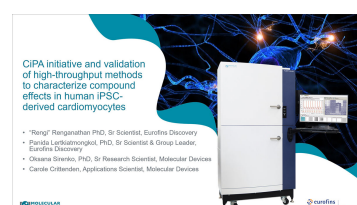


**[NEW] Product Spotlight**

**Water immersion objectives for the ImageXpress Micro Confocal system**

We're excited to announce the release of automated water immersion imaging technology for the ImageXpress® Micro Confocal High-Content Imaging System. Scientists can now capture greater physiologically-relevant phenotypic data for 3D and thick tissue samples. This technology will enable both acquisition and analysis of 3D cell model assays such as spheroids, organoids, and organ-on-a-chip biology, increasing signal up to four times, improving z-resolution, and decreasing optical aberrations for sharper, crisper images.

[Learn More](#)



**[NEW] On-Demand Webinar Spotlight**

**CiPA initiative and validation of high-throughput methods to characterize compound effects in human iPSC-derived cardiomyocytes**

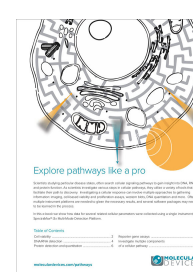
In collaboration with Eurofins, we discuss the assessment of potential cardiotoxicity of compounds with Comprehensive *in vitro* Proarrhythmia Assay (CiPA), the new cardiac safety testing paradigm that includes *in vitro* assays using human induced pluripotent stem cell derived cardiomyocytes (hiPSC-CM).

[View On-Demand Webinar](#)

**[NEW] Award Spotlight**

**Drug Discovery & Development Email of the Year!**

We were very proud to collect this year's SelectScience award for our email, 'Your guide to exploring cellular pathways'. The email provided a downloadable eBook offering guidance on how to better explore cell pathways by performing multiple techniques including cell imaging, cell-based viability and proliferation assays, western blots, DNA quantitation, and more—on a single instrument, the SpectraMax® i3x Multi-Mode Microplate Reader.



[Download eBook](#)



**Application Spotlight**

**Monitor multiple stages of apoptosis with live cell kinetic imaging**

The study of apoptosis is a critical aspect of drug discovery and development. Additionally, studying the relationship between apoptosis and other factors, such as oxidative stress, is crucial for understanding specific diseases that are associated with the dysregulation of apoptosis.

See how we ran two apoptosis assays to study the cytotoxic effects of anti-cancer compounds on HeLa cells. Long-term time-lapse imaging was performed on the ImageXpress® Pico system. On-board environmental control maintained live cell conditions by monitoring CO<sub>2</sub>, O<sub>2</sub>, temperature, and humidity.

[Learn More](#)

**[NEW] Citation Spotlight**

**Assessment of antioxidant and antidiabetic properties of *Agaricus blazei* Murill extracts**

*Agaricus blazei* Murill (ABM), a medicinal mushroom, has beneficial effects on various human metabolic diseases. The objective of this research was to evaluate the antioxidant and antidiabetic properties of ABM extracts (ethanol extract and ethyl acetate extract).

See how the SpectraMax® i3x reader was used for absorbance readings.



[Read More](#)

**Events**

**Biophysical Society**  
Booth #501  
February 15-19, 2020  
San Diego, CA USA

**PEGS**  
Booth #424  
May 4-8, 2020  
Boston, MA USA

**Society of Toxicology**  
Booth #677  
March 15-19, 2020  
Anaheim, CA USA

**World Pharma Week**  
Booth #317  
June 2-4, 2020  
Boston, MA USA

**AACR**  
Booth #3040  
April 24-29, 2020  
San Diego, CA USA

**ISSCR**  
Booth #400  
June 24-27, 2020  
Boston, MA USA



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