

Liquid Chromatography/ Mass Spectrometry



Key Features:

- Modular 3Q software with guided workflow for easy operation
- Ideal for acquiring, quantifying and reporting with flexible data viewing options
- Powerful remote diagnostics for quickly and accurately identifying system concerns

QSight Simplicity 3Q Software Suite

Software and Services to Make Lab Operations Simpler

When it comes to software, most labs need a solution that is easy-to-use and provides a streamlined workflow, from method development to results. With the QSight™ Triple Quad LC/MS/MS, that's precisely what you get. Our modular Simplicity 3Q™ software is ideal for acquiring, quantifying and reporting, has flexible data viewing options and powerful remote diagnostics. It's designed for intuitive, straightforward operation so even novice technicians find it easy-to-learn and use.

Modularity Makes It Easy to Run Samples

Our Simplicity 3Q software suite includes a collection of modules that guide users through a workflow-based experience from method development through to results processing and reporting. The suite includes Simplicity 3Q Acquire™ for data acquisition, Simplicity 3Q Quant™ for quantitation, Simplicity 3Q View™ for data viewing. Simplicity 3Q Service™ enables remote diagnostics for quick and accurate troubleshooting.

Simplicity 3Q Acquire

This data acquisition software streamlines acquisition method development and batch acquisition, with the highest levels of automation. Simplicity 3Q Acquire™ software is capable of real-time data acquisitions of more than one thousand MRMs and provides a real-time simultaneous display of large numbers of transitions. It also includes an easy concentration table editor and control options for batch acquisition.

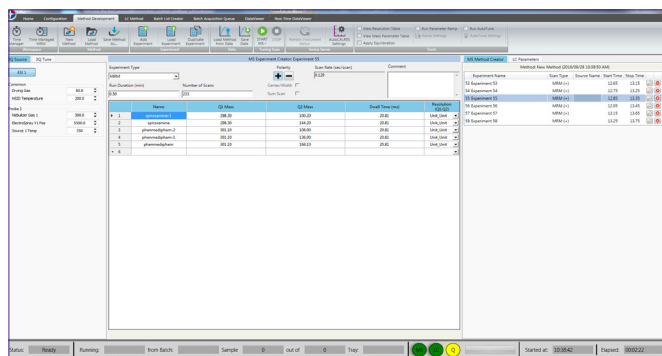


Figure 1. Simplicity 3Q Acquire enables time managed MRM method development for analyzing hundreds of pesticides in single method.

Simplicity 3Q Quant

For data analysis and reporting, Simplicity 3Q Quant™ software uses autopopulation from data acquisition files for an exceptionally fast, smooth, streamlined workflow for high throughput situations. With its full-on view of all navigation data, the software's RapidView Heads-up Display™ allows you to concentrate on your data, and your results. Multiple options are available for data processing, such as smoothing algorithms, which can be updated very easily. Adding to its flexibility, this module can be easily set to simplified views for labs that only need specific output results.

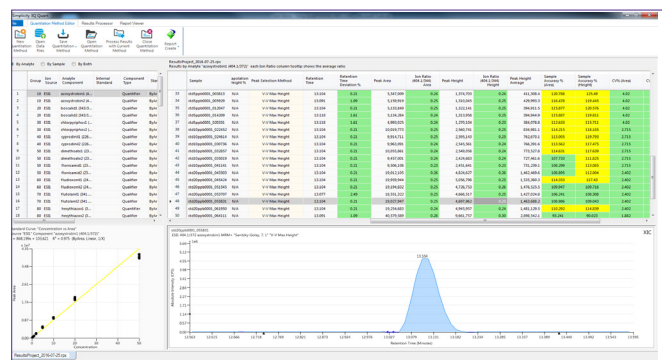


Figure 2. Simplicity 3Q Quant's intuitive result table display focuses on results needed for a quantitation workflow.

Simplicity 3Q View

Simplicity 3Q View™ provides flexible multi-analyte viewing, overlay plotting, and signal-to-noise analysis, for post-processing review and dynamic real-time display. It allows viewing of total ion chromatograms, extracted ion chromatograms, mass spectra, and various types of post-processing data results. Some of the available functions include, S/N calculations, 3D contour diagram, spectra/chromatograms overlay and background subtraction.

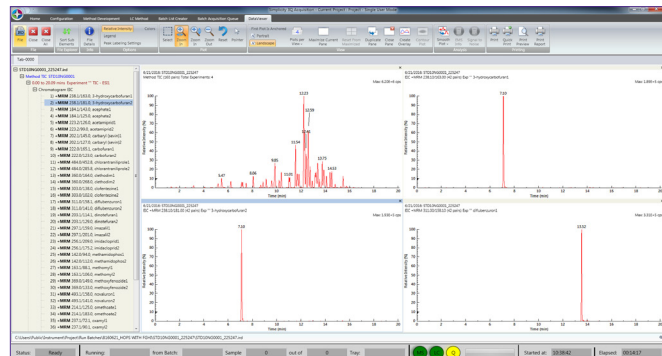


Figure 3. Simplicity 3Q View displays and processes live or post-acquired data.

