

EXTEND YOUR RANGE AND YOUR SCIENCE

Introducing the SPECTRUM 3™ FT-IR Spectrometer

Cutting-edge product development and groundbreaking innovations are built on a foundation of meaningful insights that can be trusted. From routine identification and verification of raw materials to advanced research applications, you need the sampling flexibility to quickly, confidently, and cost-effectively analyze any sample. We've distilled 75 years of innovation and leadership in Infrared spectroscopy into the Spectrum 3™ FT-IR to deliver more value, performance, and versatility than ever before.

FT-IR reimagined – with the power to empower you.

- Up to three IR ranges in one instrument for ultimate sampling flexibility
- Run more samples than ever with SmartPanel
- Better insights into reaction kinetics with faster and superior software
- Simplified hyphenation from a single provider

Simplified results, powerful collaboration

You can easily acquire, process, and share IR data with any device, anywhere, any time with Spectrum 10 software with NetPlus IR cloud connectivity. And our AssureID™ with full 21 CFR Part 11 compliance provides the most powerful solution for method development, product screening, and qualification. It's software that lets you quickly see information on the instrument and easily run routine protocols.



SPECTRUM 3™ FT-IR Spectrometer

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.