

# Agilent InfinityLab

## Quick Turn Fittings

*Engineered for efficiency and ease of use*



### [agilent.com/chem/infinitylab-supplies](https://www.agilent.com/chem/infinitylab-supplies)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2017, 2019  
Published in US, November 25, 2019  
Part Number 01200-90115  
DE.3273842593

### Procedure to make a connection with the InfinityLab Quick Turn LC fitting

1. Slide a capillary (with long socket) through the fitting.
2. Hand-tighten the fitting into the port. The connection can be tight up to 600 bar.
3. By using a 1/4 " open-wrench, tighten a maximum of 1/4 to 1/2 turn past hand-tight; the fitting tight up to 1,300 bar.

**Note:** The InfinityLab Quick Turn LC fitting can only be used with the stainless steel capillaries with long socket described in the ordering information.

### Procedure to exchange a capillary

1. Unscrew the fitting.
2. Pull the capillary out of the fitting.
3. Slide a new capillary (with long socket) into the fitting.
4. Hand-tighten the fitting into the port. The connection can be tight up to 600 bar.
5. By using a 1/4" open-wrench, tighten a maximum of 1/4 to 1/2 turn past hand-tight; the fitting is tight up to 1,300 bar.

**Note:** The InfinityLab Quick Turn LC fitting can be used multiple times with different capillaries, without loss of performance.



Agilent Technologies

## Ordering information

Agilent InfinityLab Quick Turn Parts	Part no.
<b>Fittings and ferrules</b>	
Quick Turn fitting	5067-5966
Front ferrule	5043-0924
<b>Capillaries (long socket)*</b>	
Stainless Steel 0.075 × 105 mm	5500-1198
Stainless Steel 0.075 × 150 mm	5500-1232
Stainless Steel 0.075 × 250 mm	5500-1206
Stainless Steel 0.075 × 500 mm	5500-1205
Stainless Steel 0.12 × 105 mm	5500-1188
Stainless Steel 0.12 × 150 mm	5500-1189
Stainless Steel 0.12 × 180 mm	5500-1233
Stainless Steel 0.12 × 200 mm	5500-1190
Stainless Steel 0.12 × 280 mm	5500-1191
Stainless Steel 0.12 × 500 mm	5500-1192
Stainless Steel 0.17 × 105 mm	5500-1193
Stainless Steel 0.17 × 150 mm	5500-1194
Stainless Steel 0.17 × 180 mm	5500-1234
Stainless Steel 0.17 × 200 mm	5500-1195
Stainless Steel 0.17 × 280 mm	5500-1196
Stainless Steel 0.17 × 380 mm	5500-1235
Stainless Steel 0.17 × 400 mm	5500-1236
Stainless Steel 0.17 × 500 mm	5500-1197
Stainless Steel 0.17 × 700 mm	5500-1237
Stainless Steel 0.25 × 105 mm with a female connection	5500-1261
Stainless Steel 0.25 × 150 mm	5500-1262
Stainless Steel 0.25 × 400 mm	5500-1263
Stainless Steel 0.12 × 130 mm M4	5500-1200
Stainless Steel 0.12 × 150 mm M4	5500-1288
Stainless Steel 0.17 × 150 mm M4	5500-1290

\* InfinityLab Quick Turn fittings require the capillaries specified in this table.