

CDS Solutions

Key Features:

- Easy-to-use software platform that accommodates all user needs and skill levels
- Workflow-driven interface automates both routine and complex tasks
- Provides the functionality required by modern chromatographers, such as ApexTrack™ Integration algorithm, custom calculations, and custom reports

Empower® 3 Chromatography Software: QuickStart™ Interface

Easily Harness the Powerful Capabilities of Empower® 3

The Waters® Empower® 3
QuickStart™ interface makes

it easy for users to conduct targeted tasks and workflows while harnessing the many capabilities of Empower® 3 Chromatography Software. The powerful QuickStart™ interface presents the appropriate functionality to the user – in a single window – only when they need it. For added flexibility, users may customize their own user interfaces to suit individual needs and make their work even more efficient.

A Single-Window to a Wide Range of Analytical Instruments

Through the QuickStart™ user-centric, task-oriented interface, access all the chromatographic capabilities you need as well as a wide range of analytical chromatographic and detection techniques:

- PerkinElmer Altus™ LC
- PerkinElmer Clarus® and TurboMatrix™
- Waters® HPLC, GPC, APC™, ACQUITY UPLC®, ACQUITY UPC2®, and integrated mass spectrometry systems
- Gas Chromatography (GC), Capillary Electrophoresis (CE), and Ion Chromatography (IC)
- Multi-vendor chromatographic instruments, including Agilent Technologies®, Beckman Coulter®, CTC Analytics, DANI Instruments®, Hitachi High Technologies, Shimadzu®, Thermo Scientific™/Dionex™, and others

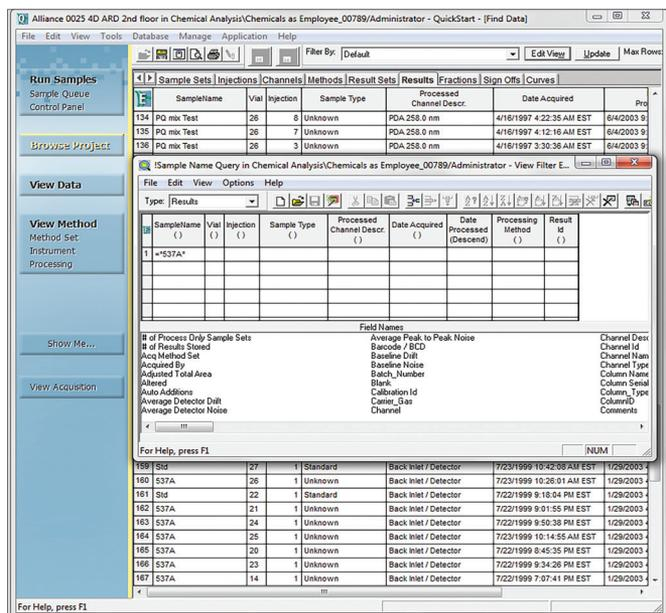


Figure 1. Quickly and easily find data using Empower® 3 View Filters that can display only what is needed and can be saved for later use.

Help Where You Need It

Setting up new acquisitions and processing chromatographic results can sometimes be challenging – especially when the task is unfamiliar to the user. The Empower® 3 QuickStart™ interface provides access to the Show Me... Help, a guided workflow wizard, to help users complete their tasks quicker and more accurately.

Data Traceability and Integrity

Empower® 3 automatically manages data through several levels of safeguards to ensure trustworthy and reliable records. Laboratories in the chemical materials, food and beverage, and environmental industries require trustworthy and reliable electronic records to protect the data integrity/authenticity of results in research and to aid troubleshooting in manufacturing. Empower® 3 QuickStart™ makes it easy to review and approve results; search for data; and verify records for authenticity.

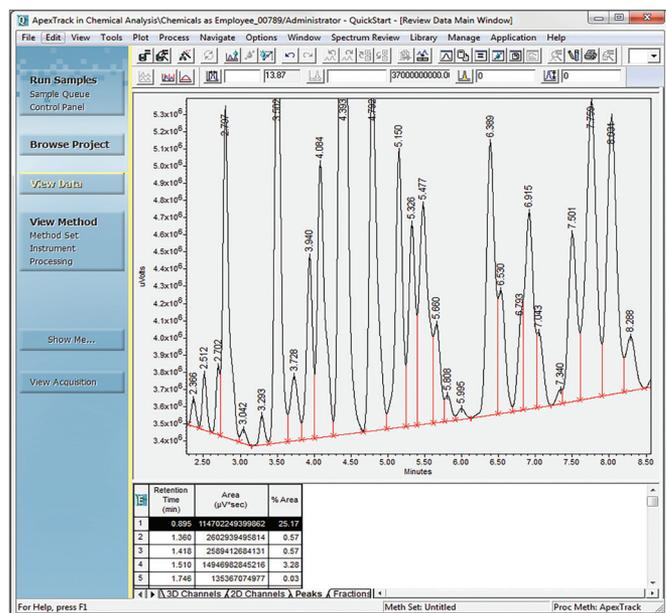


Figure 2. Focus on areas of interest by taking advantage of view settings that can be set as defaults for each type of experiment to ensure proper peak integration (automated using ApexTrack™), and quantitated values are within expected limits.

