

Agilent 9500 ICP-MS Supplies

Agilent supplies for Agilent instruments

Agilent is committed to optimizing your laboratory's productivity. This reference guide shows the most commonly ordered supplies and parts for the Agilent 9500 ICP-MS/MS. Keep this list so you can quickly find the supplies you need and minimize instrument downtime.



Sampler Cone Care Packages	
Each package includes 2 pcs of the sampler cone, 1 pc magnifier tool, 1 pk graphite gaskets, and 1 pk cotton swabs	Part Number
Ni-tip Cu base Easy-Fit sampler cone care package	M5150-67100
Ni-tip with Ni-plated base Easy-Fit sampler cone care package	M5150-67101
Pt-tip Cu base Easy-Fit sampler cone care package	M5150-67102
Pt-tip with Ni-plated base Easy-Fit sampler cone care package	M5150-67103

Sampler Cone and Supplies	Part Number
Easy-Fit Sampler cone, Ni-tip with Cu base (std for u-lens)	M5150-67000
Easy-Fit Sampler cone, Ni-tip with Ni-plated base (for sample containing > 0.5% HCl or when using UHMI aerosol dilution > 25)	M5150-67001
Easy-Fit Sampler cone, Pt-tip with Cu base (for HF digests or organic solvents)	M5150-67002
Easy-Fit Sampler cone, Pt-tip with Ni-plated base (for aggressive acid matrix, esp. HCl, HF, and HClO ₄)	M5150-67003
Graphite gasket for 9500 ICP-MS Sampler Cone, 3/pk	M5150-67029
Magnifier Tool	5190-9614
Cotton swabs, 1pk of 100 pcs	9300-2574

Skimmer Cones (Selected by Ion Lens Configuration)		
Description	9500 with u-lens	9500 with m-lens
Nickel (for common aqueous applications)	M5150-67005	G3666-67511
Platinum, with copper base (for aggressive acid digests)	M5150-67006	
Platinum, with nickel base (for organic solvent analysis with Oxygen/Argon option gas)	M5150-67007	G3666-67501

Skimmer Bases (Selection First by Ion Lens Configuration, then by the Interface Cone Tip Material (Nickel, Ni or Platinum, Pt))		
Description	9500 with u-lens	9500 with m-lens
Stainless Steel (use with Ni-tipped cones)	M5150-67018	M5150-67020
Brass (use with Pt-tipped cones)	M5150-67019	M5150-67021

Easy-Fit and Extraction-Omega Lens Assemblies	
Description	9500 with u-lens
Easy-Fit u-lens, single-piece assembly, provided in clean packaging, ready to install. For direct replacement of 9500 ICP-MS u-lens. Does not include skimmer base.	M5150-67104
Extraction-Omega Lens assembly for u-lens of 9500 ICP-MS. Includes Extraction Lens 1, Extraction Lens 2, Omega Lens, Omega Bias assembled with screws and spacers onto a stainless-steel skimmer base.	M5150-67022
Extraction-Omega Lens assembly for m-lens of 9500 ICP-MS. Includes Extraction Lens 1, Extraction Lens 2, Omega Lens, Omega Bias assembled with screws and spacers onto a brass skimmer base.	M5150-67023

Easy-Fit Torches	Part Number
Easy-Fit Torch, Quartz, 2.5 mm id for aqueous (standard)	M5150-67011
Easy-Fit Torch, Quartz, 1.5 mm id for organic solvents	M5150-67012

Grounding Manifolds	Part Number
Grounding manifold with 2 inlets for sample and internal standard online mixing (standard)	M5150-67026
Grounding manifold with 1 inlet for sample. Optional configuration recommended when using AVS MS or when internal standard online addition is not needed.	M5150-67027
Thumb screw for grounding manifold	M5150-67028

