

# Agilent Triple Quadrupole LC/MS 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 Ultivo LC/TQ	Part Number
<a href="#">Ion injector capillary</a>	
Ion injector	G6301-80004
Narrow-bore capillary cap	G1960-20310
<a href="#">Desolvation assembly, ion optics, and vacuum manifold</a>	
Canted coil spring (4)	G2571-67001
High-gain electron multiplier horn	G2571-80103
Skimmer, 2 mm	G1969-20302
AJS spray shield, small	G1958-20050
Endcap assembly, multimode	G1978-60060
<a href="#">Calibrant delivery system</a>	
50 mL CDS bottle + cap	G6301-80019
<a href="#">Rough pump (MS40+)</a>	
Nitrogen filter for main gas supply (2)	RMSN-4
AFV 60 gold oil	6040-1444
Oil mist filter, MS40+	G1960-80039
<a href="#">Gas purifiers</a>	
Nitrogen filter for main gas supply (2)	RMSN-4
Nitrogen gas filter for collision cell gas supply	RMSN-2

Agilent 6495C LC/TQ	Part Number
<a href="#">Ion injector capillary</a>	
0.8 mm id capillary cap	G1964-20028
Capillary, 6 x 0.61 mm x 90 mm, 1G	G1964-80659
<a href="#">Desolvation assembly, ion optics, and vacuum manifold</a>	
Canted coil spring (4)	G2571-67001
High-gain electron multiplier horn	G2571-80103
AJS spray shield, multi bore	G1964-20050
Endcap assembly, multimode	G1978-60060
<a href="#">Calibrant delivery system</a>	
Glass CDS bottle	9300-2576
Glass CDS bottle cap	9300-2575
<a href="#">Rough pump (MS120)</a>	
AVF 60 gold oil (2)	6040-1444
MS120 exhaust filter kit	X3702-68201
<a href="#">Gas purifiers</a>	
Nitrogen filter for main gas supply (2)	RMSN-4
Nitrogen gas filter for collision cell gas supply	RMSN-2

Agilent 6470B LC/TQ	Part Number
<a href="#">Ion injector capillary</a>	
0.6 mm id fast switch capillary	G1960-80060
Narrow-bore capillary cap	G1960-20310
Ion Injector, 180 mm, 0.6 mm id, Dielectric	G7604-60000
<a href="#">Desolvation assembly, ion optics, and vacuum manifold</a>	
Canted coil spring (4)	G2571-67001
High-gain electron multiplier horn	G2571-80103
Skimmer, 2 mm	G1969-20302
AJS spray shield, small	G1958-20050
Endcap assembly, multimode	G1978-60060
<a href="#">Calibrant delivery system</a>	
Glass CDS bottle	9300-2576
Glass CDS bottle cap	9300-2575
<a href="#">Rough pump (MS40+)</a>	
AFV 60 gold oil (2)	6040-1444
Oil mist filter, MS40+	G1960-80039
<a href="#">Gas purifiers</a>	
Nitrogen filter for main gas supply (2)	RMSN-4
Nitrogen gas filter for collision cell gas supply	RMSN-2
<b>Common-Use Parts</b>	
<a href="#">Tools</a>	
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

Common-Use Parts	Part Number
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
<a href="#">Reagents</a>	
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)	G2453-85050
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
<a href="#">Calibrant delivery system</a>	
APCI-L tuning mix	G1969-85010
ESI-L tuning mix	G1969-85000
MMI-L tuning mix	G1969-85020
Rotor seal	0101-1409
CDS PEEK tubing	0100-1919
<a href="#">Ionization source</a>	
ES nebulizer needle kit for AJS	G1958-60137
APCI nebulizer needle kit	G2428A
Corona needle	G1947-20029
Nebulizer assembly (ESI/AJS)	G1958-60098
Nebulizer assembly (APCI)	G1946-67037
<a href="#">PM kits</a>	
LC/MS PM kit, Jet Stream, 1 MS40+ RP	5190-9008
LC/MS PM kit, Jet Stream, 2 MS40+ RPs	5190-9009
LC/MS PM kit, MS120 RP	G2571-68003
Ultivo PM kit	G6301-68010



## 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](http://www.agilent.com/crosslab)

To place your order now, go to:  
[www.agilent.com/chem/lc-ms-supplies](http://www.agilent.com/chem/lc-ms-supplies)

DE.5424652778

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021  
Published in the USA, August 30, 2021  
5990-8583EN