

Agilent InfinityLab Supplies for the 1260 Infinity II Preparative LC 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 Preparative LC System.



Preparative Binary Pump supplies (G7161A)	PM*	Part number
Bottle head assembly		G7161-60042
Solvent inlet filter, stainless steel, 20 µm, 2/pk	2 each	5023-3115
Capillary kit, LC/UV 4-8 mL/min, for 9.4 or 10 mm id columns		5067-6858
Capillary kit, LC/UV 15-40 mL/min, for 21.2 mm id columns		5067-6859
Peristaltic pump, with Pharmed tubing	1 each	5065-4445
Seal wash tubing, 5 m		5065-9978
Filter frit, 2 µm	1 each	5022-2192
Piston seal (pump head, 50 mL)	4 each	5067-4299
Wash seal (pump head, 50 mL)	4 each	5067-6587
Piston (pump head, 50 mL)		G7161-60005
Check valve cartridge (pump head, 50 mL)		5067-6642
PM kit, for 1260 Infinity II preparative binary pump, 50 mL/min pump heads	Kit	G7161-68730
Preparative Autosampler supplies (G7157A)	PM*	Part number
Needle	1 each	G7157-87201
Needle seat	1 each	G7157-87101
Rotor seal injection valve	1 each	5068-0268
Peristaltic pump, with Pharmed tubing	1 each	5065-4445
Gripper finger caps, 15/pk	1 each	5063-6506
Sample loop		G7157-60515
Multi draw loop kit, 5 mL, for G7157A		G7157-68711
Drawer, for 66 x 2 mL vials		G7129-60010
Drawer, for 18 x 6 mL vials		G7129-60110
Drawer kit, for 100 x 2 mL vials		G7129-68210
PM kit, 1260 Infinity II preparative autosampler	Kit	G7157-68730

- * 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

Manual Preparative Injector (G1328D) supplies	PM*	Part number
Rotor seal	1 each	0101-1233
Stator face seal	1 each	5068-0321
PM kit, for preparative manual injector, includes stator face, rotor seal, needle port cleaner, and tools	Kit	5068-0319
Syringe, PTFE, removable Luer lock, 25 mL		5190-1544
Needle for glass syringe, Luer Lock, 16 gauge, stainless steel		9301-0407
Sample loop, 2 mL, stainless steel		0101-1244
Sample loop, 5 mL, stainless steel		0101-1243
Sample loop, 10 mL, stainless steel		RE3065023
Sample loop, 20 mL, stainless steel		RE3065025
Sample loop, 2 mL, PEEK		0101-1229
Sample loop, 5 mL, PEEK		0101-1228
Sample loop, 10 mL, PEEK		0101-1227
Sample loop, 20 mL, PEEK		0101-1226
Column Organizer (G9328A) supplies	PM*	Part number
Column holder clips (lamella type), 2/pk		G7116-68003
Column holder clamp, for columns 9.4–30 mm id		5067-6849
Column holder clamp, for columns 30-50 mm id		G7163-60002
Delay coils, for 4–8 mL/min		5067-6800
Delay coils, for 15-40 mL/min		5067-6181
Pressure relief valve kit		5067-6857
UV Detectors supplies	PM*	Part number
1260 Infinity II diode array detector (G7115A) and m wavelength detector (G7165A)	nultiple	
Long-life deuterium lamp, with RFID tag	1 each	2140-0820
Tungsten lamp		G1103-6000
Preparative flow cell, quartz, 3 mm		G7115-60001
Preparative flow cell, quartz, 0.3 mm		G7115-60002
Preparative flow cell, quartz, 0.06 mm		G7115-60003
Capillary ST flow cell, 4-8 mL/min		5500-1342
Capillary ST Flow Cell 15-40 mL/min		5500-1343
1260 Infinity II variable wavelength detector (G7114	-A)	
Long-life deuterium lamp, with RFID tag	1 each	G1314-60101
Flow cell, for VWD, 3 mm path length		G1314-60024
Flow cell, for VWD, 0.3 mm path length		G1314-60025
Flow cell, for VWD, 0.06 mm path length		G1314-60023

