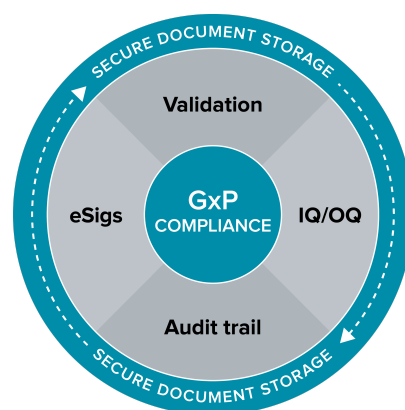


eNewsletter

November 2018

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

Software Spotlight



NEW SoftMax Pro 7.1 GxP Software

SoftMax® Pro 7.1 GxP Software is the latest, most secure software to achieve full FDA 21 CFR Part 11 compliance with streamlined workflows to ensure data integrity. Every step is optimized to simplify analysis and reporting to support our microplate readers. We will partner with you to set up single- or enterprise-level software and provide IQOQ services to establish full compliance on your microplate readers.

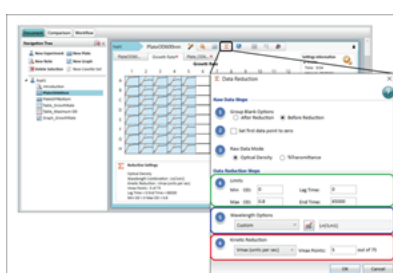
[Read More](#) ▶ [Download eBook](#) ▶

Application Spotlight

NEW Advanced kinetic analysis of a bacterial growth assay

Difficult-to-kill bacteria have become a problem in hospital-acquired infections. Identifying compounds which kill these bacteria is vital, and developing and optimizing assays to screen for the efficacy of these compounds is a challenge. Here, we show you how to set up a bacterial growth assay and easily calculate GFP signal cell density, extract growth rates, and other relevant kinetic information.

[Download Application Note](#) ▶



On-Demand Webinar

NEW Flexible monitoring of small molecules in complex matrices with Aptamer Affinity Reagents using Bio-Layer Interferometry

Are you having difficulty screening for small molecules in complex matrices? We partnered with the Aptamer Group to present data showing the versatility of our Octet BLI platforms with example data for different aptamer-based applications, with a focus on small molecule detection and quantification.

[Watch Webinar](#) ▶

On-Demand Webinar

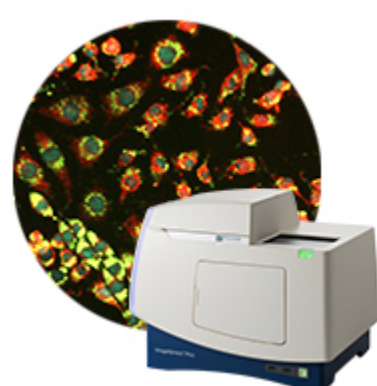
NEW FLIPR 1536-well application in high throughput screening

Did you miss our webinar last week with Pivot Park Screening Centre? Register to watch the on-demand replay of "FLIPR in Drug Discovery: Ion Channel Screening and Cardiac Safety Testing." 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 using full automation to perform complex GPCR testing, tips and tricks for 1536-well screening, and calcium kinetics in cardiac safety testing.

[Watch Webinar](#) ▶



On-Demand Webinar



NEW From cellular images to results in minutes with automated imaging

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.

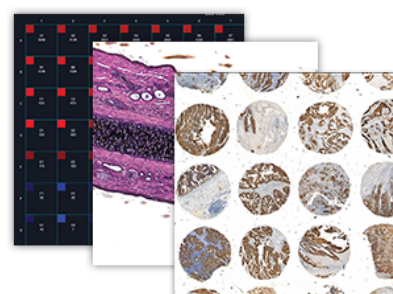
[Watch Webinar](#) ▶

Promotion Spotlight

NEW Image your cellular assay for free

Test our ImageXpress Pico Automated Cell Imaging System by letting us image your favorite cellular assay. Send us your sample and we will provide you with images and data. Get more from your images. Terms & Conditions Apply.

[Learn More](#) ▶



Upcoming Events

10th Annual PEGS Europe

November 12-16, 2018
Lisbon, Portugal
(Join ForteBio, our new Biologics business unit)

ASCB

December 8-12, 2018
San Diego, CA USA
Booth #943

French Society for Stem Cell Research

December 11, 2018
Nantes, France

SLAS

February 2-6, 2019
Washington, D.C. USA

Biophysical Society

March 2-6, 2019
Baltimore, MD USA

Follow Us:

