



SCIENCE AND INFORMATION
TECHNOLOGY ON THE SAME TEAM
IS A WINNING STRATEGY

OneSource[®]
Laboratory Services

BRIDGE THE GAP BETWEEN YOUR INSTRUMENTATION AND YOUR INFORMATION TECHNOLOGY

In today's complex lab environments, your instrumentation and your information technology need to work together seamlessly to support the critical work your scientists do. Our OneSource Information Services team is the only global life sciences organization that brings science and IT together to fully understand how your workflow works – so you can accelerate your scientific productivity and streamline all your processes.

Part of our comprehensive OneSource® Laboratory Services offering, OneSource Information Services helps labs optimize their scientific productivity at the intersection of instruments, computing, and software with effective use of information technology. These solutions combine scientific and technical expertise with well-structured project management and IT process methodologies to create predictable, measurable IT outcomes regardless of instrument technology or manufacturer.

Lab Computing: Meeting the challenge of benchtop support and management

With limited budgets and a lack of in-house expertise for creating cost-effective, reliable scientific IT processes, implementing and managing multivendor IT can be burdensome on your laboratory and IT staff. And new software initiatives often run into deployment challenges due to the unpredictability of these complex, ever-changing environments.

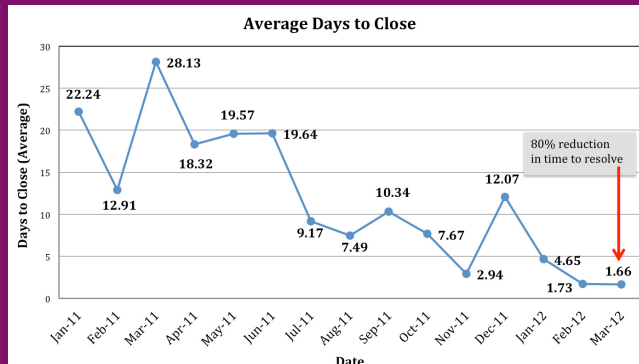
What's more, with multiple vendors, the burden of day-to-day support falls to your scientists, who can spend up to 20% their time on noncore, nonscientific issues. And you pay the cost in excessive vendor support and lost productivity.

For many organizations, a managed service is the most efficient, affordable option: So our Lab Computing offering combines deep scientific-instrumentation and IT expertise to provide a single point of accountability for a fully connected and digitally optimized laboratory environment. A Lab Computing analyst quickly gets you engaged with the right personnel, with the right skill sets in instrumentation, application software, operating systems, networking, and computing hardware, for fast problem resolution. If additional support is needed, you have access to a wealth of information via our training hubs and remote support.

Each vendor-neutral plan is tailored to your IT and compliance landscape. Scope of work can include onboarding and disposition of assets, break/fix repair, PC maintenance and software upgrades, software and instrument licensing, and security/compliance support. Plus, you'll have access to our IT Health Dashboard and custom KPI reports. The Lab Computing analyst also takes over vendor and contract management for your lab IT assets. This managed approach allows you to minimize scientific downtime and accelerate discovery.

Lab Computing's Proven Impact: Improved asset uptime and scientist productivity

On average, see an
80% drop
in time between an open and closed
Lab IT service request



Computer Systems Validation: Ensuring better compliance – and better productivity

We provide a full range of computer systems validation (CSV) services to support your laboratory, manufacturing, and enterprise systems. We take a risk-based approach according to GAMP 5 guidelines, which includes planning (Validation Master Plan), requirements design (design, configuration, function, and user) and verification (IQ/OQ/PQ, trace matrix, and summary report). This risk-based approach streamlines the validation process, helping to control costs and reduce system downtime. We can also customize our validation delivery to align with your company's policies. Each CSV project thoroughly assesses the systems' software for 21CFR Part 11 compliance – a current area of intense FDA scrutiny – and includes an evaluation of the system's data integrity, security (physical and logical), audit trail, data back-up/restore/archive, disaster recovery, as well as electronic records and signatures.

Our team of validation specialists possesses broad experience validating hundreds of systems in multiple regulated industries. What's more, we can staff your validation project with the specific expertise required for your systems and regulatory environment.

Scientific Application Services: From development to integration to deployment

Our Scientific Application Services provides end-to-end management and deployment of new and existing systems and customized application development. As your IT platforms grow over time, so does the need for proper support and management. Our scientific and laboratory applications management solutions deliver a unique blend of scientific knowledge coupled with an expertise in IT and enterprise database systems, for a wide array of needs areas: technology support, user support, production support, application enhancements and upgrades, and reporting functions through our Application Maturity Program.

For laboratory automation projects, we can create feature rich, predictable software solutions, on time and on budget. Our iterative software development framework allows for early and frequent user access to software functionality and design, enabling feedback and corrective actions early in the development lifecycle. Plus, we provide application configuration and data integration with other critical systems within the laboratory environment.

With a wide variety of third-party software applications available for science and the lab, you're faced with the challenge of selecting and ensuring the successful deployment of these technologies at a platform level. We've got the implementation experience for laboratory information management systems (LIMS), electronic laboratory notebooks (ELNs), laboratory execution systems (LES), and chemistry and biology Informatics to help ensure project success and integration.

OneSource Laboratory Services: With you every step of the way – to reach your best outcomes

Our OneSource® Laboratory Services team delivers solutions that cover all aspects of your operations and can be customized for the scientific workflows and business outcomes you're striving toward. With proven success at the world's top pharmaceutical and environmentally focused organizations, and managing more than 500,000 assets worldwide, the OneSource team provides a complete services offering with global reach and local support. Together, we're taking business results to the next level.



OneSource Laboratory Services

- Asset Services
- Analytics Platform
- Relocation Services
- Information Services
- Compliance Services
- Scientific Services

To learn more about OneSource Information Services and additional services from OneSource Laboratory Services, visit www.perkinelmer.com/InformationServices or contact your PerkinElmer OneSource representative.

Learn more at 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



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

Copyright © 2017, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

012341_03