



# QPix<sup>®</sup> Insights

Powered by Polar.ai

Optimize colony picking strategies  
with advanced AI/ML analysis

# QPix Insights Software

Microbial screening can be labor-intensive. Our QPix Microbial Colony Pickers automate colony selection and various sample preparation and plate handling processes. But how do we analyze the large volumes of multi-dimensional data generated? With QPix Insights, powered by Polar.ai, you can identify the best colonies aligned to optimal selection criteria with greater confidence in a shorter time.

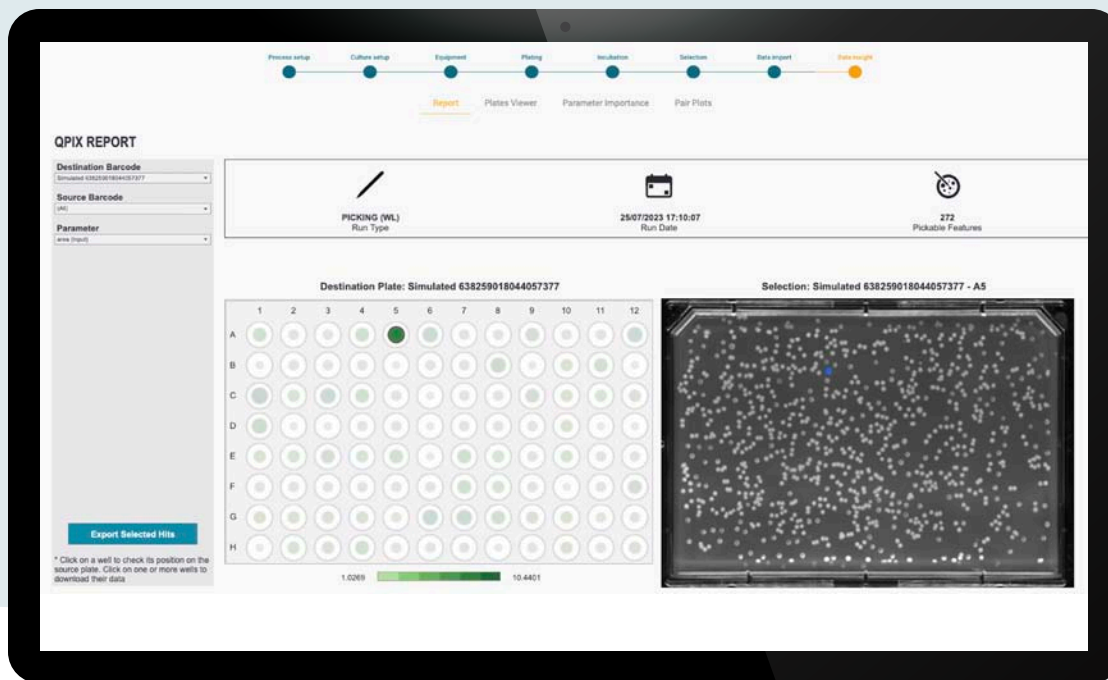


Figure 1. Plate map correlating colonies and their properties with colony on agar plate.



## AI-Powered Precision

Leveraging advanced AI and ML technologies, QPix Insights offers data interpretation, reducing errors and increasing reliability.



## Streamlined Data Integration

Say goodbye to fragmented data. QPix Insights seamlessly integrates data from various sources, providing a comprehensive overview and simplifying the analysis process.



## Optimized Colony Selection

The software's intelligent algorithms analyze colony parameters and analytical outputs to recommend the best selection criteria, improving the efficiency and quality of your selections.



## User-Friendly Interface

Designed with the user in mind, QPix Insights offers an intuitive interface that simplifies complex data analysis, making it accessible to all levels of scientific personnel.



## Cost and Time-Savings

Pick the right colony faster, leading to fewer reruns, less waste, and more time for value-adding work.

# Unlock the Full Potential of QPix

The QPix line of microbial colony pickers is known for its precision and efficiency. Now, with the integration of QPix Insights, these instruments are elevated to a whole new level of performance and functionality.



Figure 2. AI/ML model suggests refinements to selection parameters.

## Experience the Synergy

In the fast-paced world of microbial colony research, scientists often grapple with complex data analysis challenges. From time-consuming manual processes to the struggle of accurately interpreting vast data sets, these challenges can hinder the progress and reliability of your research.

Enter **QPix Insights** – a cutting-edge cloud software solution powered by Polar.ai and designed to allow easy data import and export, provide data visualization and reporting, and run AI/ML and other analytics. Co-developed by Molecular Devices and IDBS, QPix Insights is designed to address and overcome the hurdles of data analysis in microbial colony picking – allowing you to identify the best colonies aligned to optimal selection criteria with greater confidence in a shorter time.

## Key Enhancements with QPix Insights Integration



### AI/ML Technology

By utilizing AI/ML to learn from each experiment, QPix Insights suggests optimal selection parameters to apply to colonies in subsequent rounds of cloning.



### Advanced Visualization

QPix Insights provides enhanced visualization of data collected from QPix instruments. This means clearer insights into colony characteristics, leading to more informed decisions.



### Efficient Colony Selection

The software's AI-driven recommendations are tailored to the unique capabilities of QPix instruments, ensuring the selection of the best colonies with high accuracy and efficiency.



### Data Management

Manage and store data from QPix instruments in QPix Insights. This centralized data repository simplifies data retrieval and historical analysis by storing it in a single, secure location.



### Enhanced Decision-Making

With QPix Insights, you get a comprehensive view of your experiments, all in one application. From plate reader and process data to morphological and image data and analytical results, make data-driven decisions to optimize your research outcomes.

# Better Together:

## QPix Microbial Colony Pickers and QPix Insights

The combination of QPix instruments and QPix Insights represents a perfect synergy. The precision of QPix instruments, coupled with the smart analysis capabilities of QPix Insights, offers an unparalleled experience in microbial colony picking.

### Why Choose the Combined Solution?

- **Increased Productivity:** Save time and resources with streamlined processes.
- **Improved Research Quality:** Enhance the quality of your research with precise data analysis that optimizes selection criteria, yielding the highest-quality clones.
- **Future-Ready Technology:** Stay ahead in your field with the latest in technological innovation.



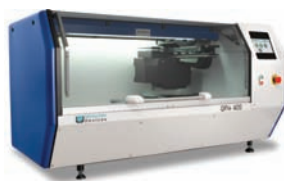
#### QPix HT system

Ready for robotics integration for maximum throughput and walkaway time.



#### QPix 450/460

From plating to picking – increase throughput with up to 210 destination plates in three stacker lanes.



#### QPix 420

Ideal system to replace manual with automated picking and allows for a flexible bed setup and labware use.



#### QPix XE\*

Ideal for small lab spaces with no compromise on efficiency, with advanced imaging and robotic capabilities.

For questions or additional support, please review QPix resources:

<https://support.moleculardevices.com/s/article/QPix-400-Series-Colony-Picking-System-Software-Download-Page>

<https://support.moleculardevices.com/s/article/Clone-Screening-software-operating-system-compatibility>

<https://support.moleculardevices.com/s/article/Molecular-Devices-Software>

For more information on IDBS Polar, please visit <https://www.idbs.com/polar/>

\*Price, time to deliver, and specifications will vary based on mutually agreed technical requirements. Solution requirements may cause adjustment to standard performance.

\*\*QPix Insights not available in China.

Custom solutions are subject to Molecular Devices Custom Products Purchase Terms available at <https://main-moleculardevices-hlxites.hlx.page/custom-products-purchase-terms>

#### Contact Us

Phone: +1.800.635.5577  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto: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

