



NEW eBook Spotlight

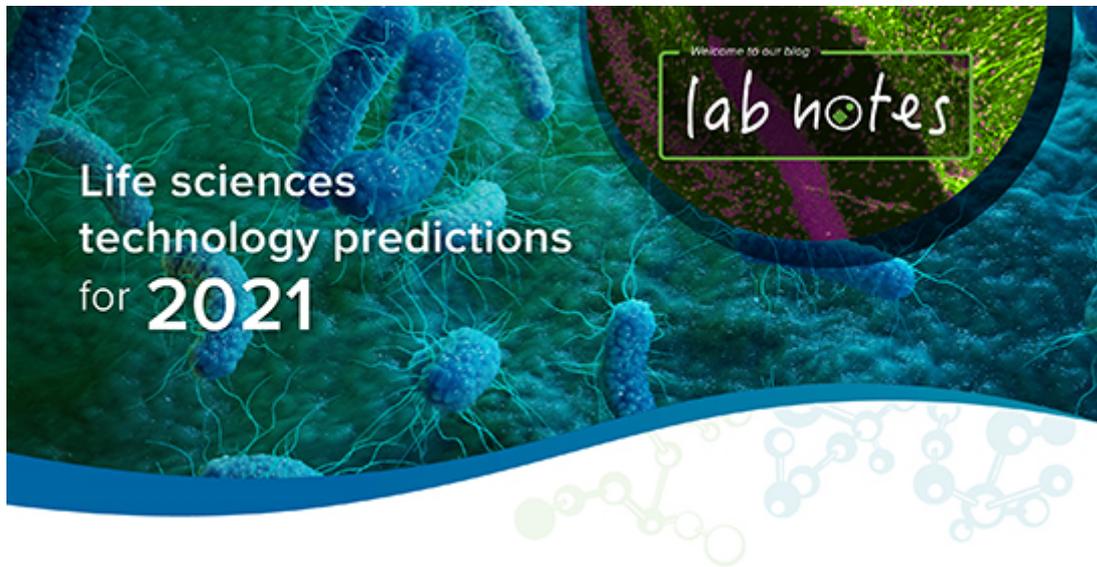
Accelerate vaccine development

Critical steps in the discovery and scale-up process can still be a challenge, but high quality monoclonal antibody producing cell lines are within your reach. In this recently published eBook, we offer 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.

Learn more about:

- Automated screening & selection of high-expressing cell lines
- Novel antibody internalization and binding assays
- Easy expression assays
- Objective verification of monoclonality

[Download eBook](#)



NEW
Lab Notes Blog

Life sciences technology predictions for 2021

In this latest blog post, our associates share their perspectives on where technology will take the life sciences industry in 2021 and beyond.

- Emergence of synthetic biology
- Increased collaborations between academia and industry
- Development of personalized medicine through gene editing and biomaterial-based drug delivery systems
- Adoption of artificial intelligence and automation platforms

Read Blog

NEW Application Spotlight

3D cell-based assays using a novel fluidics flowchip technology and high-content imaging

Here, we describe the use of high-content imaging in combination with the microfluidic-based Pu·MA System (Protein Fluidics) to perform automated assays, imaging, and image analysis with 3D cell models. We introduce a new method that's advancing the possibilities of drug discovery using microfluidics to automate 3D cell culture and cell analysis, allowing a hands-free process for media exchange and treatments while simplifying and improving workflow.

[Download Application Note](#)



NEW Application Spotlight

Determine Kd for the glucagon GLP-1 receptor with Tag-lite HTRF technology

Here, we show how the SpectraMax® i3x and SpectraMax® iD5 Multi-Mode Microplate Readers are used to perform robust, no-wash saturation binding assays using HTRF. GLP1R ligand binding was evaluated on the HTRF-certified SpectraMax readers using Tag-lite technology.

[Download Application Note](#)

NEW

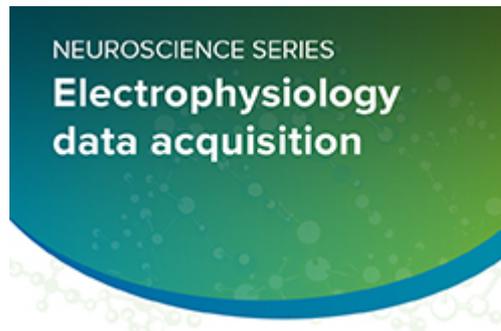
Customer Spotlight

Advancing understanding of Extracellular Vesicles (EVs)

Dr. Grégory Lavieu's team at the Université de Paris is interested in how EVs transfer their content from donor to acceptor cells, particularly the EV-delivery process. Their research could contribute to improved delivery of therapeutics. Find out how, and why, the team uses the SpectraMax® iD3 Multi-Mode Microplate Reader for sensitive luminescence and fluorescence assays.



[View Customer Story](#)



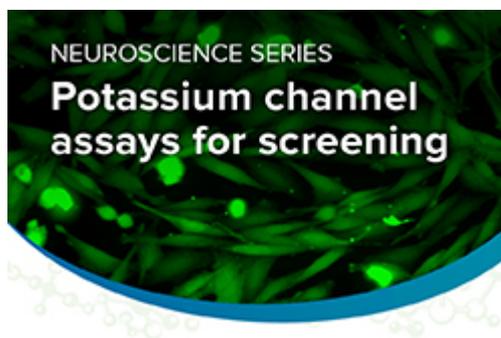
NEW

Videos Spotlight

Electrophysiology data acquisition

The Axon pCLAMP software suite provides users with greater capability in electrophysiology data acquisition and analysis. The features in Clampex 11 software enhance functionality in the Protocol Editor, increasing the user's ability to create complex experimental protocols involving many command stimulations.

[View Videos](#)



NEW

Application Spotlight

Potassium channel assays for screening

Analyze hERG liability earlier during the drug discovery process or screen for modulators of a hKCC2 cotransporter with the FLIPR® system and potassium assay kit.

[Download Application Note](#)



NEW

Video and eBook Spotlight

Eliminate line-frequency noise with HumSilencer technology

The HumSilencer™ feature within the Axon™ Digidata® system removes noise across a range of electrophysiological applications without distorting the data.

[Download Video & eBook](#)

EVENTS

[SfN Global Connectome 2021 - Virtual](#)

January 11-13, 2021 | North America

[SLAS 2021 - Virtual](#)

January 25-27, 2021 | North America

[LS² Annual e-Meeting - Virtual \(Life Sciences Switzerland\)](#)

February 17-19, 2021 | Europe Virtual

[The Biophysical Society \(BPS\) 2021 - Virtual](#)

February 22-26, 2021 | North America

[SOT Annual Meeting & ToxExpo 2021 - Virtual](#)

March 12-26, 2021 | North America

[RMS High-Content Imaging Workshop](#)

March 23-25, 2021 | UK Virtual

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