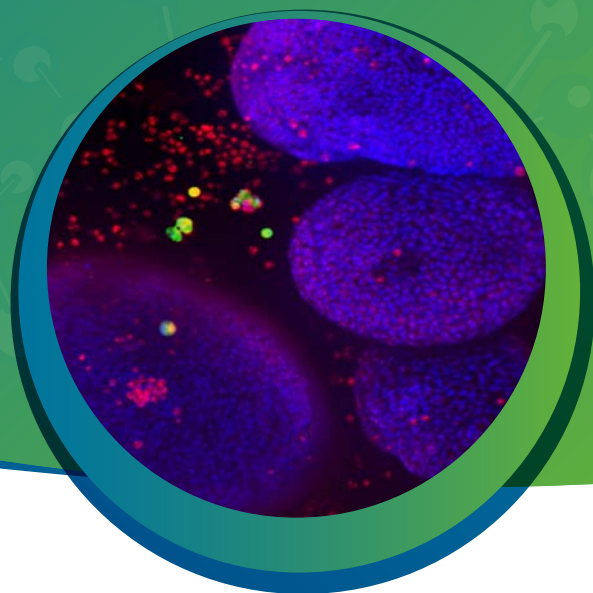


ORGANOID

INNOVATION CENTER



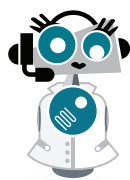
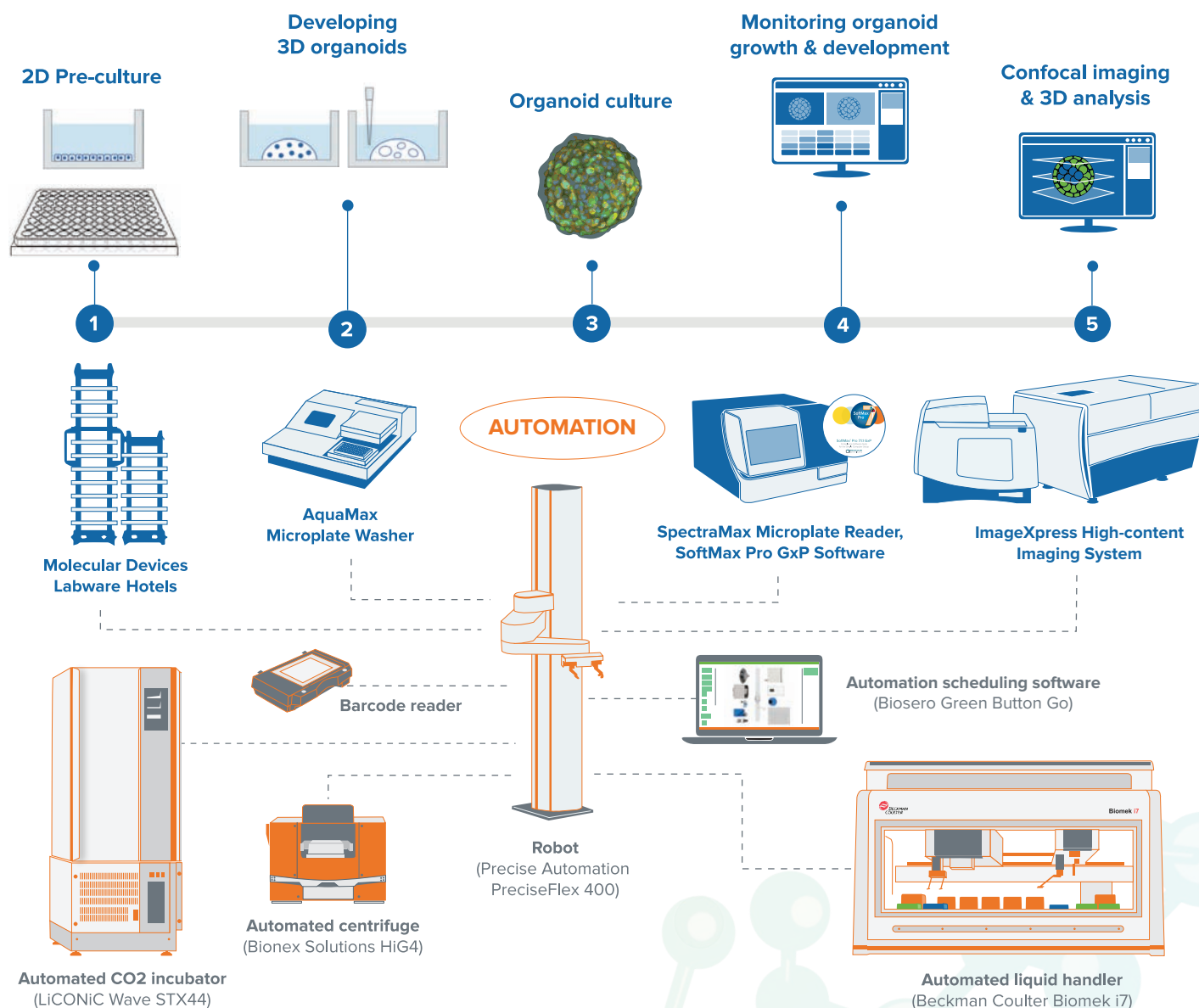
Quickly adopt innovative,
3D biological methods and
technologies for drug discovery

The Organoid Innovation Center showcases cutting-edge instruments that work harmoniously together for autonomous, long-term, 3D and 2D live cell culture growth and monitoring with intelligent imaging. This integrated workflow provides quality control alerts and readiness, 3D organoid screening, and deep learning image analysis that reveals hidden patterns other technologies miss.

The end-to-end solution standardizes the organoid development process by growing complex 3D cell cultures (incubation, treatment, passaging, and monitoring) and the downstream analysis capabilities utilizing confocal imaging, analysis, and data processing. The Organoid Innovation Center delivers consistent, unbiased, and biologically-relevant results at scale.



See the power and flexibility of an automated and customizable high-throughput screening solution



Want to know more about automating your organoid screening workflow?
Visit moleculardevices.com/organoid-innovation-center

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

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