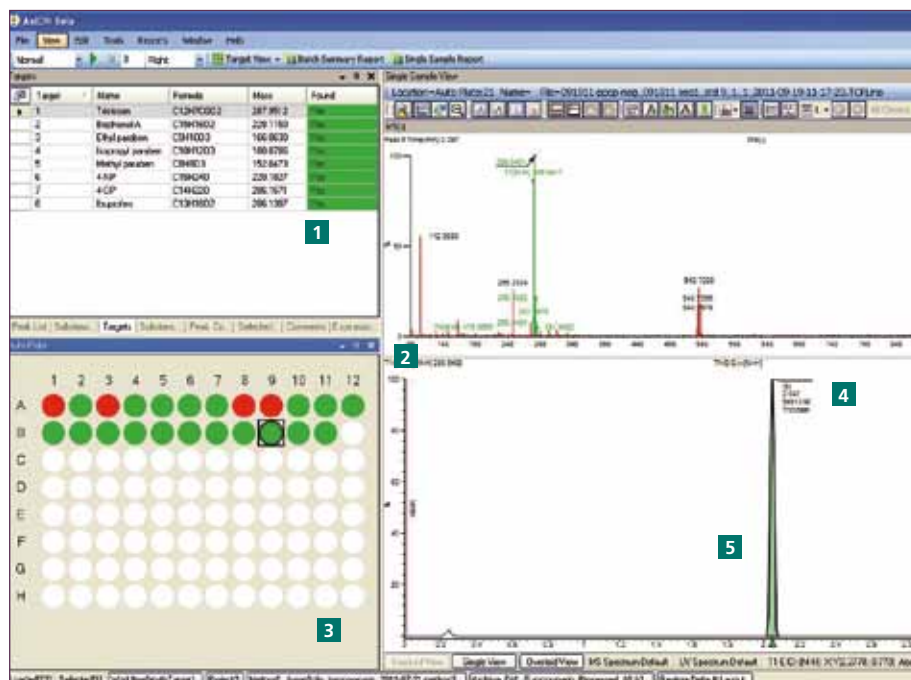


The Fast, Clear Way To Find What You're Looking For.

Designed for optimal flexibility and productivity, AxION® Solo™ streamlines the most complex LC/MS analyses by allowing you to define your own criteria and processing methods to find one or many substances by mass or formula. An intensity-driven heat mapping and graphical user interface provides superior visualization of results for quick interpretation and reporting. Whether you're searching for single or multiple targets, AxION Solo allows you to take the fast, easy way to clear, confident analytical results.

Multi-Target Analysis

- 1 Visual summary of results generates rapid decision-making.
 - Found/not found substances in all samples are displayed.
 - Quick access to unambiguous results.
- 2 Detailed data minimizes review time.
 - Supports chromatograms and spectra in the same window.
 - Outliers and anomalies quickly identified.
- 3 Heat map provides fast interpretation.
 - Signal thresholds can quickly be adjusted for each substance.
 - Multiple plates or lists of samples can be viewed at once.
- 4 Automatic peak annotation provides greater specificity.
 - Annotated substances on peaks—adducts and dimers clearly identified.
 - Substances are easily distinguished with peak labeling.
- 5 EICs displayed for greater sample insight.
 - Adjustable parameters for EIC abstraction from TIC.
 - One click to display EICs for all substances.



Speed Up And Simplify Your Multi-Target Analyses.

- Detect substances on your list of targets.
- Quickly locate a specific substance in all samples.
- Find and detect an additional substance after processing the data.

Single-Target Analysis

1 Results ranking speeds data analysis.

- Rank order rules such as purity, or found and pure.
- Multiple plates or lists can be viewed at the same time.

2 Fast found/not found/maybe found for rapid review of samples.

- Sample scoring visually shows which one has your substance.
- Visually coded to quickly understand your results.