Evaporative Light Scattering Detector (G4260B) supplies	PM*	Part number
Splitter kit, 2–10 mL/min	1 each	5023-2256
Splitter kit, 8-40 mL/min		5023-2257
MS Flow Modulator (G7170B) supplies	PM*	Part number
Stator flow modulator		0101-1393
Rebuild kit, for flow modulator, includes 3 stator face seals, 3 rotor seals, and a hex key	1 each	1535-5215
Upgrade kit, for mass-based fraction collection		5067-6861
Delay Coil Organizer (G9324A) supplies	PM*	Part number
Knitted delay kit, for 9.4 mm id columns, 4–8 mL/min flow, mass-based prep systems		5067-6800
Knitted delay kit, for 21.2 mm ID columns, 15–40 mL/min flow, mass-based prep systems		5067-6101
Preparative Fraction Collector (G1364E) supplies	PM*	Part number
Tubing kit, 4–8 mL/min, 0.5 mm id	1 each optional	G1364-68603
Tubing kit, 15–40 mL/min, 0.8 mm id	1 each optional	G1364-68604
Checkout solution	1 each	5190-6886
Exhaust tube adapter kit		G1364-68709
Tray holding 40 tubes, 30 x 100 mm, 50 mL		G1364-84543
Tray holding 60 tubes, 25 x 100 mm, 35 mL		G1364-84544
Tray holding 126 tubes, 16 x 100 mm, 14 mL		G1364-84545
Tray holding 215 tubes, 12 x 100 mm, 7 mL		G1364-84536
Tray for 40 x 2 mL vials		G1364-44612
Tray for 15 x 6 mL vials		G1364-44613
Tray for 100 x 2 mL vials		G1364-44610
Tray for 100 x 2 mL vials, thermostattable		G1364-60111
Full tray, for 4 well plates		G1364-60621
Fraction tubes, 100 x 30 mm, 45 mL, 100/pk		5042-6458
Fraction tubes, 100 x 25 mm, 35 mL, 100/pk		5042-6459
Fraction tubes, 100 x 16 mm, 14 mL, 250/pk		5022-6532
Fraction tubes, 100 x 12 mm, 7 mL, 250/pk		5022-6531

^{*} 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

Valve-Based Preparative Fraction Collector (G7166A) supplies	PM*	Part number
Back pressure regulator, 20 psi		5067-6840
Fittings, for 2.3 mm od tubing, 2/pk		5043-1215
Delay calibrant and checkout mix	1 each	5190-8223
Valve supplies		
2-position / 10-port valve head (G4730A)		
Stator head		5068-0152
Rotor seal	1 each	5068-0153
Bearing ring	1 each	1535-4045
8-position / 9-port valve head (G4731A)		
Stator head		5068-0154
Rotor seal	1 each	5068-0155
Bearing ring	1 each	1535-4045
6-position / 14-port valve head (G4734A)		
Stator head		5068-0248
Rotor seal	1 each	5068-0249
Bearing ring	1 each	1535-4045
Column selector capillary kit, 4-8 mL/min**		5067-6867
Column selector capillary kit, 15-40 mL/min**		5067-6868
Clustering Valve (G9322A) supplies	PM*	Part number
Capillary kit, for clustering up to 3 fraction collectors		5067-6811
Capillary upgrade kit, 4-8 mL/min		5043-1540

^{**}One kit per column required

Miscellaneous InfinityLab supplies for Preparative LC	PM*	Part number
System tool kit		G7120-68708
Back pressure regulator, 20 psi		5067-6840
ERI remote and leak module		5188-8053
External leak sensor		5067-6149
Tubing and fittings		
Tubing, PTFE/ESD, 1.6 mm od, 0.8 mm id, 6 m		5043-1540
Fitting, PEEK-ESD 1/4-28, for 1.6 mm od, 6/pk		5023-2898
Tubing, PTFE/ESD, 2.0 mm od, 1.2 mm id, 2 m		5023-2878
Fitting, PEEK-ESD 1/4-28, for 2.0 mm od, 6/pk		5023-2899
Tubing, PTFE/ESD, 2.5 mm od, 1.6 mm id, 6 m		5023-2882
Fitting, PEEK-ESD 1/4-28, for 2.5 mm od, 6/pk		5023-2883
Adapter, 10–32 female to 1/4-28 stainless steel male		5023-1803
Adapter, stainless steel 10–32 female to 1/4-28 female flat-bottom		5023-2875
Adapter, stainless steel 1/4–28 female flat-bottom to 10–32 male		5023-2512
Union, flat-bottom 1/4-28 ESD		5023-2508
Solvent bottles (GL45) and accessories		
Solvent bottle, clear, 125 mL, with cap		9301-6527
Solvent bottle, clear, 500 mL, with cap		9301-6523
Solvent bottle, clear, 1,000 mL, with cap		9301-6528
Solvent bottle, amber, 1,000 mL		9301-6526
Solvent bottle, clear, 2,000 mL		9301-6342
Solvent bottle, amber, 2,000 mL		9301-6341
InfinityLab identification silicone rings, 2 x 4/pk		9301-6529
Sticker for solvent bottles, removable, 100/pk		9301-6530
Vials		
Vial, screw top, amber with write-on spot, certified, 2 mL, 100/pk		5182-0716
Cap, screw, blue, pre-slit PTFE/silicone, 100/pk		5183-2076
Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk		5188-5369
Septum, pre-slit PTFE/silicone, 16 mm, 100/pk		5188-2758
Cap, screw, for 6 mL vials, 100/pk		9301-1379

^{*} 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

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

