



High Throughput Screening

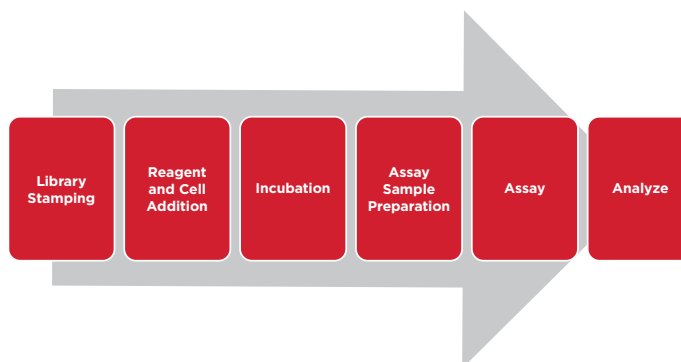
Make an impact by removing the barriers

For drug discovery and biomarker research, screening campaigns are employed to identify potential new treatments for diseases and to answer questions that remain unknown in the scientific community. From library screening, maintaining traceability of your cells and compounds, hit identification, and assay development, mining through the data collected and the required amount of hands-on time can be overwhelming. To meet the pressures associated with timelines and costs, liquid handling automation is a natural fit to remove human error and to process more scientific data in the lab which drive overall success.

Typical High Throughput Screening Workflow

Compound management, cell maintenance, and assay sample preparation for several end point analyzers all require flexibility of instrument between campaigns and to reliably manage each sample with the precision a scientist expects. The Biomek platform manages your workflow needs while keeping the flexibility in configuration as your workflows evolve.

Example workflows include small molecule screening, high content screening, phage or hybridoma screening.



Solutions That Meet Your Needs

We understand that automation is only part of your process so, based on your specific needs and challenges, our team will consult with you and engage our network of partners to provide you with access to complementary technologies and solutions in order to enhance your operation and enable you to obtain results you can repeat and rely on in an efficient manner.

Our systems answer your high throughput screening needs:

- Flexibility of the system allows you to do as little or as much as you need, whether you're doing a simple assay or complex screen
- Integration of other devices - we are continually adding to our library of over 300 device integrations
- Small to large volume pipetting mechanisms to/from low to high density plates and tubes
- Precise scheduling to provide consistent well-to-well treatment, always considering your daily throughput requirements
- Simplified data management software allows you to easily run and pull reports for analysis without the need of assistance from IT
- Confidence that the system will do its job every day, so you can do yours

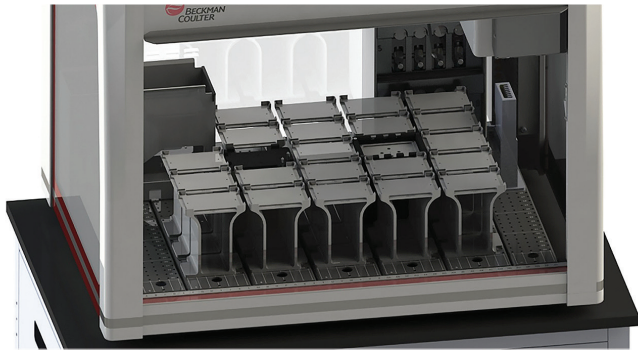


Figure 1. Have one to several plates to process on a regular basis? The i5 Span 8 or the i5 Multichannel system removes the burden of tedious manual tasks of cell-based assays. The Biomek platform aids in providing reliability for the critical step in your workflow.



Figure 2. Standardize your workflow with a complete walk-away solution. Modular design creates an adaptable system to automate your broad range of cell based assay workflows while increasing the laboratories' productivity.

Transform your Biomek liquid handler into a complete workflow solution that will optimize the efficiency, consistency, and reliability of your lab operations. From simple on-deck devices to complete robotic systems, we can develop focused, yet flexible, Biomek-based solutions by integrating Beckman Coulter and third-party instruments.

Biomek Software: Workflow Intelligence at Every Step

Every time your sample moves, data moves. Cell culture workflows are typically performed over the course of several weeks to months at a time. The Biomek Software Suite ensures data integrity to provide peace of mind and schedules activities to streamline automated workflows.

At the heart of every Biomek is the software. Biomek software provides users with the confidence to know their samples are being treated consistently every run, and with every liquid transfer, the data from your samples is being stored. The intuitive, drag and drop user interface makes setting up and maintaining your workflows easy. To meet the evolving needs of cellular workflows, we also offer:

- **SAMI EX:** provides complete automation and process control by creating planned schedules with the benefits of an optimized, predictable static schedule
- **DART 2.0 (Data Acquisition and Reporting Tool):** gathers data and synthesizes runtime information from Biomek log files to capture each manipulation of the sample during the course of the method
- **SAMI Process Management Software (SPMS):** is a calendar organizational tool that allows the addition, monitoring, and planning of SAMI EX methods and other events as part of user-defined processes
- **PowerPack:** features advanced tools to make programming data-intensive methods easy