



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 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 preventive maintenance? That's what OneSource Multivendor Service options are all about.

Reduce downtime, increase productivity, enhance compliance – and spend more time making discoveries.

CUSTOMIZED TO MATCH YOUR LAB

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

- Our highly trained and trusted service support team
- Options for regularly scheduled preventive maintenance
- One point of contact, one request, one purchase order

[Click here to simplify your lab today.](#)

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 ©2021, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.