

# TAKE YOUR ASSET MANAGEMENT PROGRAM TO THE NEXT LEVEL

**OneSource**  
Laboratory Services

## OneSource Asset Genius™ Solution: Business Intelligence Driven from Laboratory Asset Utilization

Making smarter decisions around how laboratory assets are deployed, utilized and managed throughout their lifecycle is an effective goal for laboratories as they work to drive improved overall efficiencies. Data is a powerful component in making smart business decisions, but the challenge laboratories have faced is how to quickly and easily collect and compile the data required to drive those decisions.

## OneSource Asset Genius Solution: The Data Combination You've Been Looking For

The Asset Genius solution is a cloud-based business intelligence tool that automatically collects and seamlessly combines instrument utilization, injection throughput and OneSource Asset Management service data from laboratory instruments across the enterprise, quickly turning data into insights.

- Securely capture utilization and injection throughput data from across the organization and correlate with instrument downtime, service activity, age, and analytical method use
- Scalable cloud-based architecture with encrypted communication supports data capture without the need for software installed on target instrument systems
- User-friendly interactive dashboard provides an array of visually-enhanced, multidimensional reports with easy-to-understand metrics
- 360° performance picture supports optimized resources, workloads and inventory lifecycles

## Instrument-agnostic Data Capture

The Asset Genius solution automatically captures utilization data from any PC-based instrument in the lab. Data is captured transparently without requiring any additional effort on the part of scientists or lab techs, and without workflow disruption.

## DRIVE LABORATORY OPTIMIZATION WHILE REDUCING COSTS

- Capital Expenditure Planning
- Equipment Resource Management
- Maintenance Plan Optimization

## Actionable Insight through Data Visualization

Using a pre-configured suite of reports, scientific and operations leaders can view aggregated instrument data in a single system that provides enterprise-wide insight into asset utilization metrics.

### Capital expenditure planning

Identify minimally used equipment for decommissioning or repurposing to other labs where needed; provide data driven justification for new instrument purchases.

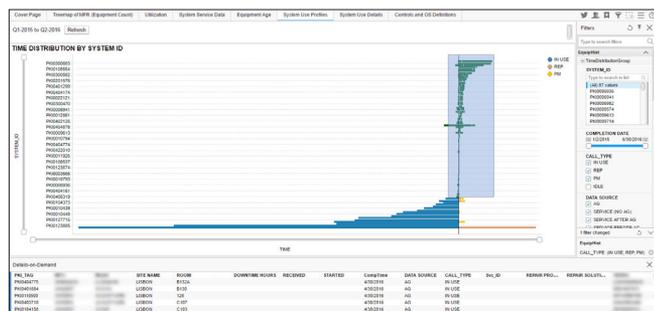


Figure 1: Quickly identify equipment with high service activity and low utilization.



Figure 2: Correlate with age at the click of a button.



Figure 3: Correlate with high service activity.

For more information about the Asset Genius solution or any of our IT solutions, please contact your PerkinElmer OneSource Laboratory Services representative, or visit [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)

### Maintenance plan optimization

Identify instruments with low utilization and high PM activity in seconds

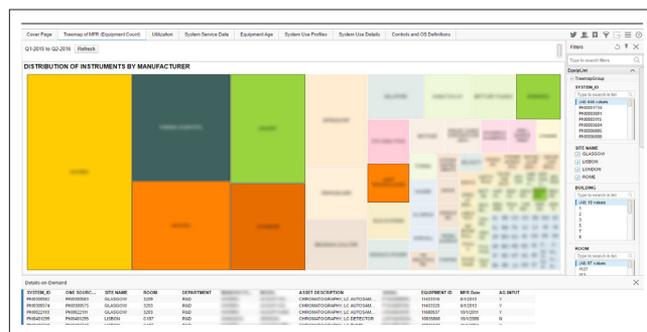


Figure 4: See the instruments with low utilization but high PM activity to identify savings opportunities.

### Increase scientific throughput and overall laboratory efficiency

The Asset Genius solution collects and reports utilization information for any PC-based laboratory asset and provides laboratory leaders with the insight they need to make data-driven decisions that capitalize on the strengths and effectiveness of their operations.

These decisions help laboratories enhance their competitive and scientific advantage by increasing scientific throughput, supporting cost certainty and increasing overall optimization of laboratory operations.



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

Copyright © 2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.