

SELECT YOUR AA HOLLOW CATHODE LAMP BY ELEMENT

| | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
|----------------------|-----------------------|-----------------------|-----------------------|--------------------------|------------------------|------------------------|-----------------------|-----------------------|------------------------|-----------------------|-------------------------|-------------------------|-----------------------|--------------------------|-----------------------|----------------------|---------------------|----------------------|---------------------|---------------------|--------------------|---------------------|----------------------|------------------|---------------------|----------------------|------------------|---------------------|----------------------|----------------------|-------------------|
| 1 H hydrogen | 2 He helium | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| 3 Li lithium | 4 Be beryllium | 5 B boron | 6 C carbon | 7 N nitrogen | 8 O oxygen | 9 F fluorine | 10 Ne neon | | | | | | | | | | | | | | | | | | | | | | | | |
| 11 Na sodium | 12 Mg magnesium | 13 Al aluminum | 14 Si silicon | 15 P phosphorus | 16 S sulfur | 17 Cl chlorine | 18 Ar argon | | | | | | | | | | | | | | | | | | | | | | | | |
| 19 K potassium | 20 Ca calcium | 21 Sc scandium | 22 Ti titanium | 23 V vanadium | 24 Cr chromium | 25 Mn manganese | 26 Fe iron | 27 Co cobalt | 28 Ni nickel | 29 Cu copper | 30 Zn zinc | 31 Ga gallium | 32 Ge germanium | 33 As arsenic | 34 Se selenium | 35 Br bromine | 36 Kr krypton | | | | | | | | | | | | | | |
| 37 Rb rubidium | 38 Sr strontium | 39 Y yttrium | 40 Zr zirconium | 41 Nb niobium | 42 Mo molybdenum | 43 Tc technetium | 44 Ru ruthenium | 45 Rh rhodium | 46 Pd palladium | 47 Ag silver | 48 Cd cadmium | 49 In indium | 50 Sn tin | 51 Sb antimony | 52 Te tellurium | 53 I iodine | 54 Xe xenon | | | | | | | | | | | | | | |
| 55 Cs cesium | 56 Ba barium | 57 La lanthanum | 58 Ce cerium | 59 Pr praseodymium | 60 Nd neodymium | 61 Pm promethium | 62 Sm samarium | 63 Eu europium | 64 Gd gadolinium | 65 Tb terbium | 66 Dy dysprosium | 67 Ho holmium | 68 Er erbium | 69 Tm thulium | 70 Yb ytterbium | 71 Lu lutetium | 72 Hf hafnium | 73 Ta tantalum | 74 W tungsten | 75 Re rhenium | 76 Os osmium | 77 Ir iridium | 78 Pt platinum | 79 Au gold | 80 Hg mercury | 81 Tl thallium | 82 Pb lead | 83 Bi bismuth | 84 Po polonium | 85 At astatine | 86 Rn radon |
| 87 Fr francium | 88 Ra radium | 89 Ac actinium | 90 Th thorium | 91 Pa protactinium | 92 U uranium | 93 Np neptunium | 94 Pu plutonium | 95 Am americium | 96 Cm curium | 97 Bk berkelium | 98 Cf californium | 99 Es einsteinium | 100 Fm fermium | 101 Md mendelevium | 102 No nobelium | | | | | | | | | | | | | | | | |

COLOR CODE LEGEND

| FLAME TYPE | SPECTROSCOPY TECHNIQUE | | | |
|----------------------------|------------------------|---|----|---------|
| Air/Acetylene | FAAS | Flame Atomic Absorption Spectroscopy | MP | MP-AES |
| N ₂ O/Acetylene | GFAA | Graphite Furnace Atomic Absorption Spectroscopy | ES | ICP-OES |
| Not detectable by AA | VGA | Vapor Generation Atomic Absorption Spectroscopy | MS | ICP-MS |

◆ Multi-Element Lamp (See ordering table on right)

ORDERING INFORMATION

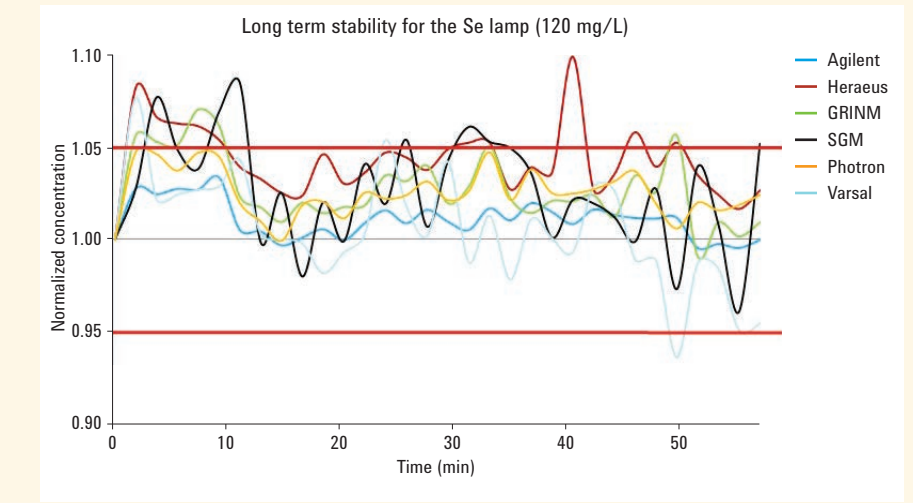
WHAT MAKES AGILENT LAMPS DIFFERENT?

