

Agilent InfinityLab LC Purification Solutions

Agilent offers tailor-made, high-efficiency LC purification solutions ranging from analytical to preparative scale. This comprehensive portfolio of state-of-the-art LC purification instruments offers you scalable LC purification solutions to meet your needs and budget. InfinityLab columns and supplies are designed to work seamlessly with Agilent InfinityLab instruments, and draw on 40+ years of technological leadership.

With Agilent OpenLab software and support from Agilent CrossLab instrument and enterprise service plans, Agilent provides a complete purification solution that will make a real difference to your lab's operational efficiency.



1290 Infinity II Preparative LC Systems – Purify with the new benchmark in efficiency

Aspire to maximum efficiency in your purification workflows. Get maximum flexibility to solve all your purification challenges with highest purity, recovery, and speed.



1260 Infinity II Preparative LC Systems – Benefit from efficient purification – every day

Achieve highest purity and recovery with the most affordable fraction collector. Get highest flexibility in fraction size while maintaining high-precision recovery.



1220/1260/1290 Infinity II Analytical-Scale LC Purification Systems – Experience the power of analytical compound isolation

(Only preparative-scale systems are covered in this guide)

Isolate your compounds of interest through high-performance separation and low-dispersion fraction collection. Choose the bio-inert option for purification of your proteins or sensitive samples

Pumps

	PM*	Part Number
Supplies for 1260 Infinity II Preparative Binary Pump (G7161A)		
Standard preventive maintenance parts		
Preventive maintenance kit, for the 1260 Infinity II preparative binary pump (G7161A), and 1290 Infinity II preparative binary pump (G7161B) equipped with 50 mL/min pump heads	PM kit	G7161-68730
Solvent bottle inlet filter, preparative, 20 µm, SST, 2/pk	2 each	5023-3115
Filter assembly, preparative, 12.7 mm id, 2 µm, SST	1 each	5022-2192
Piston Seal, for 1260 Infinity II preparative binary pump and 1290 Infinity II preparative binary pump G7161B with AtP pump head	4 each	5067-4299
Seal wash PE for 1260 Infinity II preparative binary pump G7161A and for AtP pump head, 1/pk	4 each	5067-6587
Peristaltic pump with PharMed tubing	1 each	5065-4445
Extended preventive maintenance parts		
Piston collar, 5 mm, for G7161A 1260 Infinity II preparative binary pump and G7161B 1290 Infinity II preparative binary pump with AtP pump head		G7161-60005
Check valve assembly, inlet/outlet, for G7161A/B preparative pump, with 50 mL/min pump head		5067-6642
Purge valve prep		G7161-60036
Solvent mixer		G1312-87330
Silicone tubing, 1 mm id, 3 mm od, 5 m		5065-9978
Other supplies		
Bottle head assembly for 1260/1290 Infinity II Preparative LC system		G7161-60042
InfinityLab capillary kit, for flow rate range 4 to 8 mL/min or column with 9.4 mm id		5067-7015
InfinityLab capillary kit, for flow rate range 15 to 40 mL/min or column with 21.2 mm id		5067-7016
InfinityLab capillary kit, for flow rate range 40 to 80 mL/min or column with 30 mm id		5067-7017
InfinityLab Supplies for 1290 Infinity II Preparative Binary Pump G7161B		
Standard preventive maintenance parts (200 mL/min flow pump heads)		
Preventive maintenance kit 1290 Infinity II Preparative Binary Pump	PM kit	G7161-68740
Piston seal, for 1290 Infinity II preparative binary pump G7161B with preparative pump head	4 each	5067-6589
Seal wash PE for 1290 Infinity II preparative binary pump G7161B, 1/pk	4 each	5067-6588
Filter assembly, preparative, 12.7 mm id, 2 µm, SST	1 each	5022-2192
Peristaltic pump with PharMed tubing	1 each	5065-4445