3 Unambiguous target identification for confident results.

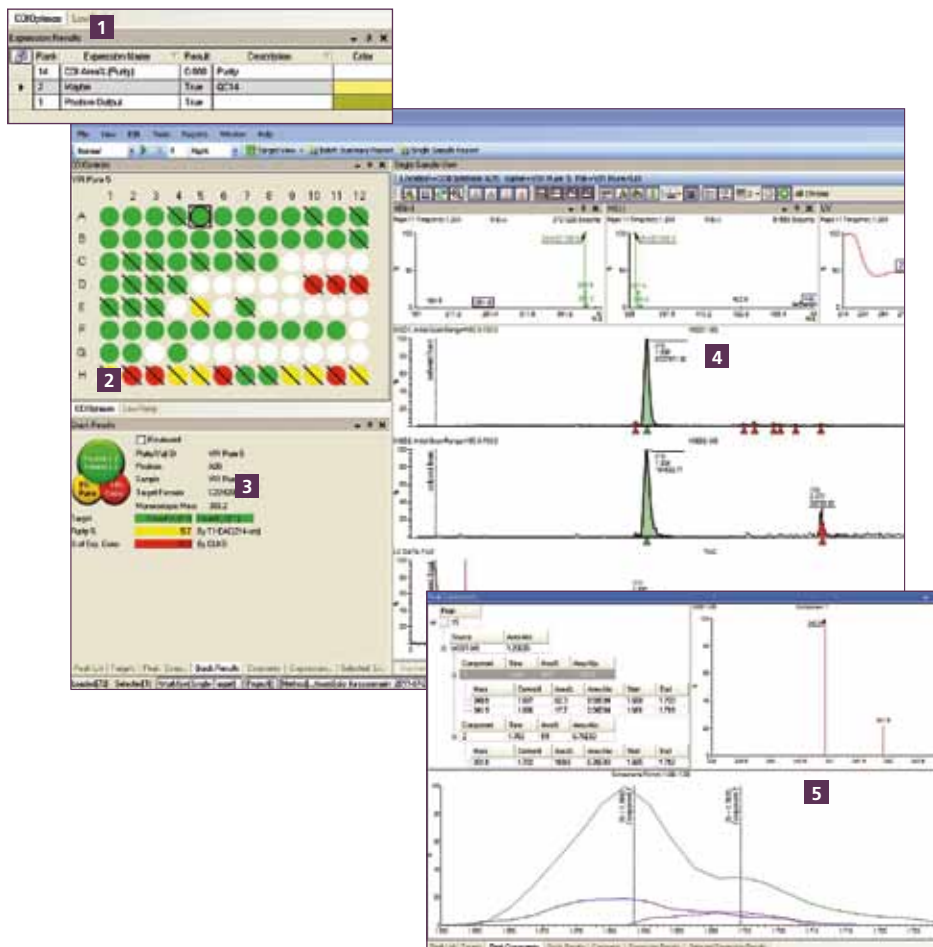
- User-defined rules such as isotope ratio, accurate mass, peak purity and peak shape.
- Multiple rules applied to substance identification at the same time.

4 Peak annotations provide specificity.

- Automatic peak annotation for substances, dimers, and adducts.
- Compare extracted peaks from each instrument mode in the same window.

5 Complete peak analysis to obtain comprehensive results.

- Easily view detailed homogeneity analysis.
- Detailed componentization for your substance.



Speed Up And Simplify Your Single-Target Analyses.

- Detect your target substance.
- Make result-driven decisions.
- Detect additional compounds.
- Find the ratio of other compounds to your target substance.
- Detect which sample contains the most of your target substance.

AxION Xpo™—An Intuitive Data Review Solution For Target Analysis.

Provides fast, powerful reviewing and reporting functionality for large quantities of LC/MS data.

Compound-centric data viewing environment offers easy access to results and allows customization to suit individual workflows.

Results are easily captured and shared using a broad array of applications.

Reports can be created and printed in customizable Microsoft® Word and Adobe® Acrobat® formats.

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

010118_01

Printed in USA