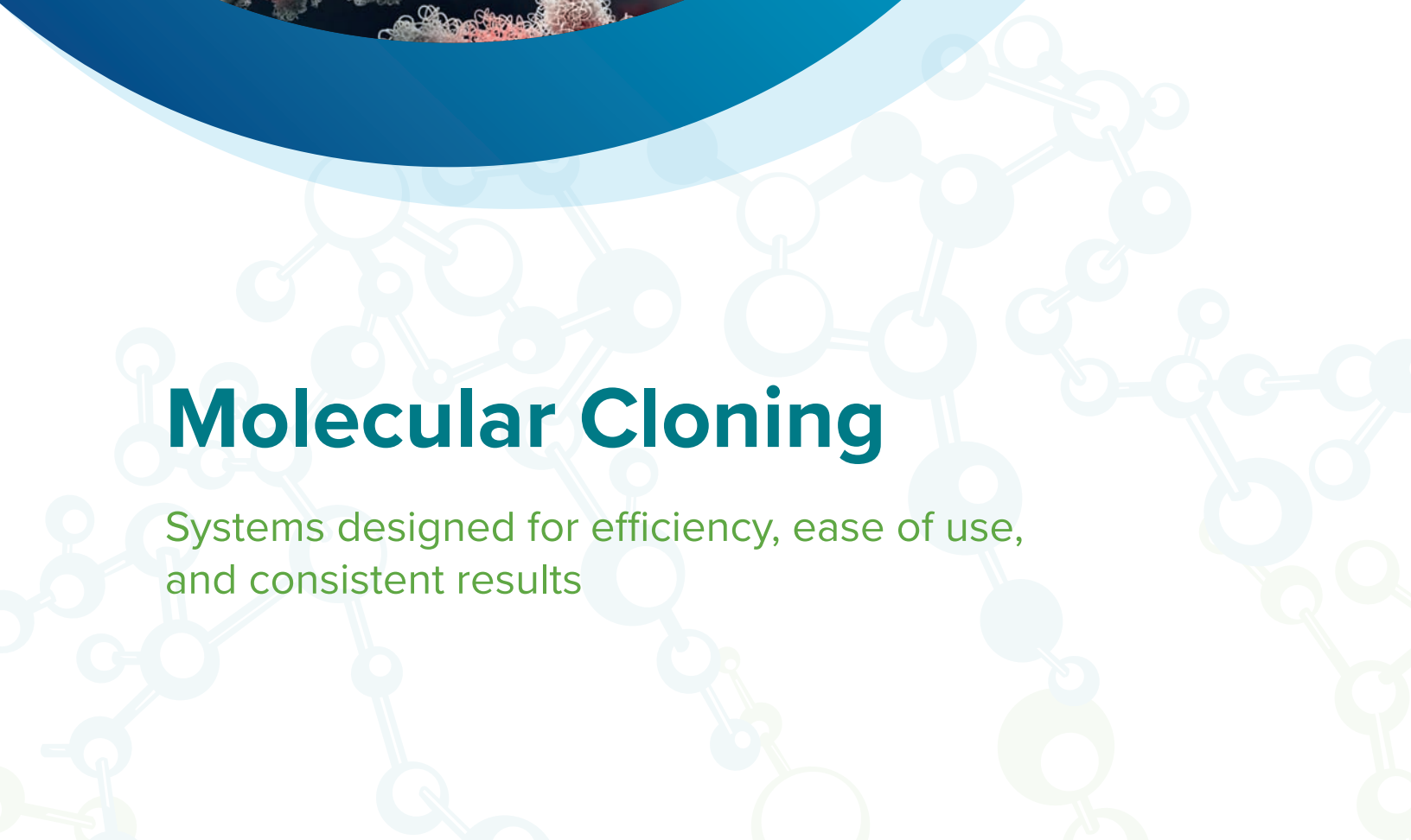
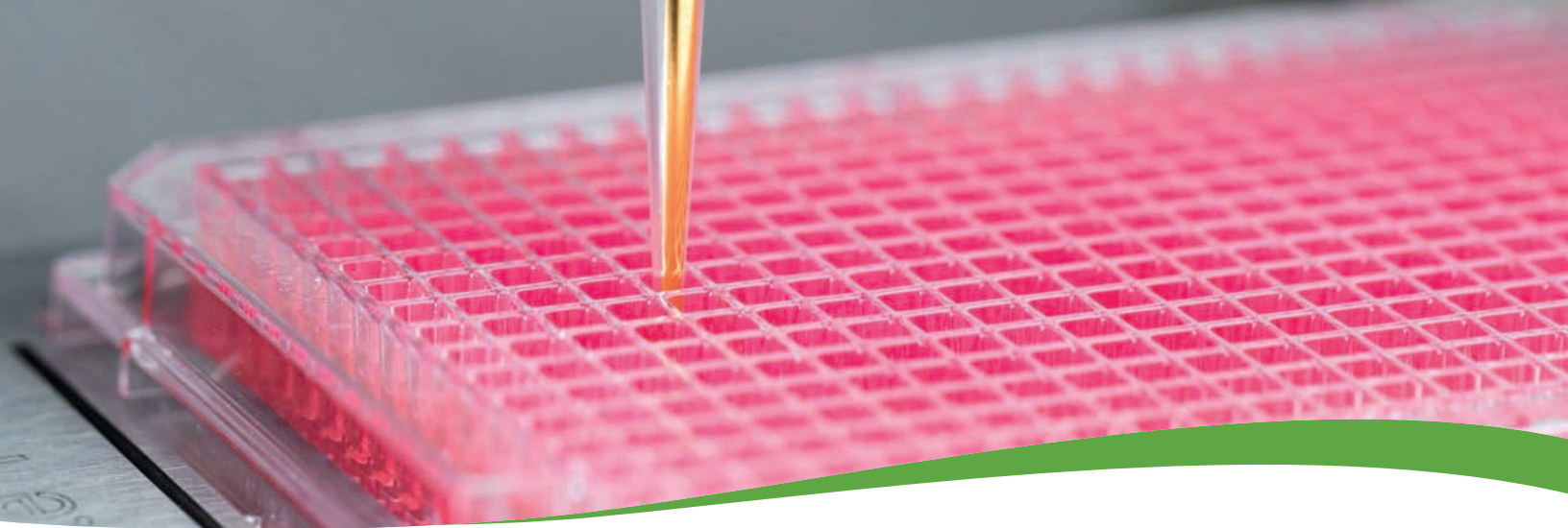