	PM*	Part Number
Standard preventive maintenance parts (optional 50 mL pump heads)		
Preventive maintenance kit, for the 1260 Infinity II preparative binary pump (G7161A), and 1290 Infinity II preparative binary pump (G7161B) equipped with 50 mL/min pump heads	PM kit	G7161-68730
Piston Seal, for 1260 Infinity II preparative binary pump and 1290 Infinity II preparative binary pump G7161B with AtP pump head	4 each/option depending	5067-4299
Seal wash PE for 1260 Infinity II preparative binary pump G7161A and for ATP pump head, 1/pk	4 each/option depending	5067-6587
Solvent bottle inlet filter, preparative, 20 µm, SST, 2/pk	2 each	5023-3115
Extended preventive maintenance parts		
Check valve assembly double seat, inlet/outlet		G7161-60052
Piston, for G7161B 1290 Infinity II preparative binary pump with preparative pump head		G7161-60028
Purge valve prep		G7161-60036
Solvent Mixer		G1312-87330
Silicone tubing, 1 mm id, 3 mm od, 5 m		5065-9978
Piston collar, 5 mm, for G7161A 1260 Infinity II preparative binary pump and G7161B 1290 Infinity II preparative binary pump with AtP pump head		G7161-60005
Check valve assembly, inlet/outlet, for G7161A/B preparative pump, with 50 mL/min pump head		5067-6642
Other supplies		
Bottle head assembly, for 1260/1290 Infinity II Preparative LC system		G7161-60042
InfinityLab capillary kit, for flow rate range 4 to 8 mL/min or column with 9.4 mm id		5067-7015
InfinityLab capillary kit, for flow rate range 15 to 40 mL/min or column with 21.2 mm id		5067-7016
InfinityLab capillary kit, for flow rate range 40 to 80 mL/min or column with 30 mm id		5067-7017
InfinityLab capillary kit, for flow rate range 80 to 200 mL/min or column with 50 mm id		5067-7018

*** PM (preventive maintenance) part information key**

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

Sample injection modules

	PM*	Part Number
InfinityLab Supplies for Preparative Open-Bed Sampler/Collector G7158B		
Parts for preventive maintenance		
InfinityLab tubing kit, 8 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60953
InfinityLab tubing kit, 50 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60952
InfinityLab tubing kit, 150 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60954
InfinityLab tubing kit, 200 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60951
Dripping adapter, for fraction valve	1 each	G9321-26600
Rotor Seal, 2 pos/14 port preparative valve head, 600 bar	1 each	5068-0323
Peristaltic pump head, for 1290 Infinity II Preparative Open-Bed Sampler/Collector G7158B	2 each	5067-6668
Needle plug assembly, for use with the 1290 Infinity II Preparative Open-Bed Sampler/Collector	1 each	G9321-60036
Tubing kit, for sampling unit	1 each	G9321-68006
Solvent distribution tubing kit, for use with 1290 Infinity II Preparative Open-Bed Sampler/Collector.	1 each	5500-1429
Preparative LC needle tube kit, 1.2 × 2.0 mm calibration, for use with the 1290 Infinity II Preparative Open-Bed Sampler/Collector	1 each	G9321-68735
Solvent bottle inlet filter, 20 µm, glass	2 each	5041-2168
Extended preventive maintenance parts		
Stator, for 2pos/14 port valve	1 each	5068-0322
Containers and vial plates		
InfinityLab drawer, ambient, with RFID tag		G9321-60085
InfinityLab tube container, 10 tubes, 30 × 150 mm, ambient		G9321-60015
InfinityLab tube container, 18 tubes, 25 × 150 mm, ambient		G9321-60025
InfinityLab tube container, 36 tubes, 16 × 150 mm, ambient		G9321-60129
InfinityLab tube container, 72 tubes, 12 × 150 mm, ambient		G9321-60131
InfinityLab tube container, 10 tubes, 30 × 100 mm, ambient		G9321-60058
InfinityLab tube container, 18 tubes, 25 × 100 mm, ambient		G9321-60035
InfinityLab tube container, 36 tubes, 16 × 100 mm, ambient		G9321-60055
InfinityLab tube container, 72 tubes, 12 × 100 mm, ambient		G9321-60045

