



ACCESS TO LATEST ASSET LOCATION AND MOVEMENT

Do you find yourself spending time on tracking and locating your lab equipment instead of spending time on science?

You're not alone. No matter what type of laboratory you work in, implementing an integrated RFID asset tracking technology makes it faster, and easier than ever, to take full inventory of your physical laboratory assets.

With PerkinElmer's passive RFID technology, you can now accurately and efficiently monitor and manage your assets.

We took the best RFID equipment available and integrated it with our state-of-the-art mobile application to deliver an end user friendly package. With the RFID packaged solution and implementation team the introduction of advanced asset location services is seamless.

Passive RFID technology from OneSource will:

- Eliminate manual data collection, analysis, and reporting for lab and assets
- Reduce manual errors, and streamline compliance reporting
- Provide accurate and up-to-date asset location information returning time spent searching for equipment

With an advanced level of inventory tracking and management through the use of a mobile app and an interactive dashboard, it is the perfect solution for labs of any size.

- Improved inventory accuracy
- Increase in service efficiency
- Reduction in resources & time to complete inventory
- Increased inventory throughput – many tags can be read at once
- Decreased instrument misplacement
- Visibility to asset chain of custody
- Correlation of asset movement and service history
- Reduction in maintenance, insurance, and depreciation costs for missing equipment

Learn more at www.perkinelmer.com/uk/product/asset-location-services-onealo

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



For a complete listing of our global offices, visit 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.