



## Which Automated Solution is Right for Your 3D Biology Research?

We're proud to offer several best-in-class automated solutions for 3D biology research, each with its own unique features and capabilities. Choose the solution that best meets the specific needs and requirements of your lab.



### CellXpress.ai™ Automated Cell Culture System

#### AI-driven, turnkey, walkaway solution

The CellXpress.ai system is an AI-driven cell culture innovation hub that gives your team total control over demanding cell culture feeding and passaging schedules—eliminating time in the lab while maintaining a 24/7 schedule for growing and scaling multiple stem cell lines, spheroids, or organoids. It's unified software environment brings together single-device initialization, protocol error protection, and machine learning-enabled decision-making to streamline the cell culture process and makes it easier than ever to navigate and monitor even the most complex workflows.



### Laboratory Automation & Customization Solutions

#### Tailored end-to-end lab solutions

Our laboratory automation and customization solutions offer the ultimate flexibility with custom-built automated workflows using cutting-edge instruments that work harmoniously for autonomous, long-term, live cell culture growth and label-free monitoring. We work with you to design and build a system that meets your exact workflow needs. Utilize flexible scheduling software to integrate all the functionalities you need to take your 3D research to the next level.



### BioAssemblyBot® 400 (BAB400)

#### More than a bioprinter

The BAB400 is a multi-functional system that brings together 3D bioprinting, liquid dispensing, and plate handling allowing you to accelerate screening of complex 3D models and workflow automation. It's 6-axis robotic arm utilizes a wide range of interchangeable tools to deliver ever-growing capabilities and reliably reproduce 3D model systems with increased throughput and precision, alleviating common issues with manual workflows.

## At-a-glance

### When you need turnkey operation for core assays:

#### CellXpress.ai Automated Cell Culture System

- 24/7/365 operation to maximize productivity.
- Protocol set-up guidance from in-house scientists.
- Onboard brightfield and fluorescent imager for day-to-day monitoring and endpoint measurement.
- Integrated software tracks the entire cell journey, including plate and well relationship after passaging events.
- Machine learning-assisted decision-making standardizes the cell growth expansion process.
- Dynamic alert system – notifies users with smart error handling
- Seamless integration with ImageXpress® high-content imager, plate centrifuge, and FLIPR available for advanced workflows.
- An approachable, truly hands-off automated solution built by scientists, for scientists.
- Protocol error protection.
- Quick start, single-device initialization.
- Compatible with a wide range of validated consumables and liquid types.
- Automated media monitoring.
- Onboard low-volume and bulk media storage.
- Select from 44 or 154 plate incubator.

### When ultimate flexibility is your goal:

#### Laboratory Automation & Customization Solutions

- 24/7/365 operation to maximize productivity.
- Integration of one or multiple high-content confocal/epifluorescence imagers available for monitoring and end-point assays.
- Protocol set-up guidance from in-house scientists.
- Dynamic alert system – notifies users with smart error handling
- Advanced, flexible scheduling software to coordinate the entire workflow.
- Designed to your specific throughput, speed, and workflow requirements.
- Several liquid handling configurations available including pipette-free dispensing.
- Unlimited expansion and configurability.
- Gives you the freedom and flexibility for assay development

### When you need multi-functional assay automation:

#### BioAssemblyBot 400

- Integration with high-content imagers available for endpoint measurements.
- Consistent dispensing, placement, and printing of organoids vs. manual workflows.
- Cost-effective solution for combining 3D printing with workflow automation.
- Compatible with custom-selected consumables.
- Small footprint.

## Need help deciding which solution best fits the needs of your lab?

Speak with an expert to find the right solution and give your lab confidence in experimental outcomes to make key decisions sooner, achieve milestones faster, and get to clinic earlier.

#### Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing of worldwide distributors.

#### Regional Offices

USA and Canada	<a href="tel:+18006355577">+1.800.635.5577</a>	Taiwan/Hong Kong	<a href="tel:+886226567585">+886.2.2656.7585</a>
United Kingdom	<a href="tel:+441189448000">+44.118.944.8000</a>	Japan	<a href="tel:+81363629109">+81.3.6362.9109</a>
Europe*	<a href="tel:0080066532860">00800.665.32860</a>	South Korea	<a href="tel:+82234719531">+82.2.3471.9531</a>
China	<a href="tel:+864008203586">+86.4008203586</a>	India	<a href="tel:+917386611198">+91.73.8661.1198</a>

\*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

