

Build more, screen more, increase your colony picking capabilities 10x!

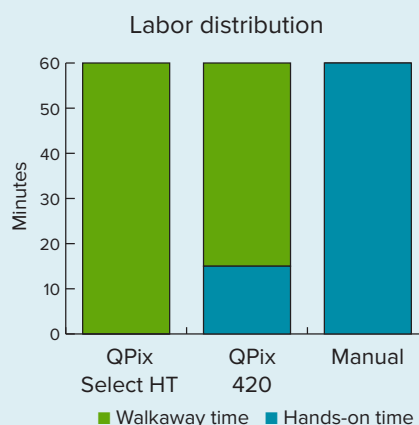
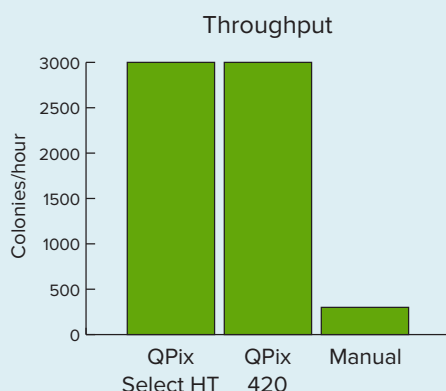
Scale is becoming an increasingly important factor in biology. The ability to scale up a process increases the chances that one is able to more readily and efficiently identify better hits earlier in the process.

Library screening is a critical step in many workflows, ranging from synthetic biology to CRISPR and gene editing. With improvements in library synthesis, it is now possible to easily and inexpensively manufacture extremely large and diverse genetic libraries. The increased quality of these genetic libraries, in theory, yields more ideal variants as researchers are able to begin with access to a larger starting pool of possibilities.

The ability to manufacture and easily engineer large libraries introduce a new bottleneck into the process—traditional colony screening and picking is slow, laborious, and time-consuming. This leaves groups with two choices:

hire additional staff for mindless colony picking sessions, or turn to proven technologies to increase the accuracy and efficiency of repetitive tasks, freeing scientists to analyze data.

The QPix™ Microbial Colony Picker leverages best-in-class colony picking technology to alleviate this bottleneck to quickly, accurately, and efficiently screen through massive genetic libraries. The easy-to-use, intuitive software guides users through setting up colony picking runs where precision robotics pick the right colonies every time. Data is automatically recorded into the machine's database, providing users with a complete audit trail and sample tracking, ensuring that no data is ever lost. Our modular, scalable series of colony pickers allows groups of all sizes to increase the accuracy and throughput of their workflow, while still allowing for future throughput growth.





QPix Select HT system

For groups looking to maximize the throughput and efficiency of their colony screening and picking process, the QPix Select HT system combines the precision and accuracy of the QPix colony picker with state-of-the-art laboratory integration that can be run 24/7 without manual intervention or supervision.

Throughput

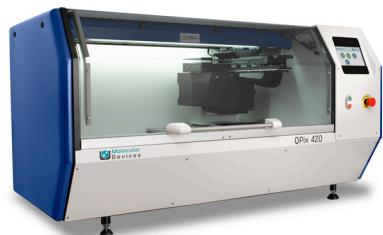
3,000 colonies/hour

Hands-on time

90 seconds to set up

Walkaway time

Entire duration of run



QPix 420 system

The QPix 420 system is a great entry level unit for groups looking to increase the throughput and efficiency of their colony screening. With the ability to pick up to 12 96-well plates before needing manual intervention, this will increase your walkaway time greatly. Additional plating and streaking and library management features with complete audit trail and sample tracking allows it to become an entire library management system—taking the guesswork out of sample tracking.

Throughput

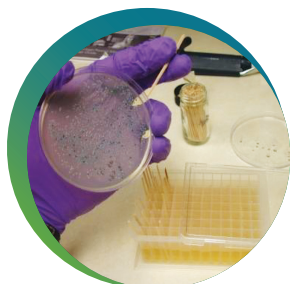
3,000 colonies/hour

Hands-on time

90 seconds to set up

Walkaway time

25 minutes at a time – only return to swap destination plates after 12 are full



Manual picking

Highly subjective and prone to human error, manual picking is a slow, laborious, and time-consuming process.

Throughput

300 colonies/hour

Hands-on time

60 minutes

Walkaway time

0 minutes

To learn more, please contact your regional Molecular Devices technical sales specialist, visit www.moleculardevices.com or call +1.800.635.5577.

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	China (Beijing)	+86.10.6410.8669	Japan	+81.3.6362.9109
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland