

Agilent InfinityLab Supplies for the Agilent Infinity II Bio LC

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 Agilent Infinity II Bio LC.

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?



Supplies for 1290/1260 Infinity II Bio Flexible Pumps (G7131A/C)	PM*	Part No.
Standard preventive maintenance parts		
PM Kit 1290 Inf II Bio Flexible Pumps	kit	G7131-68740
Pump seal	2 each	G7131-20009
Seal Wash seal	2 each	0905-1731
Peristaltic pump	1 each	5065-4445
Frit 1290 Bio Pump Outlet Filter 2/pk	1 each	5320-0048
Extended preventive maintenance parts		
1290 Passive Inlet Valve Bio		G7131-60021
Outlet Valve Universal		G7131-60028
Inlet Valve Universal		G7131-60022
Plunger Assembly		5067-5975
Rotor Seal		5068-0379
Other supplies		
Jet Weaver Mixer Kit V35 Flex Bio-C		G7131-68035
Jet Weaver Mixer Kit V100 Flex Bio-C		G7131-68100
Jet Weaver Mixer Kit V380 Flex Bio-C		G7131-68380
Jet Weaver Connection Capillary		5004-0041
Bio InLine Filter Upgrade Kit		G7131-68000

Supplies for 1290 Infinity II Bio High-Speed Pump (G7132A)	PM*	Part No.
Standard preventive maintenance parts		
PM Kit 1290 Inf II Bio High Speed Pump	kit	G7132-68740
Pump seal	4 each	G7131-20009
Peristaltic pump	1 each	5065-4445
Seal Wash seal	4 each	0905-1731
PTFE frits (pack of 5)	1 each	01018-22707
Seal cap assembly	2 each	5067-4728
Extended preventive maintenance parts		
Rotor Seal for Purge Valve (5pos/7port)		5068-0377
High pressure filter		G7132-60026
1290 Passive Inlet Valve Bio		G7131-60021
Outlet valve		G7131-60028
Plunger Assembly		5067-5975

* 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

Supplies for 1290 Infinity II Bio High-Speed Pump (G7132A) Part No.	
Other supplies	
Jet Weaver Mixer Kit V380 HS Bio-C	G7132-68380
Jet Weaver Mixer Kit V35/V100 HS Bio-C	G7132-68135
Capillary, pump head to purge valve (MP35N 0.25x250mm SI/SI)	5500-1420
Capillary, purge valve to pressure sensor (MP35N 0.25x130mm SI/SI)	5500-1421
Capillary, pump to multisampler (MP35N 0.17x500mm) SI/SI	G7132-68135

Supplies for 1290 Infinity II Bio Multisampler (G7137A) Part No.	
Parts for preventive maintenance	
PM Kit 1290 Inf II Bio Multisampler	G7137-68740
Needle Bio-compatible	G7137-87201
High-Pressure Seat Bio-Compatible 0.12	G7137-87012
Rotor Seal	5320-0010
High-Pressure Seat Bio Ultra Low Dispersion	G7137-60075

Extended preventive maintenance parts	
Metering Seal 1290 Bio 2 mm Piston	G7137-20003

Other supplies	
Bio Sample Loop 20 µL	G7137-60300
Bio Sample Loop Flex 40 µL	G7137-60400
Bio Sample Loop 100 µL	G7137-60500
Capillary, Sampler to MCT (MP35N 0.12 x 500 mm SI-SI)	5500-1279

Vials and Wellplates	
96 well plate, 1.2ml, round wells, U shape, polypropylene, 27mm high, 25/pk	5043-9308
Sealing mat, 96 wells, round, preslitted, silicone, 50/pk	5043-9317
2 mL Clear Glass Vial (deactivated)	5183-2071
2 mL PP Vial (holds 1.7 mL fill volume – similar to 2 mL glass vial)	5191-8151
250 µL PP screw vial (shape of 2 mL vial)	5190-2242
9 mm bonded screw cap with pre-slit septum (PTFE/silicone)	5190-7023
11 mm crimp cap with pre-slit septum (PTFE/silicone)	5190-9045
High Recovery glass Vial (Silanized)	5183-4497
250 uL PP insert - (needs 2 mL vial)	5190-4073
400 uL flat bottom glass insert (deactivated) - (needs 2 mL vial)	5183-2086
250 uL glass insert (deactivated) with polymer feet - (needs 2 mL vial)	5181-8872
250 µL PP insert with polymer feet - (needs 2 mL vial)	5182-0549

Multisampler Drawers and Vial plates	
Additional standard drawer kit, double-height (1/pk)	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

Supplies for Multi Column Compartment Part No.	
Quick-Connect Heat-Exchanger 1290 Bio Standard Flow (0.12 mm ID, 1.6 µL)	G7116-60071
Quick-Connect Heat-Exchanger 1290 Bio High Flow (0.12mm/3.0µL)	G7116-60081
Quick Connect Heat Exchanger 1290 Bio Ultra Low Dispersion (0.075 mm /0.6 µL)	G7116-60091
Agilent Thermal Equilibration Device	G7116-60013

Supplies for Valves Part No.	
2 Pos / 10 Port Valve Bio (G5641A)	5067-6682
Rotor Seal	5068-0205
Stator	5068-0286
6-Column Selector Valve Bio (G5640A)	5320-0025
Rotor Seal	5320-0029
Stator	5068-0019

Supplies for Detectors Part No.	
DAD supplies	
Long-life deuterium lamp, 8-pin, with RFID tag	5190-0917
Max-Light Cartridge Cell LSS 10mm	G7117-60020
Max-Light cartridge test cell	G4212-60011
VWD supplies	
Deuterium lamp, with RFID tag	G1314-60101
Bio Micro Flow Cell VWD, 3 mm, 2 µL, RFID tag	G1314-60189
Bio Standard Flow Cell VWD,10 mm, 14 µL, RFID tag	G1314-60188

