

# MEASURE PULSE PRODUCT QUALITY



DA 7250™  
Diode Array NIR Analysis System



IM 9500  
Inframatic 9500 NIR Grain Analyzer

## Maximize Performance and Quality



### Capable

Measure multiple components – moisture, protein, starch, and fat content – simultaneously in seconds. Tests whole beans, lentils, peas, flours, hulls, protein concentrates, and other sample forms with little or no sample prep. It's used at material intake, for process control, and to monitor final products.



### Accurate

Our instruments have similar accuracies to reference methods of analysis. They are designed to optimize reproducibility and transferability of calibrations. Additionally, the simple sample handling, operation, and data storage increase operational accuracy significantly.



### Reliable

The DA 7250 has no moving optical components and is housed in an IP65 case making it dust proof and water resistant. These ensure maximum instrument up time and low cost of ownership. The IM 9500 is a whole grain instrument designed for ease-of-use and rugged environments. Both come with a one-year warranty and a spare replaceable lamp.

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)

  
**PerkinElmer**  
For the Better

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

Copyright ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.