	PM*	Part Number
InfinityLab Drawer Gilson 207, ambient, contains two drawers for Gilson rack type 207, for use with 16 × 100 mm (od × length) test tubes		G9321-60201
InfinityLab Drawer Gilson 208, ambient		G9321-60208
InfinityLab SBS well plate container, ambient		G9321-60051
Vial plate, for 15 × 6 mL vials, low profile, nonstackable		5043-1826
Vial plate, for 40 × 2 mL vials		5023-2471
96-well plate, 2.0 mL, square wells, U shape, polypropylene, 41 mm, 30/pk		5043-9300
Sealing mat, 96 wells, square, preslitted, silicone, 50/pk		5043-9319
96-well plate, 0.90 mL, round wells, U shape, polypropylene, 32 mm, 50/pk		5043-9305
96-well plate, 1.0 mL, round wells, U shape, polypropylene, 27 mm, 50/pk		5043-9309
Sealing mat, 96 wells, round, preslitted, silicone, 50/pk		5043-9317
Fraction tubes and vials		
Test tubes, 30 × 150 mm, 78 mL, 100/pk		5190-9090
Test tubes, 25 × 150 mm, 55 mL, 100/pk		5190-9091
Test tubes, 16 × 150 mm, 21 mL, 250/pk		5190-9092
Test tubes, 12 × 150 mm, 11 mL, 250/pk		5190-9093
Round bottom glass vial, 30 × 100 mm 100/pk		5042-6458
Round bottom glass vial, 25 × 100 mm 100/pk		5042-6459
Test tubes, 16 × 100 mm, 250/pk		5022-6532
Test tubes, 12 × 100 mm, 250/pk		5022-6531
Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk		9301-1377
Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk		5188-5369
Cap, screw, for 6 mL vials, 100/pk		9301-1379
Septa, for 6 mL vials, 100/pk		9301-1378
Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk		5182-0716
Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk		5182-0715
Cap, screw, blue, preslit PTFE/silicone, 100/pk		5183-2076

	PM*	Part Number
Other supplies		
Sample loop, stainless steel, 400 µL, 1/16 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector (G7158B)		5068-0348
Sample loop, stainless steel, 2 mL, 1/16 in		5068-0350
Preparative sample loop, stainless steel, 5 mL		5068-0351
Sample loop, stainless steel, 5 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector G7158B		5068-0334
Sample loop, stainless steel, 10 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector		5068-0335
Sample loop, stainless steel, 20 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector		5068-0336
Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL		5190-8223
InfinityLab bottom fume hood kit, for Agilent 1290 Infinity II preparative open-bed sampler/collector		G9321-68300
Supplies for 1260 Infinity II Preparative Autosampler (G7157A)		
Parts for preventive maintenance		
Preventive maintenance kit, for the 1260 Infinity II preparative autosampler G7157A	PM kit	G7157-68730
Rotor seal, 6 port injection valve, preparative autosampler G7157A	1 each	5068-0268
Needle assembly, for G7157A preparative autosampler	1 each	G7157-87201
Needle seat, for G7157A preparative autosampler	1 each	G7157-87101
Peristaltic pump with PharMed tubing	1 each	5065-4445
Finger caps 15/pk	1 each	5063-6506
Other supplies		
Drawer, for 66 × 2 mL vials	x	G7129-60010
Drawer for 18 × 6 mL vials	x	G7129-60110
Drawer kit for 100 × 2 mL vials, classic style	x	G7129-68210
Multi draw loop kit, 5 mL, for G7157A, 1260 Infinity II preparative autosampler	x	G7157-68711
Sample loop, stainless steel, 900 µL	x	G7157-60515
Supplies for 1260 Infinity II Manual Preparative Injector (G1328D)		
Parts for preventive maintenance		
Rebuild kit, for 1260 Infinity II preparative manual injector	PM kit	5068-0319
Rotor seal for Rheodyne 3725, 3725-038, PEEK	1 each	0101-1233
Stator face seal, for 1260 Infinity II preparative manual injector	1 each	5068-0321

	PM*	Part Number
Other supplies		
Sample loop, stainless steel, 2 mL, 1.00 mm id		0101-1244
Sample loop, stainless steel, 5 mL, 1.00 mm id		0101-1243
Adapter, stainless steel, female 10 to 32 to male 5/16-24		5067-6845
Sample loop, stainless steel, 5 mL, 1/8 in od		5067-7031
Sample loop, stainless steel, 10 mL, 1/8 in od		5067-7032
Sample loop, stainless steel, 20 mL, 1/8 in od		5067-7033
Sample loop, PEEK, 2 mL, 1.60 mm id		0101-1229
Sample loop, PEEK, 5 mL, 1.60 mm id		0101-1228
Sample loop, PEEK, 10 mL, 2.00 mm id		0101-1227
Sample loop, PEEK, 20 mL, 2.00 mm id		0101-1226
Manual syringe, 25 mL, removable Luer lock with push button valve, PTFE-tip plunger		5190-1544
Glass syringe, 1 mL, Luer tip, 0.05 scale		9301-1446
Needle, sample loading, stainless steel, Luer, 16 gauge, 1/pk		RE3725086
Stator manual prep injector		5068-0320
Holder, for mounting the preparative manual injector to the prep column organizer		5067-6831
Vials and Caps for use with Autosampler and Manual Injector		
Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk		9301-1377
Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk		5188-5369
Cap, screw, for 6 mL vials, 100/pk		9301-1379
Septa, for 6 mL vials, 100/pk		9301-1378
Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk		5182-0716
Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk		5182-0715
Cap, screw, blue, preslit PTFE/silicone, 100/pk		5183-2076