Nebulizers and Supplies	Part Number
MicroMist nebulizer, borosilicate glass with PEEK gas connector. Suits Agilent 9500 ICP-MS	M5150-67024
MicroMist nebulizer gas fitting with PEEK ratchet connector for usage with M5150-67024 on Agilent 9500 ICP-MS, 750 mm length	M5150-67025
UniFit sample connector, for U-series MicroMist nebulizer, 0.5 mm id, 10/pk	G3266-80012
UniFit sample connector, for U-series MicroMist nebulizer, 0.25 mm id, 10/pk	G3266-80013
UniFit sample connector, for U-series MicroMist nebulizer, 0.75 mm id, 10/pk	G8010-80035
Nebulizer/Make-up/Dilution gas tubes for 9500 ICP-MS	M5150-67051
MicroFlow nebulizer, PFA, 200 µL/min, self-aspirating	G3139-65100
MicroFlow nebulizer, with I-AS probe, PFA, 200 µL/min, self aspirating	G3139-65102
MicroFlow nebulizer, with SPS 4 probe, PFA, 200 µL/min, self aspirating	G3139-68000
PFA-ST MicroFlow nebulizer, with demountable capillary, self aspirating. Order alternative capillary to vary flow rate from 20 to 400 µL/min available without or with autosampler probe (to suit I-AS or SPS 4/6 autosamplers).	G3139-68100
PFA-ST MicroFlow nebulizer capillary, 20 µL/min	G3139-68101
PFA-ST MicroFlow nebulizer capillary, 50 µL/min	G3139-68102
PFA-ST MicroFlow nebulizer capillary, 100 µL/min	G3139-68103
PFA-ST MicroFlow nebulizer capillary, 200 µL/min	G3139-68104
PFA-ST MicroFlow nebulizer capillary, 400 µL/min	G3139-68105
PFA-ST MicroFlow nebulizer capillary with I-AS probe, 20 µL/min	G3139-68106
PFA-ST MicroFlow nebulizer capillary with I-AS probe, 50 µL/min	G3139-68107
PFA-ST MicroFlow nebulizer capillary with I-AS probe, 100 µL/min	G3139-68108
PFA-ST MicroFlow nebulizer capillary with I-AS probe, 200 µL/min	G3139-68109
PFA-ST MicroFlow nebulizer capillary with I-AS probe, 400 µL/min	G3139-68110
PFA-ST MicroFlow nebulizer capillary with SPS 4 probe, 20 µL/min	G3139-68111
PFA-ST MicroFlow nebulizer capillary with SPS 4 probe, 50 µL/min	G3139-68112
PFA-ST MicroFlow nebulizer capillary with SPS 4 probe, 100 µL/min	G3139-68113
PFA-ST MicroFlow nebulizer capillary with SPS 4 probe, 200 µL/min	G3139-68114
PFA-ST MicroFlow nebulizer capillary with SPS 4 probe, 400 µL/min	G3139-68115

Spray Chambers and Supplies	Part Number
Spray Chamber, Quartz, for UHMI, Agilent 9500 ICP-MS (standard configuration)	M5150-67016
Dilution nozzle for UHMI	G8400-67026
Elbow gas connector for UHMI spray chamber dilution gas port	0100-2703
Connector Tube, Quartz, Straight, between spray chamber and torch	M5150-67014
Spray Chamber, Quartz, Agilent 9500 ICP-MS (Optional). Recommended if UHMI is not needed or for running organic solvents with option gas addition.	M5150-67017
Connector Tube, Quartz, with gas port for HMI aerosol dilution or Option Gas addition	M5150-67015
Straight gas connector for Connector Tube gas port	0100-2583
Endcap for Quartz Spray Chamber	G3280-60008
Clamp, torch ball joint connector, secures ball joint connector tube between torch and spray chamber, 1/pk	G8400-60327
Endcap retainer ring for spray chamber	G8400-40200
Drain tubing assembly for Quartz Spray Chamber	G3280-60555

Pump and Connecting Tubing	Part Number
Easy-Fit Peristaltic pump tubing, PVC, white/white, standard for sample uptake, recommended for aqueous, acid or alkaline matrix, 1 pack of 12 pcs	5005-0020
Easy-Fit Peristaltic pump tubing, PVC, blue/orange, standard for Internal Standard uptake, recommended for aqueous, acid or alkaline matrix, 1 pack of 12 pcs	5005-0021
Easy-Fit Peristaltic pump tubing, long-life, standard for spray chamber drain, 1 pack of 12 pcs	5005-0022
Easy-Fit Peristaltic pump tubing, PVC, black/black, optional high matrix sample and ISTD uptake, recommended for aqueous, acid or alkaline matrix, 1 pack of 12 pcs	5005-0023
Easy-Fit Peristaltic pump tubing, SolvaFlex, black/black, optional for high acid or organic solvent high matrix sample and ISTD uptake, 1 pack of 12 pcs	5005-0026
Easy-Fit Peristaltic pump tubing, SolvaFlex, blue/orange, optional for organic solvent or high acid matrix internal standard uptake, 1 pack of 12 pcs	5005-0027
Easy-Fit Peristaltic pump tubing, SolvaFlex, white/white, optional for organic solvent or high acid sample uptake, 1 pack of 12 pcs	5005-0025
Sample tubing, PFA, for sample introduction, 0.5 mm id, 1.6 mm od, 5 m	G1820-65105
Sample tubing, PFA, for internal standard introduction, 0.3 mm id, 1.6 mm od, 3 m	G1820-65478
PTFE nut for 1.6 mm od tubing, 10/pk	5064-8023
Front and back ferrule for 1.6 mm od tubing, 10/pk	5064-8024
Drain tubing, for use between peri-pump tubing and drain tank, 4.0 mm id, 2.5 m	G1820-65515
Tubing, Tygon, for drain to drain tank, 3 m	G3280-60077
Online internal standard addition kit for 9500 ICP-MS. Includes Grounding manifold with 2 inlets and connection for sample/ISTD lines	M5150-67105

