

[www.agilent.com/chem/infinitylab](http://www.agilent.com/chem/infinitylab)

DE75376504

This information is subject to change without notice.

© Agilent Technologies, Inc. 2022, 2023, 2025

Published in US, January, 2025

Part Number 01200-90114



## Agilent **InfinityLab** Quick Connect Fittings

*Engineered for efficiency and ease of use*

### Procedure to make a connection with the InfinityLab Quick Connect

1. Screw the fitting into the port until you **hear the clicks**.

2. Close the lever. The connection is tight to 1,300 bar.

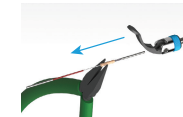
**Note:** The InfinityLab Quick Connect LC fitting can be reconnected multiple times without loss of performance. Do not turn the lever instead of the blue screw nut; this can lead to overtightening and damage of the fitting.

### Procedure to exchange a capillary

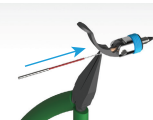
1. Disconnect the fitting from the port and grip the capillary on the PEEK adaptor with a plier.



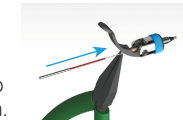
2. Pull the capillary out of the fitting.



3. Insert the new capillary into the fitting.



4. Press the PEEK adaptor into the fitting until it clicks into the final position.



# Agilent InfinityLab Quick Connect Fittings

Part	Part no.
Assemblies*	
Stainless steel 0.12 × 105 mm	5067-5957
Stainless steel 0.12 × 150 mm	5067-5958
Stainless steel 0.12 × 220 mm	5067-5959
Stainless steel 0.17 × 105 mm	5067-6166
Stainless steel 0.17 × 150 mm	5067-6167
Fittings and ferrules	
Quick Connect fitting	5067-5965
Front ferrule	5043-0924

\* Each assembly is equipped with a Quick Connect fitting, a capillary, and a Swagelok fitting, or with a female connection, where specified.

Capillaries**	
Stainless steel 0.075 × 105 mm	5500-1174
Stainless steel 0.075 × 150 mm	5500-1175
Stainless steel 0.075 × 220 mm	5500-1176
Stainless steel 0.075 × 280 mm	5500-1178
Stainless steel 0.12 × 105 mm	5500-1173
Stainless steel 0.12 × 150 mm	5500-1172
Stainless steel 0.12 × 220 mm	5500-1171
Stainless steel 0.12 × 280 mm	5500-1170
Stainless steel 0.17 × 105 mm	5500-1181
Stainless steel 0.17 × 150 mm	5500-1182
Stainless steel 0.17 × 220 mm	5500-1183

Stainless steel 0.17 × 280 mm	5500-1230
Stainless steel 0.25 × 105 mm with a female connection	5500-1258
Stainless steel 0.25 × 150 mm	5500-1259
MP35N 0.075 × 105 mm	5500-1474
MP35N 0.075 × 150 mm	5500-1475
MP35N 0.075 × 220 mm	5500-1576
MP35N 0.075 × 280 mm	5500-1577
MP35N 0.12 × 105 mm	5500-1578
MP35N 0.12 × 150 mm	5500-1579
MP35N 0.12 × 220 mm	5500-1580
MP35N 0.12 × 280 mm	5500-1581

MP35N 0.17 × 105 mm	5500-1584
MP35N 0.17 × 150 mm	5500-1585
MP35N 0.17 × 220 mm	5500-1586
MP35N 0.17 × 280 mm	5500-1587

\*\* The InfinityLab Quick Connect fitting can only be equipped with an InfinityLab capillary specified in this table. The InfinityLab capillary is designed with a spring and a holder. More dimensions are available.