

Do More in Less Time

- Simple and powerful kinetics software
- Faster data processing from complex data sets
- Greater productivity with integrated TG-hyphenation control and real time monitoring

Search Better and Deeper

- Accurate results with enhanced ATR (attenuated total reflectance) algorithm
- Search ATR spectra against more databases than ever

Interact Easily

- Interact via smart panel or PC
- Easy standard operating protocols and diagnostics
- Improved productivity for multiple runs in busy labs



Our global network of over 1500 certified service technicians can provide in-depth training, perform software and hardware upgrades, and more.

For more details on the solution, please visit <https://www.perkinelmer.com/spectrum3>

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.