Vacuum and Cooling System Supplies	Part Number
AVF Platinum oil for Agilent MS40+ foreline pump, 1 L	X3760-64004
Oil mist filter element for Agilent MS40+ (change at least annually)	G1960-80039
Cool Clear, coolant fluid, 2 Gal/pk	5799-0037

AVS MS Tubing Kits (for 9500 ICP-MS Only)	Part Number
Preconfigured kit for discrete sampling with AVS MS for 9500 ICP-MS. Includes a 1.5 mL sample loop (5005-0425).	M5171-67001
Capillary for Tune or ISTD uptake with AVS MS	M5171-67002
Tune/ISTD Uptake Connection Kit, AVS MS	M5171-67003
Carrier Uptake Connection Kit, AVS MS	M5171-67004
Carrier Grounding Connection Kit, AVS MS	M5171-67005
Nebulizer Connection Kit, AVS MS (uses UniFit type connector)	M5171-67006
AVS MS Autosampler Probe Connection Kit	M5171-67007
AVS MS Fast Pump Inlet Connection Kit	M5171-67008
AVS MS Fast Pump Outlet Connection Kit	M5171-67009

AVS MS Supplies	Part Number
Spare Valve, 7 port, for 9500 ICP-MS AVS MS	M5171-67011
AVS MS Valve Rotor Seal replacement kit for 9500 ICP-MS	M5171-67010
Stator and Rotor for AVS MS 7 port valve, 9500 ICP-MS	M5171-67012
MiraMist sample inlet connector for AVS	M5171-67013

AVS MS and ADS 2 Sample Loops	Part Number
Sample loop for AVS, 0.25 mL, 1.00 mm id, 1/pk	5005-0420
Sample loop for ADS/AVS, 0.50 mL, 1.00 mm id, 1/pk	5005-0421
Sample loop for ADS/AVS, 0.75 mL, 1.00 mm id, 1/pk	5005-0422
Sample loop for ADS/AVS, 1.00 mL, 1.00 mm id, 1/pk	5005-0423
Sample loop for ADS/AVS, 1.25 mL, 1.00 mm id, 1/pk	5005-0424
Sample loop for ADS/AVS, 1.50 mL, 1.00 mm id, 1/pk. Standard AVS MS sample loop.	5005-0425
Sample loop for ADS/AVS, 2.00 mL, 1.00 mm id, 1/pk	5005-0426
Sample loop for ADS/AVS, 2.50 mL, 1.00 mm id, 1/pk	5005-0427
Sample loop for ADS/AVS, 3.00 mL, 1.00 mm id, 1/pk	5005-0449
Sample loop for AVS, 3.50 mL, 1.00 mm id, 1/pk	5005-0450
Sample loop for AVS MS, 1.00 mL, 2.18 mm id, 1/pk	5005-0428
Sample loop for AVS MS, 1.25 mL, 2.18 mm id, 1/pk	5005-0429
Sample loop for AVS MS, 1.50 mL, 2.18 mm id, 1/pk	5005-0430
Sample loop for AVS MS, 2.00 mL, 2.18 mm id, 1/pk	5005-0431
Sample loop for AVS MS, 2.50 mL, 2.18 mm id, 1/pk	5005-0432
Sample loop for AVS MS, 3.00 mL, 2.18 mm id, 1/pk	5005-0433
Sample loop for AVS MS, 4.00 mL, 2.18 mm id, 1/pk	5005-0434
Sample tubing 1.0 mm id, 1/16" od, 5 m	G8411-67117
AVS MS CTFE nut for 1/16" od tubing, 10/pk	5043-0054
AVS MS Black flangeless fitting, 1/16" od, for 1.0 mm id sample loop, 5/pk	5023-1523
Sample tubing 2.18 mm id, 1/8" od, 5 m	G8411-67118
AVS MS CTFE nut for 1/8" od tubing, 10/pk	5043-0055
AVS MS Black flangeless fitting, 1/8" od, for 2.18 mm id sample loop, 5/pk	5023-1527

