

Optima™ 7300 V ICP-OES



Demountable Torches

Demountable design using one-piece quartz tubing for plasma and auxiliary gas flow.



Demountable Quartz Torches

# of Slots	Part No.
3	N0771500
3 ABC Fully Demountable Torch Kit	N0774075



Ceramic D-Torch

Description	Part No.
D-Torch with Quartz Outer Tube	N0777051
Quartz Outer Tube	N0777053
Base and Inner Tube	N0777054
Retaining Ring	N0777055



Torch Bonnet

Description	Part No.
Torch Bonnet	N0775289

Quick-Change Torch Modules

Description	Part No.
Quick-Change Torch Module for HF-Resistant Analysis	N0770911
Quick-Change Torch Module for Used Oil Analysis	N0770910



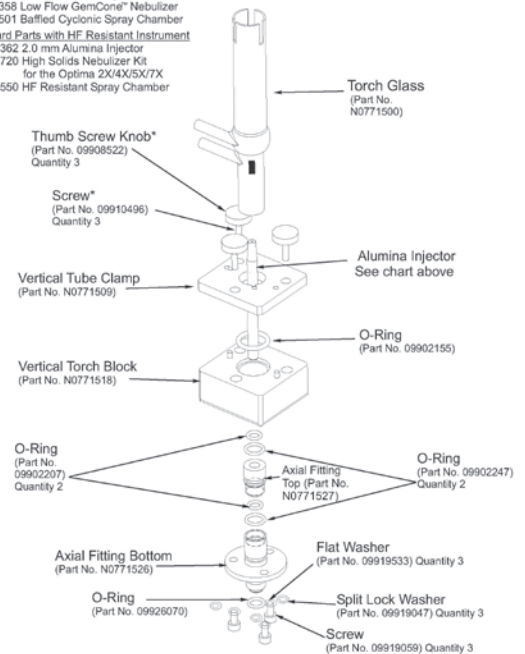
Ignitor Tape

Description	Part No.
Copper Ignitor Tape	N0775297

Torch O-Rings

Location	Dimensions	Part No.
Ball Joint	0.301 i.d., 0.070 w.d.	09926070
Complete O-Ring Kit		N0770916
Includes:	Qty	Part No.
	1	09902155
	2	09902207
	2	09902247
	1	09926070

Standard Parts with Oils Instrument
 N0771531 1.2 mm Alumina Injector
 N0770358 Low Flow GemCone® Nebulizer
 N0771501 Baffled Cyclonic Spray Chamber
 Standard Parts with HF Resistant Instrument
 N0695362 2.0 mm Alumina Injector
 N0770720 High Solids Nebulizer Kit
 for the Optima 2X/4X/5X/7X
 N0771550 HF Resistant Spray Chamber



*Thumb screw knob (Part No. 09908522) must be pressed onto screw (Part No. 09910496.)

Nebulizers

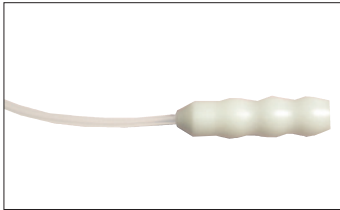
Offer high-quality sample introduction for a wide variety of applications.



MEINHARD® Nebulizers

Provide excellent sensitivity and precision for aqueous solutions and samples with few dissolved solids (less than 1%). MEINHARD® Nebulizers are self-aspirating and must not be used with solutions containing hydrofluoric acid. All require an end cap for use with Ryton® Scott spray chambers.

Type	Application	Part No.
Type A	General purpose	00472020
Type C	High dissolved solids	00472022
Type K	Reduced flow, and organics	N0681574



MEINHARD® Nebulizer Parts

Description	Part No.
Liquid-End Connector Kit, 4 mm	N0777414
Nebulizer Clip for MEINHARD® Nebulizers	N0777460

End Cap for Concentric Glass Nebulizers

Description	Part No.
End Cap	N0680504



MicroMist Nebulizers

Air Flow Rate	Uptake Rate	Part No.
0.7 L/min	0.05 mL/min	N0775349
0.7 L/min	0.1 mL/min	N0775348
0.7 L/min	0.2 mL/min	N0775347
0.7 L/min	0.6 mL/min	N0775346

Requires Quick Disconnect Part No. N0770336.



