

Quick Reference Guide

Agilent InfinityLab Supplies for the Agilent 1260 Infinity II Bio-inert LC

Agilent Technologies is committed to optimizing your laboratory's productivity, so we have produced this list of the most commonly ordered supplies and parts for the 1260 Infinity II Bio-inert LC.



Product Name Type/Group	PM Part*	Part Number
Bio-inert Pump supplies		
Bottle head assembly		G7120-60007
Glass filter, solvent inlet, 20 µm pore size		5041-2168
Purge valve		G5611-60062
Seal cap for purge valve		5067-4728
Active inlet valve		G5611-60025
Cartridge for active inlet valve		G5611-60020
Outlet ball valve		G5611-60067
Piston, sapphire		5067-4695
Piston seal	2 each	G5611-21503
Wash seal	2 each	0905-1731
Preventive maintenance kit	kit	G5611-68741
PTFE frits (5/pk)	1 each	01018-22707
Peristaltic pump (Pharmed)	1 each	5042-6401

Product Name Type/Group	PM Part*	Part Number
Bio-inert Multisampler supplies		
Needle	1 each	G5668-87200
Needle seat	1 each	G5668-87017
Rotor seal	1 each	5068-0209
Stator face seal	1 each	0100-1851
Sample loop 100 µL		G5668-60500
Drawer kit standard, double-height (1/pk)		G7167-60020
Drawer kit, single-height (1/pk)		G7167-60021
Drawer kit, triple-height (1/pk)		G7167-60022
Vial plate for 54x2 mL vials (6/pk)		G2255-68700
Multidraw kit, 400 bar		G5667-68711
Preventive maintenance kit	PM kit	G5668-68730

* Preventive maintenance part:

- kit: includes all standard parts required for PM

- part: standard part required for PM

- optional: part required for PM if referring option is installed

Product Name Type/Group	PM*	Part Number
Bio-inert manual injector supplies		
Rotor seal	1 each	5068-0082
Stator face	1 each	0100-1851
Bio-inert Multicolumn Thermostat supplies		
Bio-inert Quick Change heat exchanger (including one Quick Connect fitting, one UHP FF fitting)		G7116-60009
Bio-inert Quick Change heat exchanger, without fittings		G7116-60041
Rotor seal, bio-inert 2-position/6-port valve	1 each optional	0101-1409
Stator face, bio-inert 2-position/6-port valve	1 each optional	0100-1851
Rotor seal, bio-inert 2-position/10-port valve	1 each optional	5068-0041
Stator face, bio-inert 2-position/10-port valve	1 each optional	5068-0095
Rotor seal, bio-inert 4 column selector valve	1 each optional	5068-0045
Stator face, bio-inert 4 column selector valve	1 each optional	5068-0093
Column ID tag assembly		5067-5917
Column holder clips, soft (2/pk)		G7116-68003
Column holder clamp for larger columns (2/pk)		G7116-68004
Detector supplies		
Long-life deuterium lamp with RFID tag, for DAD and MWD	1 each	2140-0820
Tungsten lamp, for DAD and MWD		G1103-60001
Bio-inert standard flow cell with RFID, 10 mm, 13 μL , 120 bar, for DAD and MWD		G5615-60022
Titanium capillaries		
Capillary 0.17 mm id x 160 mm, with fittings		G5611-60503
Capillary 0.17 mm id x 400 mm, with fittings		G5611-60500
Capillary 0.17 mm id x 500 mm, with fittings		5500-1264
Capillary 0.17 mm id x 700 mm, with fittings		G5611-60501
Capillary 0.17 mm id x 900 mm, with fittings		G5611-60502
PEEK-lined stainless steel capillaries		
Capillary PK/ST 0.17 mm x 105 mm, incl. 2x fittings		G5667-81000
Capillary PK/ST 0.17 mm x 150 mm, incl. 2x fittings		G5667-81001
Capillary PK/ST 0.17 mm x 200 mm, incl. 2x fittings		G5667-81002
Capillary PK/ST 0.17 mm x 300 mm, incl. 2x fittings		G5667-81003
Capillary PK/ST 0.17 mm x 400 mm, incl. 2x fittings		G5667-81004
Capillary PK/ST 0.17 mm x 500 mm, incl. 2x fittings		G5667-81005
Capillary PK/ST 0.17 mm x 280 mm, for Quick Connect fitting		5500-1276
Capillary PK/ST 0.17 mm x 500 mm, for Quick Connect fitting		5500-1277
* Preventive maintenance part:		

* Preventive maintenance part:

- kit: includes all standard parts required for PM

- part: standard part required for PM

- optional: part required for PM if referring option is installed

This information is subject to change without notice.

© Agilent Technologies, Inc. 2019 Printed in the USA, August 20, 2019 5991-7669EN

Product Name Type/Group	PM*	Part Number
System supplies miscellaneous		
InfinityLab convenience kit 1260 Infinity II Bio-inert LC	InfinityLab convenience kit 1260 Infinity II Bio-inert LC	
InfinityLab capillary kit for 1260 Infinity II Bio-inert LC		5067-6621
InfinityLab Quick Connect fitting		5067-5965
InfinityLab Quick Turn fitting		5067-5966
Front ferrule for InfinityLab Quick Connect and InfinityLab Quick Turn fitting		5043-0924
Mounting tool for ultrahigh pressure front ferrule fitting and InfinityLab Quick Turn fitting		5043-0915
System tool kit		G7120-68708
Bio-inert union, 600 bar, 10-32 with PEEK insert		5067-4741
Solvent bottle, transparent 1 L		9301-6524
Solvent bottle, amber 1 L		9301-6526
Solvent bottle, transparent 1 L with cap		9301-6528
Solvent bottle, clear, 500 mL with cap		9301-6523
Solvent bottle, amber, 500 mL with cap		9301-6525
Identification rings (silicone) for solvent bottles (8pcs/4colors)		9301-6529
Stickers (removable) for solvent bottles (100/pack)		9301-6530
InfinityLab Stay Safe cap starter kit		5043-1222
InfinityLab Stay Safe cap (GL45) with one port, one venting valve, two fittings 3.2 mm		5043-1217
Venting valve for InfinityLab Stay Safe cap		5043-1190
6 L waste container including one InfinityLab Stay Safe cap GL45 with four ports		5043-1221
Charcoal filter (48 g) for waste container		5043-1193
InfinityLab Flex Bench		5043-1252
InfinityLab Flex Bench with power strip		5043-1759
InfinityLab Benchtop		5043-1711
InfinityLab Benchtop with power strip		5043-1740

Agilent columns and supplies help you get the results you need

Our wide range of parts, kits, and accessories are designed, manufactured, and tested to rigorous Agilent specifications, under a quality system registered to ISO 9001. Why risk compromising your analytical results with anything less than genuine Agilent supplies?

For a complete list of Agilent columns and supplies, visit us online at: www.agilent.com/chem/supplies

