

PinAAcle™ 500 Flame AA



Lumina Hollow Cathode and Electrodeless Discharge Lamps

Every genuine PerkinElmer Lumina™ Lamp is designed for use with and tested on our spectrometers to assure compatibility and the highest performance.



Coded Lumina Single-Element HCLs and EDLs

Element Name	Symbol	Lumina Part No.	EDL Part No.
Aluminum	Al	N3050103	
Antimony	Sb	N3050170	N3050670
Arsenic	As	N3050105	N3050605
Cadmium	Cd	N3050115	N3050615
Calcium	Ca	N3050114	
Chromium	Cr	N3050119	
Copper	Cu	N3050121	
Gold	Au	N3050107	
Iron	Fe	N3050126	
Lead	Pb	N3050157	N3050657
Magnesium	Mg	N3050144	
Manganese	Mn	N3050145	
Mercury	Hg	N3050134	N3050634
Nickel	Ni	N3050152	
Potassium	K	N3050139	
Selenium	Se	N3050172	N3050672
Sodium	Na	N3050148	
Zinc	Zn	N3050191	N3050691

Coded Lumina Multi-Element Hollow Cathode Lamps

Element Name	Symbol	Part No.
Calcium, Magnesium	Ca, Mg	N3050202
Potassium, Sodium	K, Na	N3050204
Copper, Iron, Manganese, Zinc	Cu, Fe, Mn, Zn	N3050212
Cobalt, Chromium, Copper, Iron, Manganese, Nickel	Co, Cr, Cu, Fe, Mn, Ni	N3050217
Aluminum, Calcium, Copper, Iron, Magnesium, Silicon, Zinc	Al, Ca, Cu, Fe, Mg, Si, Zn	N3050218



Burner Heads

Description	Length	Flame Type	Part No.
Single-slot	5 cm	Nitrous Oxide-Acetylene	N0400100
Single-slot*	10 cm	Air-Acetylene	N0400102*
Three-slot	10 cm	Air-Acetylene	N0400103
Single-slot	5 cm	Air-Acetylene	N0400101
Burner Head Handle			N3161040

**Standard burner head shipped with the PinAAcle 500 instrument*



Nebulizer Assemblies and Spares

Description	Type	Needle	O-ring Type	Complete Assembly Part No.	End Cap Assembly Part No.	Replacement Needle Assembly Part No.
Stainless Steel (standard)	Aqueous	Stainless	Viton	N3160143	N3160505	03030354
Plastic High Sensitivity (standard)	Aqueous	Pt/Ir	Viton	N3160144	N3160511	B3150525
Plastic High Sensitivity	Aqueous or Organics	Pt/Ir	Kalrez	N3160159	N3160189	N3160154
Plastic High Sensitivity*	Aqueous or Organics	Tantalum	Kalrez	N3160152	N3160189	
Stainless Steel	Aqueous or Organics	Stainless	Kalrez	N3160158	N3160188	
Stainless Steel	Aqueous	Plastic	Viton	N/A	N3160505	03030402
Stainless Steel	Aqueous	Platinum Alloy	Viton	N/A	N3160505	03030433

**Suitable for aqua regia in addition to organics*

Kalrez O-Ring Kits

Description No.	Type	For Stainless Nebulizers Part No.	For Plastic Nebulizers Part No.
Nebulizer O-Ring Kit	Organics	N3160192	N3160193
End Cap O-Ring Kit	Organics	N3160194	N3160194
Nebulizer, End Cap, and Burner Head O-Ring Kit	Organics	N3160156	N3160157

Autosampler Supplies

We offer a wide variety of high quality autosampler tubes and supplies. Autosampler tubes range in size from 8 mL to 50 mL and are available as conical or round bottom capped and uncapped.



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 Bottom/Freestanding	500	B0193234

Single-Element Standards

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



Single-Element Standards 1,000 mg/L in aqueous solution

Element Name	Symbol	125 mL Part No.	500 mL Part No.
Aluminum	Al	N9300184	N9300100
Arsenic	As	N9300180	N9300102
Cadmium	Cd	N9300176	N9300107
Calcium Ca		N9303763	N9300108
Chromium	Cr	N9300173	N9300112
Copper	Cu	N9300183	N9300114
Gold	Au	N9303759	N9300121
Iron	Fe	N9303771	N9300126
Lead	Pb	N9300175	N9300128
Magnesium	Mg	N9300179	N9300131
Manganese	Mn	N9303783	N9300132
Mercury	Hg	N9300174	N9300133
Nickel	Ni	N9300177	N9300136
Phosphorus	P	N9303788	N9300139
Potassium	K	N9303779	N9300141
Silver	Ag	N9300171	N9300151
Sodium	Na	N9303785	N9300152
Zinc	Zn	N9300178	N9300168

Multi-Element Standards

Quality Control Standards

Matrix	Contents	Volume	Part No.
Quality Control Standard, 21 Elements 5% HNO ₃ /tr Tartaric Acid/tr HF	100 mg/L: As, Be, Ca, Cd, Co, Cr, Cu, Fe, Li, Mg, Mn, Mo, Ni, Pb, Sb, Se, Sr, Ti, Tl, V, Zn	125 mL	N9300281
Quality Control Standard, 7A Elements 5% HNO ₃ /tr HF	1,000 mg/L: K 500 mg/L: Si 100 mg/L: Al, B, Ba, Na 50 mg/L: Ag	125 mL	N9300280

Alternate Metals I

Matrix	Concentration	Volume	Part No.
2% HNO ₃	5 mg/L: Be, Sb, Tl 10 mg/L: Co, Cu, Mn, Ni, Zn 20 mg/L: Al, Fe, V	125 mL	N9300214

Alternate Metals II

Matrix	Concentration	Volume	Part No.
2% HNO ₃	100 mg/L: K, Mg 500 mg/L: Ca, Na	125 mL	N9300215

GFAAS Mixed Standard

Matrix	Concentration	Volume	Part No.
5% HNO ₃ w/trace of HF	5 mg/L: Be, Cd 10 mg/L: Ag 20 mg/L: Cr, Fe, Mg 50 mg/L: Ba, Co, Cu, Ni 100 mg/L: Al, As, Pb, Sb, Se, Tl	125 mL	N9300244

AA Test Mix

Matrix	Concentration	Volume	Part No.
2% HCl	10 mg/L: Na, Zn 20 mg/L: K 50 mg/L: Ca, Cr, Cu, Fe, Ni	125 mL	02900540

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 AA Consumables, please visit www.perkinelmer.com/supplies

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.