

eNewsletter

June 2018

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

Product Spotlight

NEW Small but mighty absorbance readers

Bridge the gap between the affordability of filter-based readers and the flexibility of monochromator-based systems with the SpectraMax® ABS Absorbance Microplate Readers. Advanced technology provides you with the flexibility, sensitivity, and convenience for a wide range of assays such as ELISAs, microbial growth, and protein quantitation, all in an incredibly small footprint. The SpectraMax ABS readers fit in every lab and within every budget.

[Learn More >](#)



Application Spotlight

NEW Measure ROS

Reactive oxygen species (ROS) are involved in cellular processes such as cell signaling, homeostasis, and immunological defenses. However, elevated ROS levels can cause oxidative damage to cellular components. Far-red fluorescent assay kits can quickly measure ROS without interference from autofluorescence. The SpectraMax® iD5 microplate reader reduces signal background and stray light, providing wide dynamic range and sensitivity for far-red fluorescent assays.

[Download App Note >](#)

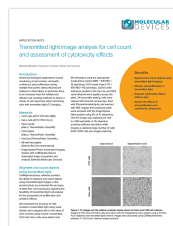


Application Spotlight

NEW Transmitted light image analysis for cell count and assessment of cytotoxicity effects

There is increasing need for reliable and efficient label-free cell counting and characterization methods to assess a variety of cell responses when monitoring cells with transmitted light imaging. Learn about the convenience, precision, and accuracy of cell characterization in transmitted light without using dyes.

[Download App Note >](#)

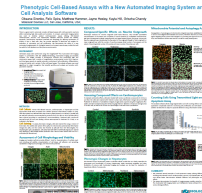


Poster Spotlight

NEW Phenotypic cell- based assays with the ImageXpress Pico

The need to automate complex cell-based assays with multi-parametric readouts while maintaining high data quality and precision is growing. In this poster presented at SLAS 2018, we demonstrated several assay models using the ImageXpress® Pico Automated Cell Imaging System.

[Download Poster >](#)



Customer Spotlight

NEW University College London

Professor Townsend-Nicholson at UCL, UK, is interested in the role cell surface receptors play in transducing extracellular signals into intracellular responses. The FLIPR Tetra® High-Throughput Cellular Screening System and SpectraMax® i3x Multi-Mode Microplate Reader are used for a range of assays, including calcium signaling, membrane potential, cAMP, hTR-FRET and dsDNA quantitation.

[Read More >](#)



Article Spotlight

NEW Rising growth of the HCS software and services market

The global high-content screening (HCS) software and services market is expected to reach approximately \$1.4 billion by 2026. Developments in informatics solutions and high-content imaging systems, increased focus on cell-based research, and rising government support of R&D are key factors driving market growth. Read more in the *Drug Target Review* article featuring Jeff McMillan from Molecular Devices.

[Read More >](#)



Upcoming Events - done

Cell Line Development & Engineering

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

ISSCR

June 20-23, 2018
Melbourne, Australia
Booth #34

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](#)