

FT-IR Spectroscopy



Spectrum Two for Polymers QA/QC

Refreshingly easy to use and maintain, Spectrum Two™ offers a readymade solution for rapid, worry-free polymer analysis. User friendly design features combined with a Polymer Application Pack ensure that with Spectrum Two, anyone can be an IR expert.

Faster Testing Implementation

In today's competitive polymer market, making efficient use of time and resources is vital. Spectrum Two can be set up in minutes to deliver the answer you need, faster than ever before.

- User-install capability allows instrument to be set up by anyone, anywhere.
- Out the box polymer reference library enables systems to go live immediately.
- Practical polymer analysis guide simplifies and speeds SOP development.
- Spectrum Two Introductory Kit allows new or inexperienced users to become familiar with the basics of IR measurements and troubleshooting common sampling issues.
- Wide range of on-site training services tailored to the needs of your laboratory, including an introduction to Spectrum Two.

Effortless Analysis

Spectrum Two is designed to be used by anyone, no matter what their level of experience. Now, operators without a scientific or technical background can achieve reliable raw material or final product verification every time, with absolute confidence:

- Configurable software allows complex analyses to be performed at the push of a single button
- Scanalyze™ technology delivers the answer you need in one single step, streamlining and accelerating operation
- Modern, robust ATR sampling minimizes sample preparation and sampling errors, decreasing analysis rework

Reduced Cost of Ownership

Minimizing instrument downtime is crucial in the polymer industry. Spectrum Two is packed with innovative technologies that ensure maximum return on investment. Patented OpticsGuard™ technology provides a unique humidity shield, protecting your instrument from detrimental environmental effects. Long life desiccant with a three year warranty reduces the maintenance burden, while the unique Dynascan™ interferometer and solid-state laser ensure reliable operation and reduce technical support calls. Plus, the unique hot-spot stabilized source increases measurement stability and extends source lifetime.

IR Ready to Go

With Spectrum Two, you can achieve accurate, dependable answers in non-traditional environments. Compact in size, battery powered and robust, the lightweight instrument can be transported to any location. You get the answer you need, wherever you are.



For more information, visit www.perkinelmer.com/spectrumtwopolymers

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 ©2011, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.