Simplicity 3Q View also allows the user to view up to 16 plots per page, which allows multiple sets of data to be opened in the same window tab and/or separate window. It also displays detailed file information specifications, which includes LC/MS method parameters and instrument information.

Simplicity 3Q Service

Troubleshoot Issues More Effectively and Efficiently

Simplicity 3Q Service™ was designed with focus on providing maximum uptime. This module uses the instrument's vast number of controls and read backs to provide accurate and timely diagnostics. At the core of Simplicity 3Q service is the AdvIO™ Electronics system, which can minimize downtime by quickly identifying or ruling out hardware concerns.

The AdvIO Electronics system enables a straightforward, powerful review and feedback for virtually every QSight instrument parameter, including voltage, temperature, pressure, and gas flow. The result: rapid diagnosis and remediation to enable greater instrument uptime.

Intuitive, Real-time Readback Facilitates Accurate Diagnostics and Resolution

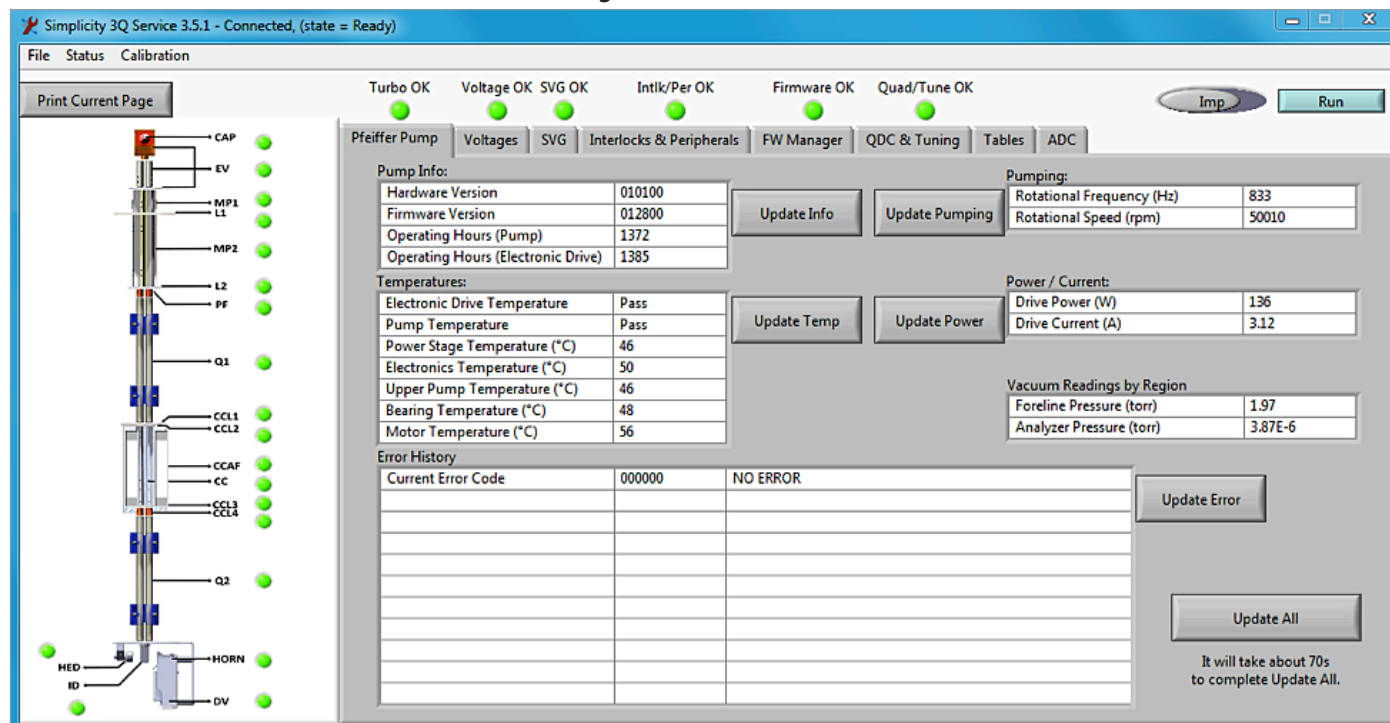


Figure 4. Each tab within the software has a specific focus on the hardware within the QSight instrument. Above each tab (after hitting the “Run” command), there will be a status LED for the tab. Status LEDs that are green, indicate that all relevant status readbacks within the tab have been satisfied. A red LED above the tab indicates that one or more indicators within the tab have not been satisfied.

Software and Systems That Are Simplicity Itself

Our QSight Triple Quadrupole LC/MS/MS with its guided workflow Simplicity 3Q software is a complete solution that delivers the functionality you need for the most efficient high-sensitivity analyses.

For more information about the QSight Triple Quadrupole LC/MS/MS, go to www.perkinelmer.com/QSight

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 ©2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.