ADS 2 Autodilutor Supplies	Part Number
Syringe, 5mL, for ADS 2 carrier on ICP-OES and ICP-MS	5299-0037
Syringe, 10 mL, for ADS 2 diluent on ICP-MS, high purity quartz	5299-0036
Spare valve, 7 ports, 2 positions for AVS 7 or ADS 2 switching valves, 1/pk	G8495-60000
Stator and rotor for 7 ports valve for AVS 7 or ADS 2 switching valve, 1/pk	G8495-60003
Rotor seal for 6-7 port valve for AVS 6/7 or ADS 2 switching valve, 1/pk	G8494-60002
ADS 2, port plug kit for valve A	5005-0473
Loop Holder, AVS/ADS 2, 1/pk	5005-0452
Tubing Clips, for ADS 2, 6/pk	5005-0453
ADS 2 tubing kit, Valve C to AVS Pump, 1/pk	5005-0107
ADS 2 tubing kit, Valve C to AVS Valve, 1/pk	5005-0182
ADS 2 tubing kit, Valve B set-up, 4/pk	5005-0102
ADS 2 tubing kit, Valve A to Valve C, 1/pk	5005-0103
ADS 2 tubing kit, Carrier/Diluent, 2/pk	5005-0105
ADS 2 tubing kit, Valve A to AVS Valve, 1/pk	G8457-68004

Carrier, Diluent, and Waste Solutions Management	Part Number
Diluent/Carrier bottle kit for ADS 2 and Autosampler, 6L HDPE with GL45 Stay Safe cap and venting valve. Suitable for all applications using an Autosampler or Switching Valve.	5005-0435
Diluent bottle kit for ADS 2 on ICP-MS, 2L PFA with GL45 Stay Safe cap and venting valve. Suitable for all applications using an Autosampler or switching valve.	5005-0436
Waste container kit, 10L HDPE with S60 Stay Safe cap and charcoal filter. Suitable for any atomic spectroscopy instrument for waste solution management.	5005-0437
InfinityLab Stay Safe cap, GL45, 3 ports	5043-1219
S60 Stay Safe cap for spectroscopy including fittings	5043-1326
Acid vapor charcoal filter with 6-months time strip for S60 Stay Safe cap	5043-1193
One-way venting valve with PTFE filter and 6-months time strip for GL45 Stay Safe cap	5043-1190

Operating Supplies Kits	Part Number
Organic Sample Introduction System for 9500 ICP-MS. Includes 1.5 mm id torch, quartz spray chamber, endcap, retaining ring, drain tubing assembly, connector tube with gas port, Black/Black SolvaFlex peri-pump tubing (12/pk), 3 x Yellow/Blue Fluran peri-pump tubing (1/pk), Tubing for drainage (4.0 mm id, 2.5 m).	M5150-67106
Standard Quartz Sample Introduction System for 9500 ICP-MS (for aqueous samples). Includes 2.5 mm id torch, UHMI quartz spray chamber, endcap, retaining ring, drain tubing assembly, straight quartz connector tube.	M5150-67107
Basic supplies kit for 9500 ICP-MS. Includes peri-pump tubing for sample, white/white, 12/pk; for ISTD, blue/orange, 12/pk; for drain, yellow/blue, 12/pk; PFA sample tubing, 0.5 mm id, 1.6 mm od, 5 m; PFA sample tubing, 0.3 mm id, 1.6 mm od, 3 m; PTFE nut for 1.6 mm od tubing, 10/pk; front and back ferrule for 1.6 mm od tubing, 10/pk; graphite gasket for 9500 ICP-MS sampler cone, 3/pk; 2.5 mm id quartz torch; rough pump oil, AVF Platinum, 1 L.	M5150-67108
Configurable supplies kit for 9500 ICP-MS. Includes a comprehensive list of supplies with mandatory options to select: 1- Sampler Cone; 2- Skimmer Cone; 3- Quartz Spray Chamber configuration; 4- Quartz Connector Tube configuration. For extra options and more information see Agilent publication: Spare Parts and Operating Supplies Kits for Agilent ICP-MS (p/n 5991-5006EN).	M5145A

