

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

Advanced MetaXpress® Software Training

April 6-8 and Oct 5-7 at Molecular Devices in Downingtown, PA
April 27-29 and Nov 9-11 at Molecular Devices in Sunnyvale, CA

This advanced course provides students with the basic tools to design and write their own custom analysis “Journals” for analyzing plate based image sets collected in MetaXpress Software. Students will be provided with a computer workstation for use during the class and will be expected to work through example journals, 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 requirements for the student to independently develop more advanced journals in their labs. Example images, journals, presentations and supplementary materials used during the training will be provided to the student upon completion.

Condensed Agenda:

Day 1 (optional) half day*

- Image Processing Fundamentals
 - Image scaling and display options
 - Image stacks and montages
 - Segmentation and thresholding
 - Filters and masks
- Measurement Fundamentals
 - Regions
 - Intensity, distance, area measurements
 - Integrated Morphometry Analysis (IMA)
 - Data logging
- Review Plate Data and Application Module Review

Day 2

- Introduction to Journaling
- Working with Journals
 - Basic journals
 - Journal editor
 - Variables
- Acquisition Journals
 - Device control
 - Program variable
 - Case studies: Image overlay display and Z stack acquisition

Day 3

- Analysis Journals - Basic Strategies
 - Working with binary masks
 - Data logging in journals
- Analysis Journals - Case Studies
- Additional Topics
 - Taskbars
 - Looping journals
 - User interaction and setup journals

*Day 1 (optional) will cover basics of image processing and image analysis and is recommended for anyone with limited experience using MetaXpress Software or MetaMorph® Software, and for those who have only used the application modules for image analysis. Day 1 or equivalent experience is a pre-requisite for the remainder of the journaling course.