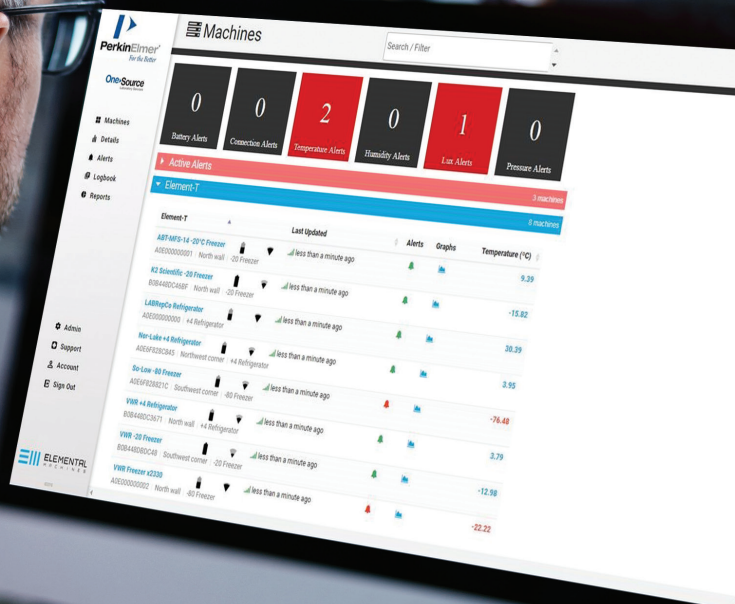


# COMPLETE CONFIDENCE WITH 24/7 LAB MONITORING

**OneSource**  
Laboratory Services



## Introducing the Asset Genius solution for wireless monitoring of assets and lab conditions.

Boost lab productivity and accelerate the pace of innovation by improving the process for monitoring laboratory instruments and environmental conditions. With the Asset Genius™ solution for laboratory monitoring, you can return time to your scientists by automating the collection, visualization, and reporting of many of your most critical lab assets. And with 24/7 monitoring and alert notifications, you can have total confidence in your assets – and your lab environment.

The Asset Genius solution includes intelligent wireless sensors that automatically measure temperature, humidity levels, and more, in a single, comprehensive solution. And by using state-of-the-art data analytics, Asset Genius can give you essential information on equipment utilization as well – all integrated into easy-to-understand interactive dashboards for deeper insights and better decision making.

With the Asset Genius solution, you're on your way to the connected lab of the future.

### Improved Productivity with Remote Monitoring

- Monitors lab and equipment via WIFI with cellular backup
- Deploy and configure sensors in minutes, with minimal IT support
- Access data anytime from your computer or mobile device
- Visualize temperature and asset data to better understand performance and trends
- View current and historical data to support compliance and exception requirements
- Manage user access to equipment and configuration parameters

For more information visit us online at [www.perkinelmer.com/onesource](http://www.perkinelmer.com/onesource)

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 ©2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.