

eNewsletter

July 2018

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

eBook Spotlight

NEW Cellular Imaging Insights

With cost efficient, high-throughput instruments, such as the [ImageXpress® imaging portfolio](#), researchers can gain insight and expedite their studies while better understanding the interaction of cells under different conditions. This eBook provides new tools to better understand complex imaging models.

[Download eBook >](#)

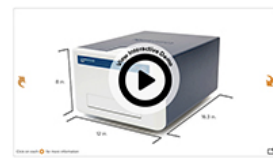


Product Spotlight

NEW Latest addition of absorbance readers to the trusted Molecular Devices family

The [SpectraMax® ABS and ABS Plus](#) are equipped with advanced technology for a wide range of assays such as ELISA, DNA, RNA, protein quantitation, and UV-VIS applications in a small footprint to fit every lab and every budget. Learn more about these feature-packed absorbance readers.

[View Interactive Demo >](#)



Infographic Spotlight

NEW Colony picking for microbiome studies

Growing and picking the right microbial colonies is an essential part of microbiome studies. Traditionally, colony picking is performed manually using sterile pipette tips or inoculation loops, which is usually slow, labor intensive, and time consuming. See how the automated QPix™ colony picker makes the entire process quicker (pick up to 3,000 clones per hour) to provide consistent and reliable results.

[Download Infographic >](#)



Infographic Spotlight

NEW Colony picking in synthetic biology

In synthetic biology, building blocks are often generated from bacteria, and the cyclical 'design-build-test' process often requires screening and picking colonies of interest. Research can therefore benefit greatly from the flexibility and throughput of an automated colony picker. See how the automated QPix colony picker makes the entire process quicker (pick up to 3,000 clones per hour) to provide consistent and reliable results.

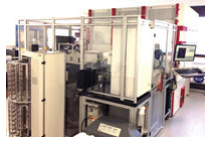
[Download Infographic >](#)



Customer Spotlight

NEW Gustave Roussy Cancer Center

The Kroemer lab at the Gustave Roussy Cancer Center in France is interested in the molecular mechanisms of cellular stress and death in cancer treatment. To enable their cutting-edge research, their Cell Biology platform needed a high-content analysis system that they could fully integrate into an automated platform. After evaluating the market, they chose ImageXpress Widefield systems.

[Read More >](#)

Citation Spotlight

The [SpectraMax® iD3 reader](#) was used to measure fluorescence intensity as shown in [Macromolecular Bioscience](#).



Citation: Synthesis and Evaluation of Chloroquine-Containing DMAEMA Copolymers as Efficient Anti-miRNA Delivery Vectors with Improved Endosomal Escape and Antimigratory Activity in Cancer Cells

[Read More >](#)

Upcoming Events

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

LabDays

September 12-13, 2018
Copenhagen, Denmark

WOTS

Oct 2-5, 2018
Utrecht, Netherlands

ELRIG Drug Discovery

October 9-10, 2018
London, UK

[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