



BACK TO WORK BACK TO COMPLIANCE

OneSource
Laboratory Services

Novel workflows, new health and safety guidelines, proper distancing in your lab – navigating the aftermath of the COVID-19 pandemic is one of the biggest challenges your lab will ever face. But you don't have to go it alone.

Our proven OneSource® Services help you understand compliance impacts from your back-to-work plan and provide you insights on what others are implementing to minimize the impact to your lab operations. Our proven track record in business continuity threats and disaster recovery will guide you in:

- Helping ensure your lab's Return To Work Plan (RTWP) is safe and successful
- Determining the impacts from new practices and finding gaps with your identified guidelines, assessing best-practice approaches, and analyzing any negative impact to regulated compliance
- Providing solutions for increased compliance, including SOP revision to capture impacted processes and workflows

Labs are our livelihood. Our customers are our passion. Whether we're assisting you remotely or on site, our experienced staff is fully devoted to getting your lab back up and running, quickly, safely and compliant.

There with You, Every Step of the Way

Resuming your lab after a major pandemic won't be easy. So we work with you – step by step – to ensure your lab is safe, compliant, and optimized for success. Our comprehensive three-phased approach includes:

Phase 1

- Defining gaps between your RTWP and identified guidelines to ensure a smooth RTW experience
- Evaluating the impact from applying new COVID-19 protocols on laboratory workflows
- Analyzing impacted and new workstreams to determine gaps in documentation and risks to regulatory compliance
- Providing suggestions for a remediation plan and proposed best practices

Phase 2 (Optional)

- We'll assist and support any remediation activities outlined in your assessment

Phase 3 (Optional)

- Periodic reassessment to be sure your lab stays on track

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 ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.