



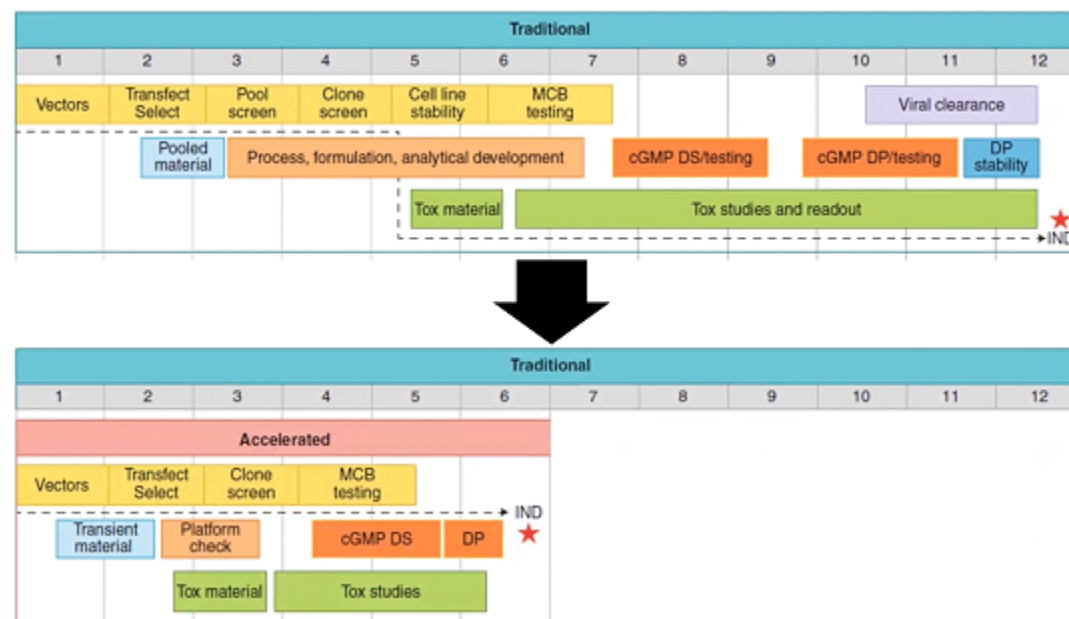
Welcome to the November edition of our newsletter!

Bringing you our latest discoveries in high-throughput screening, genomic and cellular analysis, colony selection, and microplate detection. Visit us at: moleculardevices.com

Accelerate monoclonal antibody discovery to combat COVID-19

This article reveals the steps to develop monoclonal antibodies using our clone screening instruments, from clonal selection to cell growth tracking to image-based assurance of monoclonality.

Join Rebecca Kreipke, Ph.D., as she helps scientists fast-track the FDA approval process and accelerate monoclonal antibody discovery to combat COVID-19 and other infectious diseases.



The time from discovery to proof-of-concept trials could be reduced to 5-6 months from a traditional timeline of 10-12 months.

[Read Blog Article](#)



Automated monoclonality assurance in human stem cell (hSC) engineering

Join Gargi Roy of AstraZeneca's Stem Cell Therapy team to explore high throughput human stem cell engineering workflows that automate monoclonality assurance and accelerate research timelines.

Learning Objectives

- Understand how the CloneSelect Single-Cell Printer* is an effective tool for single cell isolation of iPSCs and ESCs
- Explore how the combination of CloneSelect Single-Cell Printer and CloneSelect Imager offers confirmation of monoclonality assessments
- Discover how Clonal ESCs and iPSCs isolated by CloneSelect Single-Cell Printer maintain genomic stability and pluripotency for many generations and can be effectively programmed to differentiate into three germ layers and organ specific cell types

*Only available in North America, Mainland China, Hong Kong, Macau, and Taiwan

[Register for Webinar](#)

Rapidly quantify compound effects on mitochondrial functions of cancer cell lines

Assessment of mitochondrial function is a vital step in preclinical drug safety assessment since mitochondrial dysfunction can be triggered by certain drugs - resulting in adverse effects on cell and tissue health. Learn how the effective use of a highly sensitive cationic fluorescent dye enabled researchers to visualize and analyze mitochondrial membrane potential changes using automated imaging.



[Download the Research](#)



Use iPSC-derived cardiomyocytes for evaluation of compound effects and potential cardiotoxicity

CiPA initiative and validation of high-throughput methods to characterize compound effects in human iPSC-derived cardiomyocytes.

- Hear about the CiPA initiative
- Learn how to evaluate compound induced effects on hiPSC-derived cardiomyocytes
- See the features of the FLIPR® Penta system and ScreenWorks® Peak Pro 2 software

[Register for Webinar](#)

EVENTS

Society for Neuroscience (SfN) 50th Annual Meeting 2021

November 8-11, 2021 | Virtual

Cell Bio Virtual 2021

December 1-10, 2021 | Virtual

Antibody Engineering & Therapeutics 2021

December 12-16, 2021 | San Diego, CA

SynBioBeta: Global Synthetic Biology Conference 2022

January 25-27, 2021 | Oakland, CA

JOIN OUR TEAM

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