Kits Part No.	
Multidraw Kit for Bio Multisampler	G7137-68711
Ultra Low Dispersion Kit Bio	5004-0007
Bio Infinity II Inline Filter Kit	5720-0020
Replacement frits for Inline Filter Kit (pack of 5)	5320-0022

Fittings & Unions Part No.	
Quick Connect Fitting	5067-5965
Quick Turn Fitting	5067-5966
Blank Nut long PEEK-SST, for module diagnostic tests	5043-0277
Fitting Union MP35N (10-32 port size)	5023-2625
Restriction Capillary MP35N 0.12 x 2000mm	5005-0046

Capillaries	Part No.
Capillaries for Quick Connect Fitting	
Quick Connect Capillary MP35N 0.12 x 105	5500-1578
Quick Connect Capillary MP35N 0.12 x 150	5500-1579
Quick Connect Capillary MP35N 0.12 x 220	5500-1580
Quick Connect Capillary MP35N 0.12 x 280	5500-1581
Quick Connect Capillary MP35N 0.12 x 400	5500-1582
Quick Connect Capillary MP35N 0.12 x 500	5500-1583
Quick Connect Capillary MP35N 0.17 x 105	5500-1584
Quick Connect Capillary MP35N 0.17 x 150	5500-1585
Quick Connect Capillary MP35N 0.17 x 220	5500-1586
Quick Connect Capillary MP35N 0.17 x 280	5500-1587
Quick Connect Capillary MP35N 0.17 x 500	5500-1588
Capillaries for use with Quick Turn and other bio-compatible fitting	
Quick Turn Capillary MP35N 0.075 x 105	5500-1589
Quick Turn Capillary MP35N 0.075 x 150	5500-1590
Quick Turn Capillary MP35N 0.075 x 250	5500-1591
Quick Turn Capillary MP35N 0.075 x 500	5500-1592
Quick Turn Capillary MP35N 0.12 x 105	5500-1593
Quick Turn Capillary MP35N 0.12 x 150	5500-1594
Quick Turn Capillary MP35N 0.12 x 200	5500-1595
Quick Turn Capillary MP35N 0.12 x 280	5500-1596
Quick Turn Capillary MP35N 0.12 x 400	5500-1597
Quick Turn Capillary MP35N 0.12 x 500	5500-1598
Quick Turn Capillary MP35N 0.17 x 105	5500-1599
Quick Turn Capillary MP35N 0.17 x 150	5500-1600
Quick Turn Capillary MP35N 0.17 x 200	5500-1601
Quick Turn Capillary MP35N 0.17 x 280	5500-1602
Quick Turn Capillary MP35N 0.17 x 400	5500-1603
Quick Turn Capillary MP35N 0.17 x 500	5500-1604
Quick Turn Capillary MP35N 0.17 x 700	5500-1605

General supplies	Part No.
System tool kit	G7120-68708
InfinityLab Solvent filtration assembly	5191-6776
InfinityLab solvent filtration flask, glass, 2 L	5191-6781
Filter membrane, Regenerated Cellulose 47 mm, pore size 0.2 µm, 100/pk	5191-4340
Filter membrane, Regenerated Cellulose 47 mm, pore size 0.45 µm, 100/pk	5191-4337

Solvent Cabinet and Waste Management	Part No.
Bottle head assembly	G7120-60007
Tubing Kit, 140 mm, 2/pk, SSV to shutoff valve or degassing unit (high speed pump), degasser to MCGV (flexible pump)	G4220-60035
Bottle head assembly, ultra clean tubing quality, recommended for MS application	G7120-60017
Tubing kit, ultra clean quality, recommended for MS application, includes 4x bottle head assemblies (G7120-60017) and 3x tubing kits 140 mm (G4220-60070)	G7120-68070
Glass filter, solvent inlet, 20 µm pore size	5041-2168
Tubing for seal wash, silicone, 5 m	5065-9978
Tubing, purge valve to waste, PTFE, 5 m	5062-2462
Solvent bottle, transparent 1 L	9301-6524
Solvent bottle, amber 1 L	9301-6526
Solvent bottle, transparent 1 L, with cap	9301-6528
Solvent bottle, clear, 500 mL with cap	9301-6523
Solvent bottle, amber, 500 mL with cap	9301-6525
InfinityLab Stay Safe purging bottle	5043-1339
Identification rings (silicone) for solvent bottles (8 pcs/4 colors)	9301-6529
Sticker 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 fitting 3.2 mm	5043-1217
Venting valve for InfinityLab Stay Safe cap	5043-1190
10 L waste container and Stay Safe cap S60 with 4 ports for waste	5043-1338
InfinityLab waste canisters 5L, GL38	5043-1986
6 L waste container including one Stay Safe Cap GL45 w. 4 ports	5043-1221
Charcoal filter (48 g) for waste container	5043-1193



For a complete list of Agilent columns and supplies, visit us online at:

www.agilent.com/chem/lc-supplies

www.agilent.com/chem/biopharma-hplc-selector

InfinityLab Family: Maximize the efficiency of your LC Workflows

The InfinityLab family is an optimized portfolio of LC instruments, columns, and supplies, designed to work in perfect harmony. Combined with Agilent OpenLAB software and Agilent CrossLab Services, Agilent InfinityLab provides you with end-to-end support to make every day more productive.

Explore the Agilent InfinityLab LC Solutions

www.agilent.com/chem/infinitylab

DE44404.5460532407

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
Printed in the USA, July 26, 2021
5994-3948EN

