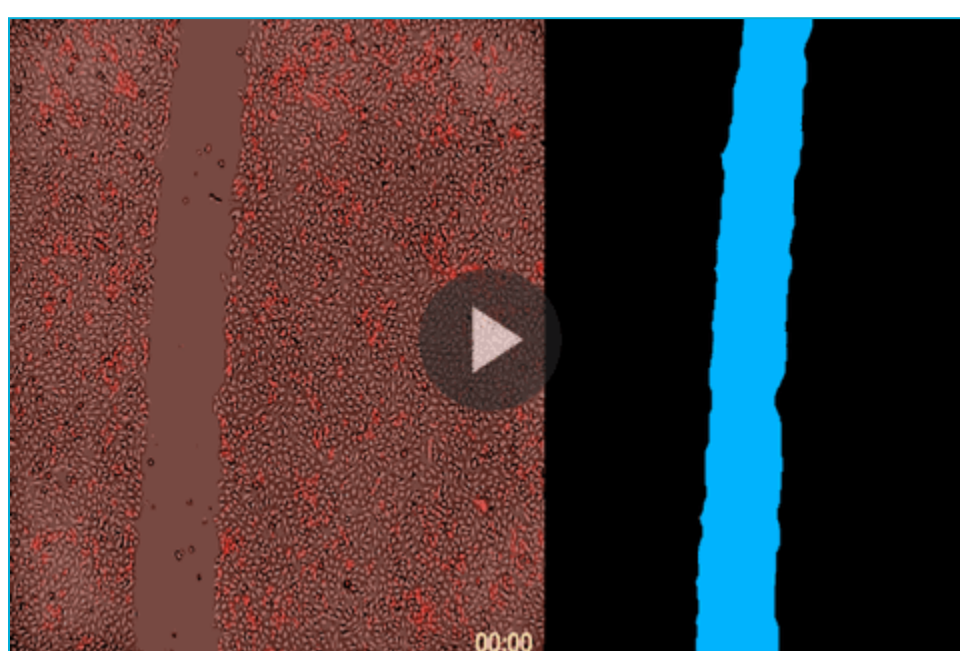


**[NEW] Video Spotlight**

The ImageXpress® Pico Automated Cell Imaging System combines high-resolution imaging with powerful data analysis. With new z-stack acquisition and environmental control functionality, the ImageXpress Pico helps you to advance your discovery even further.

To learn more, click on the video below.



**[NEW] Application Spotlight**

**Measure total protein in cell lysates**

Quantification of protein concentrations from cell lysates is a key step for many downstream applications, such as western blots and enzyme-linked immunosorbent assays (ELISAs). The Pierce Rapid Gold BCA and Pierce Detergent Compatible protein assays provide users easy methods for quantification of total protein in cell lysate samples without long incubation times or worries over the effects of detergent in lysis buffers.

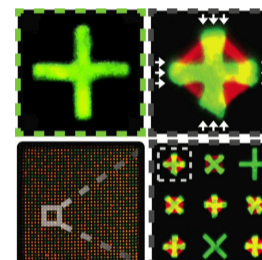
The SpectraMax® ABS Plus reader with SoftMax® Pro Software offers a complete solution for data acquisition and analysis, with preconfigured protocols that plot standard curves and calculate sample concentrations automatically. With the Speed Read setting, a 96-well plate can be read in just five seconds, further streamlining the process of protein quantification.

[Download Application Note](#) ▶

**[NEW] Poster/Application Spotlight**

**High-throughput single-cell contractility measurements using FLECS Technology**

Read about the development of an automated single-cell functional contractility assay called FLECS ([Forcyte Biotechnologies](#)) to assess both the tonic contractility of cells and the ability of test compounds to modulate the forces applied by the cells.



[Download Poster and Application Note](#) ▶



**Infographic Spotlight**

Putting together a patch-clamp rig can be intimidating. To help you get started, we've simplified the process with this quick infographic that details the importance of each component in a complete setup.

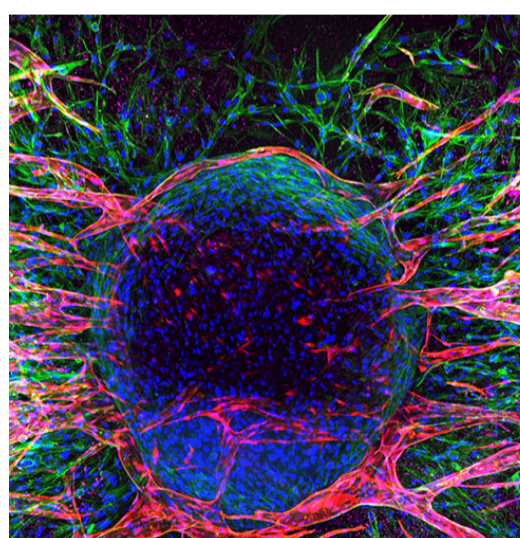
[Download Infographic](#) ▶

**Customer Spotlight**

**MIMETAS, the organ-on-a-chip company**

MIMETAS uses their OrganoPlate®, a unique 3D organ-on-a-chip platform, to develop disease, toxicology, and transport models for research, development and drug screening. They use the ImageXpress® Pico and ImageXpress® Micro Confocal systems to develop tissue models for their OrganoPlates.

[Read More](#) ▶



Vascularized liver spheroid in the OrganoPlate (Blue: DNA, Green: Actin, Red: RFP-HUVEC, Magenta: VE-Cadherin)

**Events**

**LABVOLUTION**

Booth #B72  
May 21-23, 2019  
Hannover, Germany

**ELRIG - Advances in Cell Based Screening in Drug Discovery**

May 22-24, 2019  
Gothenburg, Sweden

**3D-Culture, Organoids & Tox Screening**

Booth #E4  
June 13-14, 2019  
Rotterdam, The Netherlands

**Organ-on-a-chip & Tissue-on-a-chip**

Booth #19  
June 18-19, 2019  
Rotterdam, The Netherlands

**SLAS Europe**

June 26-28, 2019  
Barcelona, Spain

**ISSCR**

Booth #407  
June 26-29, 2019  
Los Angeles, CA, USA

Follow Molecular Devices:

