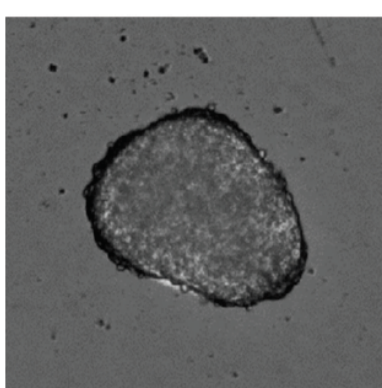


eNewsletter
September 2018

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

Application Spotlight



NEW Assay apoptosis in magnetically bioprinted HepG2 spheroids

A variety of methods for transforming 2D cultures into 3D spheroids have been devised but require a long time to form spheroids that can be difficult to handle. See how HepG2 cells can be bioprinted using the magnetic 3D cell culture system from Greiner Bio-One (m3D). The resulting spheroids are imaged on the SpectraMax® i3x reader with MiniMax™ cytometer.

[Download Application Note](#) ▶

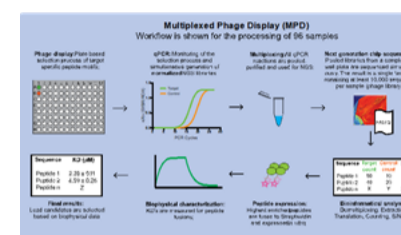
Webinar Spotlight

NEW Achieving higher throughput with multiplexed phage display

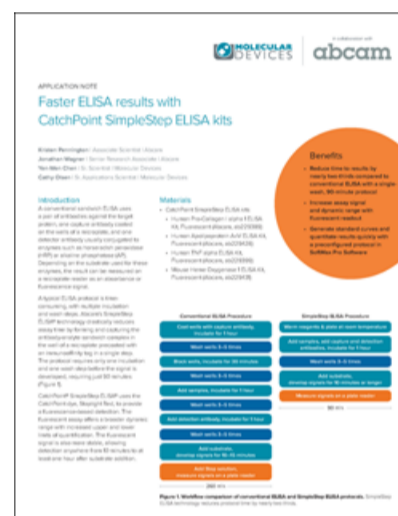
Multiplexed phage display provides a simple, robust, and seamless high-throughput-ready version of phage display that can simultaneously screen and bioinformatically analyze up to 96 peptide phage libraries in parallel. To accommodate the large number of potential lead peptides from the screening, the ForteBio Octet® System was used to characterize streptavidin peptide fusions generated by *in vitro* translation.

Webinar details: **September 27, 2018 at 8:00 AM PDT**

[Register Now](#) ▶



Application Spotlight



NEW Faster ELISA results with CatchPoint SimpleStep ELISA kits

Typical ELISA protocols are time-consuming, with multiple incubation and wash steps. CatchPoint® SimpleStep ELISA® technology drastically reduces assay time by forming and capturing the antibody-analyte sandwich complex in the well of a microplate precoated with an immunoaffinity tag in a single step.

- Reduce time to results with a single-wash, 90-minute protocol
- Generate standard curves and quantitate results quickly
- Increase assay signal and dynamic range with fluorescent readout

[Download Application Note](#) ▶

Application Spotlight

NEW Optimization of NaV1.5 channel assay

Voltage-gated ion channels are present in the excitable cell membranes of heart, skeletal muscle, brain and nerve cells. Blocking or modulating such channels can have a therapeutic effect or may interfere with normal cell function. Here, CHL cells transfected with the cardiac NaV1.5 channel are used to demonstrate differences in membrane potential assay functionality with tetrodotoxin and lidocaine & show optimization results with the FLIPR® Membrane Potential Assay Kits.

[Download Application Note](#) ▶



Application Spotlight



NEW Detection of Autophagy using ImageXpress™ Pico system

Autophagy and its dysregulation has been found to play an important role in neurodegenerative diseases and cancers, therefore the discovery of novel therapeutic targets along this process has emerged as a promising approach for drug therapies. In our application note, learn a method for detecting autophagy particles which can be used for assay development and quantitation of the compound effects on the process of autophagy.

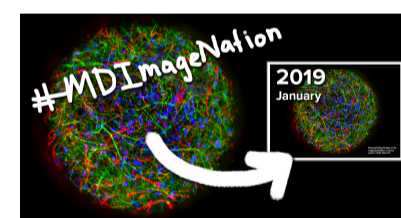
[Download Application Note](#) ▶

Contest Spotlight

NEW Imaging Contest

Get recognized for your scientific research using your ImageXpress imaging system. Submit an image (or two) to the Molecular Devices Imaging Contest. Winners will be featured in the 2019 Molecular Devices calendar. Click on the link to submit an entry or vote for your favorites.

[Submit Entry](#) ▶



Upcoming Events

ForteBio User Group Meeting

September 11, 2018
Boston, MA USA

MipTec

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

LabDays

September 12-13, 2018
Copenhagen, Denmark
Booth #19

Europhysiology

September 14-16, 2018
London, UK
Booth #21

SBI2

September 19-20, 2018
Boston, MA USA

Molecular Devices High-Content Imaging User Meeting

September 19, 2018
Boston, MA USA

Discovery on Target

September 25-28, 2018
Boston, MA USA

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)

Follow Us:

