

Agilent InfinityLab LC/MSD Series Supplies

Make sure that you don't run out of essential columns and supplies

Use this handy list to keep your shelves stocked—and keep your lab productive.



Agilent LC/MSD iQ	Part Number
Ion injector capillary	
Ultivo ion injector	G6301-80004
Narrow bore capillary cap	G1960-20310
Ionization source	
ES nebulizer needle kit for ESI/AJS	G1958-60137
Nebulizer assembly (ESI/AJS)	G1958-60098
Spray shields	
Spray shield (ESI/APCI)	G1946-20157
Desolvation assembly, ion optics, and vacuum manifold	
High gain electron multiplier horn	G2571-80103
Skimmer, 2 mm diameter	G1969-20302
Calibrant delivery system	
ESI-L tuning mix	G1969-85000
50 mL CDS bottle + cap	G6301-80019
Rough pump (MS40+)	
AFV 60 gold oil	6040-1444
Oil mist filter, MS40+	G1960-80039
Gas purifiers	
Nitrogen filter for main gas supply	RMSN-4
PM kits	
Ultivo PM kit	G6303-68010

Agilent LC/MSD	Part Number
Ion injector capillary	
0.6 mm id dielectric capillary	G7604-60000
Narrow bore capillary cap	G1960-20310
Ionization source	
ES nebulizer needle kit for ESI/AJS	G1958-60137
APCI nebulizer needle kit	G1946-68704
Corona needle	G1947-20029
Nebulizer assembly (ESI/AJS)	G1958-60098
Nebulizer assembly (APCI)	G1946-67037
Spray shields	
AJS spray shield small	G1958-20050
Spray shield (ESI/APCI)	G1946-20157
Multimode spray shield	G1978-60060
Desolvation assembly, ion optics, and vacuum manifold	
Low gain EM replacement horn, LCMS	G1960-80103
Skimmer 1	G1956-20302
Calibrant delivery system	
ESI tuning mix	G2421-60001
APCI tuning mix	G2432-60001
Glass CDS bottle	9300-2576
Glass CDS bottle cap	9300-2575

Agilent LC/MSD	Part Number
Rough pump (MS40+)	
AFV 60 gold oil	6040-1444
Oil mist filter, MS40+	G1960-80039
Gas purifiers	
Nitrogen filter for main gas supply	RMSN-4
PM kits	
LCMS PM kit, ESI, MS40+ RP	5190-9007

Agilent LC/MSD XT	Part Number
Ion injector capillary	
0.6 mm id fast switch capillary	G1960-80060
Narrow bore capillary cap	G1960-20310
Ionization source	

ES nebulizer needle kit for ESI/AJS	G1958-60137
APCI nebulizer needle kit	G1946-68704
Corona needle	G1947-20029
Nebulizer assembly (ESI/AJS)	G1958-60098
Nebulizer assembly (APCI)	G1946-67037

Spray shields	
AJS spray shield small	G1958-20050
Spray shield (ESI/APCI)	G1946-20157
Multimode spray shield	G1978-60060

Desolvation assembly, ion optics, and vacuum manifold	
Low gain EM replacement horn, LCMS	G1960-80103
Skimmer 1	G1956-20302

Calibrant delivery system	
ESI tuning mix	G2421-60001
APCI tuning mix	G2432-60001
ESI-L tuning mix (AJS)	G1969-85000
Glass CDS bottle	9300-2576
Glass CDS bottle cap	9300-2575

Rough pump (MS40+)	
AFV 60 gold oil	6040-1444
Oil mist filter, MS40+	G1960-80039

Gas purifiers	
Nitrogen filter for main gas supply	RMSN-4

PM kits	
LCMS PM kit, AJS, 1 MS40+ RP	5190-9008

Common-Use Parts	Part Number
Tools	
1/4 inch × 5/16 inch open-end wrench	8710-0510
1/2 inch × 9/16 inch open-end wrench	8710-0877
9/64 inch hex key	8710-2394
0.8 mm hex key	8710-1225
2 mm hex key	9810-0641
3 mm hex key, 12 cm long	8710-2411
10 mm hex key	8710-2612
4000 grit abrasive paper	8660-0827
Clean gloves	8650-0030
Clean chemical-resistant gloves	9300-1751
Clean lint-free cloth 15/pk	05980-60051
Cotton swab 100/pk	0580-5400
Long flat-blade screwdriver	8710-1029
Long-nose pliers	7810-0004
Nebulizer adjustment kit	G1960-67470
T10 torx screwdriver	8710-1623
T20 torx screwdriver	8710-1615
Alconox powdered precision cleaner	5190-1401

Reagents	
LC/MS-grade or de-ionized (18 MΩ/cm) water (4 L)	8500-2236
HPLC flushing solvent	G1969-85026
LC/MS-grade acetonitrile (1 L)	G2453-85060
LC/MS-grade methanol (1 L)	8500-1867
Formic acid (5 mL)	G2453-85060
InfinityLab Ultrapure LC/MS solvent 1 L: acetonitrile	5191-4496
InfinityLab Ultrapure LC/MS solvent 1 L: methanol	5191-4497
InfinityLab Ultrapure LC/MS solvent 1 L: water	5191-4498

Calibrant delivery system	
Rotor seal	0101-1409
CDS PEEK tubing	0100-1919

Desolvation assembly, ion optics, and vacuum manifold	
Nitrogen filter for main gas supply	RMSN-4
Canted coil spring (4 required)	G2571-67001
NW 20/25 clamp ring (MS40+)	0100-0549
NW 25 centering O-ring (MS40+)	0100-1551



Agilent CrossLab services. From insight to outcome.

CrossLab is an Agilent capability that integrates services and consumables to support workflow success and important outcomes like improved productivity and operational efficiency. Through CrossLab, Agilent strives to provide insight in every interaction to help you achieve your goals. CrossLab offers method optimization, flexible service plans, and training for all skill levels. We have many other products and services to help you manage your instruments and your lab for best performance.

Learn more about Agilent CrossLab, and see examples of insight that leads to great outcomes, at www.agilent.com/crosslab

To place your order now, go to:
www.agilent.com/chem/lc-ms-supplies

DE.4968171296

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020
Published in the USA, May 20, 2020
5990-8580EN