

APPLICATION NOTE

Accelerated clone selection for recombinant CHO cells using CloneSelect Single-Cell Printer f. sight fluorescence-aided sorting

[DOWNLOAD NOW](#)

NEW
Application Spotlight

Discover an effective and rapid process for generating recombinant CHO cell lines

The selection of high-producing mammalian CHO cell lines continues to represent a major bottleneck in process development for the production of biopharmaceuticals. Therefore, it is increasingly important to develop new high-throughput methods for the selection of high-expressing CHO cell lines in an efficient and cost-effective manner.

In this application note, discover an effective and rapid process for generating recombinant CHO cell lines, producing high levels of therapeutic proteins.

- Identify desirable clones at an earlier cloning stage
- Assess therapeutic protein expression label-free
- Eliminate the use of specific antibodies for screening
- Easily implement into any cell line development process

[Download Application Note](#)

EBOOK

Streamline absorbance assays for nucleic acid & protein quantitation

[DOWNLOAD NOW](#)

eBook Spotlight

Streamline absorbance assays for nucleic acid & protein quantitation

Nucleic acid and protein quantitation are essential upstream measurements of many sophisticated assays in genetics and molecular biology.

In this eBook, learn how our absorbance solutions can help you save time and effort while easily generating the data you need.

Applications covered include:

- DNA and RNA quantitation
- Measuring total protein in cell lysates
- Low-volume, high-throughput DNA and protein detection
- ELISA assays to quantitate interleukin-8 concentrations

[Download eBook](#)



Protein Quantitation and Analysis

Microplate readers with absorbance, fluorescence, and luminescence detection modes provide a versatile solution for the quantitation and study of proteins.

[Learn More](#)



Nucleic Acid (DNA/RNA) Quantitation and Analysis

Our application notes demonstrate the quantitation and analysis of nucleic acids in a microplate format, offering higher throughput compared to other methods, as well as automated calculation of results.

[Learn More](#)

Confocal Laser Illumination
High-intensity laser light source for confocal microscopy

Technology Spotlight

Reduce exposure times by up to 75% using high-output laser excitation

Using the right high-content imaging solution allows you to penetrate into deep tissue while reducing exposure times and increasing focus.

By leveraging the illumination power of an 8-channel, high-intensity laser light source, and coupled with automated workflows, you can substantially decrease exposure times and increase your imaging speed and assay throughput. In addition, you benefit from:

- Sharper images with higher signal-to-noise ratio
- A nearly two-fold boost in scan speed due to significantly reduced exposure times
- Ability to run FRET experiments using lasers for CFP and YFP with increased multiplexing requirements

[Learn More](#)

NEW
Lab Notes Blog

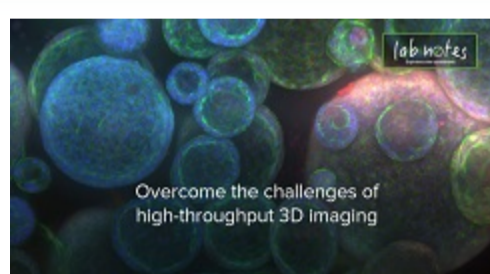
Overcome the challenges of high-throughput 3D imaging

High-throughput 3D imaging is an intricate process with many complexities, from long image scanning times to low resolution and inadequate analysis tools.

Here is a summary of the common challenges faced in 3D cell imaging and how you can overcome them with the help of Molecular Devices products.

- Out-of-focus light
- Long acquisition time
- Reliable automated focusing
- Analysis bottleneck
- Difficult light penetration

[Read Post](#)



NEW
Lab Notes Blog

FDA 21 CFR Part 11 and the importance of regulatory compliance in GMP and GLP labs

Here is a rundown of the essential components of regulatory compliance in GxP labs and our approach to achieving and maintaining these standards.

[Read Post](#)

EVENTS

[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

Ready to reach your full potential? [Click here](#) to see what life is like at Molecular Devices.