



# SIMPLIFY YOUR POLYMER ANALYSIS

## Introducing the Polymer ID Analyzer

The PerkinElmer Polymer ID Analyzer is a comprehensive Fourier Transform Infrared (FT-IR) polymer analysis system that helps you quickly and accurately identify unknown polymer samples, determine the composition of polymer blends, and verify quality of raw materials.

There is a growing demand for easy, accurate, and rapid qualitative and quantitative polymer analysis to enable increased productivity and faster product release. Whether you are a polymer producer or a buyer in food packaging, construction, or the automotive industry – accurate verification of identity, quality, and composition of materials is critical to ensure finished products hold up to your standards.

Our straightforward and easy-to-use analyzer helps reduce potential errors from your analysis, methods, and operators, enabling you to gain meaningful insights to ensure quality and optimize processes.

### Accelerate your polymer analysis

With a specialized polymer library and tools to accelerate analysis for a wide range of polymers, anyone – even those with minimal training – can easily operate the instrument.

### The Polymer ID analyzer helps you:

- Confidently analyze a wide range of polymers and complex blends
- Maximize productivity with simple and rapid verification
- Ensure quality control of raw materials, in-process samples, and finished products
- Reduce operational and training costs
- Expand revenue by analyzing at the location of your sample – don't lose time and money getting samples to the lab
- Save time and maximize revenue by analyzing at the location of your sample

## An easy to use analyzer package for maximum productivity

- Spectrum Two™ FT-IR
- Spectrum Two UATR (Attenuated Total Reflectance accessory)
- Enhanced Polymer Resource Pack – including updated starter libraries of homopolymers and blends
- Spectrum Touch Developer with Macro – a simplified, easy-to-use interface for running a pre-designed application for polymer analysis
- MultiSearch™ – for ready estimation of polymer mixtures
- User supplement document
- Workflow video



## Results in one touch

Simplify your polymer analysis workflow with easy and efficient data collection, processing, and generation of results. Create routine operating protocols and lower the likelihood of operator error with our powerful Spectrum 10™ and Spectrum Touch™ software.

Achieve easy workflow control and results in a shorter time through:

- Minimal decision points
- Intuitive one-click operation

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/polymeranalyzer>

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://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.