*** PM (preventive maintenance) part information key**

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

Fraction collectors

	PM*	Part Number
Supplies for 1260 Infinity II Preparative Fraction Collector (G1364E)		
Parts for preventive maintenance (G1364E)		
Tubing kit, for flow rate 40 to 100 mL/min, 30 mm columns, for fraction collector G1364E	1 each	G1364-68605
Tubing kit, for flow rate 4 to 8 mL/min, 9.4 mm columns, for fraction collector G1364E/F and G5664B	1 each/optional	G1364-68603
Tubing kit, for flow rate 15 to 40 mL/min, 21.2 mm columns, for fraction collector G1364E	1 each/optional	G1364-68604
Other supplies		
Tray holding 40 tubes, 30 × 100 mm, 50 mL, for G1364E fraction collectors	x	G1364-84543
Round bottom glass vial, 30 × 100 mm 100/pk	x	5042-6458
Tray holding 60 tubes, 25 × 100 mm, 35 mL, for G1364E fraction collectors	x	G1364-84544
Round bottom glass vial, 25 × 100 mm 100/pk	x	5042-6459
Tray holding 126 tubes, 16 × 100 mm, 14 mL, for G1364E fraction collectors	x	G1364-84545
Test tubes, 16 × 100 mm, 250/pk	x	5022-6532
Tray holding 215 tubes, 12 × 100 mm, 7 mL, for G1364E fraction collectors	x	G1364-84536
Test tubes, 12 × 100 mm, 250/pk	x	5022-6531
Tray holding 40 × 2 mL vials, for fraction collector G1364E/F and G5664B	x	G1364-44612
Tray, holding 15 × 6 mL vials, for fraction collector G1364E/F and G5664B	x	G1364-44613
Tray, holding 100 × 2 mL vials, for fraction collector G1364E/F and G5664B	x	G1364-44610
Tray for 100 × 2 mL vials, coolable		G1364-60111
Tray, thermostatable for 4 well plates, for use with fraction collectors G1364B/F and G5664B	x	G1364-60621
Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL	x	5190-8223
Check-out solution	x	5190-6886
Back pressure regulator, 20 psi	x	5067-6840
Exhaust tube adapter kit, for Agilent 1260 Infinity II Preparative/Analytical fraction collector, use with G1364E/F, G5664B		G1364-68709

	PM*	Part Number
InfinityLab Supplies for 1290 Infinity II Preparative OpenBed Fraction Collector (G7159B)		
Parts for preventive maintenance		
InfinityLab tubing kit, 8 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60953
InfinityLab tubing kit, 50 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60952
InfinityLab tubing kit, 150 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60954
InfinityLab tubing kit, 200 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector	1 each/option depending	G9321-60951
Dripping adapter, for fraction valve	1 each	G9321-26600
Containers and vial plates		
InfinityLab drawer, ambient, with RFID tag		G9321-60085
InfinityLab tube container, 10 tubes, 30 × 150 mm, ambient		G9321-60015
InfinityLab tube container, 18 tubes, 25 × 150 mm, ambient		G9321-60025
InfinityLab tube container, 36 tubes, 16 × 150 mm, ambient		G9321-60129
InfinityLab tube container, 72 tubes, 12 × 150 mm, ambient		G9321-60131
InfinityLab tube container, 10 tubes, 30 × 100 mm, ambient		G9321-60058
InfinityLab tube container, 18 tubes, 25 × 100 mm, ambient		G9321-60035
InfinityLab tube container, 36 tubes, 16 × 100 mm, ambient		G9321-60055
InfinityLab tube container, 72 tubes, 12 × 100 mm, ambient		G9321-60045
InfinityLab Drawer Gilson 207, ambient, contains two drawers for Gilson rack type 207, for use with 16 × 100 mm (od × length) test tubes		G9321-60201
InfinityLab Drawer Gilson 208, ambient		G9321-60208
InfinityLab SBS well plate container, ambient		G9321-60051
Vial plate, for 15 × 6 mL vials, low profile, nonstackable		5043-1826
Vial plate, for 40 × 2 mL vials		5023-2471
96-well plate, 2.0 mL, square wells, U shape, polypropylene, 41 mm, 30/pk		5043-9300
Sealing mat, 96 wells, square, preslitted, silicone, 50/pk		5043-9319
96-well plate, 0.90 mL, round wells, U shape, polypropylene, 32 mm, 50/pk		5043-9305

