

HUMAN HEALTH

ENVIRONMENTAL HEALTH



OneSource®

ALL THE  
SCIENTIFIC  
SERVICES  
YOU'VE EVER  
IMAGINED

OneSource Scientific Services

  
**PerkinElmer®**  
*For the Better*

# DISCOVERY BEGINS BY GETTING BACK TO BASICS

What does *your* ideal lab look like? In that perfect world, everything in your lab is primed and ready to deliver on your principal goal: getting first to market with innovative products of the highest quality and reliability, all while adhering to stringent regulatory and budget requirements. And getting you to that place of discovery is what we're all about.



## Efficiency-by-Design

Our design model for lab efficiency is based on these simple action items:

### Standardize



We help create a predictable lab environment that works for everyone – scientists and nonscientists alike.

### Organize



Making sure your laboratory is mapped out for maximum efficiency is one of the key components of our offering.

### Automate



We enable you to simplify laboratory tasks through automation, including robotics, informatics, and much more.

## When things work efficiently, discovery can happen

The goal of the pharmaceutical lab is to bring high-quality drug candidates to clinical trial faster than ever before. And if your lab is like most, you're being asked to do it with fewer resources. Yet, an internal Boehringer Ingelheim study concludes that up to 25% percent of research scientists' time is spent on peripheral, noncore tasks. Bottom line, it all represents resources that could be better utilized.

But what if your researchers could spend more time concentrating on their core competency – the science – and less on mundane organizational and maintenance tasks? Who knows what you could accomplish?

That's the question OneSource® Scientific Services is designed to answer.

With OneSource Scientific Services, we believe delivering higher value to your customers shouldn't mean higher costs. To that end, we deliver a "faster, smarter" approach to laboratory workflow.

This innovative suite of services – OneSource Lab Support, OneSource Scientific Support, OneSource Laboratory Efficiency Consulting, and, of course, the white-glove

OneSource Instrument Concierge™ and Instrument Check™ – delivers a simple, streamlined, integrated framework that enables you to concentrate more of your precious resources on your research, while a dedicated team of scientists takes care of noncore lab activities. This "co-sourcing" model means a more holistic approach to laboratory efficiency, allowing you to leverage the knowledge, technology, and resources of a scientific and technology leader.

The result? An engagement with OneSource Scientific Services can optimize lab operations, significantly reduce operating costs, and help maintain compliance with an ever-changing array of regulatory mandates. Best of all, it provides a comprehensive approach to lab activities – designed by scientists, for scientists – that helps your lab gain and maintain its competitive advantage.

OneSource Scientific Services. Extraordinary science is about to happen. And it starts right here.

### Specialize



Our scientific experts can complement the skill sets you already have in place – making your whole lab more productive.

### Success



Once all the pieces are in place and your scientists can concentrate on the important work ahead, you'll be amazed at what you can accomplish.



## SCIENTIFIC SERVICES DELIVERED WITH A PERSONAL TOUCH

Studies show that the highly trained scientists you employ can spend much of their time troubleshooting and maintaining instruments, tracking spare parts and warranties, and waiting for instruments or information systems to come back online. That's time that could be much better spent doing critical research.

Plus, your research itself could be compromised by poorly monitored equipment, instrumentation that's not generating reliable data, and calibration that's carried out by unqualified technicians.

To top it all off, laboratory environments are becoming more complex all the time, testing the limits of a management regimen that's already stretched to the limit. And the failure to establish a regular inspection and maintenance program can lead to inadvertent mistakes, accidents, and other equipment-related mishaps. With all this complexity, it's become a huge undertaking to ensure instrumentation is operating correctly for its intended purpose.



## OneSource Instrument Concierge: We're at your service

With our OneSource Instrument Concierge™ service, we introduce you to an even higher level of care. We start by incorporating one or more highly trained scientists into your location, dedicated to caring for your entire multivendor laboratory environment. This is the ultimate in personalized service, tailored to your lab's specific users and their requirements. Our scientific personnel are a highly skilled, onsite presence that can deliver a deep level of instrumentation expertise, enhancing your lab's daily workflow and scientific processes.

Our Instrument Concierge specialists can also provide a wide array of laboratory capabilities, tailoring instrumentation to the needs of individual users and delivering a level of knowledge and support that goes over and above what traditional services typically provide.

OneSource Instrument Concierge specialists can give your lab:

- Integrated instrument management
- Single-source accountability
- Reduced unplanned downtime
- Improved performance and productivity
- Better data integrity
- Instrument compliance
- Improved safety

## OneSource Instrument Check: We've got performance covered

With our OneSource Instrument Check™ services, you know your instrumentation is optimized to perform the way it was designed to – and will be ready to run whenever you need it. We conduct routine instrument checks that help minimize unplanned downtime, provide continuous trending of data (with control charts) to optimize instrument performance, and deliver a proactive approach to identifying problems – and solving them.

## OneSource Lab Support: Everyday science, redefined

With OneSource Lab Support services, we handle all your noncore laboratory activities, which goes a long way toward freeing your researchers to concentrate on the science at hand. We deliver the resources your lab needs to stay ahead of the competition, including managing your consumables and lab supplies, preparing your media, weighing compounds, and other basic lab activities.

## OneSource Scientific Support: We deliver change, comfortably

With OneSource Scientific Support services, we provide the people and the know-how to run key experiments to keep your lab moving forward. As valued members of your team, OneSource specialists can provide your lab with a wide variety of laboratory capabilities, including onsite flexible staffing resources and optimized occupancy and utilization of your lab space. This can mean a real competitive advantage for your organization by ensuring data confidentiality and faster turnaround times on research and development projects.

## OneSource Laboratory Efficiency Consulting: A more effective lab – by design

We start by taking our four action items – standardize, organize, automate, and specialize – and using them to help us begin a thorough assessment of all the resources you have in place (both in technology and people). From there, we classify all your activities as being either core or noncore, giving you the opportunity to see how your lab can benefit from a strategic engagement with us.

From there, we begin to design and propose a strategic improvement roadmap with lots of short-term, achievable milestones as well as longer term projects with results that are achieved over time. We make recommendations as to how laboratory tasks can be reprioritized, and how staff can be deployed to increase productivity.



# TRUST THE ONE THAT'S ALWAYS THERE FOR YOU

For nearly two decades, OneSource has been providing essential laboratory services to pharmaceuticals, contract research organizations, and other scientific enterprises. We support more than 500,000 multivendor assets in more than 120 countries. So who better to help you organize and automate your lab, making it more efficient and effective than ever before?

We take a holistic approach to your lab environment, considering a whole range of key components:

**Lab workflows** Does your lab's physical layout support the ways your scientists interact with one another? We help develop lab workflows that simply *work*.

**Instrumentation** We can design a lab environment that meets the needs of all your scientists while supporting the legacy instrumentation you've already invested in.

**Asset management** Our onsite engineers can manage and maintain instrumentation from *any* vendor – and that means your system uptime can meet the most stringent SLAs.

**Lab IT services** We combine scientific and technical expertise with information technology to create predictable, measurable IT outcomes – regardless of technology or manufacturer.

**Analytics** By bringing together data from multiple sources, our end-to-end analytics solutions help you uncover opportunities, anticipate trends, and accelerate discovery.

**Lab support** We can handle all your operational needs, from compound management and sample preparation to downstream applications for drug discovery.

**Compliance** We help you manage the complexities of the qualification process, ensuring the compliance of every instrument and computer system.

## Operating More Efficiently –For Less



At Boehringer Ingelheim, scientists were tasked with finding new ways to streamline operations and improve efficiencies, all while reducing spending. An internal operations audit revealed that scientists were spending up to one quarter of their time on non-laboratory activities.

Once Boehringer Ingelheim teamed up with PerkinElmer OneSource, scientists found they could spend more time performing vital experiments and much less time – and fewer resources – on noncore activities. In addition, with OneSource trained engineers maintaining the company's multivendor environment, downtime was further reduced as well.

## For More Information

Want to learn more about OneSource Scientific Services' holistic approach to laboratory activities and how we can help you optimize lab operations, increase lab efficiency, and maintain compliance with regulatory mandates? Our suite of OneSource Scientific Services white papers can give you a complete overview of our offerings – and how they can help you reach your scientific and business goals.

*A New Approach to Laboratory Efficiency: On-site Instrument Concierge Service*

*A Case for Co-Sourcing*

*The "New Normal" in Laboratory Efficiency: Leaner, Faster, Smarter*

*Managing Laboratory Consumables*

**Download them all at: [www.perkinelmer.com/efficiency](http://www.perkinelmer.com/efficiency)**

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)



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

Copyright ©2015, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.