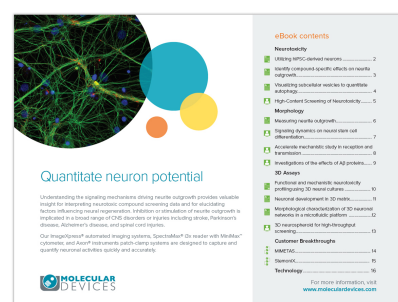


eNewsletter

October 2018

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

eBook Spotlight



NEW Quantitate Neuron Potential

Understanding the signaling mechanisms driving neurite outgrowth provides valuable insight for interpreting neurotoxic compound screening data and for elucidating factors influencing neural regeneration. In this eBook, gain valuable insights for interpreting neurobiology, from evaluation of ion channels to analysis of 3D neuron organoids.

[Download eBook](#) ▶

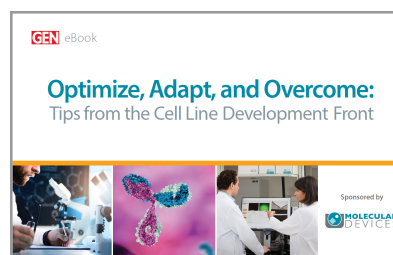
eBook Spotlight

NEW Optimize, adapt, and overcome: tips from the cell line development front

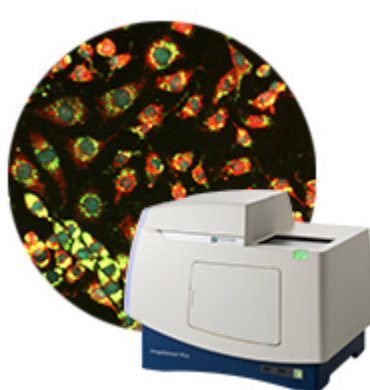
We have partnered with Genetic Engineering & Biotechnology News (GEN) to bring you this eBook, which features the following articles:

- Single-cell printing for establishing monoclonality
- Confident assurance of clonality using fluorescent dye
- How to improve cell line productivity

[Download eBook](#) ▶



Webinar Spotlight



NEW From cellular images to results in minutes with automated cellular imaging and analysis

There is a great need to automate complex cell-based assays with multi-parametric readouts while maintaining high data quality and precision. We explore the accessibility and utility of the ImageXpress™ Pico Automated Cell Imaging System for biological research and will present several assay models that are useful for both academic and bio-pharma environments.

Date and Time:

Wednesday, October 17, 2018

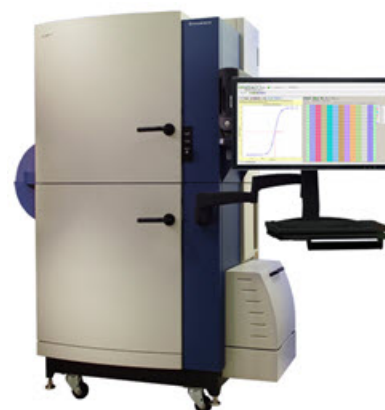
8:00 am PDT | 11:00 am EDT | 4:00 pm GMT

[Register Now](#) ▶

Webinar Spotlight

NEW Webinar: FLIPR and Pivot Park Screening Centre

Discover how Pivot Park Screening Centre leverages the FLIPR Tetra® System to perform high-throughput screening campaigns for the European Lead Factory (ELF). Also, learn about cardiomyocyte screening using full automation to perform complex GPCR testing, tips and tricks for 1536-well FLIPR screening, calcium kinetics in cardiac safety testing, and more.



Date and Time:

Tuesday, October 23, 2018

7:00 am PDT | 10:00 am EDT | 3:00 pm GMT

[Register Now](#) ▶

Application Spotlight



NEW Measure oxidative metabolism and glycolytic activity

Cells require fuel in the form of ATP to carry out most essential bioprocesses. This fuel is generated through glycolysis and mitochondrial respiration. While both produce ATP, glycolysis can function in the absence of oxygen, whereas mitochondria require oxygen for the final step in oxidative phosphorylation (OXPHOS). Here, we describe how to use the Agilent MitoXpress® Xtra Oxygen Consumption Assay and the Agilent pH-Xtra™ Glycolysis Assay on the SpectraMax® i3x reader to facilitate metabolic analysis.

[Download Application Note](#) ▶

Application Spotlight

NEW Identify compound-specific effects on neurite outgrowth

Neurite outgrowth is a commonly used assay to study neuronal development and neuronal degeneration *in vitro*. Development of neurites requires a complex interplay of both extracellular and intracellular signals. The growth of neurites can be stimulated or inhibited by neurotrophic factors. In this application note, we evaluate the efficiency of the ImageXpress Pico system to measure the ability of compounds to adversely affect the developing nervous system.

[Download Application Note](#) ▶



Customer Spotlight



NEW Phenotypic screening of physiologically relevant 3D cell models

HCS Pharma in Lille, France, is a biotechnology start-up focused on *in vitro* preclinical research and development. The team needed to implement 3D confocal imaging using powerful and reliable high-throughput systems that could be integrated within a robotic platform, allowing high-throughput imaging. They used the ImageXpress® Micro Confocal Systems for phenotypic screening of physiologically relevant 3D cell models.

[Read More](#) ▶

Contest Spotlight

NEW First ImageNation Imaging Contest Winner

Congratulations to our first ImageNation imaging contest winner, Sophia Su (Wahlin Lab, UCSD)! Her image of a 3D mini-retina organoid will be featured in the 2019 Molecular Devices calendar. Click on the link to submit an entry or vote for your favorites.

[Read More](#) ▶



Credit: Sophia Su, Wahlin Lab, UCSD

Upcoming Events

WOTS

October 2-5, 2018
Utrecht, Netherlands

ILMAC

October 3-4, 2018
Lausanne, Switzerland

BioProduction Congress

October 8-10, 2018
Dublin, Ireland
Booth #59
(Join ForteBio, our new Biologics business unit)

ELRIG Drug Discovery

October 9-10
London, UK

SLAS Europe

October 17-19, 2018
Leiden, Netherlands
Booth HCS001

SFN

November 3-7, 2018
San Diego, CA USA
Booth #3716

10th Annual PEGS Europe

November 12-16, 2018
Lisbon, Portugal

ASCB

December 8-12
San Diego, CA USA

Follow Us:

