

3D Ready™ Organoid Expansion Service

Do you have an established organoid line that grows in a gel matrix? Are you time-limited and finding it difficult to produce consistently-sized, high quality organoids for your work?

Our Organoid Expansion Service can help. Let us do the hard work and grow them for you.



We specialize in patient derived organoids



From 0.3 to over 6 million organoids per batch



Approximately 3–6 months from initial organoid receipt to delivery of vials



We grow healthy and diseased organoid types



~30–50 µm in size: optimal for imaging and biochemical assays



Assay Ready: Thaw-Plate-Test

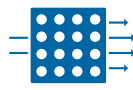
Our team of organoid experts will closely work with you throughout the process to ensure a successful expansion. Elements of the process can be customized to meet your specific needs.



1

Discussions with experts to scope out project

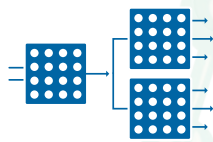
Custom requirements: Suitability, scale, QC & timelines.



2

Pre-expansion

Determining parameters for the individual line in the bioreactor



3

Bespoke bioreactor-based expansion

Harvested organoids counted, sized and frozen in cryovials



4

QC testing thawed vials:

Recovery post thaw, pathogen testing, drug validation, STR analysis, custom QC



5

Custom cryopreserved assay-ready organoids delivered to customer

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