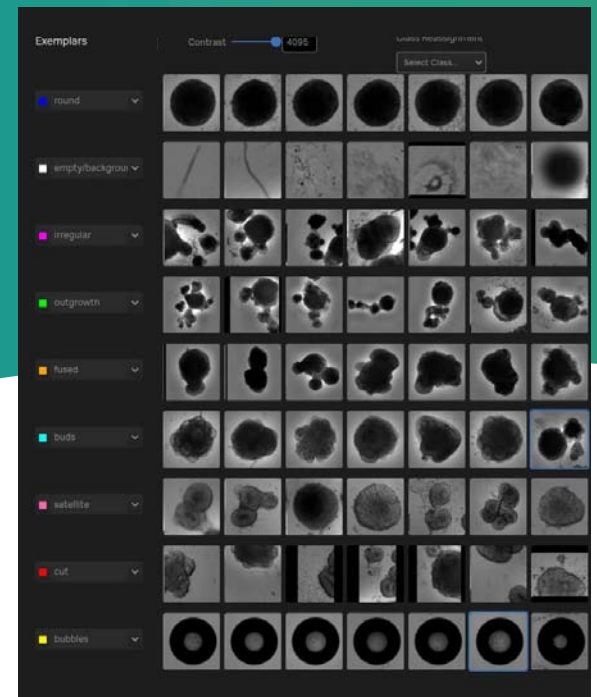


An intelligent, hands-free software environment created by scientists, for scientists.

Designed for continuous operation, the CellXpress.ai™ Automated Cell Culture System boasts a software environment that will keep you from worrying about missing a feeding or passaging step.

The CellXpress.ai system ensures workflow consistency by using automation and employs AI-based image analysis to determine the optimal time for cell culture feeding and passaging events. It even sends alerts to keep you informed of your cell culture status. All of this is backed by a comprehensive event log to confirm on-time feedings and critical task execution with complete digital microscopy records.



Preconfigured protocols and protocol elements

- Turnkey, validated protocols for organoid and cell models
- Self-contained solution eliminates the need to coordinate different instruments
- Software clearly pinpoints issues so scientists can fix problems themselves without the need for an automation engineer
- Smart media handling – offers software-controlled heating, cooling, stirring, and volume tracking in various media vessel types and volume capacities to support multiple workflow environments.

Machine learning-based classification & segmentation

- Reduce human error — label-free decision-making removes variability, maintains sterility, and increases confidence in success.
- Phenotypic classification — AI-driven classification based on an advanced set of morphological and texture measurements for a variety of biological applications, such as organoid quality control or colony formation
- Machine learning-assisted decision-making standardizes the cell growth expansion process.

Protocol failure protection

- Software detects user-entry errors and notifies user before protocols are saved
- Once a protocol is launched, users are alerted to any issues in the workflow (e.g. insufficient media)

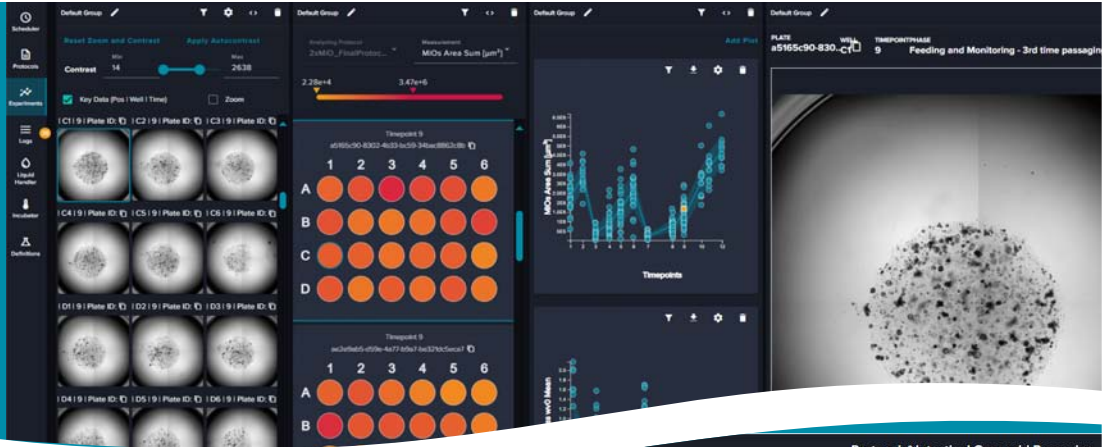
Unified software environment

- Single-source initialization — no need to initialize multiple instruments
- Remote access — receive alerts when milestones are met and access the system from anywhere to check on progress

Offering a wide range of features to make your cell culture process more efficient, reliable, and reproducible, the CellXpress.ai system brings together single-source initialization, protocol failure 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 workflow.

