



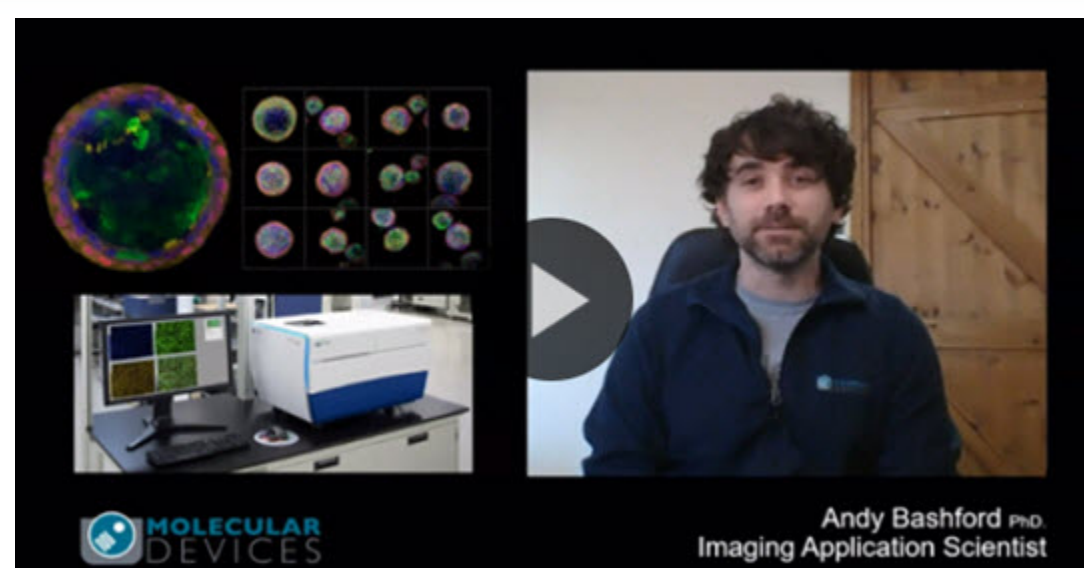
On-demand Webinar

Microplate analysis technologies for food and beverage quality testing

Absorbance microplate readers are widely used in research, drug discovery, quality control, and manufacturing processes in the pharmaceutical, biotech, food and beverage, and academic fields. They provide rapid and sensitive measurements across a wide range of concentrations for a variety of analytes that are key to assessing food or beverage quality and safety.

In this webinar, we show how SpectraMax® absorbance microplate readers and SoftMax® Pro Software are used to increase throughput and simplify analysis of data from assays used to test products including wines and beer.

[View On-demand Webinar](#)



Video Spotlight

Imaging and Analysis of a 3D Airway Organoid Model

In this 90 second video, Andy Bashford (Imaging Applications Scientist) shows a great example of imaging a 3D Airway Organoid Model using the ImageExpress® Micro Confocal High-Content Imaging System, along with some interesting ways to get the most out of this type of assay.

[View Video](#)

Application Spotlight

High-content phenotypic profiling using the Cell Painting assay

High-content phenotypic profiling is increasingly popular in research areas that span from gene function studies, drug discovery, and toxicology. The strength in this approach lies in the unbiased, multidimensional information captured at single cell resolution, which enables individual cell states to be analyzed, profiled, and compared to population-level data.



[Here](#), we show a high-content phenotypic profiling workflow based on the Cell Painting protocol by *Gustafsdottir et al**. Learn more about how this workflow utilizes easy-to-use tools without sacrificing quality.

[Download Application Note](#)



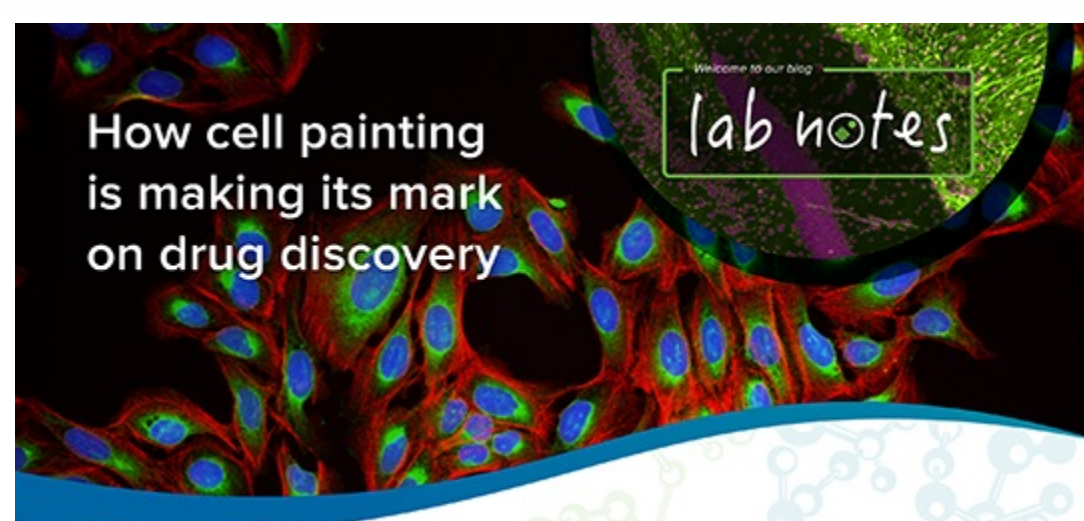
Whitepaper Spotlight

GxP regulated industry assessments of microplate readers

This document outlines references to 21 CFR Parts 58, 211 and 820 and EudraLex Annex 15 to assess the implementation of Molecular Devices microplate readers in regulated environments

- 21 CFR and EudraLex Vol. 4 overview
- Whose responsibility is it to validate the system?
- Impact of compliance vs. non-compliance

[Download Whitepaper](#)



Lab Notes Blog

How Cell Painting is making its mark on drug discovery

Cell Painting is a high-content, multiplexed image-based assay used for cytological profiling. The idea is to “paint” or stain as much of the cell as possible to get a representative image of the cell. Automated image analysis software is then used to extract hundreds to thousands of quantitative features of the cell, providing a rich phenotypic profile. Here, we shed some light on this emerging high-content image-based assay and the impact it is having on drug discovery.

[Read Blog](#)



Lab Notes Blog

Get to know our Field Applications Scientist: Cheryl Bell

In her role as Field Applications Scientist, Cheryl Bell provides scientific and sales support for our ImageExpress® Pico Automated Cell Imaging System. In this latest blog post, Cheryl tells us more about her career journey and shares some insights about making the transition from manual to automated microscopy.

[Read Blog](#)

EVENTS

[SOT Annual Meeting & ToxExpo 2021 - Virtual](#)
March 12-26, 2021 | North America

[High-Content Cellular Imaging User Meeting & Workshop - Virtual](#)
March 16-17, 2021 | Europe

[3D-Culture, Organoids and Organs-On-Chips 2021 - Hybrid](#)
March 22-23, 2021 | Boston, MA & Virtual

[RMS High-Content Imaging Workshop](#)
March 23-25, 2021 | UK

[AACR Annual Meeting 2021 - Virtual](#)
April 10-15, 2021 & May 17-21 | North America

[Organoid Discovery Symposium - Virtual](#)
April 13, 2021 | Europe

[PEGS Boston Virtual Conference & Expo - Virtual](#)
May 11-13, 2021 | North America

Join our team

Interested in a career with Molecular Devices? [Click here](#) to see our latest openings.