



## AAAnalyst™ 200/400 and HGA 900

### 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

## Graphite Components

The production of high-quality graphite components requires stringent quality control. To ensure high quality and consistency, a specific grade of high-density base graphite material has been developed for PerkinElmer's exclusive use.



### HGA Graphite Tubes

Description	Part No.
-------------	----------

#### Pyrocoated Graphite Tubes with Integrated Platforms

5-pack	B3001262
20-pack	B3001264

#### Pyrocoated Graphite Tubes

5-pack	B0105197
10-pack	B0135653
20-pack	B3000342

#### Graphite Tubes grooved for use with L'vov Platforms

5-pack	B0137111
20-pack	B3001254
50-pack	B0109322
100-pack	N3110146

#### Platforms

5-pack	B0137112
20-pack	B3001256
50-pack	B0109324
100-pack	N3110145



### HGA Contact Cylinders

Description	Part No.
-------------	----------

#### For HGA-900/850/800/700/600/300 Furnaces

1 set	B0128495
5 sets	B3130086

#### For Zeeman Furnaces

1 set	B0116823
5 sets	B0180361



### 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

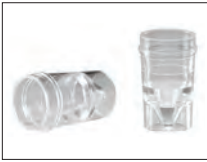


### Nebulizers

Description	Assembly Part No.	End Cap Part No.
Stainless Steel Nebulizer	03030352	00570984
High Sensitivity Nebulizer with Pt/Ir Capillary	N3150188	N0370392

## Autosampler Supplies

We offer a wide variety of high quality graphite furnace cups and autosampler tubes and supplies. Sample cups are of heavy-duty construction to prevent cracking or tipping. Their conical design allows for use of nearly the entire sample. Autosampler tubes range in size from 8 mL to 50 mL and are available as conical or round bottom capped and uncapped.



### Graphite Furnace Autosampler Cups

Volume	Composition	Quantity	Part No.
1.2 mL	Polypropylene	2,000	B0510397
2.5 mL	Polypropylene	1,000	B3001566
7.0 mL	Polypropylene	100	B3001567
2.0 mL	Polyethylene	1,000	B0087056
2.0 mL	Polystyrene	1,000	B0119079
3.5 mL	Polystyrene	100	B0129303
2.0 mL	Teflon®	30	B0087600



### Autosampler Tubes with Caps

Capacity	Size	Type	Qty	Part No.
15 mL	13 mm x 100 mm	Conical Bottom	500	B0193233
50 mL	30 mm x 115 mm	Conical Bottom	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.
<b>Quality Control Standard, 21 Elements</b> 5% HNO <sub>3</sub> /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
<b>Quality Control Standard, 7A Elements</b> 5% HNO <sub>3</sub> /tr HF	1,000 mg/L: K 1,500 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 <sub>3</sub>	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 <sub>3</sub>	100 mg/L: K, Mg 500 mg/L: Ca, Na	125 mL	N9300215

### GFAAS Mixed Standard

Matrix	Concentration	Volume	Part No.
5% HNO <sub>3</sub> 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

### Matrix Modifiers

Modifiers	Concentration	Volume	Part No.
Mg(NO <sub>3</sub> ) <sub>2</sub>	1% Mg (as nitrate)	100 mL	B0190634
Pd	1% Pd (as nitrate)	50 mL	B0109635
NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>	10% NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>	100 mL	N9303445



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](http://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](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

009547A\_01