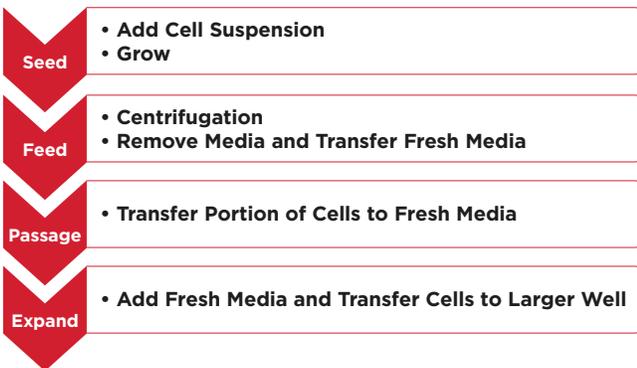


Cell Culture

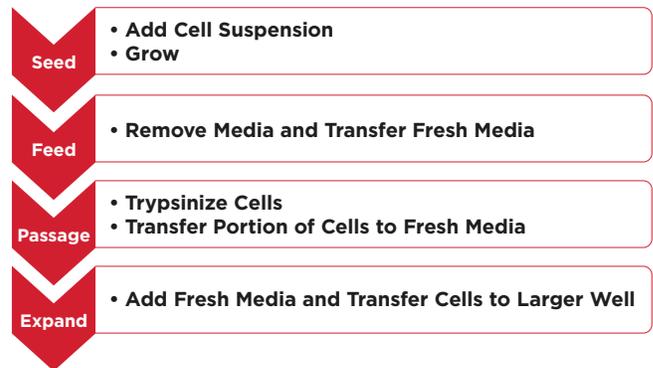
Automating Cell Maintenance

Cell culture is an essential workflow that enables studies to be performed in drug discovery and scientific research. Automation continues to aid laboratories in the time-consuming and error-prone tasks of cell maintenance. The need for automation grows as laboratories increase cell culture throughput in an effort to generate data earlier in their research studies. From seeding, feeding, and passaging cells to optimizing media conditions, automation aids in maintaining viable, healthy cells for downstream processes.

Sample Preparation for Suspension Cells



Sample Preparation for Adherent Cells



Solutions That Meet Your Needs

Minimize labor-intensive bottlenecks in your lab through automation. As your throughput increases, the Biomek automated cell culture solutions enable you to obtain results you can trace and rely on in an efficient manner, while optimizing your cell culture workflow and reducing your time at the bench.



Figure 1. Perform simple cell culture tasks on the Biomek 4000. This system is excellent for managing media exchanges and passaging cells. The Biomek 4000 fits in standard sized laminar flow hoods to maintain sterility.



Figure 2. 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 and cell culture processing. Maintain your cells in an environment that promotes efficiency and reproducibility.



Figure 3. Standardize your workflow with a complete walk-away solution. Designed to seed, monitor, passage and harvest cells based on your set criteria.

As your throughput requirements increase, the i-Series platform has integration capabilities that enable you to have a complete walk-away solution in your laboratory. From analyzers that measure confluence or cell count to centrifugation for suspension cell workflows, our team can consult with you to build a solution that will meet your needs and remain flexible as your laboratory's needs evolve.

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