



[NEW] Webinar Spotlight

Optimizing high-content screening tools for physiologically-relevant 3D *in vitro* models

Join us for this webinar with HCS Pharma, a biotechnology start-up company based in France, as we discuss the biological characterization of cells in 3D matrices for liver, brain and adipose tissues, and cancer modeling. We'll also discuss high-content screening solutions for examining 3D *in vitro* models.

- Tuesday, October 29, 2019
- 7:00 AM PDT, 2:00 PM GMT, 3:00 PM CET

[Register for Webinar ▶](#)

[NEW] User Meeting Spotlight

Axon pCLAMP 11 and Imaging User Meeting at Society for Neuroscience (SFN)

Join us at SFN in Chicago for the Axon™ pCLAMP™ 11 and Imaging User Meeting.



Speaker: Jeffrey Tang, PhD, Molecular Devices

Empower your electrophysiology studies using new Axon pCLAMP 11 software and HumSilencer adaptive noise cancellation technology

Speaker: Aaron Risinger, M.S., MBA, Molecular Devices

New Advances in Cellular Imaging using Multi-Parametric Data Acquisition and Analysis Software with the ImageXpress® Pico Automated Cell Imaging System

- Monday, October 21, 2019
- 6:30 PM – 8:00 PM CDT
- Hyatt Regency McCormick Place

[Register now ▶](#)



[NEW] eBook Spotlight

Streamline beer, wine, and food quality control and safety analyses

Absorbance microplate readers are widely used in research, quality control, and manufacturing processes in the food and beverage, pharmaceutical, and nutraceutical industries. They provide rapid and sensitive measurements of a variety of analytes for a wide range of assays including ELISAs, microbial growth, detection of key compounds and contaminants, and protein quantitation, colors, and flavors.

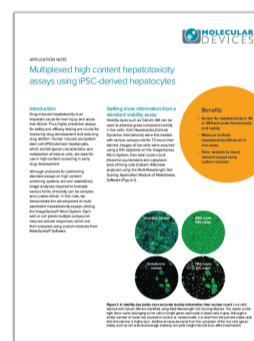
Here, you'll learn how you can use absorbance microplate readers to streamline your beer, wine, and food safety analyses using approved industry-standard methods.

[Download eBook ▶](#)

[NEW] Application Spotlight

Get more high-content tools and tips for cell painting assays

Cell painting assays are a method of using 5+ fluorescent dyes to reveal several responses of interest at once. Each well yields multiple specific and measurable cellular responses, which are then analyzed using preconfigured or custom modules from MetaXpress® software to provide a richer profile than that of multi-parametric assays.



Learn how to optimize high-content imaging capabilities and perform multiplexed assays.

[Learn more ▶](#)



[NEW] Application Spotlight

Monitor multiple stages of apoptosis with live cell kinetic imaging

The study of apoptosis is a critical aspect of drug discovery and development. Additionally, studying the relationship between apoptosis and other factors, such as oxidative stress, is crucial for understanding specific diseases that are associated with the dysregulation of apoptosis.

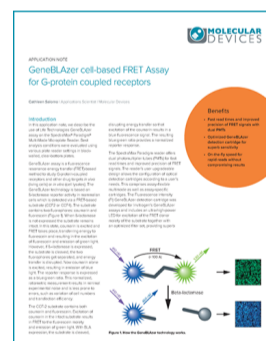
See how we ran two apoptosis assays to study the cytotoxic effects of anti-cancer compounds on HeLa cells. Long-term time-lapse imaging was performed on the ImageXpress Pico system. Live cell conditions were maintained using CellReporterXpress® software's integrated environmental control of CO₂, O₂, temperature, and humidity.

[Learn more ▶](#)

[NEW] Application Spotlight

GeneBLAzer cell-based FRET Assay for G-protein coupled receptors

GeneBLAzer assay is a fluorescence resonance energy transfer (FRET)-based method to study G-protein-coupled receptors and other drug targets *in vivo* (living cells) or *in vitro* (cell lysates).



The SpectraMax® Paradigm® Multi-Mode Microplate Reader offers dual photomultiplier tubes (PMTs) for fast read times and improved precision of FRET signals. Optical detection cartridges can be configured according to a user's needs. This comprises assay-flexible multimode as well as assay-specific cartridges. Combined with the GeneBLAzer Detection Cartridge, the reader is sensitive enough to detect even weak FRET signals, enabling miniaturization of the GeneBLAzer assay to a 1536-well microplate format with no compromise in assay performance.

[Learn more ▶](#)

Events

SynBioBeta

October 1-3, 2019
San Francisco, CA USA

Advanced Analytical Technologies for Proteins

October 2-3, 2019
Romainville, France

Stockholm LabDays

October 2-3, 2019
Stockholm, Sweden

Taiwan High Content User Meeting

5F, Palais de Chine Hotel
October 17, 2019
Taipei, Taiwan

Neuroscience

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

Cell Culture & Bioprocessing Congress

October 29-30, 2019
London, UK

BioProduction at CPHI Worldwide

November 5-7, 2019
Frankfurt, Germany

ISSCR

Booth #8
November 6-8, 2019
Toronto, Canada

American Institute of Chemical Engineers

Booth #159
November 10-15, 2019
Orlando, FL USA

Vaccines R&D

November 18-20, 2019
Newton, MA USA

PEGS Europe

November 18-22, 2019
Lisbon Portugal

Next-Generation Protein Analysis and Detection

December 2-3, 2019
Ghent, Belgium

ASCB | EMBO 2019

Booth #1103
December 7-11, 2019
Washington, DC USA

Antibody Engineering and Therapeutics

Booth #302
December 9-13, 2019
San Diego, CA USA

Follow Molecular Devices:

