



Welcome to the February edition of our newsletter!

For 40 years, we've helped scientists harness the full potential of biology with next-generation technology. Here we share the latest in automated, end-to-end solutions that span research disciplines to advance scientific discovery and improve the quality of human life worldwide.

NEW DispenCell Single-Cell Dispenser
Compact, automated cell dispenser for fast, easy single-cell isolation

Molecular Devices partners with SEED Biosciences to exclusively offer DispenCell Single-Cell Dispenser

The DispenCell™ Single-Cell Dispenser is an automated laboratory instrument designed for fast, easy, and gentle single-cell isolation. DispenCell integrates seamlessly into your laboratory workflow, with a plug-and-play approach. Flexible and effortless, DispenCell operates equally under sterile conditions in a culture hood, or on a benchtop, producing rapid, traceable results.

[Learn More](#)

APPLICATION NOTE
Sensitive RNA fluorescent quantitation with the Quant-iT RiboGreen RNA assay kit

[DOWNLOAD NOW](#)

New Application Note

Rapidly and accurately quantify RNA concentration

Accurate quantification of nucleic acid concentration is important for downstream applications including transfection, cloning, PCR, and next-generation sequencing (NGS). Often, these applications have specific target nucleic acid concentrations for optimal performance. Inaccurate quantification can increase variability in downstream assays and affect the quality of results.

[Download Now](#)

ON DEMAND
Automate your 3D biology high-throughput workflows

[WATCH NOW](#)

Aditya Kalyas, PhD
Field Applications Scientist
Molecular Devices

Jay Hoying, PhD
Chief Scientist
Advanced Solutions Life Sciences

On-Demand Webinar

Automate your 3D biology workflows with 3D bioprinting and high-content imaging

Our recent webinar with Advanced Solutions is now available to view on-demand!

In this webinar experts address the reliability and scalability challenges associated with 3D biology workflows. In addition, they share real-world examples of how you can gain crucial insights about compound efficacy, biological complexity, and physiological relevance of your 3D assay results. Key takeaways include:

- How to incorporate automation into your 3D biology workflows to generate phenotypic data from complex models in a high-throughput setting
- How to leverage bioprinting for 3D biology applications such as spheroid and organoid formation, assay performance, and vascularized human tissue
- How to monitor cell responses in real time to increase throughput and reduce variability

[Watch Now](#)

lab notes

3D organoids and automation of complex cell assays [Podcast]

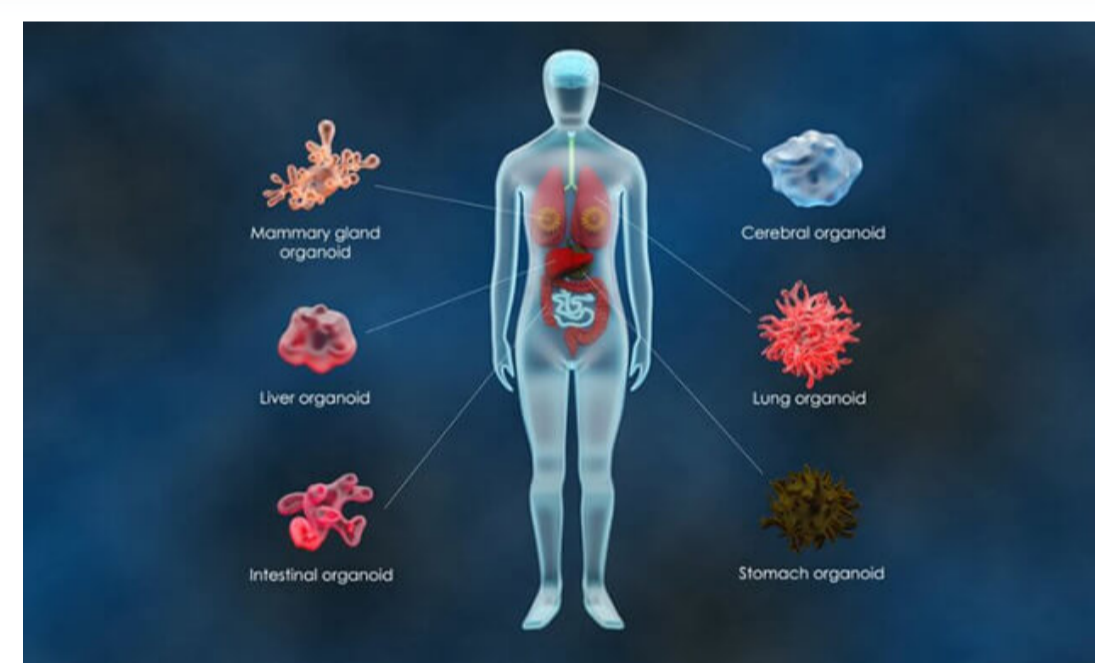
Podcast Excerpt

3D organoids and automation of complex cell assays

As we enter the era of sophisticated drug discovery with gene therapy and personalized medicine, we need to be prepared to study complex diseases, assess the therapeutic effect of drugs and identify adverse effects that can pose risks to patient health.

In this podcast excerpt, Senior scientist Oksana Sirenko discusses the advantages of 3D cell models while addressing challenges in 3D cell imaging, such as image quality, high throughput, automation, and analysis.

[Learn More](#)



Drug Discovery World Article

Gene editing in organoids: accounting for complexity in drug discovery

More researchers are using gene editing to build disease models that better represent human tissues' complex biology, signaling a shift away from 2D cell culture or animal models to organoids.

In this Drug Discovery World article, Dr. Oksana Sirenko explains how this trend has the potential to improve drug discovery.

[Read Now](#)

Imaging User Meeting 2023
Copenhagen | 9–10 May

Save the Date!

Join us in Copenhagen for our 2023 Imaging User Meeting!

This is an excellent opportunity to discover how your peers are utilizing cutting-edge imaging technologies to advance their own research, have the opportunity to share your own work, and meet our expert team of imaging scientists.

We're also pleased to be able to offer a tour of the Core Facility for High-Content CRISPR Screens at the University of Copenhagen's Biotech Research and Innovation Center.

[Learn More](#)

SCALE UP ELECTROPHYSIOLOGY
Fast track from patch-clamp to publication

[DOWNLOAD NOW](#)

New Brochure

Fast track from patch-clamp to publication with the Axon Instruments portfolio

Turn your idea into a publication with our comprehensive patch-clamping solutions, which enable the full range of patch-clamp technique experiments, from the smallest single-channel recordings to the largest macroscopic recordings.

The addition of Axon pCLAMP 11 Software Suite creates a streamlined workflow, allowing for sophisticated and efficient experiments, and higher quality data generation so you can make a scientific breakthrough even faster.

[Download Brochure](#)



Millions in FDA fines and thousands of warning letters:

How GxP Compliance Software can help avoid them

Since 2015, the number of warning letters issued by the FDA has doubled and more than 50 percent of observations impact data integrity, posing hefty fines of up to \$1 million per incident. Is your regulated lab audit ready?

[Learn More](#)

EVENTS

BPS (Biophysical Society)
February 18-22, 2023 | San Diego, CA

SLAS (Society for Laboratory Automation & Screening)
February 25 - March 1, 2023 | San Diego, CA

BPI West (BioProcess International - US West)
February 27 - March 2, 2023 | San Diego, CA

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. ©2022 Molecular Devices, LLC 3880 N First Street, San Jose, CA 95134 USA