

WEBINAR

## Expand your assay applications with Multi-Mode Microplate Readers – From ELISA to AlphaScreen

Jenna Bedard, M.S.  
Field Applications Scientist, BioResearch  
Molecular Devices



**NEW**  
On-demand Webinar

**Expand your assay applications with multi-mode microplate readers—from ELISA to AlphaScreen**

Discover the multitude of avenues you can pursue with multi-mode microplate readers. In this webinar you will learn:

- What is absorbance, fluorescence, luminescence, TRF, FP, and AlphaScreen
- What are common assays for each read mode
- How to detect varying levels of Insulin in Type 1 Diabetic Patients' pancreatic beta cells
- How each multi-mode microplate reader can fit your lab's specific needs
- How SoftMax® Pro GxP Software can help you achieve data compliance in GMP/GLP labs

[View On-demand Webinar](#)



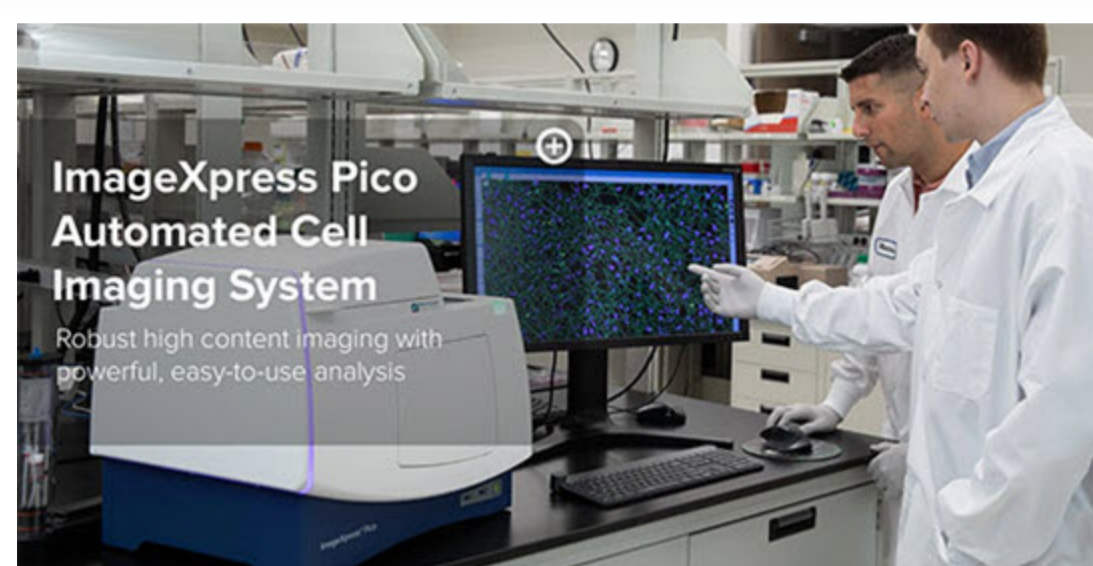
**NEW**  
Application Spotlight

**Maximize your data using multiplexed high-content screening assays**

There has been growing interest in the use of multiplexed high-content screening assays for drug discovery and biological research, as it enables scientists to gather more meaningful data in a single experiment—saving both time and money.

Here, we present two application notes that describe the development of multiplexed assays for toxicity assessment using the ImageXpress® Pico Automated Cell Imaging System and ImageXpress® Micro 4 High-Content Imaging System.

[Download Application Notes](#)



**NEW**  
Interactive Experience

**ImageXpress Pico Automated Cell Imaging System**

If you're looking to image your cells, learn more about the ImageXpress Pico system with our new interactive tools! Watch a demo video, learn more about available objectives and filters, and view the various available configurations. The system can be configured to fit your assay needs!

[Take Interactive Tour](#)



**NEW**  
eBook Spotlight

**Get a front row view of your cells**

Our SpectraMax® i3x Multi-Mode Microplate Reader with MiniMax™ 300 Imaging Cytometer enables rapid cell imaging and analysis, giving you a better view of cell proliferation, marker expression, and cytotoxicity.

- Image cells, run cell-based assays, and detect western blots, all on one platform
- Count cells and estimate confluence without fluorescent dyes
- Improve accuracy of cell counting for difficult-to-image cell types

[Download eBook](#)