	PM*	Part Number
96-well plate, 1.0 mL, round wells, U shape, polypropylene, 27 mm, 50/pk		5043-9309
Sealing mat, 96 wells, round, preslitted, silicone, 50/pk		5043-9317
Fraction tubes and vials		
Test tubes, 30 × 150 mm, 78 mL, 100/pk		5190-9090
Test tubes, 25 × 150 mm, 55 mL, 100/pk		5190-9091
Test tubes, 16 × 150 mm, 21 mL, 250/pk		5190-9092
Test tubes, 12 × 150 mm, 11 mL, 250/pk		5190-9093
Round bottom glass vial, 30 × 100 mm 100/pk		5042-6458
Round bottom glass vial, 25 × 100 mm 100/pk		5042-6459
Test tubes, 16 × 100 mm, 250/pk		5022-6532
Test tubes, 12 × 100 mm, 250/pk		5022-6531
Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk		9301-1377
Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk		5188-5369
Cap, screw, for 6 mL vials, 100/pk		9301-1379
Septa, for 6 mL vials, 100/pk		9301-1378
Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk		5182-0716
Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk		5182-0715
Cap, screw, blue, preslit PTFE/silicone, 100/pk		5183-2076
Other supplies		
20 psi Backpressure regulator		5067-6840
Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL		5190-8223
Supplies for 1260 Infinity II Preparative Valve-Based Fraction Collector (G7166A)		
Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL		5190-8223
InfinityLab solvent bottle, clear 125 mL, with cap		9301-6527
InfinityLab solvent bottle, clear, 500 mL, with cap		9301-6523
InfinityLab solvent bottle, clear, 1,000 mL, with cap		9301-6528
InfinityLab identification silicone rings, 4 colors, 8/pk		9301-6529
InfinityLab stickers, for solvent bottles, 100/pk		9301-6530
InfinityLab fitting, for 2.3 mm od tubing, 2/pk		5043-1215

UV detectors

	PM*	Part Number
1260 Infinity II Diode Array & Multi Wavelength Detector (G7115A/G7165A)		
InfinityLab long-life deuterium DAD/MWD lamp, with RFID tag	1 each	2140-0820
Tungsten VIS lamp		G1103-60001
InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A), 3 mm path length		G7115-60001
InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A), 0.3 mm path length		G7115-60002
InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A) 0.06 mm path length		G7115-60003
Capillary, stainless steel, for preparative flow cells (DAD/MWD), use for flow rates 4 to 8 mL/min. Two capillaries required per flow cell.		5500-1342
Capillary, stainless steel, for preparative flow cells (DAD/MWD), use for flow rates 15 to 40 mL/min. Two capillaries required per flow cell.		5500-1343
Tubing, PTFE-ESD, 80 cm length, 0.5 mm id, 1.6 mm od, with fittings for preparative flow cell		G1315-67306
1260 Infinity II Variable Wavelength Detector (G7114A)		
InfinityLab long-life deuterium VWD lamp, with RFID tag	1 each	G1314-60101
Preparative flow cell for VWD, 3 mm path length, 2 µl cell volume, with RFID tag		G1314-60024
Preparative flow cell for VWD, 0.3 mm path length, with RFID tag		G1314-60025
Preparative flow cell for VWD, 0.06 mm path length, with RFID tag		G1314-60023
Preparative flow cell, Quartz, 3 mm, for VWD		G7114-60024
Supplies for 1260 Infinity II Evaporative Light Scattering Detector		
Splitter kit, 2 to 10 mL/min, analytical, for Agilent G4260B evaporative light scattering detector		5023-2256
Splitter kit, 8 to 40 mL/min, preparative, for Agilent G4260B evaporative light scattering detector		5023-2257

