



Exploring absorbance based assay applications: from virus to cannabis research

May 7, 2020
8:00 AM PDT | 11:00 AM EDT
4:00 PM BST | 5:00 PM CEST

NEW Webinar Spotlight

Exploring absorbance based assay applications: from virus to cannabis research

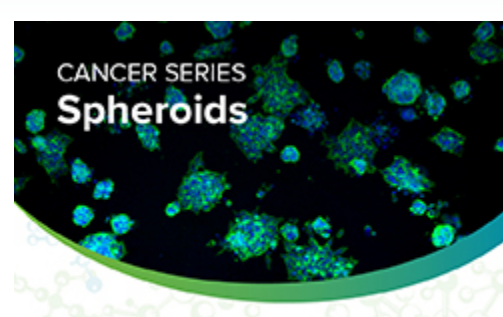
Microplate readers with absorbance detection capability are widely used in biological and chemical research, drug discovery, and quality control. There are many absorbance-based assay applications that can be utilized for research such as:

- Detecting immunoglobulins (Ig) with ELISA
- Streamlining analysis of beer and wine measuring absorbance
- Determining total aflatoxin in cannabis using an ELISA

Join our webinar to learn more:

Speaker: YoungMee Yoon, PhD, Field Applications Scientist, Molecular Devices
Date: May 7, 2020
Time: 8:00 AM PDT | 11:00 AM EDT | 4:00 PM BST | 5:00 PM CEST

[Register for Webinar](#)



NEW Application Spotlight

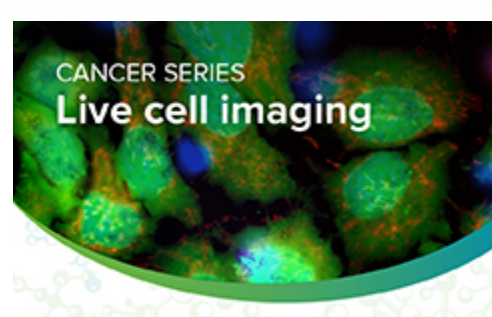
Spheroids in cancer research

Cancer is a group of related diseases where abnormal cells grow uncontrolled and can spread throughout the body. With over 200 different types, cancer is a leading cause of death worldwide.

The rapid progress of cancer research over the past few decades has seen the rise of spheroids—3D aggregates of cells in culture—as a valuable research tool that provides more physiological relevance than traditional 2D cell culture.

Learn how spheroids formed using a variety of cancer cell types are analyzed using methods featured in our application notes.

[Download Application Notes](#)



NEW Application Spotlight

Live cell imaging in cancer research

Live cell imaging is the study of cellular structure and function in living cells via microscopy. It enables the visualization and quantitation of dynamic cellular processes in real time.

Live cell imaging encompasses a broad range of biological applications, from long-term kinetic assays to fluorescently labeling live cells.

Learn how cellular processes are analyzed using methods featured in our application notes.

[Download Application Notes](#)



NEW On-demand Webinar

Automated Imaging and You—Quantitative microscopy for every lab, powerful data for all

In this webinar, after presenting the ImageXpress® Pico system and CellReporterXpress® software, we will introduce the benefits that automated microscopy and imaging can offer your laboratory, with reference to a number of case studies—from basic cell scoring and analysis to more complex applications such as neurite tracing.

[View On-demand Webinar](#)

NEW Application Spotlight

Optimized settings for Transcreener TR-FRET assays

Transcreener® HTS is a universal, high throughput biochemical assay platform based on the detection of nucleotides, which are formed by thousands of cellular enzymes—many of which catalyze the covalent regulatory reactions that are central to cell signaling and are of great value as targets in drug discovery.

Here we demonstrate the validation of SpectraMax® iD5 and i3x microplate readers for use with the Transcreener ADP2 TR-FRET Red Assay.



[Download Application Note](#)



NEW Customer Spotlight

University of Rouen

The Cell Imaging Platform of the University of Rouen in Haute-Normandie, France, chose the SpectraMax® iD3 microplate reader for ease-of-use in its teaching laboratories, and the FlexStation® 3 microplate reader for high-performance measurement of intracellular calcium mobilization studies.

[Read More](#)

NEW Customer Spotlight

Bioneer

Based in Denmark, Bioneer A/S is revolutionizing early drug discovery for immuno-oncology and neurodegenerative disease modeling with high-throughput imaging of 3D disease models using the ImageXpress® Micro Confocal High-Content Imaging System.



[Read More](#)

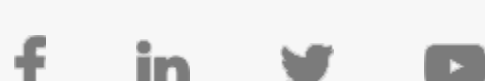
EVENTS

ISSCR (Virtual)
June 23-27, 2020 | Boston, MA USA

FENS (Virtual)
July 11-15, 2020 | Glasgow, United Kingdom

ECS2020
August 23-27, 2020 | Cork, Ireland

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



[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