



Luciferase Assays

NEW

COVID-19 Research

Luciferase assays to aid in your COVID-19 vaccine research

Luciferases are enzymes that use a substrate called luciferin, along with oxygen and ATP, in an energetic process that produces light—like the yellow glow of fireflies. The power of luciferase has been harnessed by scientists to devise reactions whose light output is used to monitor biological processes including gene expression, biomolecular binding, and cell viability.

- Measuring luciferase expression using the SpectraMax® Glo Steady-Luc™ Reporter Assay Kit
- Dual-Luciferase Reporter (DLR) Assay
- Monitor NF-κB activation with a sensitive dual reporter assay
- Detect dual luciferase expression
- Highly sensitive dual luciferase detection with the SpectraMax® DuoLuc™ Reporter Assay Kit

To view application notes related to COVID-19 research? Visit our COVID-19 research site.

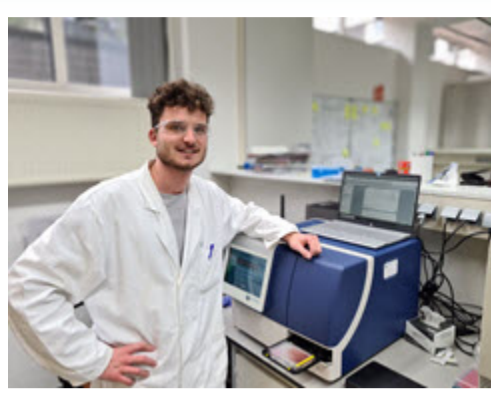
[Learn More](#)

NEW

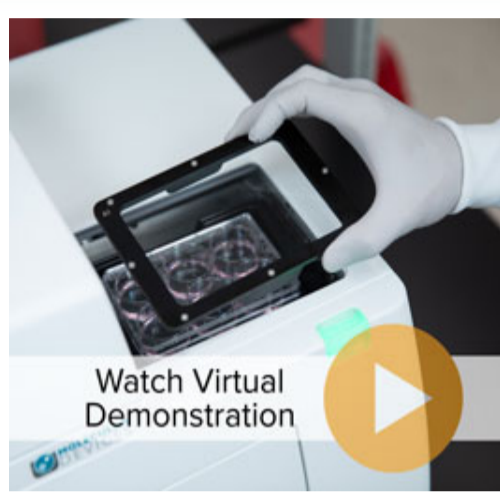
Customer Spotlight

Advancing research into life-changing mRNA therapeutics

Based in The Netherlands, the RiboPro team design and produce high-quality messenger RNA (mRNA) for academic and industrial researchers. Discover how they are using the SpectraMax® iD3 Multi-Mode Microplate Reader to advance research into life-changing mRNA therapeutics.



[Read Customer Story](#)



Watch Virtual Demonstration

NEW

Virtual Demo

ImageXpress Pico Environmental Control System

In this video, cellular imaging application scientist, Matthew Hammer demonstrates on the ImageXpress® Pico Automated Cell Imaging System, how to install and use the Environmental Control System, which is fully integrated to control temperature, humidity, CO₂ and O₂ allowing users to perform multi-day, live-cell, time-lapse experiments.

[Watch Virtual Demo](#)



NEW

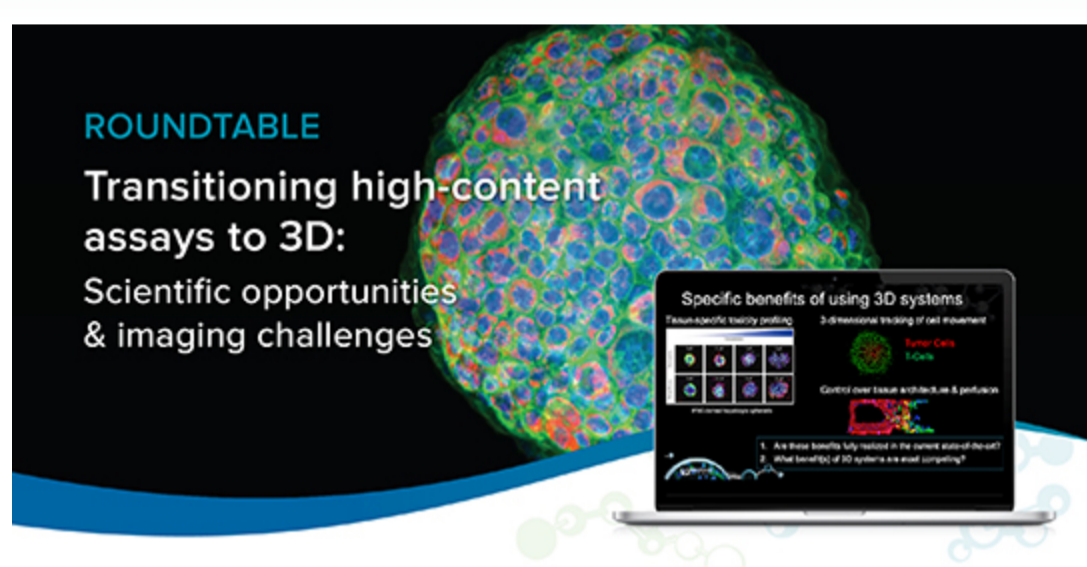
Application Spotlight

High-throughput screening of 3D cell cultures with multiple high density scaffold-free spheroids for cancer toxicity studies

