



PRODUCT NOTE

FT-IR and NIR Spectroscopy

Spectrum Quant Quantitative Analysis Software

Unleash the potential of your PerkinElmer FT-IR and NIR spectrometers with Spectrum® Quant software. Designed for novice and advanced spectroscopists who need to integrate accurate quantitative analysis into their IR or NIR analysis. Spectrum Quant combines very simple method development with a range of

data processing options to ensure optimum quantitative analysis that can be achieved quickly and reliably. Spectrum Quant is designed by spectroscopists for spectroscopists. Like the highly popular Spectrum 10 Spectroscopy software for the PerkinElmer spectrometers, Spectrum Quant manages your data to enable you to easily work with multiple spectra, methods and results to maximize your productivity. The common design approach with Spectrum 10 enables quantitative methods to be easily incorporated into Spectrum 10, macros and Spectrum Touch methods, to provide simple turnkey analyzer operation with PerkinElmer IR and NIR systems.

Smart, simple quantitative analysis

Spectrum Quant is ideal for laboratories that need to deliver results efficiently, without the need for extensive software training or methods development. If you use Spectrum 10 to collect your IR or NIR data, Spectrum Quant requires virtually no additional training to get up and running (Figure 1). Method development can be fast-tracked thanks to the close linkage between the two packages. For example, when calibration spectra are collected using Spectrum 10, concentrations can be entered with sample information and all data automatically incorporated into a quantitative method using the smart method

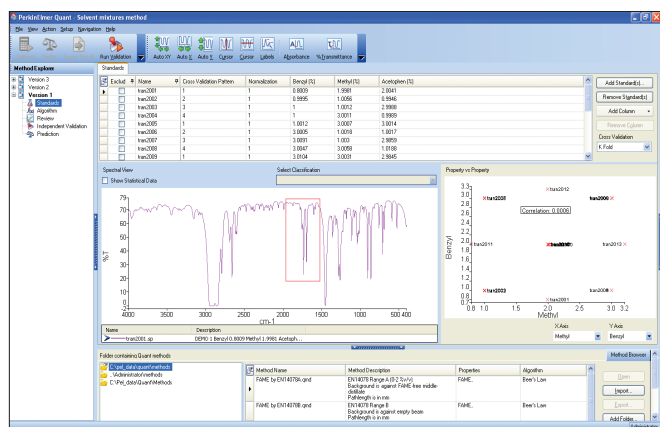


Figure 1. Simple layouts allow you to manage Quant methods, data and results.

Wizard (Figure 2) or custom data fields settings in Spectrum 10 – minimizing the possibilities of transcription error. As the method is developed and refined using the powerful interactive Quant tools, the project is managed with the Method Explorer feature of the software. Method versions allow changes to be tracked and stored enabling you to converge on optimum choice of method parameters as quickly as possible. With Spectrum Quant you can get methods implemented quickly and easily.

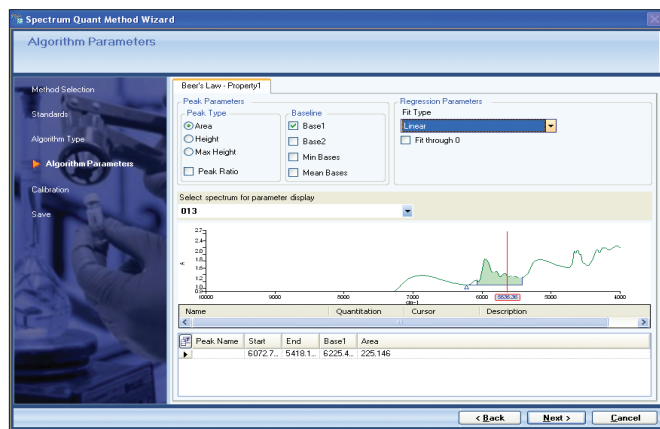


Figure 2. Quant method Wizard for simple method building from Spectrum 10.

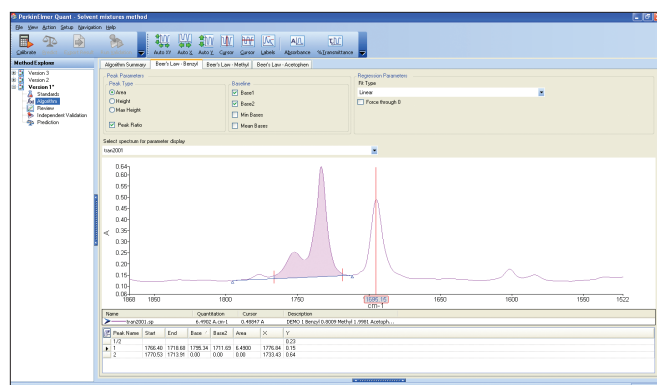


Figure 3. Interactive parameter setup from spectral displays.

