



AAAnalyst™ 200/400 and HGA 900

Lumina Hollow Cathode and EDL Lamps

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



Lumina Coded Single-Element Lamps/EDL Lamps

| 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 |

Multi-Element Coded 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 µg/mL 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 µg/mL: 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 µg/mL: K 1,500 µg/mL: Si 100 µg/mL: Al, B, Ba, Na 50 µg/mL: Ag | 125 mL | N9300280 |

Alternate Metals I

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

Alternate Metals II

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

GFAAS Mixed Standard

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

AA Test Mix

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

Matrix Modifiers

| Modifiers | Concentration | Volume | Part No. |
|--|--|--------|----------|
| Mg(NO ₃) ₂ | 1% Mg (as nitrite) | 100 mL | B0190634 |
| Pd | 1% Pd (as nitrite) | 50 mL | B0109635 |
| NH ₄ H ₂ PO ₄ | 10% NH ₄ H ₂ PO ₄ | 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

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

009547_01