

ADVANCING DISCOVERIES TOGETHER



Welcome to the August 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

NEW

Lab Automation Solutions

Explore lab automation-ready workflows for an end-to-end, high-throughput screening solution

No matter the scientific question or application at hand, an automated workflow can bring answers to light, faster. By integrating leading technology from across the industry into a flexible, customizable workcell, researchers can greatly optimize their time in the lab and free up resources for other critical tasks, speeding discovery.



- **Lab automation for high-content screening (HCS):** Our automated, high-content screening workcell provides an end-to-end solution that helps standardize live 2D/3D cellular development process with cell culture, treatment, and incubation, through to imaging, analysis, and data processing, delivering consistent, unbiased, and biologically-relevant results at scale.
- **Lab automation for high-throughput ELISA workflow:** Enzyme-linked Immunosorbent Assay is one of the most popular quantitative methods to detect a target antigen such as toxin or foreign substance within a sample. The assay is easy to set up and the range of potential analytes is vast, but the assay procedure is time consuming and labor intensive.
- **Lab automation for high-throughput clone screening:** Scalable and simple-to-use products establish clonal populations for antibody discovery and cell line development. Our automated clone screening workflows integrate laboratory devices to increase your throughput and efficiency while reducing human interaction.

[Explore Lab Automation](#)

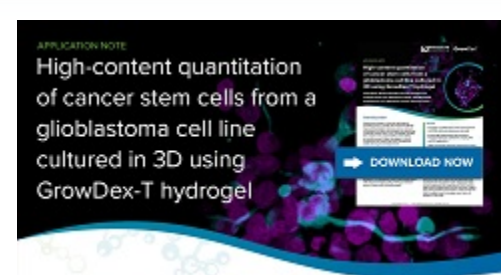


Application Spotlight

Cellular signaling and cellular responses

Many tools have been developed to measure cellular responses occurring through a wide range of signaling pathways. Here, we present a variety of cell signaling applications that can be run using microplate readers with fluorescence, TR-FRET, fluorescence polarization, and luminescence detection modes.

[Learn More](#)



Application Spotlight

Develop scalable, high-content 3D assays for cancer stem cells

Developing 3D high-content assays that are suitable for high-throughput screening requires careful consideration. Read how GrowDex-T[®] hydrogel provides well-defined support for cells that are easy to work with and provides a reliable means of assay development.

[Download Application Note](#)



Transform your cell imaging

Simplify your complex cell imaging and analysis with robust, easy-to-use workflows

Maintaining sample integrity during the course of your cell imaging experiments, while acquiring images with sufficient resolution to generate meaningful and reproducible data, is critical for revealing complex cellular and biological pathways. This eBook highlights how you can perform a wide variety of complex cellular imaging assays easily.

[Download eBook](#)

Product Spotlight

Advance your microbial and mammalian clone screening with proven, automated technologies

Check out our updated offerings to automate and expand your bioproduction processes. Our proven solutions provide dedicated, scalable, and easy-to-use products for establishing clonal populations.



Clone screening solutions
Automated, high-throughput clone screening solutions

[View Brochure](#)

EVENTS

SIMB (Society of Industrial Microbiology & Biotechnology) Annual Meeting
August 8-11, 2021 | Austin, TX

The 13th Annual Bioprocessing Summit - Hybrid
August 16-19, 2021 | Boston, MA

LabDays Aarhus 2021
September 15-16, 2021 | Aarhus, Denmark

ELRIG DE Digital Forum - Virtual
September 16, 2021

19th Annual Discovery on Target Conference - Hybrid
September 27-30, 2021 | Boston, MA

SynBioBeta 2021
October 4-6, 2021 | Oakland, CA

SB² (Society of Biomolecular Imaging and Informatics) 2021 - Virtual
October 5-6, 2021

SLAS 2021 Building Biology in 3D Symposium - Hybrid
October 26-27, 2021 | Jupiter, FL

JOIN OUR TEAM

Ready to reach your full potential? Click [here](#) to see what life is like at Molecular Devices.



[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 3860 N First Street, San Jose, CA 95134 USA