

Agilent Supplies for the Agilent 1260 Infinity II SFC 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 SFC System.



SFC control module	PM*	Part number
Seal	1 each	G4301-20200
Booster pump upgrade kit		G4301-68500
Low dispersion nozzle assembly		G4301-60501
Pump supplies	PM*	Part number
Bottle head assembly		G7120-60007
Bottle head assembly, ultraclean tubing quality, recommended for MS applications		G7120-60017
Tubing kit, ultraclean quality, including four bottle head assemblies (G7120-60017) and three tubing kits 140 mm (G4220-60070)		G7120-68070
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), for channel B	1 each	5063-6589
Piston seal PE, 2/pk (normal phase), for channel A	1 each	0905-1420
Piston, sapphire		5067-4695
Inlet valve (channel B)		G1312-60066
SFC inlet valve (channel A)		G4302-60066
Outlet valve (channel B)		G1312-60067
Outlet valve N-type (channel A)		G1312-60167
Purge valve		G1312-60071
PTFE frits for purge valve (5/pk)	1 each	01018-22707
Seal cap, for purge valve	1 each	5067-4728
Wash seal PE (normal phase)	2 each	0905-1718
Gasket, seal wash (6/pk)	1 each	5062-2484
Peristaltic pump for seal wash (Pharmed tubing)	1 each	5065-4445

* Preventive maintenance options:

- kit: includes all standard parts required for PM
- part: standard part required for PM
- optional: part required for PM 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	1 each	5068-0279
Additional standard drawer kit, double-height (1/pack)		G7167-60020
Additional drawer kit, single-height (1/pk)		G7167-60021
Additional drawer kit, triple-height (1/pk)		G7167-60022
Vial plate for 54 x 2 mL vials (6/pk)		G2255-68700
Sample loop flex 100 µL		G4267-60500
Peristaltic pump for seal wash (Pharmed tubing)	1 each	5065-4445
Multi column thermostat supplies	PM*	Part number
InfinityLab Quick Connect heat exchanger (0.12 mm id, 1.6 µL volume)		G7116-60015
InfinityLab Quick Connect heat exchanger (0.12 mm id, 3.0 µL volume)		G7116-60031
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 (PEEK) for 4-column switching valve	1 each optional	5068-0264
Rotor seal (PEEK) for 2 position/6 port valve	1 each optional	0101-1409
Rotor seal (PEEK) for 2 position/10 port valve	1 each optional	0101-1415

* Preventive maintenance options:

- kit: includes all standard parts required for PM

- part: standard part required for PM

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

Capillaries and fittings	PM*	Part number
Capillary ST 0.17 mm id x 500 mm length, pump to sampler & sampler to MCT		5500-1246
Capillary ST 0.17 mm id x 380 mm length, column to detector		G1315-87311
Capillary ST 0.12 mm id x 280 mm length, column to detector		5500-1191
InfinityLab Quick Connect fitting assembly, 0.17 mm id x 105 mm length		5067-6166
InfinityLab Quick Turn LC fitting		5067-5966
Supplies and kits	PM*	Part number
SFC/MS splitter kit		G4309-68715
InfinityLab capillary kit for 1260 Infinity II LC		5067-6614
InfinityLab capillary kit for 1290 Infinity II LC		5067-6615
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 1L, 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
Sticker (removable)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 fittings 3.2 mm		5043-1217
Venting valve for InfinityLab Stay Safe Cap		5043-1190
6 L waste container including with one InfinityLab Stay Safe cap GL45 with four ports		5043-1221
Charcoal Filter (48g) 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/supplies

This information is subject to change without notice.