

**[NEW] Product Spotlight**

**Introducing Digital Confocal, Live Preview and other new features for ImageXpress Pico system**

We are proud to announce several new features for the ImageXpress® Pico Automated Cell Imaging System. Features such as Digital Confocal, 2D on-the-fly deconvolution, Live Preview, and multi-wavelength cell scoring offer you the ability to advance your discoveries in a small, affordable imager. More than a digital microscope, the system combines high-resolution imaging with powerful analysis.

[Learn More](#)



**[NEW] Application Spotlight**

**Measure cancer cell viability using a homogeneous, stable luminescence assay**

Luminescent cell viability assays use firefly luciferase reaction as a way to determine the relative numbers of living cells under different treatments or experimental conditions. Metabolically active cells produce ATP, which is required by the luciferase reaction. When ATP is the limiting component in the reaction, the amount of light produced is proportional to and serves as the readout for the number of viable cells.

Here, we validated performance of Promega's CellTiter-Glo 2.0 Cell Viability Assay on the SpectraMax® iD5 Multi-Mode Microplate Reader. Both ATP and viable cells were used to generate standard curves demonstrating the sensitivity and linearity of the assay.

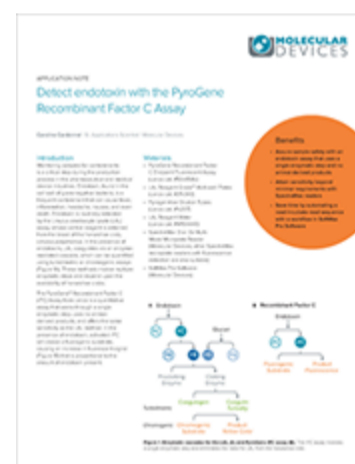
[Download Application Note](#)

**[NEW] Application Spotlight**

**Detect endotoxin with the PyroGene Recombinant Factor C Assay**

Monitoring samples for contaminants is a critical step during the production process in the pharmaceutical and medical device industries. Endotoxin, found in the cell wall of gram-negative bacteria, is a frequent contaminant that can cause fever, inflammation, headache, nausea, and even death.

Assure sample safety with the PyroGene™ Recombinant Factor C (rFC) Assay that uses a single enzymatic step and no animal-derived products. Plus, attain sensitivity beyond minimal requirements with the SpectraMax® readers.



[Download Application Note](#)



**[NEW] Application Spotlight**

**Scaling up automated tube formation assay for targeting tumor angiogenesis**

Tube formation assay a model that is representative of an *in vivo* microenvironment which allows researchers to rapidly evaluate anti-angiogenic effects by quantifying capillary-like tubular structures derived from endothelial cells grown in an extracellular matrix.

Here, we describe the development of an automated tube formation assay that streamlines image acquisition and analysis, which is a major bottleneck for scaling up similar assays.

[Download Application Note](#)

**[NEW] Application Spotlight**

**Simplify 3D cell imaging and analysis**

Utilization of 3D cellular assays adds value to research and screening campaigns, spanning the translational gap between 2D cell cultures and whole-animal models. By reproducing important parameters of the *in vivo* environment, 3D models can provide unique insight into the behavior of stem cells and developing tissues *in vitro*.

Learn how to overcome challenges of 3D workflows by visiting our new application page highlighting 3D imaging and analysis techniques, tips, and resources.



[Learn More](#)

**Events**

**BioProduction at CPHI Worldwide**  
November 5-7, 2019  
Frankfurt, Germany

**PEGS Europe**  
November 18-22, 2019  
Lisbon Portugal

**ISSCR**  
Booth #8  
November 6-8, 2019  
Toronto, Canada

**Next-Generation Protein Analysis and Detection**  
December 2-3, 2019  
Ghent, Belgium

**American Institute of Chemical Engineers**  
Booth #159  
November 10-15, 2019  
Orlando, FL USA

**ASCB | EMBO 2019**  
Booth #1103  
December 7-11, 2019  
Washington, DC USA

**Vaccines R&D**  
November 18-20, 2019  
Newton, MA USA

**Antibody Engineering and Therapeutics**  
Booth #302  
December 9-13, 2019  
San Diego, CA USA

