

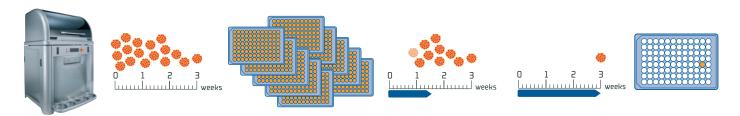
Accelerate cell line development

Get to your high value hits faster

Transform your workflows to elevate productivity

Fast screening and selection of secretory cell lines with the ClonePix[™] 2 Mammalian Colony Picker

Screen 10X more clones in weeks, not months. Productivity comparison: ClonePix: 10,000 clones in three weeks, Limiting dilution: 1000 clones in two months.

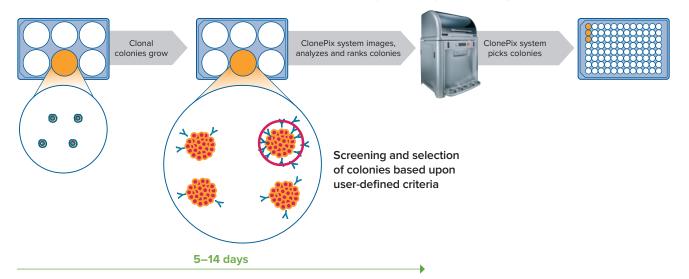


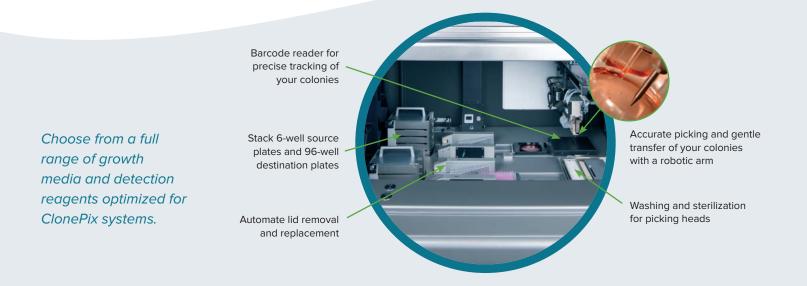
User selects colonies based on

system's analysis and ranking

Select and pick with more accuracy and confidence

Cells plated into semi-solid medium

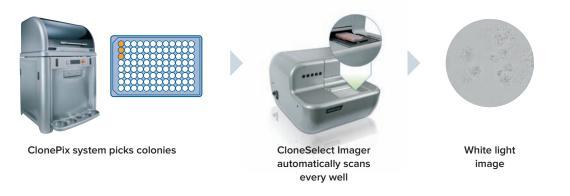




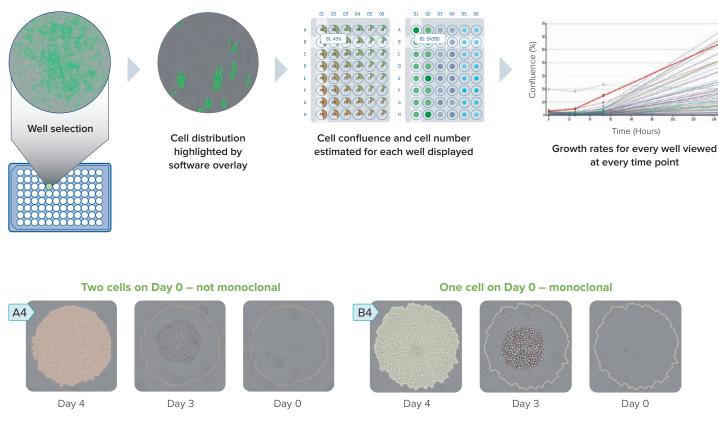
Advance to automated screening with intelligent imaging and quantitative analysis

Assessment of colony outgrowth with CloneSelect Imager[™] replaces subjective, time-consuming manual screening

Following screening and selection of clones, the ultra-fast CloneSelect Imager provides objective assessment of cell growth. Track and quantify cell growth over time automatically.



Analyze cell confluence and cell number in every well to verify monoclonality



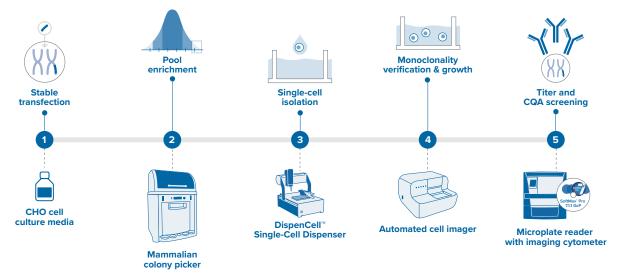
Verification of monoclonality - viewing the origin of a colony

The growth (image) history of each well can be tracked back to its starting point-providing evidence of monoclonality.

Develop stable cell lines

Typical cell line workflow

Cell line development is the process of establishing a clonally-derived cell population which has been genetically engineered to express a desirable phenotype (such as producing large amounts of recombinant protein) for a stable period of time. Single cells proliferate to form colonies that can then be assessed for the desirable characteristic.



Accelerate cell line development with a range of Molecular Devices platforms



ClonePix 2 Mammalian Colony Picker

Automatically screen more clones in less time than conventional techniques, select cells with optimal expression levels, and pick colonies with accuracy. ClonePix systems are now used in over 100 laboratories around the world to increase workflow productivity, leaving more time to better characterize target proteins and run new projects.



DispenCell[™] Single-Cell Dispenser

The DispenCell[™] Single-Cell Dispenser is an automated laboratory instrument developed for fast, easy and gentle single-cell isolation. Designed by scientists for scientists, DispenCell has been designed to integrate seamlessly into your lab routine, with a plug-and-play approach.



CloneSelect Imager – standard and fluorescence

With high quality imaging and intelligent image analysis, CloneSelect Imager allows you to assess cell confluence objectively and quantitatively. Cell growth is viewed and tracked in every well in every plate.

Phone: +1.800.635.5577 Web: www.moleculardevices.com Email: info@moldev.com Check our website for a current listing of worldwide distributors.

USA and Canada +1.800.635.5577 United Kingdom +44.118.944.8000 00800.665.32860 Europe* China +86.4008203586

Taiwan/Hong Kong +886.2.2656.7585 Japan South Korea India

+81.3.6362.9109 +82.2.3471.9531 +91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners. DispenCell is a trademark of SEED Biosciences SA. Specifications subject to change without notice. Patents: www.moleculardevices.com/productpatents FOR RESEARCH USE ONLY, NOT FOR USE IN DIAGNOSTIC PROCEDURES.

Regional Offices

©2023 Molecular Devices, LLC 2/23 1966D Printed in USA