



## OPTIMIZE YOUR NEW INSTRUMENT WITH PREVENTATIVE MAINTENANCE (PM) VISITS

**OneSource**  
Laboratory Services

### OneSource Warranty Plus Option for New Instruments

During your first year of new instrument ownership, you can rest assured you have the support you need to keep it up and running through your instrument warranty. But what if you need to have a documented Preventative Maintenance (PM) completed, too? With the OneSource Services Warranty Plus option, you can add a PM to your first-year warranty, allowing you to meet regulatory compliance requirements. Warranty Plus is only available during the first year warranty. Take advantage now.

#### With Warranty Plus, you will benefit from:

- Maximizing instrument performance by reducing downtime risk
- Proactively ensuring accurate and reproducible results
- Detecting and correcting instrument issues before failure occurs
- Optimizing both operating and capital expenses

To learn more about Warranty Plus option, visit: [www.perkinelmer.com/warrantyplus](http://www.perkinelmer.com/warrantyplus) and ask your Sales Representative to add it to your purchase.

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.