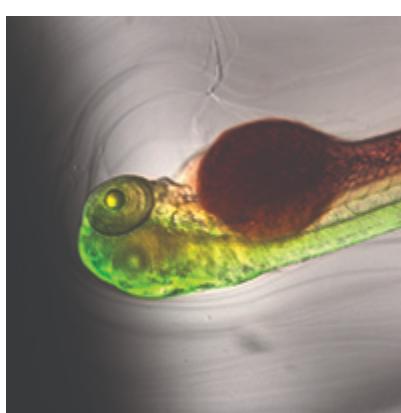


eNewsletter
December 2018

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

Product Spotlight



NEW Product introduction at ASCB | EMBO 2018
Join us December 8-12, Booth 943 to learn more about our new ImageXpress® Pico Automated Cell Imaging System and the SpectraMax® ID3 and ID5 Multi-Mode Microplate Readers.

New ImageXpress Pico system features:
-Environmental control
-Z-stack acquisition
-Double marker expression analysis
[Read More](#) ▶

eBook Spotlight

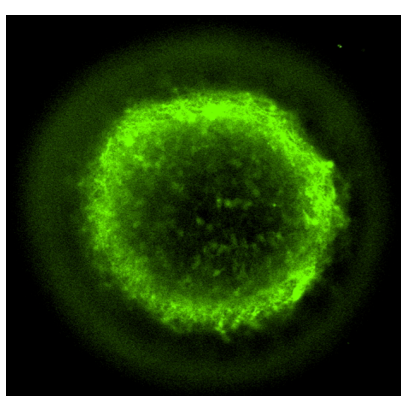
NEW Phenotypic screening with iPSC-derived cardiomyocytes and neurons

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](#) ▶



Article Spotlight



NEW Neurotoxicity profiling using 3D human iPSC-derived neural spheroids

Learn how we used an iPSC-derived 3D neurospheroid assay platform to develop more complex, biologically relevant, and predictive *in vitro* assays to screen larger sets of compounds with potential for neurotoxicity. Written in collaboration with [Stemonix](#) and the National Institute of Environmental Health Sciences. Read the peer-reviewed article online.

[Read Article](#) ▶

Application Spotlight

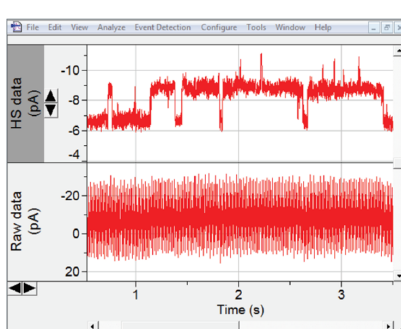
Normalize HTRF cytokine assays to cell viability

Pro- and anti-inflammatory cytokines play a central role in autoimmunity and inflammatory and infectious diseases. They are also key players in metabolic disorders and oncology, particularly the anti-tumor immune response. Homogeneous Time Resolved Fluorescence (HTRF) from Cisbio Bioassays provides a versatile platform for the quantitation of various cytokines & chemokines.

[Download App Note](#) ▶



Article Spotlight



NEW Combating electrical hum with HumSilencer technology

Electrical hum is the most common source of background noise in electrophysiology experiments. It is unwanted electrical activity detected by your recording equipment, which interferes with or hides the activity you are trying to measure. The HumSilencer™ Adaptive Noise Cancellation system provides a cleaner signal with less interference from electrical noise and better results.

[Read Article](#) ▶

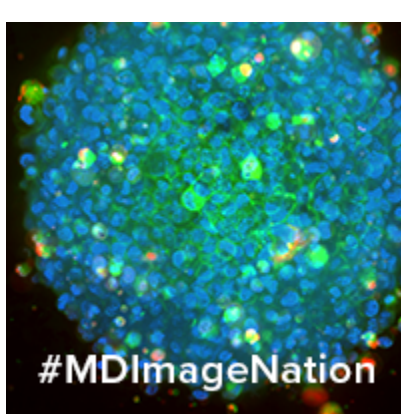
Automation Spotlight

Automated workflow solutions

The Octet HTX and RED384 systems are the industry-leading platforms for protein quantitation and kinetic characterization, performing assays at unmatched speeds with the ability to read up to 96 samples in parallel. Measure titers or perform kinetic screens in minutes, not hours. Increase the throughput and unattended run time of Octet HTX and RED384 systems with integrated automation bundles. Pick and choose a variety of possible configurations—each with varying amount of components like robotic arms, plate hotels, incubators, and more—that best fit your unique research needs.

[Download Brochure](#) ▶

Contest Spotlight



NEW ImageNation Imaging Contest Winners

Congratulations to our ImageNation imaging contest winners! Winners will be included in the 2019 Molecular Devices calendar. Download wallpapers of the winning images from the ImageNation landing page.

[Download Wallpapers](#) ▶

Upcoming Events

ASCB | EMBO
December 8-12, 2018
San Diego, CA USA
Booth #943

UK Light Microscopy Facility Managers Meeting
January 3-4, 2019
Liverpool, UK

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

An Introduction to High Content Imaging (RMS)
February 12, 2019
Nottingham, UK

3D Cell Culture
February 20-21, 2019
London, UK

Biophysical Society
March 2-6, 2019
Baltimore, MD USA

Society of Toxicology
March 10-14, 2019
Baltimore, MD USA

Forum Labo
March 26-28, 2019
Paris, France

Follow Us:

