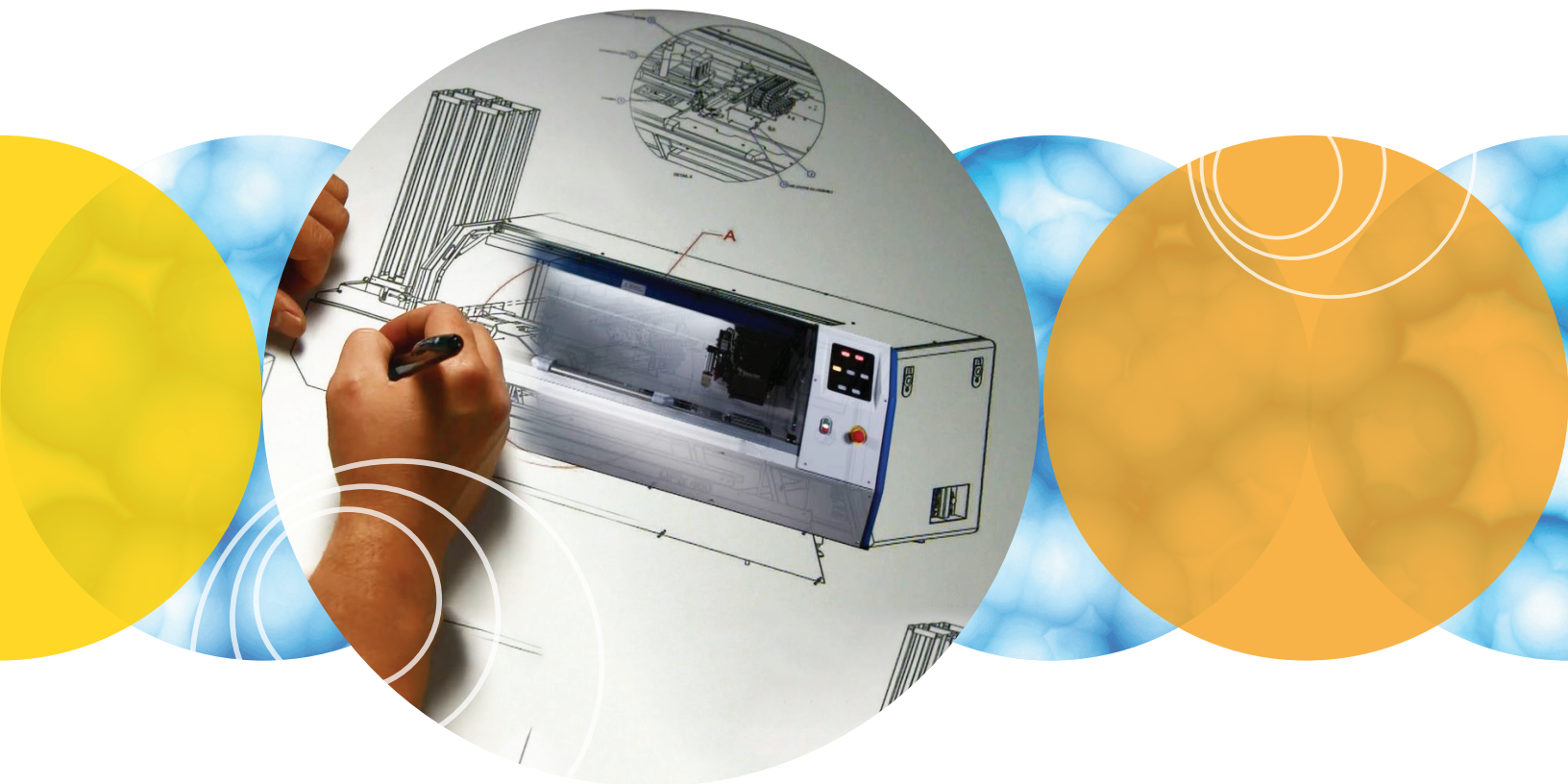


Advanced Workflow Engineering Solutions



Tailor our instruments. Transform your research.

**WHEN STANDARD DOESN'T FIT,
CUSTOMIZE YOUR QPIX SYSTEM,
NOT YOUR RESEARCH**

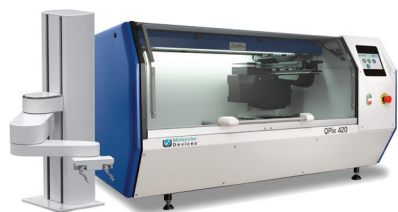
KEY STRENGTHS

- Turnkey automation solutions
- Services from consultation to implementation
- Hardware/software customization and support
- Validation of workflows with actual biological samples



Customized Colony Picking & Plating

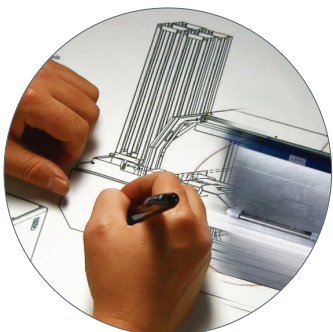
Our expert team of engineers and application scientists tailors our best-in-class QPix™ systems to meet your workflow demands for automated colony picking and plating. Proven applications include synthetic biology, antibody production & screening, strain evolution, and metagenomic research.



Customization example: integration of QPix Microbial Colony Picker with a robotic arm.

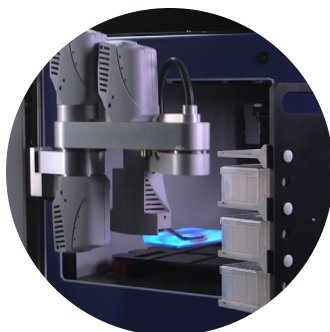
Our Advanced Workflow Engineering Solutions (AWES) team provides:

- Comprehensive package of services from consultation through implementation
- Complete setup and support for your fully integrated workflow
- Validation testing with actual biological samples
- In-house expertise for both application needs and instrument design
- Global service and support with over 30 years of experience in life science research



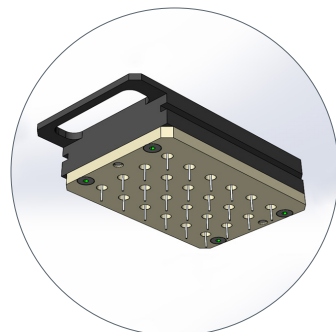
Turnkey service

Our AWES team works with you to understand your needs and challenges. We share our application and instrumentation expertise and propose solutions. We then test the chosen solution in-house to ensure it works for your application. Upon validation, we install the system in your lab and follow through on-site to support you as you advance your research.



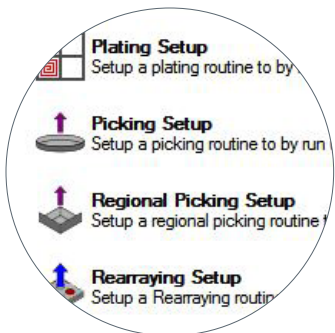
Scale-up automation

We adapt the QPix system to seamlessly integrate into larger automation workstations. AWES can design and test the whole workstation from scratch, incorporate it into an existing automated solution, or deliver a customized, automation-ready system for future integration.



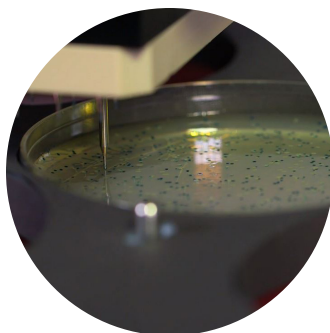
Hardware customization

From custom labware, to unique transfer pins, to special handling requirements, we'll work to find a solution and prove it through validation.



Software customization

We can modify the software to match specific Application Programming Interface (API) needs, enable unique workflow capability, integrate into LIMS or data management systems, customize a User Interface (UI), and of course control any custom designed hardware.



Biological validation

Our application scientists work with you to understand your process and biological requirements, so that we can reproduce and test the actual performance in-house prior to shipment.



Global service and support

With over 30 years of experience in the life science industry, you can count on Molecular Devices to deliver quality products and worldwide support.

Check out our customer case studies at
www.moleculardevices.com/custom-solutions

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

Regional Offices

USA and Canada
[+1.800.635.5577](tel:+1800.635.5577)
United Kingdom
[+44.118.944.8000](tel:+44.118.944.8000)
Europe*
[00800.665.32860](tel:00800.665.32860)

China (Beijing)
[+86.10.6410.8669](tel:+86.10.6410.8669)
China (Shanghai)
[+86.21.3372.1088](tel:+86.21.3372.1088)
Hong Kong
[+852.2248.6000](tel:+852.2248.6000)

Japan (Osaka)
[+81.6.7174.8331](tel:+81.6.7174.8331)
Japan (Tokyo)
[+81.3.6362.5260](tel:+81.3.6362.5260)
South Korea
[+82.2.3471.9531](tel:+82.2.3471.9531)

Brazil
[+55.11.3616.6607](tel:+55.11.3616.6607)

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