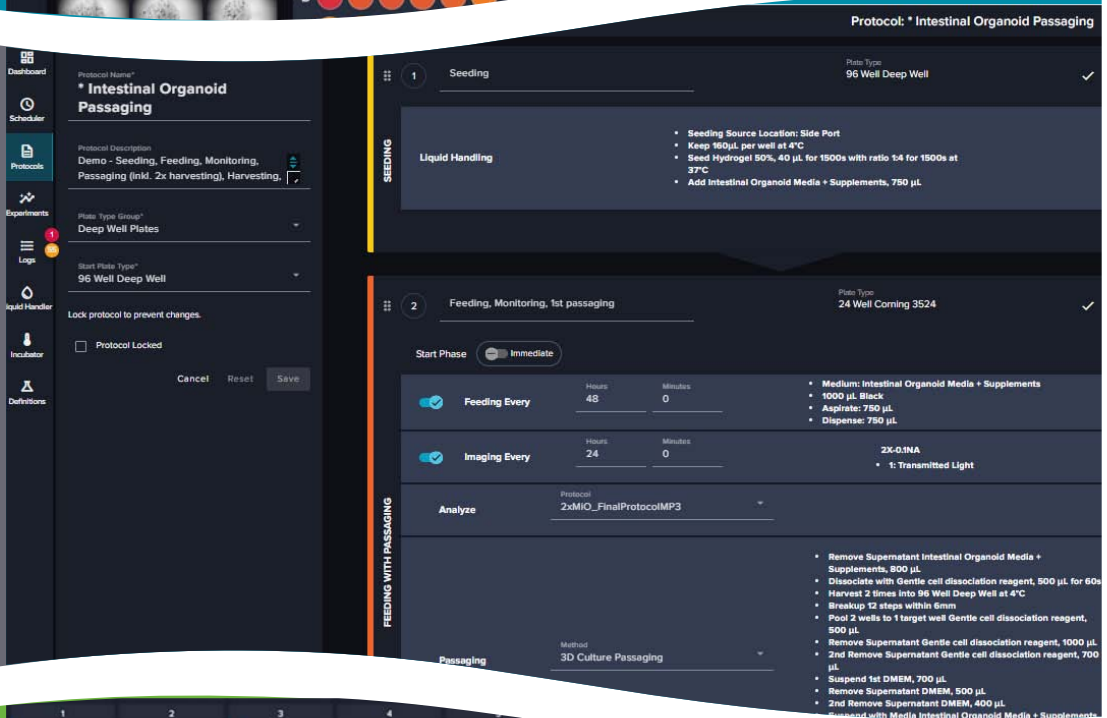
EXPERIMENTAL RESULTS PANEL

Flexible 'widget space' allows the user to construct their own data view appropriate for their experiment.



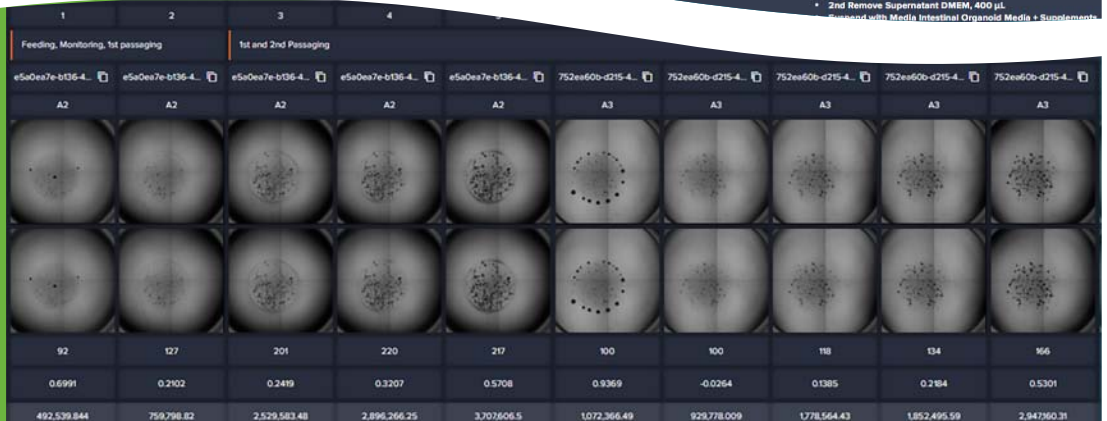
PASSAGING PROTOCOL

Easily create protocols with an intuitive interface designed for biologists. Add building blocks for each phase of the workflow, including seeding, feeding, incubation, harvesting, passaging, imaging, and analysis.



CELL JOURNEY

Track your cell culture experiments over time and across passage events.



Contact Us

Phone: +1.800.635.5577
 Web: www.moleculardevices.com
 Email: info@moldev.com
 Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1.800.635.5577
 United Kingdom +44.118.944.8000
 Europe* 00800.665.32860
 China +86.4008203586
 Taiwan/Hong Kong +886.2.2656.7585
 Japan +81.3.6362.9109
 South Korea +82.2.3471.9531
 India +91.73.8661.1198

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