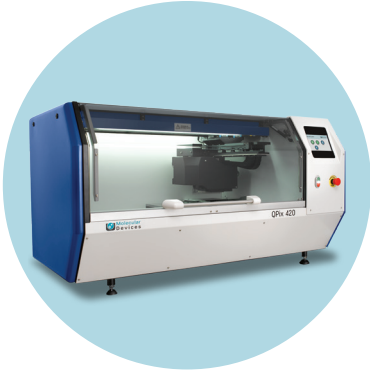
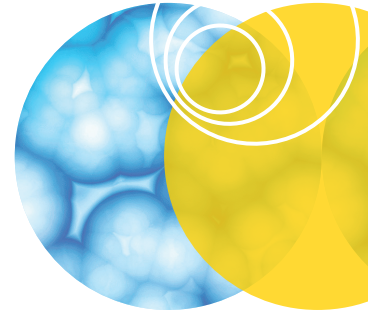


Molecular Devices

Advancing your cell line development workflow through automated and proven technologies



QPix™ 400 Systems

Empowering you to select the best colonies faster

QPix™ systems earned a well-deserved reputation for performance and reliability at major sequencing centers throughout the Human Genome Project. Get ready to discover why research institutes, sequencing facilities, biotech and pharmaceutical companies worldwide use over 600 QPix 400 systems every day.

Key features

- Scalable automation to suit your throughput needs
- True walk-away operation with easy setup
- Electronic data tracking for well-documented data control
- Sterile environment with customizable HEPA filtration options



ClonePix 2

Transform development workflows: screen more clones in less time

Cut cell line and antibody development times. Screen more clones in less time with the ClonePix system's fully automated workflow. Our users see major improvements both in productivity and overall costs compared to conventional techniques. Now you can reduce your timelines for monoclonal antibody generation by screening 10x more clones in weeks, not months!

Key features

- Screen more clones in less time
- Accurate, automatic colony picking avoids errors associated with limiting dilution
- Excellent image quality allows for screening of stable, high-producing clones
- Increased productivity of a cell line development workflow



CloneSelect Single-Cell Printer

Demonstrate monoclonality with greater confidence

The CloneSelect™ Single-Cell Printer™ (SCP) by Cytena and Molecular Devices is a fully automated system that utilizes proprietary microfluidics-based technology and real-time image analysis to sort and deposit single cells into standard microplates—while simultaneously providing assurance of monoclonality through image documentation.

Key features

- Confidently isolate single cells and provide consistent, high-quality image evidence of monoclonality to the FDA
- Highly efficient single-cell sorting with an average printing efficiency >80%
- Achieve an average of 75% cell viability with gentle handling of cells — comparable to limiting dilution method
- Minimize cross contamination through disposable cell printing cartridges
- Fast turnaround time—prints a 96-well plate within 5-10 minutes



CloneSelect Imager FL

Verify monoclonality confidently

Single cell identification with the aid of fluorescence offers a more conclusive method to meet regulatory demands in cell line development. The CloneSelect™ Imager FL can help you meet those demands with its automated analysis of cells in both white light and fluorescence, and can be used for concurrent confluency and monoclonality studies.

Key features

- 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 & fluorescence imaging
- Deliver real-time results with on-the-fly analysis
- Automation and integration ready

Cell Biology Reagents



XP Media and CloneMedia for Mouse Hybridoma Generation

The XP Media and CloneMedia for Mouse Hybridoma Generation is a complete solution that supports all stages of hybridoma cell line development from fusion to scale up.



CHO Growth A

Complete semi-solid and liquid CHO media optimized for increased antibody production.



CloneMedia

Ensures formation of discrete, clonal colonies.



CloneMatrix

Ensures formation of discrete, clonal colonies.



Recombinant CloneDetect

Recombinant CloneDetect is a unique, anti-human IgG (Fc-region) detection agent.



CloneDetect

Detection agents for visualizing secreted antibodies. All products are animal-derived, sterile, and azide free.



Cleaning And Sterilizing Solutions

Ideally suited to clean instruments and accessories.



Calibeads

Fluorescent microspheres for the verification of imaging and picking functions on ClonePix systems.

Contact Us

Phone: [800.635.5577](tel:800.635.5577)
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing of worldwide distributors