SPS 4/6 Autosampler Consumables	Part Number
Racks, Tubes, and Storage Bottles	
Standards rack, 5-position, (standard configuration). Provides 5 locations for 250 mL 60 mm od narrow neck sample containers.	G8410-40001
Standards rack, 34-position (optional configuration). Provides 12 locations for 50 mL 30 mm od tubes and 22 locations for 18 mm od tubes. (Only supported with MassHunter software version 4.5 or greater).	G8410-40000
Sample rack for 13 mm od tubes, 90 positions, 1/pk	6610026400
Sample rack for 16 mm od tubes, 60 positions, 1/pk	6610025400
Sample rack for 20 mm od tubes, 40 positions, 1/pk	6610025500
Sample rack for 25 mm od tubes, 24 positions, 1/pk	6610026500
Sample rack for 30 mm od tubes (50 mL centrifuge tubes), 21 positions, 1/pk	6610026600
Rack overlay for 60-position sample rack (2 per rack recommended), for Agilent SPS 4/6 autosampler. Allows 13 mm and some 16 mm od tubes to be used in 60-position rack.	6610026100
Rack overlay for 40-position sample rack (2 per rack recommended), for Agilent SPS 4/6 autosampler. Allows 18 mm od tubes to be used in 40-position rack.	6610026000
Bottle 250 mL for 5-position standards rack, 1/pk	5043-0064
Centrifuge tubes, polypropylene, 5 to 50 mL at 2.5 mL intervals, 500/pk, screw cap, 29 mm od	190065200
Polypropylene tubes, 17 mm od, 16.5 mL, 1,000/pk	00000001600L
PFA Lab Bottle, 31.5 mm od, 20 mL with 33 mm closure	G8410-68011
PFA Lab Bottle, 38 mm od, 50 mL with 33 mm closure	G8410-68012
PFA Lab Bottle, 57.5 mm od, 100 mL with GL45 closure	G8410-68013
PFA Lab Bottle, 72 mm od, 200 mL with GL45 closure	G8410-68014
PFA Lab Bottle, 87 mm od, 500 mL with GL45 closure	G8410-68015
SPS sample rack 21 bottles × 31.5 mm od	G8410-68001
SPS sample rack 8 bottles × 38 mm od	G8410-68002
SPS sample rack 4 bottles × 57.5 mm od	G8410-68003
SPS sample rack 2 bottles × 72 mm od	G8410-68004
SPS sample rack 2 bottles × 87 mm od	G8410-68005
Pump Tubing for the Rinse Pump	
Peristaltic pump tubing, PVC SolvaFlex, 3-bridged, gray/gray tabs, for Agilent SPS 4/6 when pumping acidic or aqueous rinses, 12/pk	3710049000
Peristaltic pump tubing, Santoprene, 3-bridged, gray/gray tabs, for ketone-based solvents on AA, MP-AES, and ICP-OES instruments, 6/pk	3710057500
Peristaltic pump tubing, FPM, 3-bridged, black/black tabs with flared ends, for use with Agilent SPS 4/6 autosampler when pumping aromatic organic samples, 12/pk	3710069100

Spare Probes	
SPS 4 probe, 0.25 mm id, red marker, inert, FEP sleeved, 1/pk	G8410-80100
SPS 4 probe, 0.5 mm id, yellow marker, inert, FEP sleeved, 1/pk (standard probe for ICP-MS)	G8410-80101
SPS 4 probe, 0.8 mm id, green marker, inert, FEP sleeved, 1/pk	G8410-80102
SPS 4 sample probe with red marker, 1/pk. Required when using the Agilent SPS 4 autosampler with microtiter tray for sampling from a 96 position well plate. Requires, but does not include, the well plate tray adaptor kit or any compatible well plates.	G8410-80105
SPS 4 probe 1.00 mm id, blue marker, inert, FEP sleeved, 1/pk (recommended for AVS/ADS 2)	G8410-80103
SPS 6 probe 0.25 mm id, red marker, inert, FEP sleeved, 1/pk	G8009-64800
SPS 6 probe 0.5 mm id, yellow marker, inert, FEP sleeved, 1/pk (standard probe for ICP-MS)	G8009-64801
SPS 6 probe 0.8 mm id, green marker, inert, FEP sleeved, 1/pk	G8009-64802
SPS 6 probe 1.00 mm id, blue marker, inert, FEP sleeved, 1/pk (recommended for AVS/ADS 2)	G8009-64803
SPS 6 sample probe with red marker, 1/pk. Required when using the Agilent SPS 6 autosampler with microtiter tray for sampling from a 96 position well plate. Requires, but does not include, the well plate tray adaptor kit or any compatible well plates.	G8009-64805

