

Agilent InfinityLab Supplies for the Agilent 1260 Infinity II Prime System

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 Prime System.



Flexible pump supplies	PM*	Part number
Preventive maintenance kit, includes piston seals (2x), seal-wash seals (2x), peristaltic pump, outlet filter frit (1x)	kit	G7104-68741
Solvent cabinet		5067-5760
Bottle head assembly		G7120-60007
Solvent inlet filter, glass, pore size 20 µm		5041-2168
Solvent inlet filter, stainless steel, pore size 10 µm		01018-60025
Inlet valve		G4204-60022
Outlet valve		G4220-60028
Piston assembly		5067-5975
Frit for outlet filter (2/pack)	1 each	5067-5716
Peristaltic pump with Pharmed tubing	1 each	5065-4445
Inline filter upgrade kit		G7104-68000
Frit for inline filter, 0.3 µm (5/pack)	1 each optional	5023-0271
Jet Weaver mixer kit V35 µL, includes mixer and connection capillaries		G4204-68035
Jet Weaver mixer kit V100 µL, includes mixer and connection capillaries		G4204-68100
Jet Weaver mixer kit V380 µL, includes mixer and connection capillaries		G4204-68380
Inlet valve		G4220-60022
Outlet valve		G4220-60028
Pump seal PE (2/pk)	1 each	0905-1420
Piston assembly		5067-5938
Wash seal PE (1/pk)	2 each	0905-1718
Gaskets seal wash (6/pk)	1 each	5062-2484

* 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 (G7167A) supplies	PM*	Part number
Preventive maintenance kit, includes needle, needle-seat assembly, rotor seal	Kit	G7167-68730
Needle assembly	1 each	G4267-87201
High-pressure needle seat	1 each	G4267-87012
Rotor seal for injection valve	1 each	5068-0209
Metering seal for 100 µL analytical head		0905-1719
Sample loop, flexible, 100 µL, for single-needle option		G4267-60500
Drawer kit, standard, double-height (1/pack)		G7167-60020
Drawer kit, single-height (1/pack)		G7167-60021
Drawer kit, triple-height (1/pack)		G7167-60022
Vial plate for 54× 2-ml vials (6/pack)		G2255-68700
Vial plate for 40× 2-mL vials		5023-2471
Multidraw kit, includes 500 and 1500 µL seat-capillary extensions		G7167-68711
Large-volume injection kit, includes 80 µL seat-capillary extension for multidraw operation		G4216-68711
Supplies for Multisampler with Dual-Needle Option		
Rotor seal (PEEK) for peripheral valve	1 each optional	5068-0229
Rotor seal (PEEK) for injection valve	1 each optional	5068-0232
Sample loop, flexible, right, 40 µL		G4267-60411
Sample loop, flexible, right, 100 µL		G4267-60511
Sample loop, flexible, left, 40 µL		G4267-60401
Vial Sampler (7129C) supplies	PM*	Part number
Preventive maintenance kit	kit	G7116-60051
Finger caps for gripper fingers (15/pk)	1 each	6563-6506
Drawer for 66× 2-mL vials		G7129-60010
Drawer for 18× 6-mL vials		G7129-60110
Drawer kit 100× 2-mL vials, classic style		G7129-68210
Needle assembly	1 each	G7129-87200
Needle-seat assembly	1 each	G7129-87017
Rotor seal for injection valve	1 each	0101-1416
Sample loop 100 µL		G7129-60500
Multicolumn Thermostat (G7116A) supplies	PM*	Part number
InfinityLab Quick Connect heat exchanger, 0.17 mm id, 3.0 µL volume		G7116-60051
InfinityLab Quick Connect heat exchanger, 0.17 mm id, 6.0 µL volume		G7116-60061
Rotor seal for 4-column switching valve	1 each	5068-0264
Rotor seal for 2-position/6-port valve	1 each	0101-1409
Rotor seal for 2-position/10-port valve	1 each	0101-1415
InfinityLab Column ID tag assembly		5067-5917
InfinityLab Column ID tag assembly		G7116-68003
Column holder clips soft for small columns		G7116-68004
DAD and MWD (G7115A/G7165A) supplies	PM*	Part number
Long-life deuterium lamp with RFID tag	1 each	2140-0820
Standard flow cell with RFID tag		G1315-60022
Semimicro flow cell with RFID tag		G1315-60025

* 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

VWD (G7114A) 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
Semimicro flow cell, 6 mm, 5 µL, with RFID tag		G1314-60183
High-pressure flow cell, 10 mm, 14 µL, with RFID tag		G1314-60182
Inlet tubing assembly, PEEK, 600 × 0.2 mm id		5062-8522
DAD (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
Inline pressure relief valve kit		G4212-68001
FLD supplies	PM*	Part number
Detector lamp		2140-0600
Flow cell, 8 µL, 20 bar		G1321-60005
Flow cell, 4 µL, 20 bar		G1321-60015
Wavelength calibration kit, includes syringe, calibration sample, filters, fittings		G7121-68001
Capillaries and fittings	PM*	Part number
InfinityLab Quick Connect fitting assembly with capillary, 0.17 mm id, 105 mm length		5067-6166
InfinityLab Quick Connect capillary, 0.17 mm id, 105 mm length		5500-1181
InfinityLab Quick Turn LC fitting		5067-5966
Replacement ferrule for InfinityLab Quick Connect and Quick Turn fitting		5043-0924
Capillaries and fittings	PM*	Part number
Convenience kit for 1260 Infinity II LC, includes capillary kit, vials, solvent bottles		5067-6617
Capillary kit for 1260 Infinity II LC		5067-6614
System tool kit for Infinity II LC instruments		G7120-68708
InfinityLab Flex Bench		5043-1252
InfinityLab Flex Bench with power strip		5043-1759
InfinityLab Benchtop		5043-1711
InfinityLab Benchtop with power strip		5043-1740
Inline filter kit, 0.3 µm		5067-6189
Frit, 0.3 µm, for inline filter (5/pack)		5023-0271
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
InfinityLab Stay Safe cap (GL45) with 1 port, 1 venting valve, and 2 fittings, 3.2 mm		5043-1217
Venting valve for InfinityLab Stay Safe cap		5043-1190
Waste container, 6 L, including InfinityLab Stay Safe cap with 4 ports		5043-1221
Charcoal filter for waste container, 48 g		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.