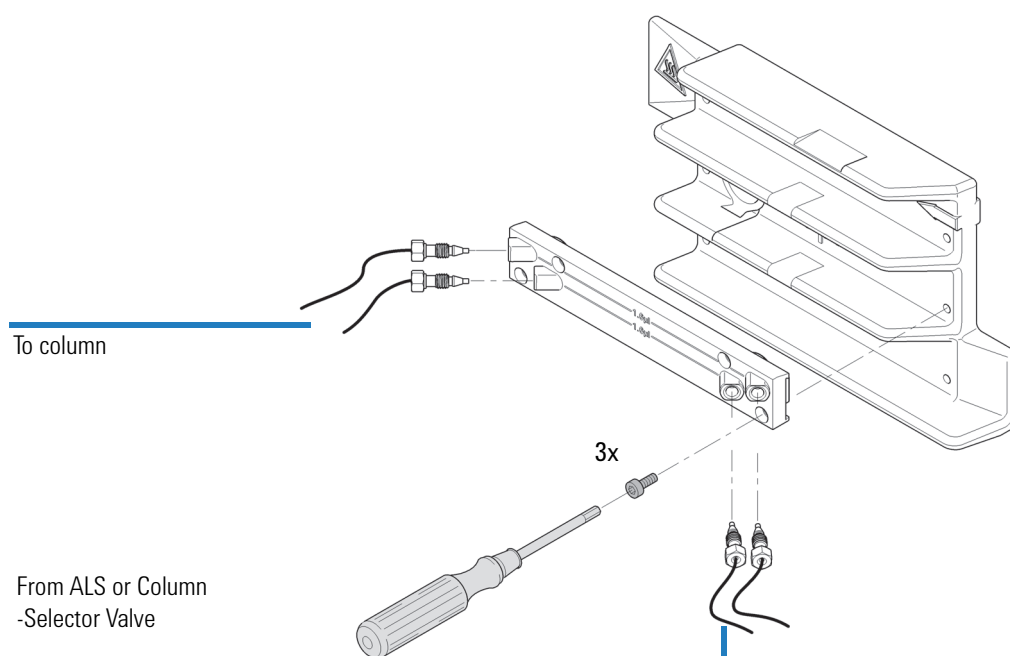
Agilent Technologies

Table 1 Delivery Checklist

Amount	Part	Comment
1	Low Dispersion Heat Exchanger Double Assembly	Incl. 3 screws for mounting
2	Capillary 0.12mm ID, 105mm length (5500-1188)	Connection Heat Exchanger to Column
	Capillary Fitting kit with long screw (G1314-68703)	Heat Exchanger Outlet Port and Column Inlet Port (alternative fittings shown below)

Table 2 Tools Required

Part	Comment
Hex Key 2.5mm 15cm long straight handle (p/n 8710-2412)	Required for fixing the Low Dispersion Heat Exchanger to the Heater Assembly
Hex driver, 1/4in, slitted (p/n 5023-0240)	Recommended for tightening the fitting nuts

**Figure 2** Fixing Heater or Cooling Devices (G1316C)