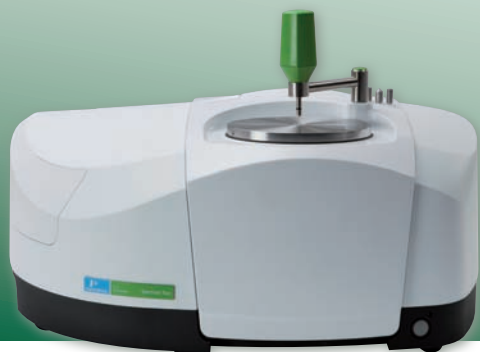


FT-IR Spectroscopy



Spectrum Two for Pharmaceuticals and Regulated Industries

Spectrum Two™ delivers everything required to perform IR analyses within regulated environments with complete confidence. From rapid raw material identification through to sensitive analysis of formulated products, Spectrum Two offers the best user experience without any compromise in analytical performance.

PerkinElmer works with you to assure absolute compliance of your IR testing, every time. Our in-depth knowledge of the pharmaceutical market has been distilled into a dedicated Pharmacopoeia Compliance Resource Kit compatible with the seven major global pharmacopoeias. As well as removing the need to source detailed pharmacopoeia testing information, it allows you to focus on the implementation of your regulated IR analyses.

Choose from a range of Pharmacopoeia Compliance Resource Kits:

USP: United States Pharmacopoeia

EP: European Pharmacopoeia

IP: Indian Pharmacopoeia

JP: Japanese Pharmacopoeia

CP: Chinese Pharmacopoeia

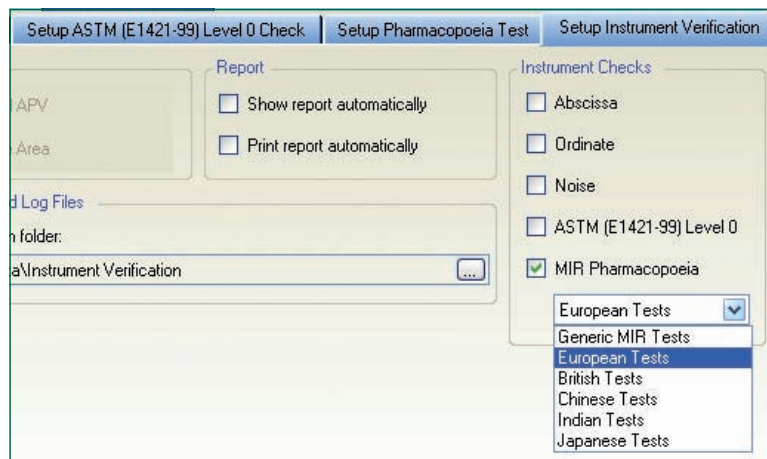
IntPhr: International Pharmacopoeia

Stress-free IR System Commissioning

With Spectrum Two, implementing an IR testing solution in all pharmacopoeial environments is made easy. State-of-the art accessories, configurable software and our Pharmacopoeia Compliance Resource Kit make learning and

setting up compliant testing protocols faster than ever before. Meanwhile, a range of instrument-specific qualification services ensure maximum uptime and productivity.

To facilitate SOP writing, the resource pack includes detailed operating instructions and troubleshooting guides covering IR sampling, testing and materials identification. Straightforward documentation enables operators to understand 21 CFR part 11 technical compliance and reduce the validation burden. For the set-up and implementation of training programs, you can access PerkinElmer's knowledge and experience via the Educational Resource and Introductory Kits also available.



On-demand facility for user specific pharmacopoeial testing.

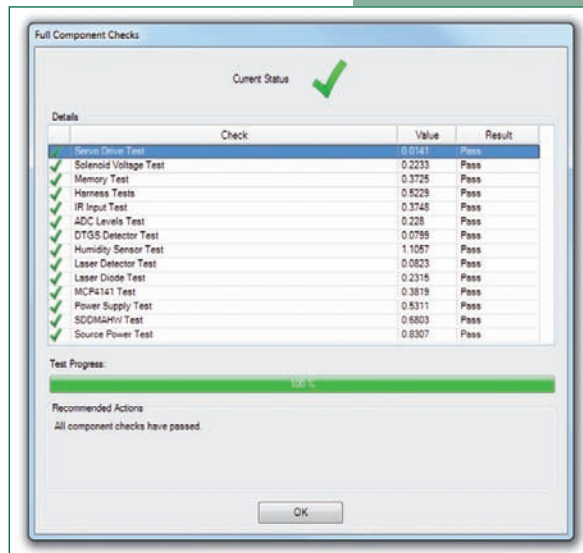
Trusted Results

Whether you are in research, product development, or excipient, API or final formulated product testing, compliance is mandatory and quality essential. Spectrum Two's patented AVI and AVC technologies use high resolution gas-phase spectra to reduce spectral interferences and inconsistencies, without slowing your day-to-day operations. Additionally, a multitude of automated system checks assure reliable answers, every time.

Driven by Spectrum Enhanced Security (ES) software, the Spectrum Two has 21 CFR part 11 technical compliance built in from the ground up. With extensive and configurable access controls, audit trailing, system validations tests, reporting and electronic signatures, you can be secure without delay.

Multi-Level System Validation Tests

- Instrument component checks certify the instrument module is operating correctly.
- Automated instrument performance testing ensures regulatory IR criteria are met.



Automated component checks help to ensure regulatory compliance.

- Contamination checks test for sample carry over.
- Data quality assessments monitor for sampling or operator error.
- Result evaluation limits guarantee unequivocal answers.

IR Confidence for Everybody

Thanks to its unique design features, Spectrum Two can be used by everybody, no matter what their level of experience. Spectrum software is configurable, enabling protocols to be started at the click of a tool-bar button, and operators are guided to the result with a clear and thorough system-user dialog.

If a problem does occur, illustrated troubleshooting tips provide advice on common sampling or operator errors.

Measurements in Non-Laboratory Environments

Spectrum Two is designed with a number of features to bring laboratory performance to non-laboratory situations. Wherever you are, you can rely on Spectrum Two to deliver the compliant results you need – fast.

- OpticsGuard™ technology provides a unique humidity shield to protect vital optical components.
- The battery-powered option ensures protection against power interruptions and enables measurements in the absence of access to mains power.
- Its compact size allows movement between measurement points.
- Wireless connectivity enables operation in environments without the inconvenience of network cabling.

A Commitment to Service

PerkinElmer offers a range of support services designed to help you get the best from Spectrum Two and drive the productivity of your IR testing processes.

With a global team of certified, factory-trained customer support engineers in more than 150 countries around the world, OneSource Laboratory Services from PerkinElmer lives up to its name by being the single source for all your instrument maintenance and repair needs. Flexible and scalable, OneSource programs are tailored to the specific needs and goals of individual customers. From complete, all-inclusive long-term contracts to an à-la-carte offering of individual services, we have the people, experience and resources to deliver the ideal solution to labs of all types and sizes.

OneSource offers the most comprehensive portfolio of professional laboratory services in the industry, including complete care programs for virtually every technology and manufacturer. By allowing you to consolidate all your service contracts under a single supplier, and by providing responsive, expert technical advice and support at a moment's notice, we ensure your instrumentation – and your lab – is running at optimum levels at all times. Whether it's care and repair, validation and compliance, asset management and laboratory relocation, software and hardware upgrades, or education and training, OneSource is the one you can count on.



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

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.

009524_07

Printed in USA