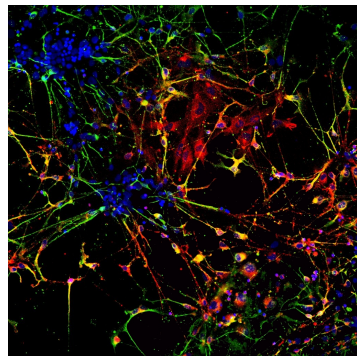


eNewsletter

January 2019

cellular imaging • high-throughput screening • colony selection
microplate detection • analysis

Tutorial Spotlight



NEW Phenotypic screening in 3D culture including ECM: Advantages and challenges

Join us at our tutorial during SLAS 2019 in Washington, D.C. with Nathalie Maubon and Grégory Maubon, PhD from HCS Pharma. During this talk, we will discuss how the cellular microenvironment impacts the proliferation and/or differentiation of cells. A few examples will be presented in oncology, CNS, and metabolic diseases.

Date: Monday, February 4, 2019

Time: 12:30 – 1:15 PM

Room: 143C

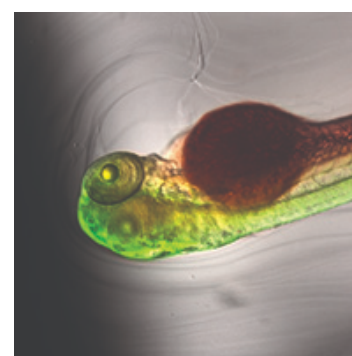
[Register Now](#) ▶

Product Spotlight

NEW Live cell imaging and 3D acquisition using the ImageXpress Pico system

We unveiled last month new features for the ImageXpress® Pico Automated Cell Imaging System including environmental control, z-stack acquisition, LEDs and filters, and double marker expression analysis. These new features enable live cell imaging and 3D acquisition in an easy to use system to better address phenotypic imaging.

[Learn More](#) ▶



eBook Spotlight



NEW Phenotypic screening with iPSC-derived cardiomyocytes and neurons

Here, we use both imaging and calcium oscillation analysis to develop profiles of compounds in iPSC-derived cardiomyocytes such as hERG blockers, β -adrenergic agonists, and environmental toxins. iPSC-derived neuronal cultures were evaluated with neuromodulators as well as environmental toxins. It includes data acquired with the FLIPR Tetra® System, ImageXpress® Micro Confocal system, FlexStation® 3 reader, and more.

[Download eBook](#) ▶

Application Spotlight

NEW Streamline BCA-based protein quantitation

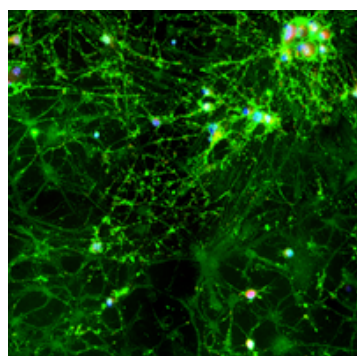
The BCA assay is a two-step colorimetric assay used to quantitate the total protein in a sample. Learn how to quantitate a cellular protein sample with BCA protein assay kits on the SpectraMax® iD5 Multi-Mode Microplate Reader.

- Quantitate proteins with a 5-minute, room-temperature incubation
- Measure assay absorbance values quickly with minimal background noise
- Graph standard curves & calculate sample protein concentrations easily

[Download Application Note](#) ▶



Article Spotlight



NEW Forming 3D Neuronal Models of the Brain

The ability to acquire and analyze in 3D provides a more physiologically relevant model to study. With 3D imaging and analysis, researchers can study physiology and pathophysiology in the 3D context, gain a greater understanding of neurodegenerative disease, and discover potential therapeutics.

[Read More](#) ▶

Citation Spotlight

The **SpectraMax® iD3** microplate reader was used to evaluate optical density as shown in the International Journal of Molecular Medicine.

Citation: Genipin protects against H₂O₂-induced oxidative damage in retinal pigment epithelial cells by promoting Nrf2 signaling

[Read More](#) ▶



Upcoming Events

SLAS

February 2-6, 2019
Washington, D.C. USA

Biophysical Society

March 2-6, 2019
Baltimore, MD USA

An Introduction to High Content Imaging (RMS)

February 12, 2019
Nottingham, UK

Society of Toxicology

March 10-14, 2019
Baltimore, MD USA

3D Cell Culture

February 20-21, 2019
London, UK

Forum Labo

March 26-28, 2019
Paris, France



[Privacy Policy](#) | [Terms & Conditions](#) | [Trademarks & Logos](#)

For Research Use Only. Not for use in diagnostic procedures. The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners.

©2019 Molecular Devices, LLC 3860 N First Street, San Jose, CA 95134 USA