Accelerated method optimization

Using Spectrum Quant means you can get the best results out of your valuable IR and NIR data. A range of quantitative analysis and data visualization options are available in a modular structure – from simple peak height/area analysis using the Beers Law module to powerful full spectrum multivariate methods such as partial least squares (PLS) or principal components regression (PCR) (Table 1).

Table 1. Spectrum 10 Quantitative Analysis modules.

| Quantitative Technique | Software Module |
|--|---|
| Univariate analysis including peak height, area regression methods | Spectrum 10 with Beers Law Quant (LX108873) |
| Multivariate analysis with partial least squares and principal components regression (PLS1 and PCR+) with model diagnostics and validation | Spectrum Quant Algorithm Pack L1101024 |

A suite of preprocessing algorithms are included to assist with common sampling variability issues and improve method robustness. Once calibration spectra are defined, the powerful graphical interaction allows calculation ranges (Figure 3), baselines and other parameters to be setup graphically using the spectra and cursors, and scatter plots can quickly reveal potential problems with spectral or concentration data in the calibration.

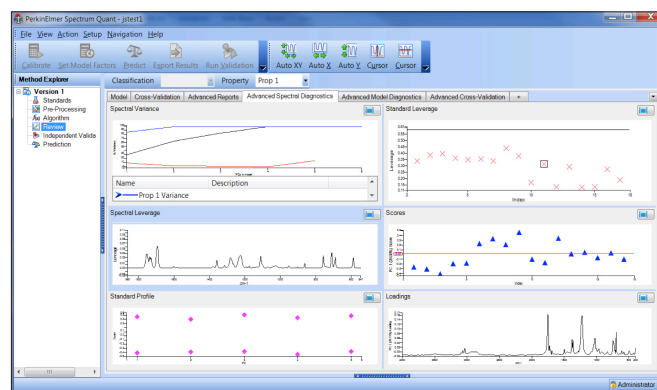


Figure 4. Review plots for easier evaluation of statistical data.

Put your Quant methods into action

Medida de la muestra

◀

Muestra 1 de 1

▶

effluent sample

Editar



Limpie la cubeta con la muestra dos veces. Llene la cubeta con la muestra.

Coloque la cubeta en el soporte para-cubetas del portamuestras.

Presione "seguir" para empezar el escaneo de la muestra.

▶

Barrido

Macro Settings

Name: Macro 1

Description:

Group Name: Macros

Process

1 Build a Mark

2 Kerning

3 Diacritics

4 Search and Compare

5 Quantization

6 Equalization

7 Resampling and Output

Name: Equation_1, Description: Formulas Area (M1, 1750, 1700). Use for next step table. 1. Generate input speech table.

If condition: Equation_1

Operator: >

Criterion: 3

Criterion 2:

Then conditional formatting:

Add text to result: [Red]

Apply to text: [X]

Else conditional formatting:

No formatting: []

Help OK Cancel

Using the unique Spectrum Touch interface allows simple touch/tap navigation through your analysis rather than the traditional point and click methods, to fully exploit modern touchscreen PC functionality (Figure 6). Performing multi-component analysis, even simultaneously on multiple spectra in Spectrum 10 is a single-click operation and exploring the results tables couldn't be easier with the sophisticated tabular/graphics display in Spectrum 10 and Quant. You can also specify control limits on calculated concentrations so that out-of-specification results can be visibly highlighted in results tables (Figure 7). With Spectrum Quant method integration you can concentrate on the results rather than the tasks.

With a familiar Microsoft® Outlook®-style interface, Spectrum Quant allows you to adjust your views to suit your preferences. Customizable Quant action and graphics toolbars can show your most-used tasks – while methods information, setups and navigators can be positioned in panes side by side – or pushed aside to avoid clutter as you would a normal work area.

You can send selected elements of the methods or results to other applications such as Microsoft® Word® for incorporating into your reports (Figure 8), or use the built-in Report Editor to customize Quant reports – or transfer results tables into third party spreadsheet and trending packages for specific views. Access control into the software is also provided via a separate customizable login screen.

Local language selection and Microsoft® support

Spectrum Quant is compatible with Windows® 7/8 with 64 bit support. Local language support is automatically selected from the Windows® locale setting, including English, Spanish, Chinese, Japanese, German, French and Portuguese versions.

Microsoft Gold Certified Partner

The Spectrum Quant software development site is a Microsoft® Gold certified partner. Spectrum team members have demonstrated the highest levels of technical excellence, marketplace impact and satisfaction of customers using Microsoft® products and services.

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 ©2012-2014, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

010556A_01