* PM (preventive maintenance) part information key

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

Column organizers

	PM*	Part Number
Supplies for Preparative Column Compartment (G7163B) /Column Organizer (G9328A)		
Column holder clips (lamella type), 2/pk		G7116-68003
Column holder clamp for columns 9.4 to 30 mm id		5067-6849
Column holder clamp, for columns 30 to 50 mm id		G7163-60002

Supporting modules

	PM*	Part Number
Supplies for 1290 Infinity II MS Flow Modulator (G7170B)		
Rebuild Kit, for flow modulator	1 each	1535-5215
InfinityLab capillary kit for mass-based fraction collection systems		5067-7023
Supplies for 1260 Infinity II Delay Coil Organizer (G9324A)		
InfinityLab knitted delay coils kit, for use in mass-based prep systems, for flow rate 4 to 8 mL/min or 9.4 mm id columns		5067-6800
InfinityLab knitted delay coils kit, for use in mass-based prep systems, for flow rates 15 to 40 mL/min or 21.2 mm id columns		5067-6801
InfinityLab knitted delay coils kit, for use in mass-based preparative systems, for flow rate 40 to 80 mL/min		5067-6802
InfinityLab knitted delay coils kit, for use in mass-based preparative systems		5067-6804

* PM (preventive maintenance) part information key

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

Valve supplies

	PM*	Part Number
Supplies for 2Pos/10Port Valve Head Prep Scale (G4730A)		
Rotor seal, five-groove, 600 bar, for 2pos/10 port preparative scale valve head (p/n 5067-4193)	1 each	5068-0153
Supplies for 8Pos/9Port Valve Head Prep Scale (G4731A)		
Rotor seal, single-groove, 600 bar, 2 pos/10 port preparative scale valve head (p/n 5067-4193)	1 each	5068-0155
Supplies for 6-Column Selection Valve Head Prep Scale (G4734A), obsolete		
Rotor seal, for the 6-column selector valve head, Preparative scale, 600 bar	1 each	5068-0249

	PM*	Part Number
Supplies for 6-Column Selection Valve Head Prep Scale (G4734B)		
Rotor, for use with 6-column selector valve head, preparative scale (p/n 5067-6722)	1 each	5068-0331
Bearing ring, for 2D-LC valve head, 1,300 bar		1535-4045
Capillary kit, for Preparative column selection valve, flow rate 4 to 8 mL/min		5067-6867
Capillary kit, for Preparative column selection valve, flow rate 15 to 40 mL/min		5067-6868
Capillary kit, for Preparative column selection valve, flow rates 40 to 80 mL/min		5067-6869
Capillary kit, for Preparative column selection valve, flow rates 80 to 200 mL/min		5067-6870
Supplies for 1260 Infinity II Clustering Valve (G9322A)		
Tubing, PTFE (ESD), 0.8 mm id, 1.6 mm od, 6 m		5043-1540
Capillary kit, for clustering up to 3 fraction collectors		5067-6811
Rail system, for fixation of clustering valve G9322A or recovery valve G7166A, on Agilent 1290 Infinity II preparative open-bed fraction collector G7159A/B		G9321-60013
Tubing guide assembly, for Agilent 1290 Infinity II preparative open-bed fraction collector G7159A/B		G9321-60026
Supplies for 2Pos/14Port Valve Head Prep (G4738A)		
Rotor Seal, 2 pos/14 port preparative valve head, 600 bar	1 each	5068-0323
InfinityLab upgrade capillary kit, for Autoscale Fraction Collection		5067-7019
Miscellaneous InfinityLab supplies		
System tool kit		G7120-68708
Preparative pressure relief valve kit		5067-6857
Back Pressure Regulator, 100 psi		5042-6443
Back pressure regulator, 20 psi		5067-6840
Stay Safe Caps and Waste Management		
InfinityLab Stay Safe cap GL45, 1 port for preparative LC		5043-1333
InfinityLab Stay Safe cap GL45, 2 ports for preparative LC		5043-1334
InfinityLab thread adapter, GL45(M) to GPI 38(F), PTFE		5043-1192
Fitting for 4.7 mm tubing, for prep LC Stay Safe caps, 2/pk		5043-1335
2 Screw plugs for Stay Safe caps prep LC		5005-2970
Kit, including InfinityLab Stay Safe cap (S60 thread, 4 ports) and 10 L waste can		5043-1338
Waste can, 10 L, with S60 thread		5043-1337
InfinityLab Stay Safe cap, S60 thread, 4 ports for waste can, including 2 x 3.2 mm od fitting, 2.3 mm od fitting, and 1.6 mm od fitting		5043-1336
InfinityLab charcoal filter with time strip, 58 g		5043-1193