SPS 4/6 Autosampler Options	Part Number
High uptake volume rinse kit for Agilent SPS 4/6 autosampler to connect to ISIS 3. Allows rinse uptake rate to be increased to 25 mL/min.	G8410-68260
SPS 4/6 rinse tube kit, for low volatile organic solvents in rinse station. Includes 1 m capillary, 3-bridged gray/gray PVC Solvaflex (4 pump tubes), barb connector, 2.5–1.5 mm, solvent-resistant tubing for inlet + outlet.	G8410-68261
Pump drain kit for Agilent SPS 4/6 autosampler. Allows the liquid overflow to be pumped from the rinse station. Includes wide-bore tubing, purple/white tabs (4 pump tubes), barb connectors, 2.5 m PVC Solvaflex to connect to a remote pump (not supplied).	G8410-68262
Well plate startup kit for Agilent SPS 4/6 autosampler. Allows the SPS 4/6 to be adapted to sample from 96-well plates. Includes well plates (10/pk), well plate adapter, and sample probe. Only supported with the ICP-MS MassHunter software.	G8410-68270

ICP-MS Stock Tuning Solutions	Part Number
Li, Y, Ce, Tl, and Co; 100 mL, 10 mg/L; matrix 2% HNO ₃	5188-6564
ICP-MS stock tuning solution: Li, Mg, Y, Ce, Tl, Co; 100 mL, 10 mg/L; matrix 2% HNO ₃	5190-0465
Tuning solution: Li, Y, Ce, Tl, and Co; 2 × 500 mL, 10 µg/L; matrix 2% HNO ₃	5184-3566
PA tuning solution kit: 2 × 100 mL bottles	
Tuning 1:	
20 ppm each of Zn, Be, Cd, As;	
10 ppm each of Ni, Pb, Mg; 5 ppm each of Tl, Na Al, U, Cu, Th, Ba, Co, Sr, V, Cr, Mn, 6Li, Sc, In, Lu, Bi;	
2.5 ppm each of Y, Yb; matrix 2 to 5% HNO ₃	5188-6524
Tuning 2:	
10 ppm each of Mo, Sb, Sn, Ge, Ru, Pd; 5 ppm of Ti, Ir; matrix 10% HCl/1% HNO ₃ /trace HF	

10 µg/mL Single Element ICP-MS Standards Supplied in 100 mL bottles	Part Number
Aluminum (Al) in 2% HNO ₃	5190-8561
Antimony (Sb) in 1% HNO ₃ , trace tartaric acid	5190-8562
Arsenic (As) in 2% HNO ₃	5190-8563
Barium (Ba) in 2% HNO ₃	5190-8564
Beryllium (Be) in 2% HNO ₃	5190-8565
Boron (B) in H ₂ O	5190-8566
Cadmium (Cd) in 2% HNO ₃	5190-8567
Chromium (Cr) in 2% HNO ₃	5190-8568
Copper (Cu) in 2% HNO ₃	5190-8569
Iridium (Ir) in 2% HCl	5190-8570
Lead (Pb) in 2% HNO ₃	5190-8571
Lithium (Li) in 2% HNO ₃	5190-8572
Lutetium (Lu) in 2% HNO ₃	5190-8573
Manganese (Mn) in 2% HNO ₃	5190-8574
Mercury (Hg) in 5% HNO ₃	5190-8575
Nickel (Ni) in 2% HNO ₃	5190-8576
Platinum (Pt) in 5% HCl	5190-8577
Scandium (Sc) in 2% HNO ₃	5190-8578
Selenium (Se) in 2% HNO ₃	5190-8579
Silver (Ag) in 2% HNO ₃	5190-8580
Strontium (Sr) in 2% HNO ₃	5190-8581
Terbium (Tb) in 2% HNO ₃	5190-8582
Tin (Sn) in 5% HCl	5190-8583
Uranium (U) in 2% HNO ₃	5190-8584
Vanadium (V) in 5% HNO ₃	5190-8585
Yttrium (Y) in 2% HNO ₃	5190-8586
Zinc (Zn) in 2% HNO ₃	5190-8587