- Optimum performance:** Proprietary cathode composition and unique lamp processing procedures ensure best long-term intensity, sensitivity, and stability.
- Longer service life:** Typical lifetime for Agilent lamps exceeds 5,000 mA hours for a lower cost of ownership.
- Better stability:** Agilent lamps feature an active "getter" spot that prolongs lamp life and improves stability.
- High sensitivity:** Our proprietary cathode composition and optimized operating parameters deliver the best S/N performance extending detection capabilities, even at trace levels.
- Lamps work right out of the box:** Prior to shipment, every Agilent lamp is analytically tested and conditioned to ensure that it meets demanding standards for intensity, noise, and stability.

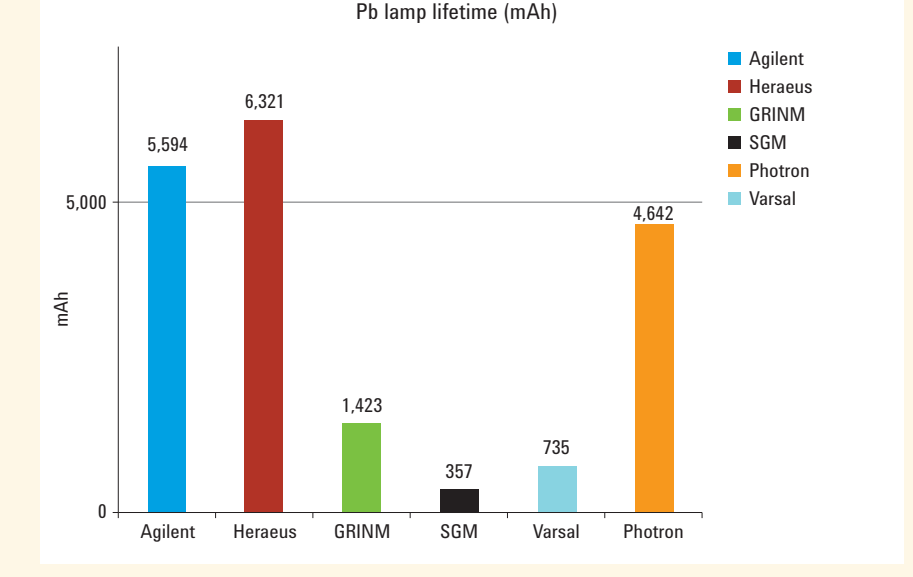
Lamp selection guidelines

- Coded lamps**
- Element coded for automatic recognition by the Agilent AA instrument
 - Instrument automatically selects the right lamp for analysis—even if you've moved it
 - Reduces operator error when working with multiple lamps
- Uncoded lamps**
- The most economical choice
 - Compatible with all Agilent and most other AA systems (except PerkinElmer and Shimadzu systems using SR correction)
 - Same excellent performance provided by Agilent coded single-element lamps

Proof of long-term stability
Agilent lamps provide stable output after 10 minutes of warm-up, eliminating drift and reducing measurement errors.



Comparison of (Se) lamps for a 120 mg/L calibration standard. The average precision for the Agilent lamp was < 1% RSD over this 1-hour period, compared with < 3% RSD for the Agilent lamp.



Comparison of lead (Pb) lamp lifetimes. The Agilent lamp gave the second longest lifetime, even with this volatile element.

- ### AGILENT MULTI-ELEMENT HOLLOW CATHODE LAMPS
- Unique element combinations extend the versatility of any AA instrument
 - Economical operation, since a single lamp covers multiple elements
 - Save time by eliminating lamp warm-up when switching elements
 - Same performance as Agilent single-element lamps (at suggested conditions)

| Coded Multi-Element Lamps | Part No. |
|--|------------|
| Aluminum/Calcium/Magnesium – Al/Ca/Mg | 5610108900 |
| Calcium/Magnesium – Ca/Mg | 5610107100 |
| Cobalt/Chromium/Copper/Iron/Manganese/Nickel – Co/Cr/Cu/Fe/Mn/Ni | 5610107600 |
| Copper/Iron/Manganese/Zinc – Cu/Fe/Mn/Zn | 5610109600 |
| Copper/Iron/Silicon/Zinc – Cu/Fe/Si/Zn | 5610109700 |
| Copper/Zinc – Cu/Zn | 5610119200 |
| Silver/Cadmium/Lead/Zinc – Ag/Cd/Pb/Zn | 5610108700 |
| Silver/Chromium/Copper/Iron/Nickel – Ag/Cr/Cu/Fe/Ni | 5610109500 |
| Sodium/Potassium – Na/K | 5610107000 |

