

Agilent InfinityLab Supplies for the Agilent 1260 Infinity II LC with Multisampler

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 LC with Multisampler.



Pump supplies	PM*	Part number
Bottle head assembly		G7120-60007
Glass filter, solvent inlet, 20 µm pore size		5041-2168
Stainless steel filter, solvent inlet, 10 µm pore size		01018-60025
Piston seal PTFE, 2/Pk (reversed phase)	1 each	5063-6589
Piston seal PE, 2/Pk (normal phase)	1 each optional	0905-1420
Sapphire piston		5063-6586
Passive inlet valve		G1312-60066
Passive inlet valve (normal phase)		G1312-60166
Cartridge for active inlet valve, 600 bar		G1312-60020
Outlet valve		G1312-60067
Outlet valve (normal phase)		G1312-60167
Purge valve		G1312-60071
PTFE frits, for purge valve (pack of 5)	1 each	01018-22707
Seal cap for purge valve	1 each	5067-4728
Wash seal PTFE (reversed phase)	2 each	0905-1175
Wash seal PE (normal phase)	2 each optional	0905-1718
Gasket seal wash (pack of 6)	1 each	5062-2484
Peristaltic pump for seal wash (pharmed tubing)	1 each	5065-4445
PM kit for isocratic / quaternary pump	kit	G1310-68741
PM kit for binary pump	kit	G1312-68741
Seal wash preventive maintenance kit for quaternary and binary pumps		G1310-68742

* PM (preventive maintenance) relevant part information:
 - x each: amount of part needed per PM
 - kit: preventive maintenance kit that includes all standard parts for PM
 - optional: part required for PM only if referring option is installed

Multisampler supplies	PM*	Part number
Needle assembly standard	1 each	G4267-87201
High pressure needle seat 0.12 mm PEEK	1 each	G4267-87012
Metering seal PE for 100 µL analytical head		0905-1719
Piston assembly for 100 µL analytical head		5067-5678
Rotor seal (PEEK) for injection valve (600 bar)	1 each	5068-0209
Additional standard drawer kit, double-height (1/pk)		G7167-60020
Additional drawer kit, single-height (2/pk)		G7167-60021
Additional drawer kit, triple-height (2/pk)		G7167-60022
Vial plate for 54 x 2 mL vials, 6/pk		G2255-68700
Multidraw kit Infinity II, includes 500 and 1500 µL seat capillary extensions		G7167-68711
Large volume injection kit, includes 80 µL seat capillary extension for multi-draw operation		G4216-68711
Sample loop flex 100 µL, for single needle option		G4267-60500
Preventive maintenance kit 1260 Infinity II multisampler	kit	G7167-68730

Supplies for dual needle option

Rotor seal (PEEK) peripheral valve	1 each	5068-0229
Rotor seal (PEEK) injection valve	1 each	5068-0232
Sample Loop Flex 100 µL, right		G4267-60511
Sample Loop Flex 100 µL, left		G4267-60501
Preventive maintenance kit Multisampler dual needle	kit	G7167-68720

Multicolumn Thermostat supplies	PM*	Part number
InfinityLab Quick-Connect heat exchanger (0.17 mm id, 1.6 µL volume)		G7116-60051
InfinityLab Quick-Connect heat exchanger (0.17 mm id, 3.0 µL volume)		G7116-60061
Rotor seal (PEEK) for 4-column switching valve, 600 bar		5068-0264
Rotor seal (PEEK) for 2-position/6-port valve, 600 bar		0101-1409
Rotor seal (PEEK) for 2-position/10-port valve, 600 bar		0101-1415
Column ID Tag assembly		5067-5917
Column holder clips, soft (2/pk)		G7116-68003

VWD detector supplies	PM*	Part number
Deuterium lamp with RFID Tag	1 each	G1314-60101
Standard flow cell 10 mm, 14 µL, with RFID tag		G1314-60186
Semi-micro flow cell 6 mm, 5 µL, with RFID tag		G1314-60183
High pressure flow cell 10 mm, 14 µL, 400 bar, with RFID tag		G1314-60182
Inlet tubing assembly, PEEK 600 mm x 0.2 mm id		5062-8522

DAD/MWD detectors (G7115A, G7165A) supplies	PM*	Part number
Standard flow cell with RFID tag		G1315-60022
Semi-micro flow cell with RFID tag		G1315-60025
Long-life deuterium lamp with RFID tag	1 each	2140-0820
Tungsten lamp		G1103-60001

* PM (preventive maintenance) relevant part information:

- x each: amount of part needed per PM

- kit: preventive maintenance kit that includes all standard parts for PM

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

DAD detector (G7117A) supplies	PM*	Part number
Long-life deuterium lamp (8-pin) with RFID tag	1 each	5190-0917
Max-Light cartridge cell (10 mm, V(σ) 1.0 µL)		G4212-60008
Max-Light cartridge cell (60 mm, V(σ) 4.0 µL)		G4212-60007
FLD detector supplies	PM*	Part number
Flow cell 8 µL, 20 bar		G1321-60005
Flow cell 4 µL, 20 bar		G1321-60015
FLD wavelength calibration kit		G7121-68001
Capillaries and fittings	PM*	Part number
Capillary ST 0.17 mm x 500 mm, pump to sampler and sampler to MCT		5500-1246
Capillary ST 0.17 mm x 380 mm, column to detector		G1315-87311
Capillary ST 0.12 mm x 280 mm, column to detector		5500-1191
InfinityLab Quick Connect fitting assembly, 0.17 mm id x 105 mm length		5067-6166
Supplies and kits	PM*	Part number
1260 Infinity II convenience kit		5067-6617
1260 Infinity II capillary Kit		5067-6614
Inline filter kit (0.3 µm)		5067-6189
System tool kit		G7120-68708
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 (8 pcs/4 colors)		9301-6529
Sticker for solvent bottles (100/pk)		9301-6530
InfinityLab Stay Safe cap starter kit		5043-1222
InfinityLab Stay Safe cap (GL45) with one port, one venting valve, two fitting 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 4 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/lc-supplies