Multi-Element Calibration Standards	Part Number
Multi-element calibration standard 1, 100 mL: 10 mg/L of Ce, Dy, Er, Eu, Gd, Ho, La, Lu, Nd, Pr, Sc, Sm, Tb, Th, Tm, Y, Yb; matrix 5% HNO ₃	8500-6944
Multi-element calibration standard 2A, 100 mL: 10 mg/L of Ag, Al, As, Ba, Be, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Ti, U, V, Zn; matrix 5% HNO ₃	8500-6940
Multi-element calibration standard 3, 100 mL: 10 mg/L of Sb, Au, Hf, Ir, Pd, Pt, Rh, Ru, Te, Sn; matrix 10% HCl/1% HNO ₃	8500-6948
Multi-element calibration standard 4B, 100 mL: 10 mg/L of B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr; matrix H ₂ O/trace HF	8500-6942

Semiquantitative Standards Supplied in 100 mL bottles	Part Number
ICP-MS semiquantitative calibration standard 1. Contains: 10 µg/mL Ag, Al, As, Ba, Bi, Ca, Cd, Ce, Dy, Er, Eu, Ga, Gd, Ho, La, Lu, Mg, Na, Nd, P, Pb, Pr, Rb, Sc, Se, Sm, Sr, Tb, Th, Tl, Tm, U, Y, Yb; matrix 40% aqua regia	5190-8594
ICP-MS semiquantitative calibration standard 2. Contains: 10 µg/mL Au, B, Be, Co, Cr, Cu, Fe, Ge, Hf, Ir, K, Li, Mn, Mo, Nb, Ni, Os, Pd, Pt, Re, Rh, Ru, Sb, Si, Sn, Ta, Te, Ti, V, W, Zn, Zr; matrix 40% aqua regia, trace HF	5190-8595

Environmental Standards	Part Number
Environmental calibration standard, 100 mL: 1,000 mg/L of Fe, K, Ca, Na, Mg; 10 mg/L of Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Ti, V, Zn, Th, U; matrix 10% HNO ₃	5183-4688
Initial calibration verification standard, 100 mL: 1,000 mg/L of Fe, K, Ca, Na, Mg, Sr; 10 mg/L of Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Ti, V, Zn, Th, U; matrix 5% HNO ₃	5183-4682
Internal standard mix, 100 mL: 100 ppm of Be, Sc, Ge, Rh, In, Tb, Lu, Bi; matrix 10% HNO ₃	5188-6525
Internal standard mix, 100 mL: 10 mg/L of 6Li, Sc, Ge, Y, In, Tb, Bi; matrix 5 to 10% HNO ₃	5183-4681
Environmental spike mix, 100 mL: 1,000 mg/L Fe, K, Ca, Na, Mg, 100 mg/L Ag, Al, As, Ba, Be, Cd, Co, Cr, Cu, Mn, Mo, Ni, Pb, Sb, Se, Ti, V, Zn, U; matrix 5% HNO ₃	5183-4687
6020 interference check solution A, 100 mL: 20,000 ppm of Cl; 3,000 ppm of Ca; 2,500 ppm each of Fe, Na; 2,000 ppm of C; 1,000 ppm each of Al, Mg, P, K, S; 20 ppm each of Ti, Mo; matrix 5% HNO ₃ /trace HF	5188-6526
6020 interference check solution B, 100 mL: 20 ppm each of Cr, Co, Cu, Mn, Ni, V; 10 ppm each of As, Cd, Se, Zn; 5 ppm of Ag; matrix 5% HNO ₃	5188-6527

Agilent supplies to help you get the results you need

Agilent supplies are designed, manufactured, and tested to rigorous specifications under a quality system registered to ISO 9001. Why risk compromising your analytical results with anything less than genuine Agilent supplies?

For a complete list of Agilent atomic spectroscopy supplies, including AA lamps and supplies for AA, ICP-OES, ICP-MS and MP-AES products, visit: www.agilent.com/chem/atomic

DE-014150

This information is subject to change without notice.

© Agilent Technologies, Inc. 2026
Printed in the USA, April 22, 2026
5994-9096EN