| Uncoded Multi-Element Lamps | Part No. |
|--|------------|
| Calcium/Magnesium – Ca/Mg | 5610129100 |
| Cobalt/Chromium/Copper/Iron/Manganese/Nickel – Co/Cr/Cu/Fe/Mn/Ni | 5610129200 |
| Copper/Zinc – Cu/Zn | 5610129300 |
| Sodium/Potassium – Na/K | 5610129000 |

AGILENT HIGH INTENSITY ULTRAA LAMPS

UltraAA lamps are high-intensity, boosted discharge hollow cathode lamps for Agilent AA instruments that provide increased emission intensity, enhance sensitivity by up to 40%, and improve detection limits, allowing determinations at even lower levels.

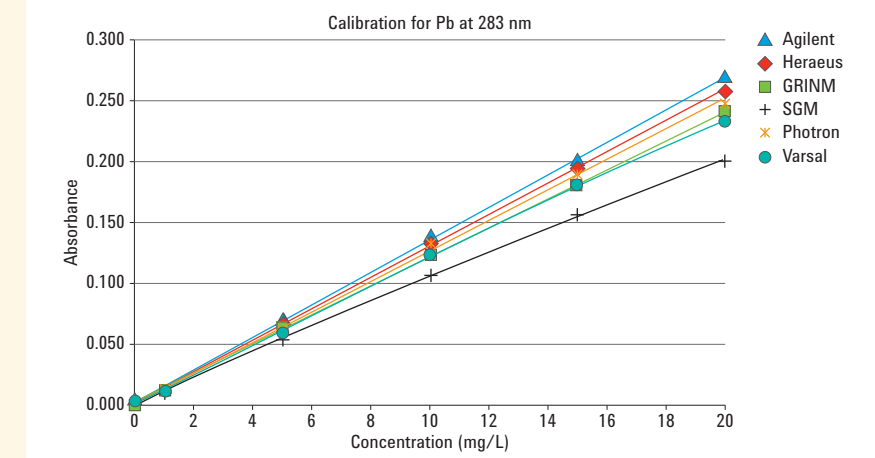
| Coded Multi-Element UltraAA Lamps | Part No. |
|--|------------|
| Aluminum/Calcium/Magnesium – Al/Ca/Mg | 5610133600 |
| Cobalt/Chromium/Copper/Iron/Manganese/Nickel – Co/Cr/Cu/Fe/Mn/Ni | 5610134500 |
| Cobalt/Molybdenum/Lead/Zinc – Co/Mo/Pb/Zn | 5610135200 |
| Copper/Iron/Manganese/Zinc – Cu/Fe/Mn/Zn | 5610135000 |
| Copper/Iron/Silicon/Zinc – Cu/Fe/Si/Zn | 5610135100 |
| Copper/Zinc – Cu/Zn | 5610134600 |
| Silver/Cadmium/Lead/Zinc – Ag/Cd/Pb/Zn | 5610108900 |
| Silver/Chromium/Copper/Iron/Nickel – Ag/Cr/Cu/Fe/Ni | 5610134900 |

| Uncoded Multi-Element UltraAA Lamps | Part No. |
|-------------------------------------|------------|
| Arsenic/Copper/Iron – As/Cu/Fe | 5610135300 |
| Nickel/Selenium – Ni/Se | 5610135400 |

| Background correction lamps | Part No. |
|---------------------------------------|--------------|
| Deuterium lamp for Agilent AA systems | 68431-800000 |

Lamps for PerkinElmer instruments
Agilent offers a range of single-element lamps that are equivalent in performance—and a suitable alternative—to PerkinElmer Lumina lamps used with PerkinElmer AA systems.
For more details, see: www.agilent.com/chem/pelamps

Reduce troubleshooting by choosing Agilent AA lamps
High-quality Agilent lamps generate a narrow, interference-free emission line, with good intensity and signal-to-noise to ensure proper calibration. Lamps from alternate suppliers can show inconsistent performance, which affects the accuracy and reliability of your results.



Calibration curves for lead lamps: The Agilent lamp provides the best sensitivity and linearity.

Get more from your elemental analysis:
www.agilent.com/chem/productivityspectro

To learn more about the advantages of Agilent lamps and testing methodologies, see: www.agilent.com/chem/aalampcomparison
Protect our environment by recycling your used lamps. Learn more at: www.agilent.com/chem/aarecycle



To order your AA hollow cathode lamps, visit www.agilent.com/chem/aalamps
To find a local Agilent representative, go to www.agilent.com/chem/contactus

