

1311 Orleans Drive
Sunnyvale, CA
94089-1136
Ph: +1-800-635-5577
Fx: +1-408-548-6461

In-Depth AcuityXpress™ Software Training

May 12-13 at Molecular Devices in Sunnyvale, CA
July 21-22 at Molecular Devices in Downingtown, PA

This advanced course provides students with an introductory/intermediate level exposure to many of the features in AcuityXpress Software. Students will be provided with a computer workstation for use during the class and will be expected to work through example data sets, with the instructor and teaching assistants, which clearly illustrate the topics being covered. These hands-on examples will be representative of actual customer data and will progress from the basic to the intermediate level illustrating the key concepts required for the student to work independently to analyze HCS data in their labs. Example data sets, presentations and supplementary materials used during the training will be provided to the student upon completion.

Condensed Agenda:

Day 1

- Introduction to AcuityXpress Software
 - Overview
 - Demo dataset
 - Supporting Materials
- The AcuityXpress Software Interface
 - Data table
 - Datasets folder
 - Common tasks
 - Analysis visualizations
 - MDCEarth configuration
- Multi-Parametric Analysis
 - Simple Match
 - PCA and SOM
 - Selection and reverse selection
- Creating a Dataset - Screening Example
 - Plate annotation
 - Plate normalization
 - Dataset query
 - Data table configuration

Day 2

- Data Analysis - Screening Example
 - Applying multiparametric tools (SOM, PCA) to screening
 - Creating and working with quick lists
- Preparing a Dataset - Dose Response Example
 - Plate annotation
 - Plate normalization
 - Dataset query
 - Data table configuration
- Data Analysis - Dose Response Example
 - Curve fitting
 - Using PCA and SOM with dose-response data
- Additional Topics
 - Reports
 - Web Links
 - Importing and exporting
 - Q&A