

Agilent 1260 Infinity III Max Uptime Kit Bio

Technical Note

The Agilent 1260 Infinity III Max Uptime Kit Bio is for Agilent 1260 Infinity III Bio-Inert Systems.

Max Uptime Kit Bio

1260 Infinity III Max Uptime Kit Bio (G5654-68707):

#	p/n	Description
1	0890-1763	PEEK tubing 0.18 mm x 1500 mm
3	5067-4741	ZDV union (Bio-inert)
1	5065-4426	Colored finger-tight PEEK fittings, 10/PK
1	5971-5210	Assortment Box (not orderable separately)
1	8710-1930	Plastic tubing cutter
1	8710-2391	Rheotool socket wrench ¼ inch
2	5041-2168	Solvent inlet filter, 20 µm pore size
1	01018-22707	PTFE frits (pack of 5)
1	5043-0915	Fitting mounting tool
1	G5611-60502	Capillary Ti 0.17 mm x 900 mm, L (Bio-inert) Connection from Pump to Sampler
1	5500-1264	Capillary Ti 0.17 mm x 500 mm, SL/SLV Connection from Pump to Sampler
1	G5667-81005	Capillary PK/ST 0.17 mm x 500 mm, RLO/RLO (Bio-inert) Connection from Sampler to Bio Heat Exchanger

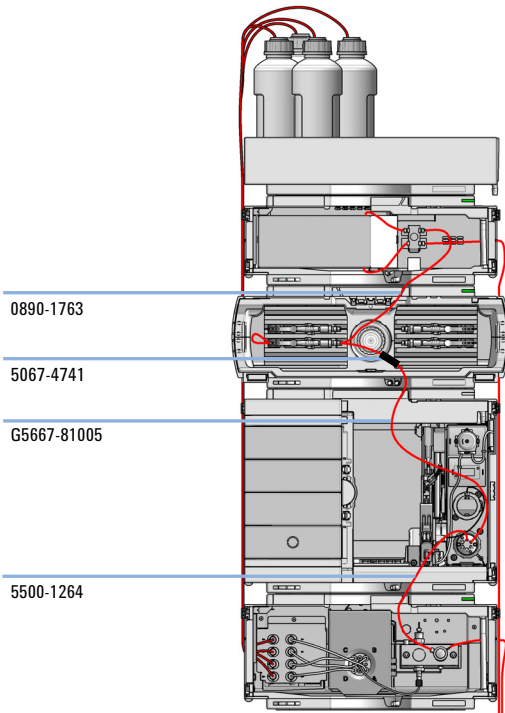


Figure 1 Capillary Connections

The information in this document also applies to Infinity II modules.

www.agilent.com

©Agilent Technologies Inc. 2016-2024

Printed in Germany
Edition: 10/2024



Part No: G5654-90130 Rev. C
Document No: D0026609 Rev. C