Here, we demonstrate the use of the Corning® Elplasia® 96-well plates with a 3D culture workflow that includes spheroid generation, compound treatment, cytotoxicity assay, 3D imaging on the ImageXpress® Micro Confocal High-Content Imaging System, and 3D image analysis using MetaXpress® High-Content Image Acquisition and Analysis Software.

- Easily increase the number of spheroids per experimental condition
- Grow, stain, and image large numbers of spheroids simultaneously
- Use high-content imaging for simultaneous 3D analysis of multiple spheroids or organoids

[Download Application Note](#)



NEW

Roundtable Discussion

Transitioning high-content assays to 3D: Scientific opportunities and imaging challenges

In this roundtable, our panel of experts discuss the benefits and inherent challenges of making the transition from traditional 2D high-content assays to more complex 3D biology.

Key highlights:

- Benefits of using 3D cellular models for high-content imaging
- Main barriers to entry into using 3D cell models for high-content imaging and analysis
- Features that will be essential to accelerating high-content imaging of 3D biology

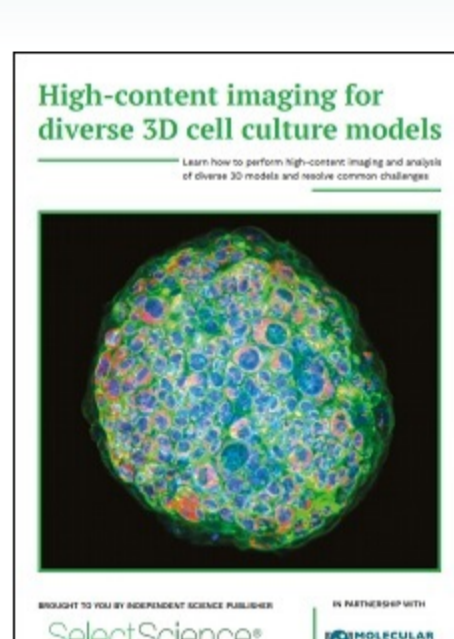
[View Roundtable](#)

NEW

eBook Spotlight

High-content imaging for diverse 3D cell culture models

In this recently published eBook with SelectScience, we present a collection of applications and case studies using the ImageXpress Micro Confocal system and MetaXpress software to investigate diverse 3D models and resolve common challenges experienced in 3D cell culture assays.



[Download eBook](#)



NEW

Video Spotlight

GxP compliance solutions for GMP/GLP labs

Our comprehensive suite of proven compliance solutions can advance your compliance journey, giving your team more time to focus on research.

- IQ/OQ/PM Services preserve microplate reader and washer documentation in a digital and compliant format
- Validation plates provide automated and traceable validation of reader performance
- SoftMax® Pro GxP Software helps you achieve full FDA 21 CFR Part 11 compliance
- Software installation and on-site software validation services support regulatory guidelines and are conducted by our certified specialists

[View Video](#)



NEW

Application Spotlight

Neurotoxicity assessment

The investigation of neuronal function and neurotoxicity has benefited from the development of assays that use a variety of physiologically relevant readouts: intracellular calcium flux, analysis of neuronal 3D co-cultures, and cell-based phenotypic assays.

[Download App Notes](#)



NEW

Application Spotlight

Neurite outgrowth

Neurite outgrowth is a commonly used assay to study neuronal development and neuronal degeneration *in vitro*. It is an efficient and effective high-throughput screening experiment suitable for evaluation of neurotoxicity across large libraries of chemical compounds.

[Download App Notes](#)

EVENTS

[Future Labs Live - Virtual](#)
November 17-18, 2020 | North America

[Cell Bio - Virtual](#)
December 2-16, 2020 | North America

[Antibody Engineering - Virtual](#)
December 13-17, 2020 | North America

[SfN Global Connectome 2021 - Virtual](#)
January 11-13, 2021 | North America

[SLAS 2021 - Virtual](#)
January 25-27, 2021 | North America

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. ©2020 Molecular Devices, LLC 3860 N First Street, San Jose, CA 95134 USA