

SUSTAINABLE SCIENCE

A Greener Lab with QPix®

Transforming Lab Sustainability

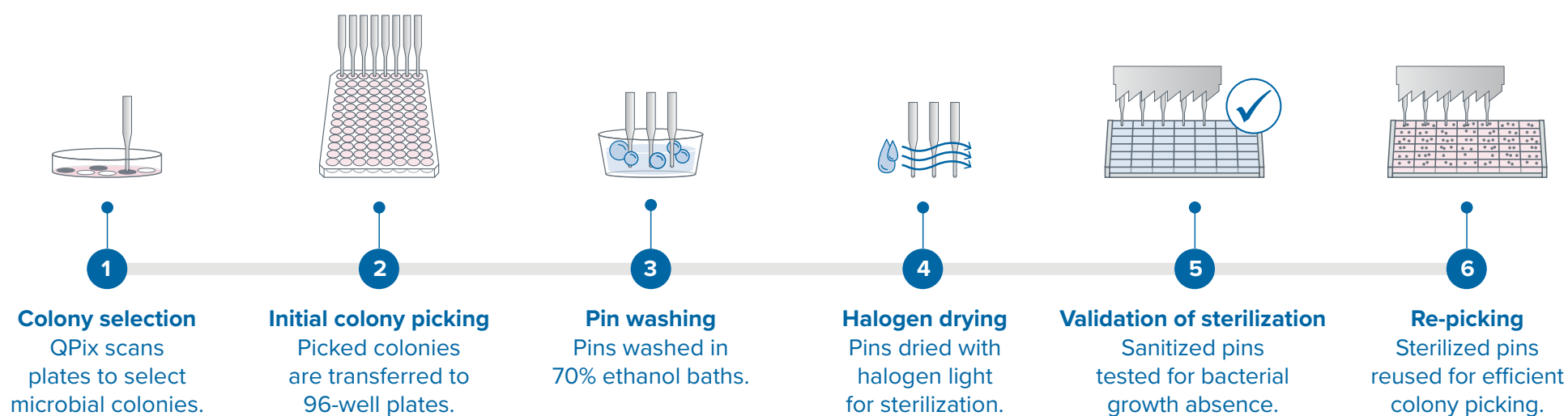
Harnessing QPix technology for an environmentally friendly laboratory



In today's laboratories, finding the right balance between efficiency and environmental responsibility is crucial. The QPix Microbial Colony Picker offers a game-changing solution that boosts productivity while promoting sustainability. With advanced technology and eco-friendly methods, QPix helps reduce waste, lower energy consumption, and cut operational costs. This innovative tool sets a new standard for green lab practices, allowing labs to operate efficiently without compromising their commitment to the environment.

Step-by-Step: The QPix Pin Process

Discover how QPix ensures sterility and efficiency through its innovative reusable pin process.



Impact of QPix in Colony Picking



Enhanced Sterility

Ensures contamination-free colony picking with rigorous sterilization protocols.



Cost Savings

Reusable pins significantly cut down on consumable costs.



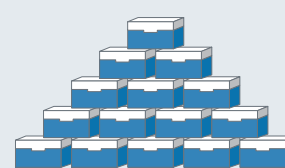
Increased Efficiency

Automates the picking process, reducing manual labor and errors.



Environmental Sustainability

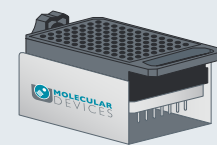
Reduces waste and energy consumption, promoting eco-friendly lab practices.



Disposable tips

96 tips/box
→ 80 boxes/month
*assuming to pick 384 clones/working day

VS



QPix re-usable pins

96 pins/picking head
→ Re-usable saving

Find out more moleculardevices.com/qpix

www.moleculardevices.com