

NEW PRODUCT
Discover More

Introducing the ImageExpress Confocal HT.ai High-Content Imaging System

The ImageExpress® Confocal HT.ai High-Content Imaging System combines premium hardware technology with a self-learning software analysis workflow for increased signal and ultra-fast 3D organoid screening.

Built on the success of our flagship model, the ImageExpress Confocal HT.ai system utilizes a seven-channel laser light source with eight imaging channels to enable highly multiplexed assays while maintaining high-throughput by using shortened exposure times. The powerful combination of MetaXpress® software and IN Carta™ software simplifies workflows for advanced phenotypic classification and 3D image analysis with machine learning capabilities and an intuitive user interface so that you can quickly uncover new insights that other technologies miss.

See the exponential power of the ImageExpress Confocal HT.ai system.

[Learn More](#)



NEW
Application Spotlight

Assessment of angiogenesis with automated time-lapse imaging

Here, we demonstrate the capability of the ImageExpress® Pico Automated Cell Imaging System to accurately capture the effects of pro-angiogenic and inhibitory compounds while maintaining optimal live cell environmental conditions.

Benefits:

- Visualize angiogenic network formation over time with live cell, time-lapse imaging
- Conduct quantitative assessment of HUVEC cells throughout the various stages of angiogenesis with preconfigured analysis modules
- Generate multiple readouts to evaluate angiogenic networks including tube length and area, as well as numbers of branches, segments, and connected sets

[Download Application Note](#)



NEW
Application Spotlight

Sensitive fluorescent quantitation of DNA with the Quant-iT Pico Green dsDNA Assay Kit

Here, we demonstrate how you can reliably measure concentrations as low as 50 pg/mL of double-stranded DNA (dsDNA) with our SpectraMax® Multi-Mode Microplate Readers and the Quant-iT Pico Green assay kit.

Key benefits include:

- Sensitive fluorescent quantitation of DNA down to 50 pg/mL
- Linear dynamic range spanning over four orders of magnitude
- Easy analysis of results with preconfigured protocols in SoftMax® Pro Software

[Download Application Note](#)



NEW
Lab Notes Blog

Get to know our Field Applications Scientist: Dwayne Carter

Dwayne Carter is a cell biologist and educator who joined Molecular Devices in November 2020. In this latest blog post, we asked Dwayne to tell us more about his background, career milestones, current role, and the future he envisions for clone screening.

[Read Blog](#)

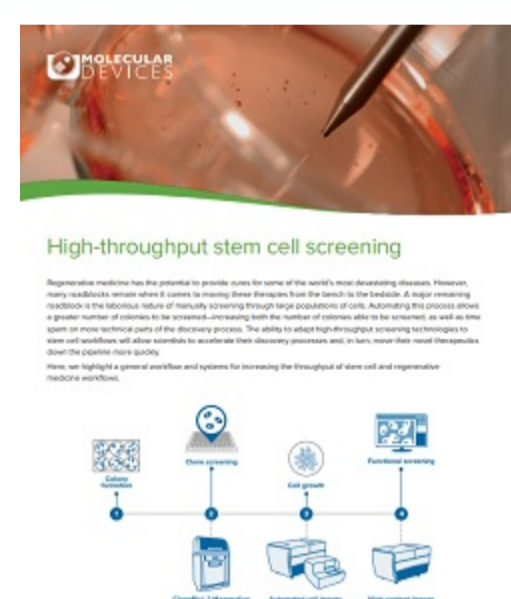
NEW
Customer Spotlight

A UK centre of excellence for the bioscience industry

The National Horizons Centre (NHC) in the UK has a major focus on training the next generation of scientists and biotechnologists. Discover how they use the SpectraMax® iD5 Multi-Mode Microplate Reader and SoftMax Pro Software to give students hands-on experience of biochemical and *in vitro* assays, and analysis, for challenging bioscience projects.



[Read Customer Story](#)



NEW
Product Spotlight

High-throughput stem cell screening workflow

Here, we highlight a general workflow and systems for increasing the throughput of stem cell and regenerative medicine workflows.

The ability to adapt high-throughput screening technologies to stem cell workflows will allow scientists to accelerate their discovery processes and, in turn, move their novel therapeutics down the pipeline more quickly.

[Download Flyer](#)

NEW
Product Spotlight

Build more, screen more, increase your colony picking capabilities 10X

Library screening is a critical step in many workflows, ranging from synthetic biology to CRISPR and gene editing. The ability to manufacture and easily engineer large libraries introduces a new bottleneck into the process—traditional colony screening and picking are slow, laborious, and time-consuming.

The QPix™ Microbial Colony Picker leverages best-in-class colony picking technology to alleviate this bottleneck to quickly, accurately, and efficiently screen through massive genetic libraries.



[Download Flyer](#)

EVENTS

LS⁺ (Life Sciences Switzerland) Annual e-Meeting - Virtual
February 17-19, 2021 | Switzerland

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

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. ©2021 Molecular Devices, LLC 3860 N First Street, San Jose, CA 95134 USA