**NEW**  
Application Spotlight

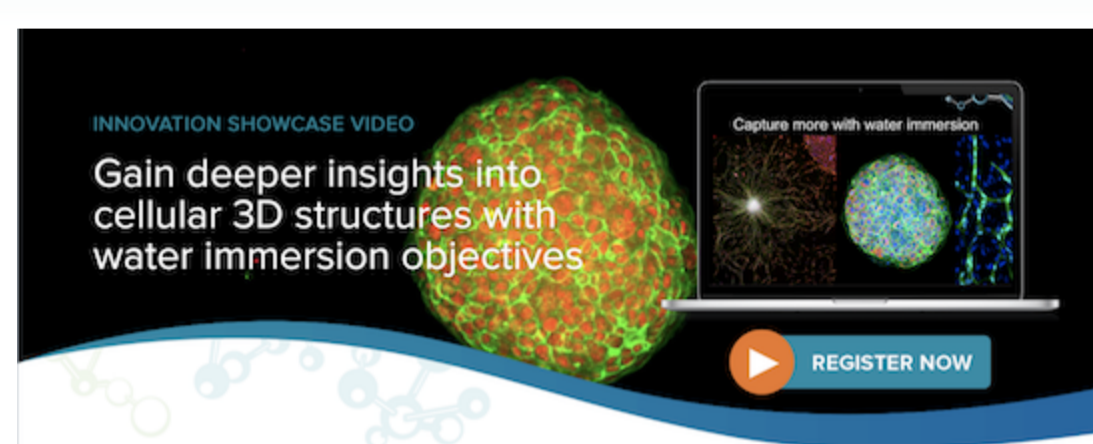
**Spheroids**

Cancer is a leading cause of death worldwide, but researchers are making progress towards understanding its underlying mechanisms.

Spheroids, 3D aggregates of cells in culture, provide more physiological relevance than traditional 2D cell culture. Learn how spheroids formed using a variety of cancer cell types.



[Download Application Notes](#)



**NEW**  
Innovation Spotlight

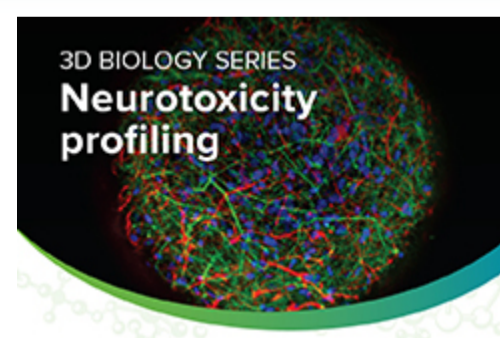
**Gain deeper insights into cellular 3D structures with water immersion objectives for high-content imaging**

With the increasing interest in using 3D cultures for assay development and phenotypic screening for a range of cellular models, water immersion objectives are essential for capturing more data at greater depths in 3D structures, such as spheroids and thick tissues.

This presentation details how high-performance water immersion objectives on the ImageXpress® Micro Confocal High-Content Imaging System enhance the resolution, sensitivity and throughput for a variety of complex 3D cell-based assays.

As a supporter of, and participant in, ISSCR's 2020 Virtual, we share with you our Innovation Showcase presentation. This presentation does not represent an endorsement from or support of the ISSCR.

[View Video](#)

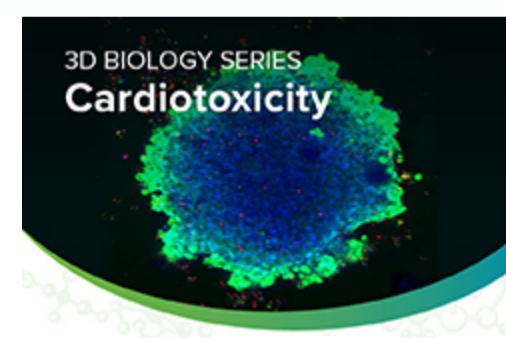


**NEW**  
Application Spotlight

**Neurotoxicity profiling**

We show that functional and morphological assays using 3D neural spheroids formed with human iPSC-derived cells can be used for evaluation of drug candidates and neurotoxicity assessment.

[Download Application Notes](#)



**NEW**  
Application Spotlight

**Cardiotoxicity**

Live-cell imaging and 3D models provide greater predictivity for compound efficacy and toxicity. We present methods for the formation of 3D spheroids derived from human iPSC-derived cells.

[Download Application Notes](#)

**EVENTS**

**LabDays Copenhagen**  
September 2-3, 2020 | Copenhagen, Denmark

**Organ-on-a-Chip, Tissue-on-a-Chip & Organoids**  
September 9-10, 2020 | Rotterdam, The Netherlands

**SB12 - Virtual**  
September 16-17, 2020 | North America

**Discovery on Target - Virtual**  
September 16-18, 2020 | North America

**SynBioBeta - Virtual**  
September 29-October 1, 2020 | North America

**Future Labs Live - Virtual**  
November 17-18, 2020 | North America

**Cell Bio - Virtual**  
December 2-16, 2020 | North America

**Antibody Engineering - Virtual**  
December 13-17, 2020 | North America



[Privacy Policy](#) | [Terms & Conditions](#) | [Trademarks & Logos](#)

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. ©2020 Molecular Devices, LLC 3800 N First Street, San Jose, CA 95134 USA