

ONESOURCE MULTIVENDOR SERVICE OPTIONS

OneSource
Laboratory Services

Are you getting the experienced service you need for all of your instruments?

With so many different vendors' instruments in your lab, it can be challenging to ensure everything is being maintained properly. Some labs struggle to get the most productivity and efficiency from all their instruments, while others need to streamline and simplify workflows to maintain regulatory compliance – and reduce the risk of noncompliance. Either way, you're always scrambling to figure out who to call for service, as quickly as possible, before you lose too much time...and money.

What if there was a one-stop service contract option for your lab – from a company with decades of deep-seated multivendor experience – that repaired all your instruments, offered state-of-the-art validation and compliance services, and provided reliable preventative maintenance? That's what OneSource Multivendor Service options is all about.

Ready to reduce downtime, increase productivity, enhance compliance – and spend more time making discoveries?

Visit www.perkinelmer.com/onesourcemultivendor for more information.

BUT “ONE-STOP” DOESN'T MEAN “ONE SIZE FITS ALL”:

We offer multiple service contract options that ensure your instruments – even those that aren't manufactured by PerkinElmer – are well maintained, keeping your lab running at peak performance. Other benefits include:

- Access to a dedicated multivendor support team
- A continued relationship with your trusted PerkinElmer service engineer
- One point of contact, one request

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


PerkinElmer
For the Better

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.

014731_01

PKI