Comprehensive Data Processing and Reporting

Accurate and consistent chromatographic results are important to universities and corporations alike. Empower® 3 QuickStart™ streamlines the collection, processing, reviewing, and reporting of chromatographic results. Capabilities accessed through the QuickStart™ interface include:

- Instrument acquisition
- Peak detection (including the ApexTrack™ Integration algorithm for consistent results)
- Quantitation
- GPC processing and calculations
- Data review
- Searching for data
- Mass spectrometry
- Custom reporting

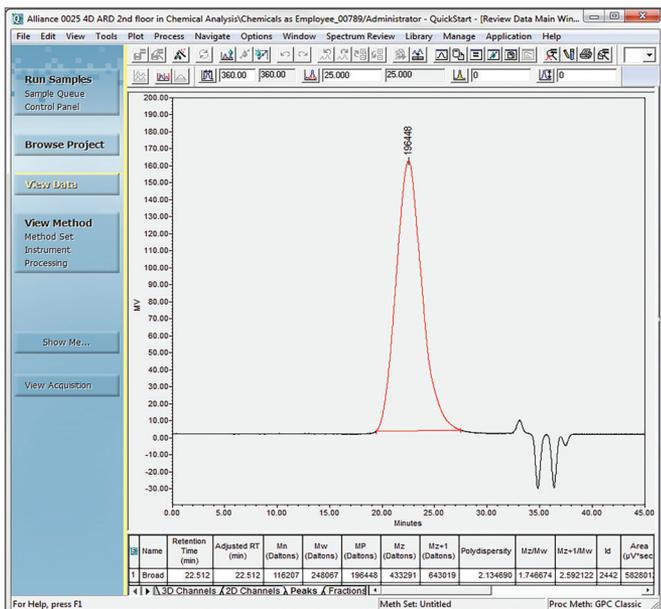


Figure 3. Have fast access to all areas of the software using the QuickStart™ interface, including GPC calculations.

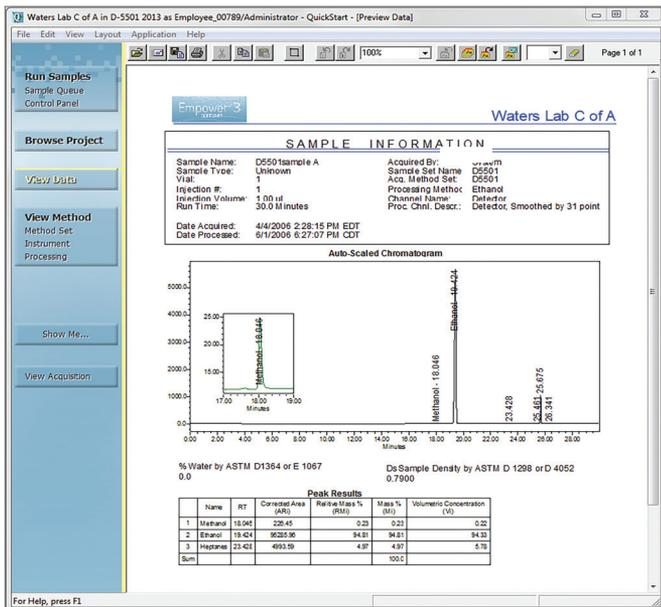


Figure 4. Empower® 3's customized reporting capabilities can show exactly what you need in the format you want with easy creation of templates using a drag-and-drop approach.

Benefits of Using the Empower® Quickstart

The QuickStart™ interface provides users with an easy-to-use, single-window software interface for acquiring, processing, and reporting chromatographic results. Benefits of using Empower® 3 QuickStart™ include:

- Users presented only with the functionality they need – greatly improving speed and accuracy of laboratory results
- A reduction in the time and cost associated with training
- Harmonized control of and streamlined access to a wide variety of analytical chromatographic and detection techniques
- Easily confirm trustworthiness and reliability of chromatographic records
- Show Me... Help leads users through unfamiliar tasks thereby boosting productivity

Comprehensive Support for All of Your Chromatography Needs

PerkinElmer has more than 60 years of knowledge and experience in chromatography. Our team of specialists serve dynamic and diverse industries and understand your application needs. We offer the most comprehensive portfolio of professional services for virtually every technology and manufacturer. More importantly, we are responsive and help ensure your lab is running optimally, at all times. Whether you need validation and compliance, software and hardware upgrades or education and training and more, PerkinElmer can empower your lab.