

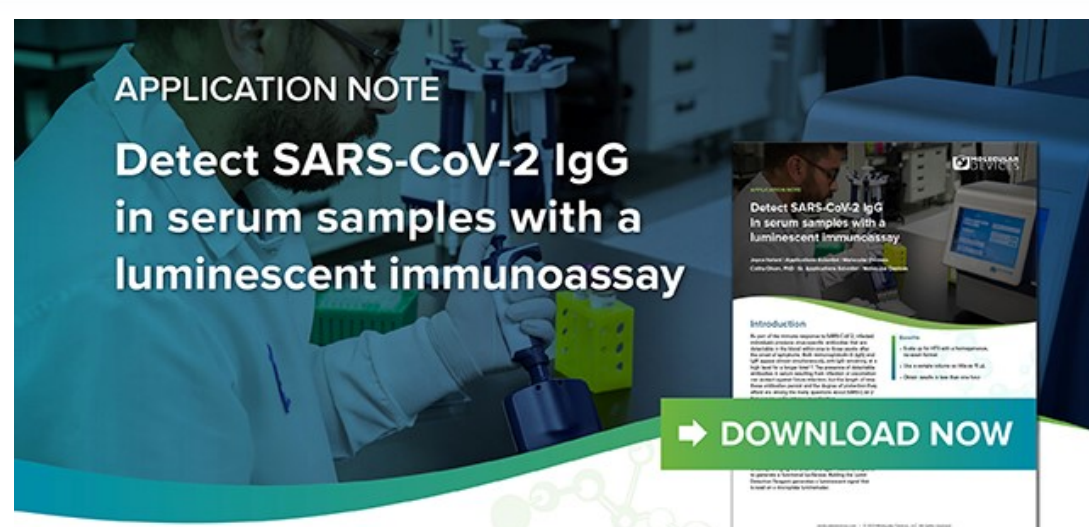
NEW

We proudly announce the opening of our Organoid Innovation Center

Modular, automated 3D cell culture and image analysis lab enables customers to streamline and scale organoid interrogation, advancing drug discovery research

The Organoid Innovation Center at our global headquarters expands beyond imaging to demonstrate an end-to-end solution that addresses challenges associated with every step in the sample prep-to-report pipeline for assays performed on complex 3D biological models. On display are state-of-the-art methods and technologies to test automated workflows for organoid culturing to deliver consistent, unbiased, biologically-relevant results at scale.

[Explore the Organoids Innovation Center](#)



Application Spotlight

Detect SARS-CoV-2 IgG in serum samples with a luminescent immunoassay

Discover how the Promega Lumit™ Dx SARS-CoV-2 Immunoassay, used in conjunction with the SpectraMax® iD5 Multi-Mode Microplate Reader and SoftMax® Pro Software, provides a simple workflow for faster results and enables higher throughput with less hands-on time.

- Scale up for HTS with a homogeneous, no-wash format
- Use a sample volume as little as 15 µL
- Obtain results in less than one hour

[Download Application Note](#)

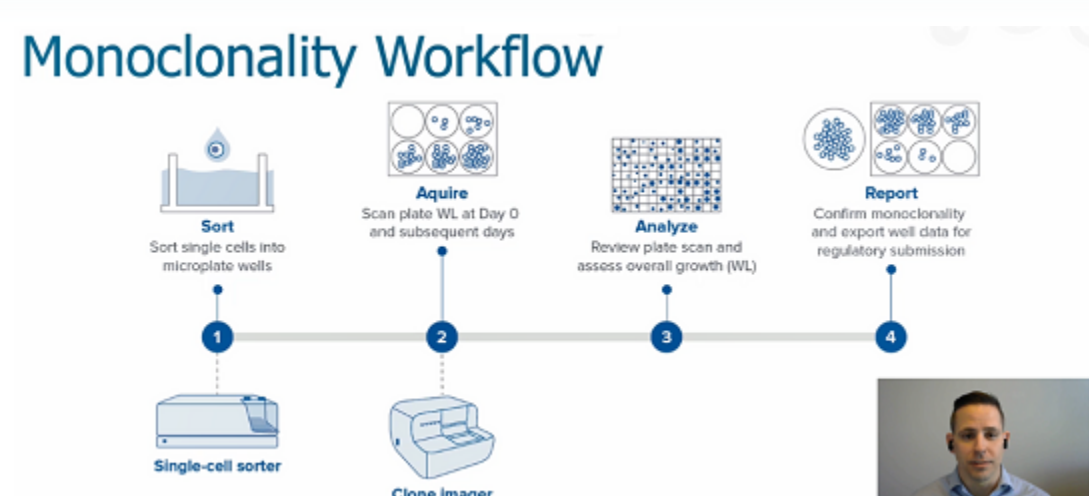


Application Page

Streamline beer, wine, and food quality control and safety

To ensure quality, authenticity and safety, manufacturers of foods and beverages need to perform a wide range of tests on their products. Many of these tests are based around ELISA, which are used in areas as diverse as checking for the presence of bacterial endotoxins, mycotoxins, various allergens, growth promoter hormones and antibiotics, and for meat species identification. Discover how you can streamline beer, wine, and food quality control and safety analyses on our new food and beverage page.

[View Application Page](#)



Video Spotlight

Monoclonality Workflow Overview

Achieve high assurance of monoclonality with immediate image-based evidence in a single workflow!

In this quick video, discover how our [monoclonality workflow solution](#) can help you confidently verify clonality with proven, automated technologies for a wide range of cell types.

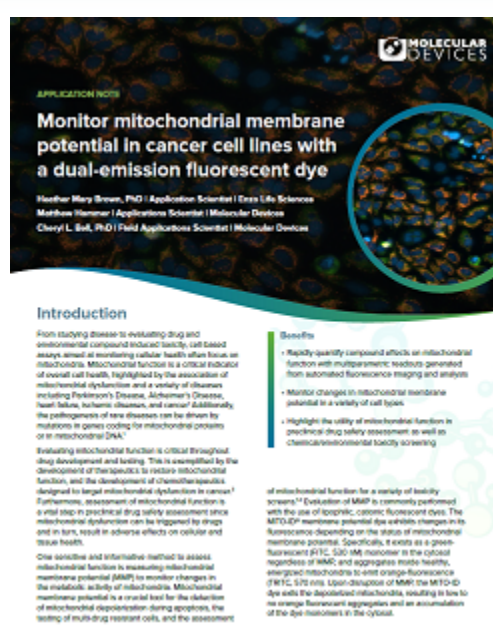
[Watch Video](#)

Application Spotlight

Monitor mitochondrial membrane potential changes using automated imaging

Assessment of mitochondrial function is a vital step in preclinical drug safety assessment since mitochondrial dysfunction can be triggered by drugs, and in turn result in adverse effects on cellular and tissue health.

We demonstrate the effective use of a highly sensitive cationic fluorescent dye (MITO-ID) to visualize and analyze mitochondrial membrane potential changes using automated imaging.



[Download Application Note](#)

EVENTS

PEGS Boston Virtual Conference & Expo
May 11-13, 2021 | North America

ISSCR 2021 Virtual Conference - Virtual
June 21-26, 2021 | North America

SLAS Europe 2021 Digital - Virtual Conference & Exhibition
June 23-25, 2021

3D-Culture, Organoids & Organ-on-a-Chip Europe 2021
June 28-30, 2021
Hybrid event - Rotterdam, The Netherlands

Join our team

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