	PM*	Part Number
Tubing and Adapters		
Tubing, PTFE (ESD), 1.6 mm od, 0.8 mm id, 6 m		5043-1540
Tubing, PTFE (ESD), 1.6 mm od, 0.5 mm id, 80 cm length, with fittings for preparative flow cell		G1315-67306
Tubing, PTFE, ESD-striped, 2.0 mm od, 1.2 mm id, 2 m length		5023-2878
Tubing, PTFE, ESD-striped, 2.5 mm od, 1.6 mm id, 6 m length		5023-2882
Fitting, 1/4-28, for ESD-PEEK tubing with 1.6 mm OD, 6/pk		5023-2898
Fitting, 1/4-28, for ESD-PEEK tubing with 2.0 mm OD, 6/pk		5023-2899
Fitting, 1/4-28, for ESD-PEEK tubing with 2.5 mm OD, 6/pk		5023-2883
Fitting, for ESD-PEEK tubing, 5/16 to 3/16 inch, 12/pk		5043-1847
Union, flat-bottom, 1/4-28 ESD		5023-2508
Fitting adapter, 1-4 28 to 10-32 coned, ESD protective material		5500-1410
Adapter nut, flat bottom, 1/4-28-M, 5/16-24-F, stainless steel		5023-3093
Adapter, stainless steel, female 10-32 to male 5/16-24		5067-6845
Adapter, stainless steel, female to male 10-32 to 1/4-28		5023-1803
Adapter, stainless steel, 10-32 female to 1/4-28 female, for prep LC		5023-2875
Adapter, stainless steel, female 1/4-28 to male 10-32, for prep LC		5023-2512
System Capillaries		
Capillary, stainless steel, 0.12 × 70 mm, SI/SI		5500-1329
Capillary, stainless steel, 0.12 × 300 mm, NX/NX		5500-1432
Capillary, stainless steel, 12 × 350 mm, SX/U6		5500-1433
Capillary, stainless steel, 0.12 × 2000 mm, SX/U6		5500-1434
Capillary, stainless steel, 0.17 × 800 mm, NX/NX		5500-1431
Capillary, stainless steel, 0.17 × 800 mm, SX		5500-1430
Capillary, stainless steel, 0.17 × 1150 mm, SX		5500-1423
Capillary, stainless steel, 0.17 × 1200 mm, NX/NX		5500-1462
Capillary, stainless steel, 0.25 × 220 mm, NX/NX		5500-1425
Capillary, stainless steel, 0.25 × 700 mm, SI/SI		5500-1330
Capillary, stainless steel, 0.25 × 800 mm, NX/NX		5500-1461
Capillary, stainless steel, 0.25 × 1100 mm, NX/NX		5500-1422
Capillary, stainless steel, 0.25 × 1200 mm, NX/NX		5500-1424
Capillary, stainless steel, 1.0 × 1900 mm, SX/SI		5500-1415

