

# HIGH QUALITY AND PERFORMANCE IN ATOMIC ABSORPTION GRAIN TESTING

## How can the PinAAcle 900 aid in my grain analysis?

Consumers are exposed to low levels of heavy metals on a daily and long term exposure can have negative health impacts. Since the elements themselves are distributed unevenly throughout, for example, cereal grain, with the germ and the outer layers having the highest concentrations, analyzing these grains to detect low analyte levels with accuracy and controlled reproducibility is a challenge.

The atomic absorption analysis technique provides a high performance option with features like a closed-furnace design that is sealed at both ends with easily removable bayonet-mount windows. In addition, independently controlled external and internal gas streams provide maximum flexibility, tube life, and sensitivity.

Systems like the PinAAcle™ 900 AA spectrometer make it faster and easier to get from sample to results by reducing your grain method development time, while PerkinElmer consumables and superior services will keep your lab at peak performance.



PinAAcle 900 Series  
Atomic Absorption Spectrometers



Lumina Hollow Cathode Lamp



THGA Graphite Tubes



# What consumables are needed for my grain applications?

Invest in the consumables that are tested and validated to fit your analytical needs. For repeatable and accurate results that are on-time throughout the lifetime of your PinAAcle 900, trust the consumables engineered for your instrument. From Lumina hollow cathode lamps that allow automatic setup to the breadth of certified standards to yield reliable results, you can relax knowing you have the best.

Our specialized consumables have the ability to control operating costs and increase your laboratory uptime. Try them in your lab today.

## Confidently select your PinAAcle 900 consumables here:

- ▶ [AA Lamps](#)
- ▶ [Flow Injection Supplies](#)
- ▶ [Graphite Tubes](#)
- ▶ [Nebulizers](#)
- ▶ [Sample Cups](#)

[www.perkinelmer.com/category/aa-consumables](http://www.perkinelmer.com/category/aa-consumables)



Get the complete solution by getting the most out of your instruments, systems, processes and data. OneSource instrument services, including comprehensive instrument service and repair, asset management, and analytics services, you have the information you need, in a format you need it, to maximize your lab's productivity.

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 ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014791\_01

PKI

# Some Grain Analysis Applications to Suit Your Lab



[Fast Digestion Analysis of Lead and Cadmium in Rice Using GFAAS with Longitudinal Zeeman Background Correction](#)



[Fast Digestion Analysis of Lead and Cadmium in Rice Using GFAAS with Deuterium Background Correction](#)



[Analysis of Micronutrients in Fortified Breakfast Cereal by Flame AA Using Microwave Digestion and FAST Flame Sample Automation](#)

