

Let's talk about: Secondary Screening

Biomek Workstation from Beckman Coulter
- A solution as flexible as the biology you study



Liquid handling automation can turn weeks of work into a hands-free, walk-away project, but only if you choose the right platform and software package for the job. Biomek liquid handlers from Beckman Coulter offer a wide selection of integrations and the most straightforward user experience.

• Intuitive, user-friendly software



Biomek software, including **SAMI EX** and **DART 2.0**, is consistently hailed by end users for being easy to learn and even easier to use.

*"I find Biomek software to be one of the most reliable and user-friendly software packages, compared to competitive alternatives."**

Walt Dewolf
Researcher, **Array BioPharma**

*"[The Biomek] allows us to use all volume ranges through the use of pin tools in addition to the pipetting head."**

Martin Neuenschwander
Researcher, **FMP-Berlin**

• Flexible hardware for optimal throughput



Expand your Biomek from a single head to a double head, or add components to create a complete assay system.

• Set it and forget it



State-of-the-art automation enables you to program your entire screen, from cell seeding and feeding, to dose-response curves and viability assays, so you can keep your evening plans. Come back tomorrow for lunchtime analysis.

*"With my Biomek, I can now start my screenings at night and analyze my data after lunch."**

Lab Technician
Large Enterprise
Consumer Products Company

• D.I.Y. - Be IT free



Pull reports on your scans at any time, without needing to call IT for support. Biomek makes data analysis painless.

DID YOU KNOW?

73% of Biomek users never have to consult IT to pull data reports.

• Save your thumbs



Choosing a Biomek means your thumbs won't tire from the repetitive stress of pipetting an entire screening campaign. Your thumbs will thank you!

*"The Biomek workstations provide reproducible and efficient assay set up with significant reduction in ergonomic stress for scientists."**

Researcher
Global 500 Pharmaceuticals Company

• Take your pick: pipette tip or pin tool



With the option for two pods, you can pick hits and scale up, without either exhausting your sample or touching the machine.

*"With my Biomek, I can now pin tool compounds like there's no tomorrow."**

Operations Manager
Educational Institution

*"Beckman made it much easier for us to perform hit pickings - because of the software.. [I now can] leave students to perform hit-picking unsupervised."**

Susanne Hasse, Researcher
Max Planck Institute of Molecular Cell Biology and Genetics

• Point-and-click programming



Design your screening program with Biomek's simple drop-down menus, double-check your program's logic, and review your program's Estimated Time to Completion (ETC).

• Faster screens for faster results



Get your data faster, without reducing the rigor or quality of your screens.

*"With my Biomek workstation, I can now screen multiple samples in a matter of minutes."**

Researcher
Small Business
Health Care Company

*Research by TechValidate