Molecular Cloning

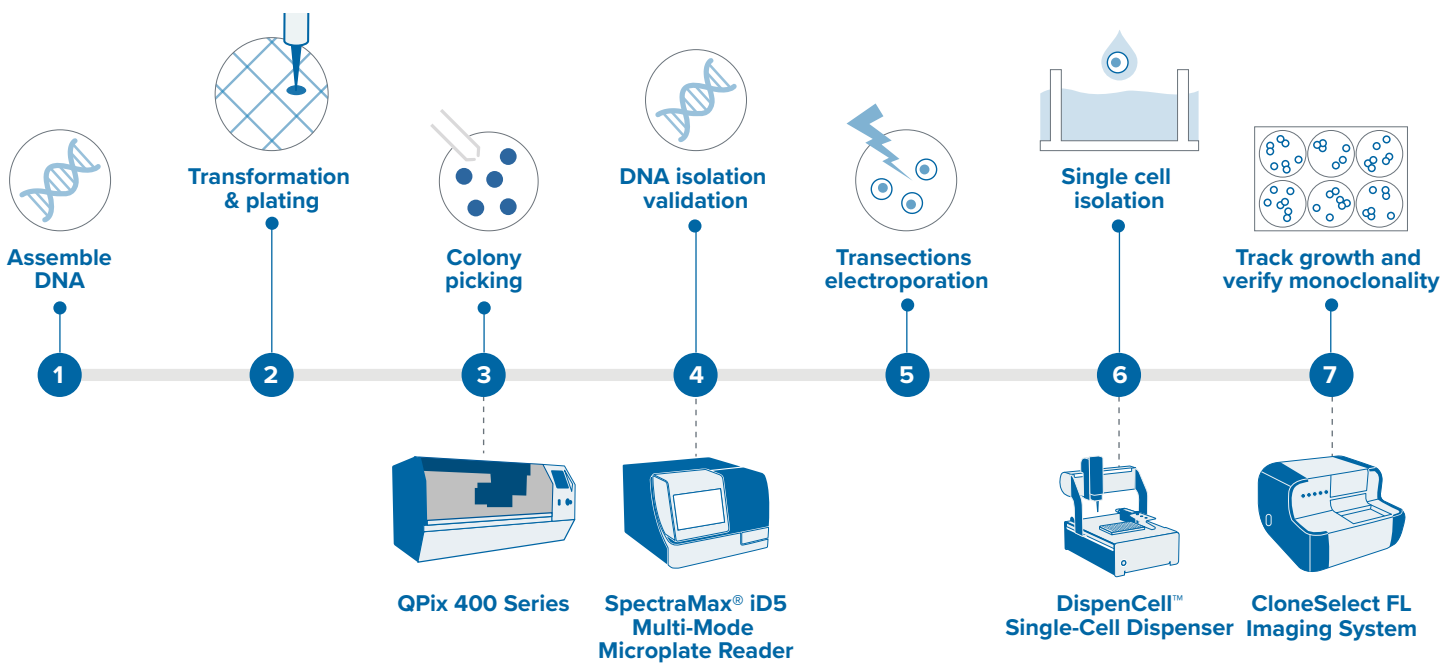
Systems designed for efficiency, ease of use,
and consistent results



