

CUSTOM SOLUTIONS

QPix 400 Series Colony Pickers



ACCELERATE *YOUR* SUCCESS

Achieve your scientific, operational, and financial goals with a QPix™ Colony Picker perfectly tailored to your specific workflow

When you can lay claim to the world's first and most respected colony-picking instrument, how do you set about making it even better? What can you do to improve on a technology that already offers superior speed, robustness, and accuracy? At Molecular Devices, we believe the answer lies in tailoring it to your lab, your applications, and your goals.

That's why our Custom Solutions group does more than simply integrate components from different manufacturers. We can provide a customized QPix platform specifically designed around your workflow, a truly one-of-a-kind solution developed to help you achieve faster, better results.

When you start with a foundation of the world's best technology and set no limits on customization possibilities, the sky's the limit.

QPix Highlights

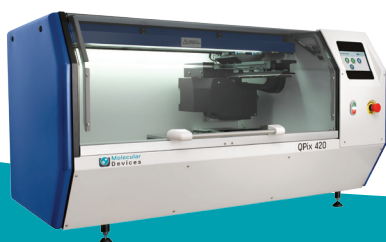
The instrument entrusted to map the human genome offers:

- Superior picking speed & accuracy
- Unique liquid handling & spreading capabilities
- Compatibility with a broad range of labware
- A wide range of fully pneumatic pin picking tools and biology-specific pin designs
- Exclusive, application-specific software algorithms for colony selection
- Complete data-tracking software

We Can Do That

A partial list of customizations that have enabled customers to translate time-consuming manual assays into robust, reliable automated solutions:

- Robotic integration from almost any manufacturer
- Specialized plate transport systems
- Filtration options
- Modified colony-picking heads
- Tailored software operating system modules



Contact your sales representative for more information:
Accelerate.My.Success@moldev.com
+1-800-635-5577

www.moleculardevices.com

Most Accurate Technology

1

+

Most Flexible Customization

1

=

Most Amazing Results

3

LET'S CHANGE THE WORLD TOGETHER

No matter what your area of focus, a custom QPix system can make a world of difference to the efficiency and productivity of your lab:

- Pharmaceuticals
- Agrochemicals
- Bioequivalents

CUSTOMIZING QPIX TO AUTOMATE AGAR-TO-AGAR MICROBIAL SAMPLE TRANSFERS

The QPix colony picker is now capable of performing agar-to-agar transfers of microbial samples. From single colonies to complex biology to dense microbial grids, QPix can be customized to automate the entire process, eliminating the hands-on time and potential errors of manual transfers.

By performing agar-to-agar transfers on an instrument instead of by hand, you can free up virtually 8 hours of time for other tasks, making it easy to streamline your workflow, increase sample throughput, and optimize lab productivity. No more manual gridding or stamping of colonies. No more repetitive stress. Just an efficient, ergonomic, walkaway solution.

Flexible and customizable, the QPix agar-to-agar module can be tailored to a variety of specific needs, from sample re-arranging and gridding to replicating and cherry picking.

Simple Steps To Greater Efficiency

STEP 1

Capture sample biology grown on solid agar or in liquid culture.

STEP 2

Transfer biological samples to a destination plate filled with solid agar.

STEP 3

Create dense arrays by pinning or re-arranging specific colonies to facilitate high throughput requirements.

Our Unique Pins. Your Unique Biology.

Specially engineered QPix pins ensure the reliable, efficient transfer of microbial samples for enhanced downstream screening capabilities. Designed for specific organisms, the pins optimize the transfer of everything from e-coli and yeast to algae and streptomyces—whatever your unique biology may be.

Simple, efficient, fully automated agar-to-agar microbe transfers on the QPix colony picker. Just one more way Molecular Devices can accelerate your success.

For more information or to see how QPix can enhance lab efficiency by automating your agar-to-agar applications, please contact us at: Accelerate.My.Success@moldev.com or +1-800-635-5577



POTENTIAL USES

- Library construction, screening and management
- Microbial strain evolution
- Antibiotic resistance screening
- Generation of microbial colony arrays for downstream screening
- Protein-protein interaction studies
- Phage display
- Culturing organisms which do not tolerate liquid media