SeaSpray Nebulizers

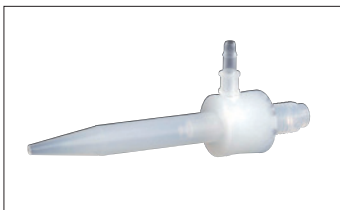
Air Flow Rate	Uptake Rate	Part No.
0.7 L/min	2 mL/min	N0775345



PEEK MiraMist® Nebulizer

The MiraMist® fits regular spray chambers as a direct replacement for glass concentric nebulizers. Its features include standard pressure (45 psi), low sample flow and enhanced parallel path design.

Description	Uptake Rate	Part No.
PEEK MiraMist®	0.2 to 2.5 mL/min	N0775330



MicroFlow Nebulizers

MicroFlow nebulizers are constructed entirely from chemically resistant fluoropolymers, ideal for strong acids, alkalis, organics, high salt solutions and high solid concentrations. They are resistant to clogging, are reliably self-aspirated or pumped and produce a fine aerosol for high transport efficiency and high sensitivity.

Description	Part No.
PolyPro ST Nebulizer	N0777219
PFA-ST Nebulizer	N0777049
PFA-ST3 Microflow Nebulizer	N0777676

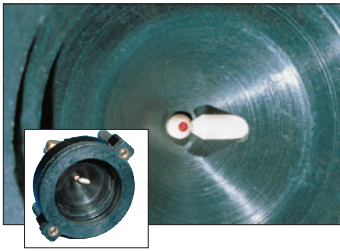


GemCone™ Nebulizers

Low-flow GemCones™ permit lower nebulizer gas flow rates, useful for spectral lines with high excitation energies and for providing a more robust plasma. GemCone™ Nebulizers require an end cap for use with Ryton® Scott Spray Chambers.

Description	Part No.
GemCone™ Low-Flow Nebulizer	N0770358
GemCone™ High Dissolved Solids	N0690670
GemCone™ Sample Capillary Adapter	N0371505
End Cap for GemCone™ Nebulizers	N0680504
Quick Disconnect Tubing	N0770336

Nebulizers, continued



GemTip Cross-Flow II Nebulizers

Excellent general purpose nebulizer for the analysis of strong mineral acids (including HF) and samples with less than 5% dissolved solids.

Description	Part No.
GemTip Cross-Flow II Nebulizer	N0770546
GemTip Cross-Flow II Tip Kits	N0690676
O-Ring Kit for GemTip Cross-Flow Nebulizers	N9300067

Spray Chambers

The spray chamber is an integral part of the sample introduction system responsible for filtering the sample mist to permit the appropriate droplet size distribution to reach the plasma.

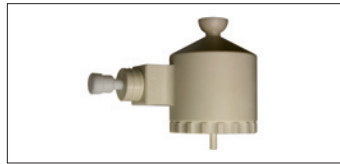


Baffled Cyclonic Spray Chambers

Type	Part No.
Standard	N0771501
O-Ring Free	N0775356
HF Resistant	N0771550

Unifit Drain Connectors with Tubing for Baffled Cyclonic Spray Chambers

Sample Tube Size	Qty	Part No.
1/16 in. o.d. x 0.75 mm i.d. x 700 mm	3	N0774080
1.3 mm o.d. x 0.75 mm i.d. x 700 mm	3	N0774077



Tracey TC44 Spray Chambers

Type	Part No.
TC-44 Cyclonic Spray Chamber	N0771550
Tracey TFE Spray Chamber with Helix	N0777497

Injectors and Injector Support Adapters



Alumina Injectors

Alumina injectors are corrosion-resistant to all mineral acids including HF and aqua regia. They are also used with less volatile organic solvents, e.g., xylene.

Dimensions	Part No.
2.0 mm i.d.	N0695362
1.2 mm i.d.	N0771531



Sapphire Injectors

Sapphire injectors are ideal for silicon applications.

Dimensions	Part No.
2.0 mm i.d.	N0695494



Quartz Injectors

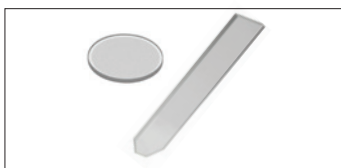
Quartz injectors are designed for organic solvents and non-HF matrices.

