



NEW
Complete compliance solutions for GMP/GLP labs

Molecular Devices is a leader in comprehensive compliance solutions with microplate detection systems and software. Combined with validation services and support, our solutions assure data integrity and compliance in GMP/GLP labs.

- IQ/OQ/PM services preserve instrument documentation in a digital and compliant format
- Software installation services verify and document that required components are installed to operational specifications
- Software validation service supports FDA 21 CFR Part 11 guidelines

[Learn More](#)



NEW
COVID-19 Spotlight

Applications and resources for COVID-19

See how our technology can help support your immunology and vaccine development research in our new COVID-19 applications resource hub. Learn about common applications used in infectious disease research:

- Cell count/cell viability
- Cell line development
- Protein/antigen binding affinity
- Protein/nucleic acid quantification
- Viral neutralization/ pathogenesis

[Learn more](#)

NEW
Cover Spotlight

Feature in Drug Target Review

Our Sr. Imaging Applications Scientist, Jayne Hesley, discusses how cellular imaging can be brought within the reach of every lab—from 3D high-throughput to simple automated imaging—in the latest edition of Drug Target Review. Check out the stunning image of co-cultured Normal Human Dermal Fibroblast (NHDF) and HUVEC cells, acquired using the ImageXpress® Micro Confocal system, as shown on the front cover.



[Read more](#)



NEW
Virtual Tour

ImageXpress Pico Automated Cell Imaging System

Watch our Biolmaging Specialist, Shawn Alvarado, PhD, as she takes you on a virtual tour of the ImageXpress® Pico system—a powerful, high-content imaging and analysis system which includes environmental control, objectives and filters, and powerful analysis tools.

[Take Virtual Tour](#)



NEW
On-demand Webinar

Exploring absorbance-based assay applications: from virus to cannabis research

Watch this webinar and learn:

- How absorbance detection works
- How to detect immunoglobulins (Ig) with ELISA—more ways to investigate infectious diseases and immune responses
- How to streamline analysis of beer and wine measuring absorbance
- How to determine total aflatoxin in cannabis using an ELISA

[View On-demand Webinar](#)



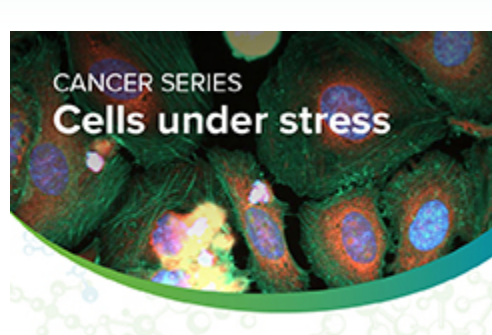
NEW
Application Spotlight

Cell migration

Cell migration is broadly defined as the movement of cells from one location to another. It is an essential process required for many biological events including embryonic development, wound healing, and immunological responses. The invasion of tumor cells into surrounding tissues, as well as metastasis, are areas of cancer that can be studied using *in vitro* cell migration methods.

Learn how to measure cellular migration over time and perform real-time analysis.

[Download Application Notes](#)



NEW
Application Spotlight

Cells under stress

Discovery and evaluation of anti-cancer therapies include development of cell-based models, screening for novel drugs, and understanding the relevant mechanisms of action. The cellular effects of processes like DNA damage and autophagy, a regulated process of degrading and recycling damaged proteins and organelles in response to cellular stress, can be analyzed efficiently using automated cell imaging.

Learn how automated cellular imaging provides an efficient method for analyzing the cellular effects of anti-cancer compounds.

[Download Application Notes](#)

EVENTS

AACR (Virtual)
June 22-24, 2020 | North America

ISSCR (Virtual)
June 23-27, 2020 | North America

FENS (Virtual)
July 11-15, 2020 | Europe

ECS2020
August 23-27, 2020 | Cork, Ireland

PEGS
August 31-September 4, 2020 | Boston, MA USA