

5 TIPS to Improve Your Molecular Cloning Process with Synthetic Biology Automation

1 EMBRACE AUTOMATED LIQUID HANDLING

BENEFIT Accurate & reproducible pipetting.
IMPACT Minimize errors & increase throughput.

2 UTILIZE ROBOTIC WORKSTATIONS FOR EFFICIENCY

BENEFIT Increased overall efficiency.
IMPACT Perform multiple tasks, reduce errors, & increase throughput.

3 OPT FOR HIGH-QUALITY AUTOMATED IMAGING SYSTEMS

BENEFIT Fast & reliable colony picking.
IMPACT Precise identification, elimination of false positives, & accelerated screening process.

4 IMPLEMENT AUTOMATED DATA ANALYSIS & MANAGEMENT

BENEFIT Streamlined workflow & enhanced collaboration.
IMPACT Eliminate errors, enhance accuracy, & improve reporting.

5 EMPHASIZE CONNECTIVITY AND INTEGRATION

BENEFIT Streamlined workflow & improved productivity.
IMPACT Reduced manual interventions, cohesive workflow, & increased efficiency.

Automate your synthetic biology processes with us.
 Visit moleculardevices.com/syntheticbiology

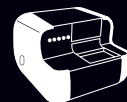
Accelerate your research with our products.



QPix



DispenCell™



CloneSelect™
Imager FL



ClonePix 2

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