Dimensions	Part No.
0.8 mm i.d.	N0695461
1.2 mm i.d.	N0695460
1.6 mm i.d.	N0695449
2.0 mm i.d.	N0695442

Injector Support Adapters

Spray Chamber	Part No.
(bottom) Cyclonic	N0771526
(top) Cyclonic	N0771527

Purge Windows and Load Coil



Purge Extension Windows

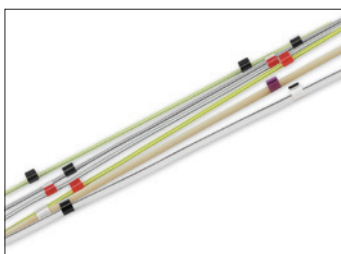
View	Part No.
Radial	N0581497



Load Coil

Description	Part No.
Load Coil	N0771536

Autosampler Pump Tubing



i.d. (mm)	Stop Colors	Standard PVC (Pkg 12)	Flared PVC (Pkg 12)	Solvent Flex (Pkg 12)	Viton® (Pkg 12)	Santoprene® (Pkg 12)	Silicone (Pkg 6)
0.19	orange red	N0695476	N0773111				
0.25	orange blue	N0773117	N0773112				
0.38	orange green	N0777110	N0777042				
0.44	green yellow		N0773113				
0.51	orange yellow		N0777476		N0777446		
0.76	black black	09908587	N0777043	00473550	N0773118		00473552
1.02	white white				N0777442		N0777441
1.14	red red	09908585		09923037	N0773115	N0773119	N0691595
1.30	gray gray				N0777445	N0777444	N0777443
1.42	yellow yellow	N0777521					
1.52	yellow blue	09923536					
1.65	blue blue				N0777447		
1.85	green green					N0773116	
2.79	purple white			02508080			09923448
3.18	orange white	N8122012					

Autosampler Tubes with Caps



Capacity	Size	Type	Qty	Part No.
15 mL	17 mm x 120 mm	Conical Bottom	500	B0193233
50 mL	28 mm x 115 mm	Conical Skirted Bottom	500	B0193234

Autosampler Tubes

Capacity	Size	Type	Qty	Part No.
8 mL	13 mm x 100 mm	Round Bottom	1,200	N0777156
8 mL	13 mm x 100 mm	Round Bottom	250	N0777159
15 mL	17 mm x 100 mm	Round Bottom	1,200	N0777167
15 mL	17 mm x 100 mm	Round Bottom	250	N0777168
16 mL	17 mm x 100 mm	Round Bottom	1,000 (packs of 125)	N9301205
50 mL	30 mm x 115 mm	Round Bottom	500 bulk	N0777158
50 mL	30 mm x 115 mm	Round Bottom	500 (50 tubes per bag)	N0777164
	Plug for 15 mL N0777167 Vials		1,200	N0777599



Single-Element Standards – 1,000 mg/L

Whatever your application, you can depend on PerkinElmer Pure Standards to yield reliable, accurate results. PerkinElmer offers a wide selection of atomic spectroscopy calibration standards. Each solution is supplied with a comprehensive Certificate of Analysis that documents the quality and reliability.

Element	Symbol	Matrix	Pure Grade 125 mL	Pure Grade 500 mL
			Part No.	Part No.
Aluminum	Al	2% HNO ₃	N9300184	N9300100
Cadmium	Cd	2% HNO ₃	N9300176	N9300107
Calcium	Ca	2% HNO ₃	N9303763	N9300108
Chromium	Cr	2% HNO ₃	N9300173	N9300112
Copper	Cu	2% HNO ₃	N9300183	N9300114
Iron	Fe	2% HNO ₃	N9303771	N9300126
Lead	Pb	2% HNO ₃	N9300175	N9300128
Mercury	Hg	10% HNO ₃	N9300174	N9300133
Sodium	Na	2% HNO ₃	N9303785	N9300152
Zinc	Zn	2% HNO ₃	N9300178	N9300168