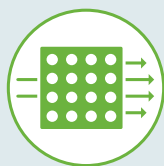


Introduction

Our solutions for antibody discovery and cell line development provide dedicated, scalable, and easy-to-use products for establishing clonal populations. The systems feature a selection of options and models to address your specific research including multiple imaging modes, biology-specific pins, fluidics, and environmental control. These solutions combine intelligent imaging with analysis and automation to increase throughput, consistency, and provide image-based documentation.



Verify monoclonality easily



Sort viable single cells efficiently



Streamline workflows



Improved probability of finding rare clones of interest

QPix® 400 Series Microbial Colony Pickers

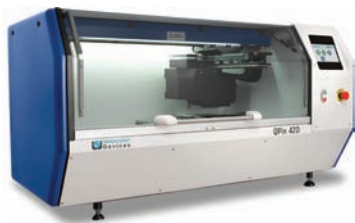
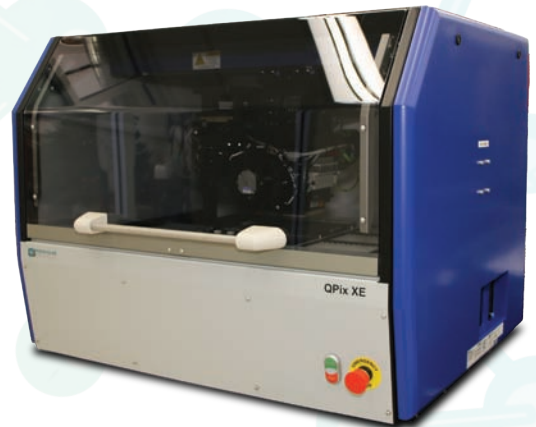
Featuring the all-new QPix XE

The QPix XE is a compact microbial colony picker ideal for labs with reduced space, but need efficient colony picking through put. The QPix XE is a great entry level system to replace highly subjective and error prone manual picking with automated colony picking.

Automated microbial clone screening and library management system

The QPix 400 Series Microbial Colony Pickers combine intelligent image analysis with precise automation for fast and efficient screening of large libraries. With a variety of data tracking and assay tools, the QPix Software streamlines the control and management of complex and iterative processes.

- Use the QPix system for synthetic biology, DNA assembly, antibody discovery, protein engineering, and phage display workflows
- Streamline your workflow with scalable-automation – pick up to 30,000 colonies per day
- Electronic data tracking for well-documented data control



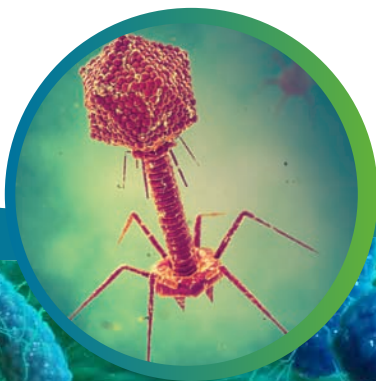
QPix 420 system



QPix 450/460 systems



QPix HT system



SpectraMax® iD5 Multi-Mode Microplate Readers

Microplate readers are commonly used to measure the quantity and quality of DNA during DNA isolation and purification protocols. By using a microplate reader to measure DNA concentration and purity, scientists can ensure that their DNA isolation and purification protocols are working correctly, and that they are obtaining high-quality DNA for downstream applications.



