

HOW ARE YOU MEASURING UP?

OneSource
Laboratory Services

The OneSource® Analytics Platform you have come to rely on just got better.

These capabilities allow you to analyze not only your performance across your own laboratories and sites, but also across the OneSource global customer base. These capabilities allow for peer group benchmarking comparisons to identify opportunities to improve your lab efficiency and performance.

Global Peer Benchmarking provides an overall program review across key metrics such as downtime, service response time and first time fix rates. Quickly drill into details relating to these metrics to understand the underlying factors that are driving the performance. These comparisons across peers groups can be as detailed as by manufacturer and model. With the interactive dashboards, analysis such as summary of your assets with the highest service events, by age and with the most downtime is quickly available enabling faster decisions to improve lab performance.

Get the 360 degree view of lab performance with the OneSource Analytics platform now with Global Peer Benchmarking to optimize your lab performance. Understand your competitive positioning relative to your peers to focus and improve your capabilities to gain a strategic advantage.



Global Peer Benchmarking is the newest dashboard within the OneSource analytics platform.

With Global Peer Benchmarking you can now:



Benchmark against your peers across regions and asset types to identify improvement areas such as service levels



Isolate best-in-class performance for regions and sites to adapt as best practices



Learn about key metrics and trends down to specific instrument models



Identify competitive strengths and weakness relative to your peers



Make data driven decisions for key initiative such as capital planning

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com


PerkinElmer
For the Better

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® and OneSource® are registered trademarks of PerkinElmer, Inc. All other trademarks are the property of their respective owners.
014312_01