

[NEW] Webinar Spotlight

Empower electrophysiology data analysis with the new Clampfit 11 Advanced Event Detection



October 2, 2019 | 9:00 AM PDT

The Clampfit software module is a useful tool to manipulate and analyze electrophysiological data acquired by pCLAMP™ software. Recently, Molecular Devices released Clampfit 11 software with Advanced Event Detection functions to automate data analysis and improve accurate measurement. In this webinar, Dr. Jin Yan will provide an overview of new event detection features in Clampfit 11. She will also demonstrate the use of Template search, Population Spike search, and Action Potential search.

- Wednesday, October 2, 2019
- 9:00 AM PDT

[Register for Webinar ▶](#)

[NEW] Product Spotlight



Water immersion objectives for ImageExpress Micro Confocal system

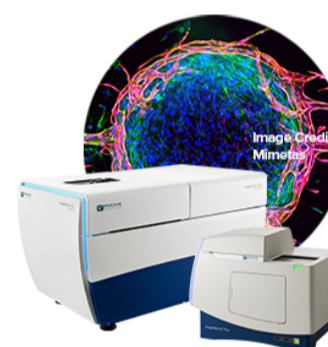
Water immersion objectives can offer signal increases of up to 4X which can help scientists see deeper into 3D and thick tissue samples at lower exposure times. Learn more about our newest high-performance customization.

[Learn More ▶](#)

[NEW] User Meeting Spotlight

High-Content Imaging

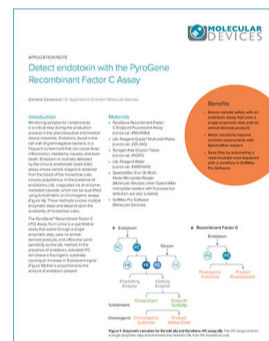
Advances in 3D imaging allow researchers to gain more insights than ever before. Join us to network with colleagues and hear key academic and industry leaders share their research and workflow solutions.



- Wednesday, September 11, 2019
- 9:30 AM – 3:00 PM PDT
- Farmer & The Seahorse – The Alexandria at Torrey Pines
- San Diego, California

[Register for User Meeting ▶](#)

[NEW] Application Spotlight



Detect endotoxin with the PyroGene Recombinant Factor C Assay

Monitoring samples for contaminants is a critical step during the production process in the pharmaceutical and medical device industries. Endotoxin, found in the cell wall of gram-negative bacteria, is a frequent contaminant that can cause fever, inflammation, headache, nausea, and even death.

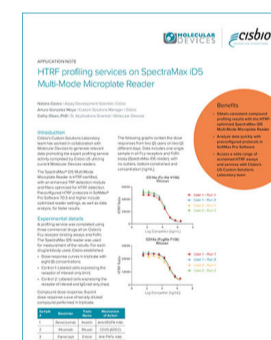
Assure sample safety with the PyroGene™ Recombinant Factor C (rFC) Assay that uses a single enzymatic step and no animal-derived products. Plus, attain sensitivity beyond minimal requirements with the SpectraMax® readers.

[Download Application Note ▶](#)

[NEW] Application Spotlight

HTRF profiling services on SpectraMax iD5 reader

Due to the collaboration with Molecular Devices and the SpectraMax® iD5 Multi-Mode Microplate Reader, Cisbio can now offer custom HTRF profiling services as a reliable, fast and reproducible approach to drug discovery using technology suitable for antibody affinity and the potential of biosimilarity studies using relevant drugs available on the market.



The SpectraMax iD5 reader is HTRF-certified, with an enhanced TRF detection module and filters optimized for HTRF detection. Preconfigured HTRF protocols in SoftMax® Pro Software include optimized reader settings, as well as data analysis, for faster results.

[Download Application Note ▶](#)

Events

Organ-on-a-Chip Workshop (MIMETAS)
September 5-6, 2019
Reading, UK

SynBioBeta
October 1-3, 2019
San Francisco, CA USA

High-Content Imaging User Meeting
September 11, 2019
San Diego, CA USA

Advanced Analytical Technologies for Proteins
October 2-3, 2019
Romainville, France

LabDays Aarhus
September 11-12, 2019
Aarhus, Denmark

Stockholm LabDays
October 2-3, 2019
Stockholm, Sweden

Organ-on-a-chip Workshop (MIMETAS)
September 12-13, 2019
San Jose, CA USA

Neuroscience
Booth #1216
October 19-23, 2019
Chicago, Illinois USA

SBI2
Booth #9
September 17-19, 2019
Boston, MA USA

Cell Culture & Bioprocessing Congress
October 29-30, 2019
London, UK

Ilmac
Booth #E152
September 24-27, 2019
Basel, Switzerland

CRI-CIMT-EATI-AACR International Cancer Immunotherapy
Booth #16
September 25-28, 2019
Paris, France

Follow Molecular Devices:

