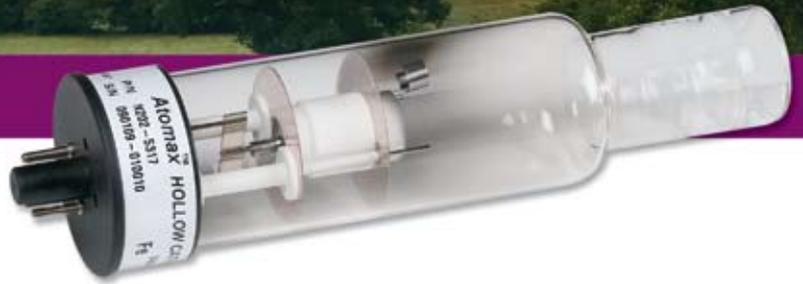




Atomax Hollow Cathode Lamps



High Quality 1.5" Hollow Cathode Lamps – Designed for Any AA Instrument

Atomax 1.5" AA Lamps from PerkinElmer, the World Leader in Atomic Absorption

New Atomax™ 1.5" hollow cathode lamps are crafted to the same high standards that have made PerkinElmer 2" Lumina™ hollow cathode lamps the best in the industry. Drawing on more than 40 years of HCL design and manufacturing expertise, Atomax lamps now offer laboratories around the world a dependable, top-quality lighting source, regardless of the brand of AA instruments used.

Value-Priced and Fully Compatible with 1.5" Format AA Instruments

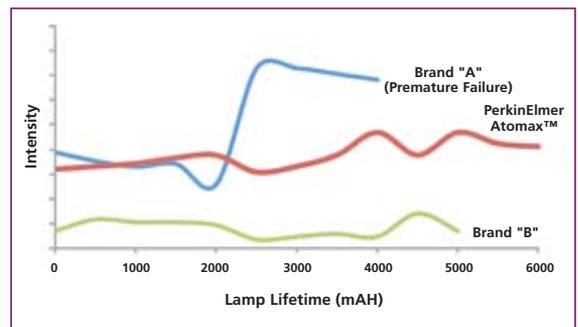
PerkinElmer Atomax lamps are designed to fit most commercially available atomic absorption instruments including those manufactured by Varian, Inc., Thermo Fisher Scientific Inc., Shimadzu Scientific Instruments, GBC Scientific Equipment Pty Ltd., Analytik Jena AG and Hitachi, Ltd. Combining affordability and industry-wide compatibility allows users of other brands of AAs to benefit from PerkinElmer's proven lamp performance. Each lamp is thoroughly tested to ensure that it will stand up to the most demanding requirements. Atomax lamps represent an exceptional value, and our money-back guarantee means you can order with complete confidence.



Every Atomax lamp includes a certificate of quality, our promise of outstanding performance levels from the leader in AA.

Tested and Validated for Highest Quality Performance

All 1.5" Atomax hollow cathode lamps are thoroughly tested to ensure maximum light output and spectral purity. Comparative results show that Atomax lamps provide equal or better performance than other AA lamps in the market. Precision engineered for high sensitivity, stable light output, low noise, and long life, you can rely on Atomax lamps for the accurate results you need.



Atomax lamps demonstrate high initial intensity and minimal variation to the lamp warranty period (5,000 mAH) and beyond, comparing favorably with other AA lamps tested under the same conditions.

PerkinElmer Offers a Complete Selection of Atomax 1.5" Single-Element Atomic Absorption Lamps

Element Name	Symbol	Part Number
Aluminum	Al	N2025301
Antimony	Sb	N2025347
Arsenic	As	N2025302
Barium	Ba	N2025305
Beryllium	Be	N2025306
Bismuth	Bi	N2025307
Boron	B	N2025304
Cadmium	Cd	N2025309
Calcium	Ca	N2025308
Cesium	Cs	N2025312
Chromium	Cr	N2025311
Cobalt	Co	N2025310
Copper	Cu	N2025313
Dysprosium	Dy	N2025314
Erbium	Er	N2025315
Europium	Eu	N2025316
Gadolinium	Gd	N2025319
Gallium	Ga	N2025318
Germanium	Ge	N2025320
Gold	Au	N2025303
Hafnium	Hf	N2025321
Holmium	Ho	N2025323
Indium	In	N2025324
Iridium	Ir	N2025325
Iron	Fe	N2025317
Lanthanum	La	N2025327
Lead	Pb	N2025339
Lithium	Li	N2025328
Lutetium	Lu	N2025329
Magnesium	Mg	N2025330
Manganese	Mn	N2025331
Mercury	Hg	N2025322
Molybdenum	Mo	N2025332
Neodymium	Nd	N2025335

Element Name	Symbol	Part Number
Nickel	Ni	N2025336
Niobium	Nb	N2025334
Osmium	Os	N2025337
Palladium	Pd	N2025340
Phosphorus	P	N2025338
Platinum	Pt	N2025342
Potassium	K	N2025326
Praseodymium	Pr	N2025341
Rhenium	Re	N2025344
Ruthenium	Ru	N2025346
Rhodium	Rh	N2025345
Rubidium	Rb	N2025343
Samarium	Sm	N2025351
Scandium	Sc	N2025348
Selenium	Se	N2025349
Silicon	Si	N2025350
Silver	Ag	N2025300
Sodium	Na	N2025333
Strontium	Sr	N2025353
Tantalum	Ta	N2025354
Terbium	Tb	N2025355
Tellurium	Te	N2025356
Thallium	Tl	N2025359
Thorium	Th	N2025357
Thulium	Tm	N2025360
Tin	Sn	N2025352
Titanium	Ti	N2025358
Tungsten	W	N2025362
Vanadium	V	N2025361
Ytterbium	Yb	N2025364
Yttrium	Y	N2025363
Zinc	Zn	N2025365
Zirconium	Zr	N2025366

For secure and easy-to-use online ordering, please go to www.perkinelmer.com/atomax.

With our precision-designed products, simple ordering process and best-in-class customer service, you can enjoy the peace of mind that comes with PerkinElmer's reliable support. For more information or to discuss your specific needs, contact us today.

PerkinElmer Warranty

Every Atomax lamp is warranted for 5,000 milliampere-hours.*

*See Atomax warranty card for complete details.

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 © 2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

008991_01

Printed in USA