



Welcome to the September edition of our newsletter!

Bringing you our latest discoveries in high-throughput screening, genomic and cellular analysis, colony selection, and microplate detection. Visit us at: moleculardevices.com



From customer feedback to workflow improvements

The path to understanding complex biological processes and diseases is paved with a lot of challenges. As the desired level of insight increased, so did the requirements from state-of-art laboratory solutions. That's why solutions that worked last year may be inadequate in solving today's research questions.

Paying close attention to the complex research goals and methods of collaborators, we've implemented several updates to our high-content imaging and analysis solutions. This article summarizes customer feedback and the recent product updates.

[Let's Innovate](#)

Interested in the research shared at ISSCR 2021?

This year's ISSCR (The International Society for Stem Cell Research) Annual Meeting brought leaders together from across the globe to discuss new technologies, share research insights, and explore the newest breakthroughs in stem cell science and regenerative medicine.

If you didn't get a chance to attend our Innovation Showcase or poster presentations during the show—we've got you covered. We've gathered all of our poster sessions including the latest research on organoid automation, machine learning analysis, cancer tumoroids, brain organoids, and cardiac assay—all in a handy, one-stop-shop just for you.



[Download the Research](#)



Proven antibody discovery solutions from target validation to scale up

Accelerate vaccine development with a complete solution that addresses all stages of antibody discovery research—from target validation, screening, and clone selection to characterization and process scale up. Our comprehensive portfolio suite is designed to optimize your productivity to get safe and efficacious vaccines and therapeutics to market faster.

[Download eBook](#)

Analyzing 3D spheroids for high-throughput screening

Imaging and analysis of spheroids are crucial to understanding the formation and behavior of tumor microenvironments. It can also aid in obtaining more insightful and realistic data from anti-cancer drug screening. Here we developed and optimized methods for the formation of 3D spheroids, as well as confocal imaging and analysis methods for high-throughput screening.



[Learn More](#)



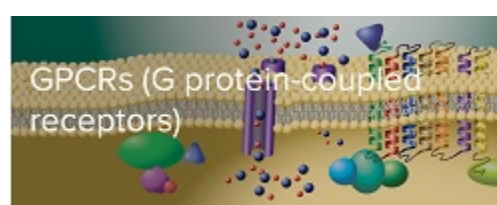
Highly robust homogeneous (no-wash) assay with simple and fast workflow

Discover how the cost-effective THUNDER Phospho-ERK1/2 (T202/Y204) assay can be used to measure EGF-induced phosphorylation of ERK1/2 in HEK293 cells with the SpectraMax® iD5, i3x, and M5e Multi-Mode Microplate Readers and SoftMax® Pro Software to generate robust assay data with streamlined data analysis and graphing.

[Download Application Note](#)

High-throughput cellular screening solutions

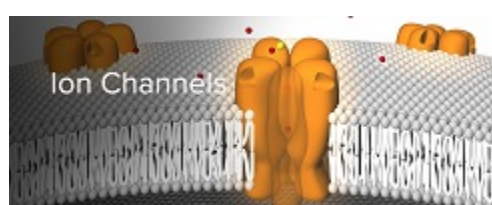
FLIPR® Penta High-Throughput Cellular Screening System is an industry-renowned instrument for monitoring GPCRs and ion channels. It supports a wide range of assay kits such as GPCRs using Calcium kits and Ion Channel assays using optogenetics, Membrane Potential, Calcium, and Potassium kits. Check out our latest research on:



Solutions for identifying early leads against GPCRs

Here we focus on applications to provide a high-throughput kinetic screening solution for toxicology and lead compound identification.

[Learn More](#)



High-throughput screening of ion channel in drug discovery

Discover solutions to support studies of ion channel function including assay kits, cellular screening and imaging systems, and microplate readers.

[Learn More](#)

EVENTS

LabDays Aarhus 2021

September 15-16, 2021 | Aarhus, Denmark

ELRIG DE Digital Forum - Virtual

September 16, 2021

Ecole thématique sur les organoïdes

September 13-17, 2021

19th Annual Discovery on Target Conference - Hybrid

September 27-30, 2021 | Boston, MA

SynBioBeta 2021

October 4-6, 2021 | Oakland, CA

SBI® (Society of Biomolecular Imaging and Informatics) 2021 - Virtual

October 5-6, 2021

TEDD Annual Meeting 2021

October 14 | Wädenswil, Switzerland

ILMAC

October 19-21 | Basel, Switzerland

SLAS 2021 Building Biology in 3D Symposium - Hybrid

October 26-27, 2021 | Jupiter, FL

Discovery UK

October 26-27 | London, UK

JOIN OUR TEAM

Interested in a career with Molecular Devices? Click [here](#) to see our latest openings.



[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. ©2021 Molecular Devices, LLC 3860 N First Street, San Jose, CA 95134 USA