To use a microplate reader for DNA isolation validation, the following steps can be taken:

1. First, the DNA samples are prepared and added to a UV transparent 96-well plate in appropriate dilutions or concentrations.
2. Next, the microplate reader is programmed to measure the absorbance of the DNA samples at specific wavelengths, usually at 260 nm and 280 nm. The absorbance values at these wavelengths can be used to calculate the concentration and purity of the DNA.
3. The software in the microplate reader can then be used to generate a standard curve to convert the absorbance values to DNA concentration.
4. Once the standard curve is generated, the absorbance values of the DNA samples can be measured and the DNA concentration can be calculated using the standard curve.
5. The 260/280 absorbance ratio is also calculated to determine the purity of the DNA samples. A ratio of 1.8 is considered to be indicative of pure DNA.
6. Plate reader also can be used to run spectrum scan (230–320nm). This will show if there are any contaminants.



DispenCell™ Single-Cell Dispenser

Automation for fast, easy and gentle single-cell isolation

DispenCell integrates seamlessly into your laboratory workflow, with a plug-and-play approach. Flexible and effortless, DispenCell operates equally under sterile conditions in a culture hood, or on a simple benchtop.

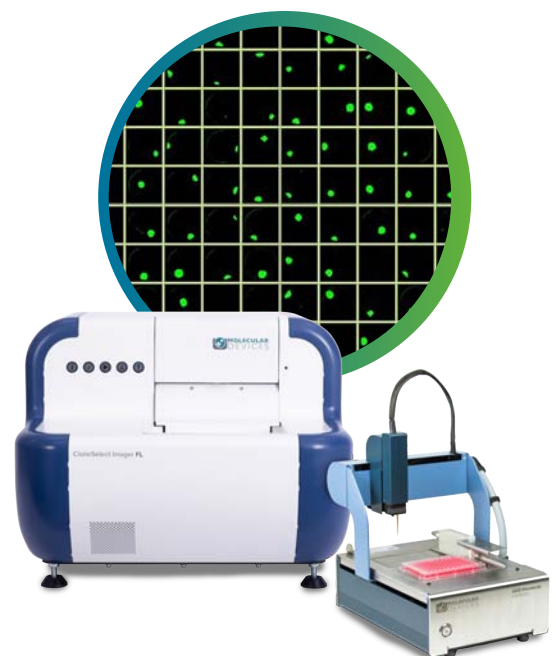
- Confidently isolate single cells and provide consistent, high-quality evidence of monoclonality to the FDA
- Highly efficient single-cell sorting with an average monoclonality reliability of > 90% observed
- Extremely gentle handling of the cell sample (less than 0.1 si), while preserving cell viability and outgrowth
- Minimize cross contamination through disposable cell dispensing cartridges
- Fast turnaround time—prints a 96-well plate in less than 10 minutes (384-well plate in less than 40 minutes)
- Sort wide range of cell lines: CHO, HEK, SF9, iPSC, primary cells
- Compact footprint takes up a small amount of valuable bench space

Verify monoclonality confidently at day zero with a CloneSelect Imager FL and DispenCell bundle

Streamline your workflow when you pair the DispenCell with the CloneSelect Imager FL

The all-new CloneSelect Imager FL adds high contrast multichannel fluorescent technology in addition to the standard white light imaging that allows for accurate single-cell detection and proof of monoclonality at day 0. Identify and verify gene edits with comparative confluence assays.

- Document evidence of single cells and confluency digitally for auditing and submission to regulatory authorities
- Image cells non invasively at multiple time points to monitor colony formation
- Screen using high resolution white light imaging
- Deliver real-time results with on-the-fly analysis
- Automation and integration ready





Get started. Explore single-cell dispensing with us.

Contact Us

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

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

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