

eNewsletter

May 2018

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

Application Spotlight

NEW Improved cell counting method using transmitted light

Hematoxylin is a commonly used dye for staining protocols in histology. It reacts like a basic dye with a purplish-blue color and stains acidic or basophilic structures, including the cell nucleus and organelles that contain RNA. See how you can increase the contrast in transmitted light images to improve cell counting for cell lines with relatively flat morphology or indistinguishable intercellular space.

[Download Application Note >](#)

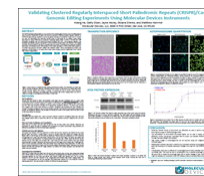


Poster Spotlight

NEW Validating CRISPR/Cas9 Genomic Editing Experiments

The CRISPR/Cas9 gene editing system is a popular tool for studying gene function, and validating it is necessary to ensure that genes of interest are successfully knocked down or knocked out. See how the SpectraMax® i3x microplate reader and the ImageXpress® Micro Confocal imaging system can be utilized in gene editing experiments by using CRISPR/Cas9 to knock down autophagy-related protein 5 (ATG5) in HEK293 cells.

[Download Poster >](#)



Demo Spotlight

NEW Capture and analyze cellular images

Compact and affordable, the ImageXpress® Pico Automated Cell Imaging System does more than imaging—it offers intelligent analysis capabilities for various cell-based assays ranging from simple cell counting to sophisticated neurite tracing.

[View Demo >](#)

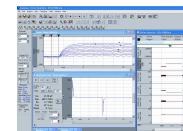


Software Spotlight

Advanced software for flexible data acquisition and analysis

The pCLAMP™ 11 Software for patch-clamp electrophysiology controls and records voltage-clamp, current-clamp, and patch-clamp experiments. It allows you to design advanced protocols easily, streamline workflows, and perform analysis faster.

[Read More >](#)



Customer Spotlight

University of Edinburgh

The Edinburgh Genome Foundry (EGF) manufactures genetic material for their customers using a fully automated robotic platform. To meet the needs of the EGF, the Molecular Devices Advanced Workflow Engineering Solutions (AWES) Team modified a QPix™ 420 system to offer hardware for liquid handling and modified structural design to allow for robotic arm access.

[Read More >](#)



Citation Spotlight

SoftMax® Pro 7 Software preconfigured protocols and custom assay workflows simplify microplate data acquisition and analysis. Read more in the citation below.

Citation: Validation of the ELISA Method for Quantitative Detection of TNF- α in Patients with Intracerebral Hemorrhage

[Read More >](#)



Upcoming Events

PEGS

April 30-May 4, 2018
Boston, MA USA
Booth #212

Cell Line Development & Engineering

June 12-14, 2018
Park Central Hotel
San Francisco, CA

SLAS Europe

June 27-29, 2018
Brussels, Belgium
Booth #401

FENS

July 7-11, 2018
Berlin, Germany
Booth #132

SIMB

August 12-16, 2018
Chicago, IL USA

MipTec

September 11-14, 2018
Basel, Switzerland
Booth #A48



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