



HUMAN HEALTH

ENVIRONMENTAL HEALTH

WE'RE TURNING
GOOD STRATEGIES
INTO GREAT SCIENCE

OneSource Strategic Lab Solutions


PerkinElmer[®]
For the Better

TO SOLVE PROCESS CHALLENGES, START HERE

We hear it every day: “We *know* the direction we want to go. We just don’t have the resources” – or the people, or know-how, or buy-in – “to get there.” Like most businesses, laboratories are being pressured on all sides to reduce costs by automating processes, improving data management and data sharing, and moving toward a more streamlined, paperless approach to lab operations. And with data generated by scientific instrumentation and sample analysis growing exponentially, all the time, it’s a situation that can overwhelm even the best-run labs.

IT'S WHAT GETTING BACK TO SCIENCE IS ALL ABOUT

Most labs don't maintain the in-house expertise needed to plan and implement sweeping projects to transform the everyday chaos of most lab operations into the sleek, efficient, effective lab of the future.



After all, you should be concentrating on your core competency – your *science* – and leaving the infrastructure, operations, and even lab management to the professionals.

That's why so many smart labs are looking to outsourcing to help them plan, design, deploy, and manage their lab environments. Problem is, few companies possess the grounding in both scientific instrumentation, lab-specific business processes, *and* information technology – skills that are absolutely necessary when undertaking major lab projects.

But one organization has those skills in abundance. OneSource® Laboratory Services.

OneSource Strategic Lab Solutions: We deliver *change*

As an integral part of our OneSource Laboratory Services, OneSource Strategic Lab Solutions addresses your need to continuously improve your lab's efficiency, while reducing its costs. So we take into account not only your instruments, but also your IT systems, staff, and processes. We then go to work identifying and delivering on change opportunities that make sense for your environment – without sacrificing your ability to meet your lab's scientific goals.

The result? You'll see increased productivity, with less risk. More effective data management and streamlined and efficient workflows. Improved access to critical information, with data going where it needs to go, right on time. Better value from your investments in people, processes, and technologies. And the knowledge that you're in compliance with all statutes and regulations that apply to your industry – and your lab. Best of all, you have a single vendor for all your lab's requirements and a single point of contact who's with you throughout your project, from inception to final signoff.

You have a lot invested in instruments, systems, processes – and people. Let us help you make the most of them.

GREAT SCIENCE STARTS WITH A SINGLE ENGAGEMENT

With OneSource Strategic Lab Solutions, our expertise in lab processes and information technology is totally focused on helping you meet your lab's specific business and scientific goals. We take a holistic approach to improving your lab environment, delivering a comprehensive suite of services that can reduce costs, improve operations, and mitigate the risks that stand in the way of taking your lab in the direction you want it to go.

Strategic Prioritization Services

Which project is the right one for your lab to invest in? It's a simple question, but a critical one. We've helped scores of labs identify, prioritize, select, and plan projects that map closely to organizational strategies and goals – and that deliver the quick ROI these labs were counting on.

Efficiency and Productivity Services

Working closely with your key stakeholders, we map out the hardware, software, services, and people power you'll need to make the transition from your current lab environment to an efficient, effective, highly automated scientific enterprise. (See sidebar, next page.)

Scientific Knowledge Management Services

Ensuring that all your stakeholders have the information they need to make sound business and scientific decisions should be a key priority for your lab. Strategic Lab Solutions can deliver a sound Scientific Knowledge Management strategy in which data can be collected, analyzed, stored, reported – and used to support the key decisions you make every day.

Partnership and Outsourcing Services

Few labs have the resources necessary to maintain in-house expertise in both lab instrumentation and information technology – which makes the development of an outsourcing strategy and partner relationships more important than ever. OneSource Strategic Lab Solutions applies technical and financial analysis along with scientific and operations expertise to help you select the right outsourcing partners for your lab – and manage them effectively over the long haul.

Personnel Services

Every lab needs to develop its people – and to train and develop effectively, you need a plan. OneSource Strategic Lab Solutions can construct a Master Training Plan to make sure that all your personnel – both scientists and support staff – are trained in existing and newly developed procedures and that SOPs are clearly understood and followed.

Compliance Services

Compliance with the statutes and regulations that govern your industry can be a real burden on your resources – and your people. Strategic Lab Solutions provides a Master Validation Plan that identifies the best path to regulatory compliance for your lab and lays out the scope and schedule of validation activities you need to accomplish to stay fully compliant.



Strategic Lab Solutions takes a holistic approach to all your lab assets.



Digging Deeper

Success Begins with a Good Plan

If you need it, we work with you on a personalized **Lab Automation Plan** that looks at your lab as it exists today – and shows you the path to the lab you've always envisioned.

This is a tailored plan that does as little or as much as you need it to. But a complete plan describes the hardware, software, processes, and effort it will take to move to a seamless, fully automated, and paperless lab environment.

The keyword here is *integration*. The Lab Automation Plan shows you the path to integrating all your lab's information technology systems and specialized instrumentation, enabling you to protect and enhance the value of the copious amounts of data you generate every day.

Plus, the plan can assist with:

- Developing methods that speed research and protect data quality
- Providing recommendations for lab automation, incorporating systems such as ELN, SDMS, LIMS, and DMS.
- Leveraging your existing investments in systems such as SAP and Sharepoint by integrating them into your laboratory workflows
- Prioritizing the procurement activities that can help you realize your lab's business and scientific goals
- Laying out the regulatory requirements and metrics you'll need to understand to maintain compliance

What's more, a **Lab Dashboard and Analytics Plan** delivers reporting tools that improve capacity planning, productivity, resource utilization, and situational awareness – all key components of any well-functioning lab.



SAY HELLO TO THE LAB YOU ALWAYS IMAGINED

No two labs are entirely alike. And neither are any two engagements with OneSource Strategic Lab Solutions.

We've built and managed some of the most complex, sophisticated laboratories in the world. And we've helped startup labs – labs with more ideas than resources – identify and prioritize the projects that can provide a quick return on investment in the short term and support long-term growth when the time comes. No engagement is too small or too overwhelming.

What's more, we can deliver the expertise necessary to give an objective review of existing projects that may be in need of some recalibration. Maybe you've lost some key resources, missed some milestones, or are over budget. We can deliver unbiased, as-is analysis of where your project stands today, and then reset expectations with expert recommendations, risk analysis, and, if necessary, a revised plan.

Most of all, we can deliver confidence that your lab can meet its goals – and that all you need is a little support where and when it's really needed.

You do the science, we'll do the rest

Strategic Lab Solutions is part of the OneSource suite of services. OneSource is built around a team of more than 1,600 highly skilled service professionals with more than 400,000 assets under their care, in more than 120 countries around the world – all dedicated to helping you reach your financial, operational, and scientific goals. How? By becoming an integral part of your business, providing the highest level of technical support and scientific expertise.

Strategic Lab Solutions can actually serve as a front end to other OneSource services, including Analytical Method Services, Asset Procurement and Disposition, Business Intelligence Solutions, Instrument Service and Repair, Laboratory Relocation, Qualification and Validation, and Scientific IT Services.

So no matter what scientific discipline you work in or how complex a lab you run, there will come a time when a change in operations is essential to help your lab meet its short- and long-term challenges. But nothing should take your focus from where it needs to be – on your science. That's when you need a little outside expertise.

That's when you need OneSource Strategic Lab Solutions.

stra•te•gic [*struh-tee-jik*], *adjective*

1. Of, relating to, or marked by strategy <a *strategic* plan>
2. Necessary to or important in the initiation, conduct, or completion of a strategic plan
3. Of great importance within an integrated whole or to a planned effect <emphasized *strategic* points>

so•lu•tion [*suh-loo-shuhn*], *noun*

1. An action or process of solving a problem
2. An answer to a problem :EXPLANATION; *specifically*: a set of values of the variables that satisfies an equation
3. Strategic Lab Solutions : A trusted partner helping you achieve your business and scientific goals

con•sul•tant [*kuhn-suhl-tnt*], *noun*

1. One who consults another
2. One who gives professional advice or services :EXPERT

For more information, contact us at StrategicLabSolutions@perkinelmer.com



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 ©2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.