

eNewsletter

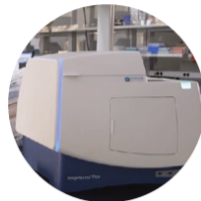
March 2018

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

Product Spotlight

NEW ImageXpress® Pico Automated Cell Imaging System

More than a digital microscope, the ImageXpress® Pico system combines cellular imaging and powerful analysis for individual labs that need an easy automated solution. Capable of brightfield, fluorescence, colorimetric, and live cell imaging, the system can analyze data on the fly and allows users to view results in various formats including image montages, scatter plots, videos, and more.



[Read More >](#)

Application Spotlight

NEW Calcium flux assay for *in vitro* neurotoxicity studies and drug screening

See the feasibility of measuring intracellular calcium flux in primary rat cortical neurons in a microplate format using the EarlyTox™ Cardiotoxicity Kit with the FlexStation® 3 Multi-Mode Microplate Reader, which enable compound handling for real-time monitoring of calcium kinetics.

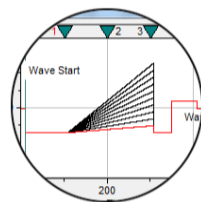


[Download App Note >](#)

Webinar Spotlight

New Speed up data analysis with the new Clampfit 11 Advanced Analysis Module

There are many new and enhanced features in the pCLAMP™ 11 Clampfit Advanced Analysis Module including population spike, action potential, and batch data analyses. Register for our webinar and learn how to use those tools.



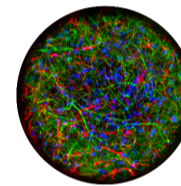
Date: March 20, 2018
Time: 11:00AM PDT, 2:00PM EDT

[Register Now >](#)

Webinar Spotlight

New StemoniX microBrain 3D Assay Ready Plates as a Human iPSC-based Neural Spheroid Platform for High Throughput Drug Screening

Learn about StemoniX® microBrain® 3D Assay Ready, a highly homogenous human induced pluripotent stem cell (hiPSC)-derived cortical spheroid screening platform.



Date: April 17, 2018
Time: 8:00AM PDT, 11:00AM EDT

[Register Now >](#)

Video Spotlight

HCS Pharma uses augmented reality to validate its industrial platform

In this video interview, Nathalie Maubon from HCS Pharma (Lille, France) and Cedric Point from Molecular Devices show how they used augmented reality to validate the HCS Pharma automated imaging platform, which includes two ImageXpress® Micro Confocal systems.



[Watch Video >](#)

See HAPix, our automation platform from inside

This short immersive video from HCS Pharma in Lille, France gives you a whole new perspective on a cellular imaging lab—from the inside of an automated imaging platform! See the HCS Pharma HAPix platform, a HCS automation platform for imaging that features two ImageXpress Micro Confocal systems.



[Watch Video >](#)

Upcoming Events

Society of Toxicology

March 11-15, 2018
San Antonio, TX USA

5th Microbiome R&D and Probiotics Europe

March 21-22, 2018
Rotterdam, The Netherlands

Forum LABO 2018

March 28-29, 2018
Lyon, France

Analytica 2018

April 10-13, 2018
Munich, Germany

11th Annual Proteins & Antibodies Congress

April 16-17, 2018
London, UK

PEGS

April 30-May 4, 2018
Boston, MA USA



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

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. ©2018 Molecular Devices, LLC

3860 N First Street, San Jose, CA 95134 USA