

# OPTIMIZE YOUR LABORATORY'S PRODUCTIVITY WITH AGILENT 1260 INFINITY PURIFICATION LC SUPPLIES

The Measure of Confidence

## Quick reference guide

### Agilent supplies for Agilent instruments

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 Purification LC.

Keep this list handy so you can quickly find the supplies you need and minimize instrument downtime.

Pump supplies	Part No.
Piston seal, for G1361A preparative pump	5022-2188
Sapphire plunger	G1361-22402
Preventative maintenance kit for 1100/1200 Infinity Series and 1260 Infinity LC prep pump. Includes filter cup (3150-0942), 4 seal prep flanges (5022-2188), filter assembly (5022-2192), peristaltic pump (0890-1764) tubing, 4 prep pump seals	G1361-68710
Accessory kit, prep pump/gradient, for G1361A. Includes stainless steel connecting capillaries, solvent mixer, 2 L solvent bottle, bottle head assembly, filter, glass stop valve, stainless steel union, tubing, and other parts	G1361-68707

Prep autosampler supplies	Part No.
Needle assembly, for G2260A	G2260-87201
Needle seat, prep autosampler, 0.5 mm id, 20 µL	G2260-87101
Valve assembly, inlet/outlet single seat	G1361-60012
Valve assembly, inlet/outlet double seat	G1361-60052

Dual loop autosampler supplies	Part No.
Peristaltic pump for 1100/1200 Infinity Series LC	5065-4445
Rotor seal 10-port valve, high flow (Vespel)	0100-2415

Fraction collector supplies	Part No.
Preparative tubing kit, PTFE, 0.8 mm id	G1364-68711
Analytical tubing kit, PTFE, 0.25 mm id	G1364-68712
Low-dispersion tubing kit, PTFE, 0.15 mm id	G1364-68723



The Measure of Confidence



**Agilent Technologies**

Valves	Part No.
External switching valve, 12-position/13-port preparative solvent selection valve, for G1160A	0101-1288 (RheBuild Kit)
Stator, for 0101-1288	0101-1365
Manual injection valve with position sensing switches, 3725i, stainless steel prep valve	0101-1232
Rotor seal, for 0101-1232	0101-1233
Manual prep injection valve kit, stainless steel. Includes position sensing, 10 mL loop, 25 mL syringe, ring mounting bracket, start cable, connecting capillaries, 0.5 mm id, 40 cm and 60 cm	5065-9922

Detector supplies	Part No.
DAD preparative flow cell, for G1315A/B, G1365A/B, 3 mm, 120 bar, stainless steel	G1315-60016
Repair kit, for G1315-60016	G1315-68712
DAD preparative flow cell, 0.3 mm, 20 bar, quartz	G1315-60017
DAD preparative flow cell, 0.06 mm, 20 bar, quartz	G1315-60018

Connectors	Part No.
Capillary, stainless steel, from injector valve to prep head, 0.5 mm id, 160 mm	G2258-87301
High-flow union, no fittings	5022-2133

Filters	Part No.
Bottle head assembly	G1361-60022
Glass filter, solvent inlet, 40 µm pore	3150-0944
Frit adapter, PTFE, 4 mm	G1361-23204
Semi-prep inline filter, 0.5 µm pore, 12.7 mm id, 1 to 5 mL/min	5064-8273
Replacement frits, for 5064-8273	5022-2185
High pressure semi-prep filter, 10 µm pore, 19 mm id, 5 to 10 mL/min	5022-2165
Replacement frits for 5022-2165, 10/pk	5022-2166
Prep filter, 10 µm pore, 10 to 100 mL/min	5065-4500
Replacement glass cartridge, for 5065-4500	5065-9901

## Agilent supplies - to help you get the results you need

For greater confidence in your chromatography results, you can optimize your total system with quality Agilent columns and supplies. Our wide range of columns, supplies, sample prep, kits and accessories is 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 Agilent's complete listing of LC supplies, including our own range of genuine Agilent parts and supplies for our LC product range, visit us online at:

[www.agilent.com/chem/supplies](http://www.agilent.com/chem/supplies)

This information is subject to change without notice.

© Agilent Technologies, Inc. 2014  
Published in UK, February 13, 2014  
5991-3965EN



**Agilent Technologies**