	PM*	Part Number
0.3 mm id capillaries		
Capillary, stainless steel, 0.3 × 500 mm, SX/-		5500-1316
Capillary, stainless steel, 0.3 × 500 mm, long socket, no fittings		5500-1448
Capillary, stainless steel, 0.3 × 550 mm, SX/-		5500-1449
Capillary, stainless steel, 0.3 × 550 mm, SX/-/SI		5500-1408
Capillary, stainless steel, 0.3 × 650 mm, SX/-		5500-1314
Capillary, stainless steel, 0.3 × 650 mm, SX/-/SX		5500-1308
Capillary, stainless steel, 0.3 × 750 mm, SX/SI		5500-1437
Capillary, stainless steel, 0.3 × 800 mm, SX/-		5500-1406
Capillary, stainless steel, 0.3 × 900 mm, no fittings		5500-1450
Capillary, stainless steel, 0.3 × 1000 mm, SX/-		5500-1407
Capillary, stainless steel, 0.3 × 1150 mm, SX		5500-1435
Capillary, stainless steel, 0.3 × 1450 mm, NX/NX		5500-1436
Capillary, stainless steel, 0.3 × 1900 mm, SX/SI		5500-1412
0.5 mm id capillaries		
Capillary, stainless steel, 0.5 × 150 mm, no fittings		5022-6509
Capillary, stainless steel, 0.5 × 280 mm, no fittings		5500-1336
Capillary, stainless steel, 0.5 × 400 mm, no fittings		5500-1334
Capillary, stainless steel, 0.5 × 500 mm, no fittings		5500-1451
Capillary, stainless steel, 0.5 × 500 mm, SX/-		5500-1317
Capillary, stainless steel, 0.5 × 550 mm, SX/-/SI		5500-1386
Capillary, stainless steel, 0.5 × 550 mm, SX/-		5500-1452
Capillary, stainless steel, 0.5 × 600 mm, no fittings		5500-1333
Capillary, stainless steel, 0.5 × 650 mm, SX/-		5500-1315
Capillary, stainless steel, 0.5 × 650 mm, SX/-/SX		5500-1309
Capillary, stainless steel, 0.5 × 750 mm, SX/SI		5500-1440
Capillary, stainless steel, 0.5 × 800 mm, no fittings		5500-1335
Capillary, stainless steel, 0.5 × 800 mm, SX/-		5500-1384
Capillary, stainless steel, 0.5 × 900 mm, no fittings		5500-1453
Capillary, stainless steel, 0.5 × 1000 mm, SX/-		5500-1385
Capillary, stainless steel, 0.5 × 1150 mm, SX/-		5500-1438
Capillary, stainless steel, 0.5 × 1450 mm, NX/NX		5500-1439
0.6 mm id capillaries		
Capillary, stainless steel, 0.6 × 100 mm, SX/SX		5500-1427
Capillary, stainless steel, 0.6 × 500 mm, SX/-		5500-1320
Capillary, stainless steel, 0.6 × 500 mm, no fittings		5500-1454
Capillary, stainless steel, 0.6 × 550 mm, SX/-		5500-1455
Capillary, stainless steel, 0.6 × 650 mm, SX/-/SX		5500-1318
Capillary, stainless steel, 0.6 × 650 mm, SX/-		5500-1319
Capillary, stainless steel, 0.6 × 800 mm, SX/-		5500-1387
Capillary, stainless steel, 0.6 × 900 mm, no fittings		5500-1456
Capillary, stainless steel, 0.6 × 1000 mm, SX/-		5500-1388
Capillary, stainless steel, 0.6 × 1450 mm, NX/NX		5500-1442
Capillary, stainless steel, 0.6 × 1900 mm, SX/SI		5500-1413

	PM*	Part Number
0.7 mm id capillaries		
Capillary, stainless steel, 0.7 × 550 mm, SX/-/SI		5500-1389
Capillary, stainless steel, 0.7 × 750 mm, SX/SI		5500-1443
Capillary, stainless steel, 0.7 × 1900 mm, SX/SI		5500-1414
0.94 mm id capillaries		
Capillary, stainless steel, 0.94 × 60 mm, SI/-		5500-1447
Capillary, stainless steel, 0.94 × 500 mm, no fittings		5500-1457
Capillary, stainless steel, 0.94 × 500 mm, SX/-		5500-1391
Capillary, stainless steel, 0.94 × 550 mm, SX/-		5500-1458
Capillary, stainless steel, 0.94 × 550 mm, SX/-/SI		5500-1395
Capillary, stainless steel, 0.94 × 650 mm, SX/-		5500-1393
Capillary, stainless steel, 0.94 × 650 mm, SX/-/SX		5500-1390
Capillary, stainless steel, 0.94 × 750 mm, SX/SI		5500-1446
Capillary, stainless steel, 0.94 × 800 mm, SX/-		5500-1392
Capillary, stainless steel, 0.94 × 900 mm, no fittings		5500-1459
Capillary, stainless steel, 0.94 × 1000 mm, SX/-		5500-1394
Capillary, stainless steel, 0.94 × 1150 mm, SX/-		5500-1444
Capillary, stainless steel, 0.94 × 1450 mm, NX/NX		5500-1445

Capillary fitting codes key

SI fitting = Stainless steel fitting with short screw

SX fitting = Stainless steel fitting with extra long screw

U6 = Stainless steel female 10-32 port thread size connector

NX = Stainless steel fitting with long screw with SW5 hex head

www.agilent.com/chem

DE.445

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
 Printed in the USA, December 21, 2020
 5994-2810EN