Universal Data Acquisition Standards Kit

Matrix	Contents	Vol.	Part No.
Solution Kit (Includes items listed below – 1 bottle of each)			N9306225
5% HNO ₃	10 mg/L: Ce, Dy, Er, Eu, Gd, Ho, La, Lu, Nd, Pr, Sm, Sc, Tb, Th, Tm, Y, Yb	125 mL	N9300232
5% HNO ₃	10 mg/L: Al, Ag, As, Ba, Be, Bi, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, In, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Ti, V, U, Zn	125 mL	N9300233
10% HCl/ 1% HNO ₃	10 mg/L: Au, Hf, Ir, Pd, Pt, Rh Ru, Te, Sb, Sn	125 mL	N9300234
5% HNO ₃	10 mg/L: Hg	125 mL	N9300253
H ₂ O/tr HF	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235

Multi-Element Standards

Matrix	Contents	Vol.	Part No.
Multi-Element Solution 1			
2% HNO ₃	10 mg/L: Be, Bi, Ce, Co, In, Mg, Ni, Pb, U	125 mL	N9300231
Multi-Element Solution 2			
5% HNO ₃	10 mg/L: Ce, Dy, Er, Eu, Gd, Ho, La, Lu, Nd, Pr, Sc, Sm, Tb, Th, Tm, Y, Yb	125 mL	N9300232
Multi-Element Solution 3			
5% HNO ₃	10 mg/L: Ag, Al, As, Ba, Be, Bi, Ca, Cd, Co, Cr, Cs, Cu, Fe, Ga, Hg*, In, K, Li, Mg, Mn, Na, Ni, Pb, Rb, Se, Sr, Tl, U, V, Zn *Supplied in separate bottle.	125 mL	N9300233
Multi-Element Solution 3 without Mercury			N9301720
Multi-Element Solution 4			
10% HCl/1% HNO ₃	10 mg/L: Au, Hf, Ir, Pd, Pt, Rh, Ru, Sb, Sn, Te	125 mL	N9300234
Multi-Element Solution 5			
H ₂ O/tr HF	10 mg/L: B, Ge, Mo, Nb, P, Re, S, Si, Ta, Ti, W, Zr	125 mL	N9300235

Interference Check Standards

Matrix	Contents	Vol.	Part No.
Interferents A			
5% HNO ₃	5,000 mg/L: Al, Ca, Mg 2,000 mg/L: Fe	500 mL	N9300226
Alternate Interferents A			
5% HNO ₃	1,000 mg/L: Cr, Cu, Mn, Ni, Ti, V	500 mL	N9300228
Analytes B			
5% HNO ₃ /trTartaric Acid/tr HF	100 mg/L: Cd, Ni, Zn 60 mg/L: Sb 50 mg/L: Ba, Be, Co, Cr, Cu, Mn, V 20 mg/L: Ag 10 mg/L: As, Tl	125 mL	N9300227
Alternate Analytes B			
5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: Al, As, B, Mo, Na, Sb, Se, Tl 10 mg/L: Ca, Fe, Mg, Si	125 mL	N9300229

Instrument Setup Solutions

Matrix	Contents	Vol.	Part No.
Vis Wavecal Solution			
2% HNO ₃	50 mg/L: K 10 mg/L: La, Li, Mn, Na, Sr 1 mg/L: Ba, Ca	250 mL	N9302946
UV Wavecal Solution			
5% HCl	100 mg/L: K, P, S 20 mg/L: As, La, Li, Mn, Mo, Na, Ni, Sc	125 mL	N0681470
5% HCl	100 mg/L: K, P, S 20 mg/L: As, La, Li, Mn, Mo, Na, Ni, Sc	500 mL	N0582152
Low UV Standard			
2% HNO ₃	10 mg/L: Al, P, S	250 mL	N0691580
Calcium Stray Light Standard			
H ₂ O	10,000 mg/L: Ca	125 mL	N0691581

Every day you count on PerkinElmer to provide you with solutions that deliver reliable performance, control operating costs and maximize operational time. Our complete portfolio of consumables, parts, supplies, training and service helps you meet both routine and demanding measurement challenges. We invest heavily in testing and validating our products to ensure you receive guaranteed compatibility and performance – on-time, for every instrument in your laboratory.

For a complete listing of consumables and supplies, please visit www.perkinelmer.com/supplies