

Agilent Supplies for the Agilent 1260 Infinity Isocratic, Binary and Quarternary LC Systems

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 Binary, Quarternary and Isocratic LC Systems.



Pump supplies	PM*	Part number
Purge valve assembly		G1312-60071
PTFE frits, 5/pk	1 each	01018-22707
Seal cap gold, for purge valve	1 each	5067-4728
Passive inlet valve for G1310B/G1311B/C		G4280-60036
Active inlet valve for G1312B/C		G1312-60025
Cartridge for active inlet valve, 600 bar		G1312-60020
Outlet ball valve		G1312-60067
Outlet valve (normal phase)		G1312-60167
Sapphire piston		5063-6586
Pump seal (reversed-phase) 2/pk	1 each	5063-6589
Pump seal (normal-phase) 2/pk	1 each	0905-1420
Wash seal (normal-phase)	1 each	0905-1718
Wash seal gasket (reversed phase) 6/pk	kit	5062-2484
PM kit for binary pump, includes pump seals (reverse phase), PTFE frits, seal cap	kit	G1312-68741
PM kit for isocratic and quaternary pump, includes pump seals (reverse phase), PTFE frits, seal cap	kit	G1310-68741
Seal wash PM kit (2x required for binary pump)	kit	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

High Performance Autosampler supplies	PM*	Part number
Needle assembly for G1367E	1 each	G4226-87201
Needle seat assembly	1 each	G1367-87012
Peristaltic pump	1 each	5065-4445
Rotor seal, injection valve	1 each	0101-1416
100 µL loop capillary for G1367E		5067-4710
Piston seal, PTFE, 2/pk	1 each	5063-6589
Vial plate, for 54 x 2 mL vials, 6/pk		G2255-68700
PM kit for G1367E, includes needle assembly, needle seat, peristaltic pump, rotor seal, metering seal	kit	G1367-68741
Autosampler supplies	PM*	Part number
Needle assembly for G1329B	1 each	G1313-87201
Seat assembly, 0.17 mm id	1 each	G1329-87017
Seat assembly, 0.12 mm id		G1329-87012
Rotor seal, injection valve	1 each	0101-1416
100 µL loop capillary for G1329B		01078-87302
Piston seal, PTFE, 2/pk	1 each	5063-6589
Multidraw upgrade kit		G7167-68711
100 position tray, for 2 mL vials		G1364-60111
PM kit for G1329B, includes rotor seal, needle, seat assembly 0.17mm, seal for metering piston	kit	G1313-68719

* 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

System supplies	PM*	Part number
InfinityLab Quick Connect fitting assembly, 0.17 mm id, 105 mm length		5067-6166
LC tool kit		G7120-68708
Restriction capillary, 0.12 mm id, 2 m length		5022-2159
Blank nut, long, 10-32, PEEK with stainless steel core, finger-tight, for system diagnostic tests		5043-0277
Inline Filter 0.3µm		5067-6189
Replacement Frits (for Inline Filter), 5/pk		5023-0271
InfinityLab Flex Bench		5043-1252
InfinityLab Flex Bench with power strip		5043-1759
InfinityLab Benchtop		5043-1711
InfinityLab Benchtop with power strip		5043-1740
Solvent bottle, transparent 1 L		9301-6524
Solvent bottle, amber 1 L		9301-6526
Solvent bottle, transparent 1 L, with cap		9301-6528
Identification rings (silicone) for solvent bottles (8 pcs/4 colors)		9301-6529
Stickers (removable) for solvent bottles (100/pack)		9301-6530
InfinityLab Stay Safe cap starter kit		5043-1222
6 L waste container including InfinityLab Stay Safe cap GL45 with 4 ports		5043-1221
Charcoal filter, 48 g, for waste container		5043-1193

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

This information is subject to change without notice.