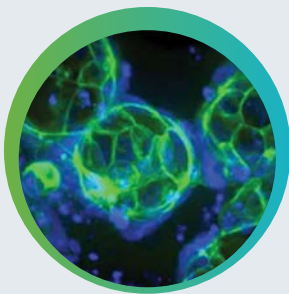


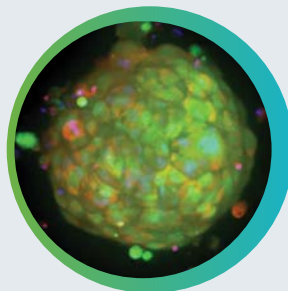
Gain human-relevant insights with patient-derived assay-ready organoids

Cryopreserved, highly reproducible patient-derived organoids ready for immediate use!

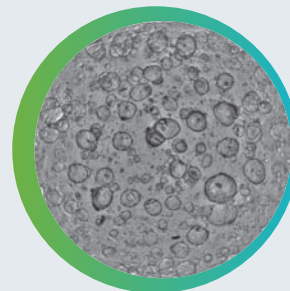
Our 3D Ready™ organoids are patient-derived models offering more physiological relevance and are ready to use straight from cryopreserved vials. Manufactured using our proprietary bioprocess and provided with full protocols and comprehensive scientific support, our organoids empower you to achieve reliable, reproducible research quickly and easily – with low upfront investment in time and resources.



Colorectal Cancer (CRC) Organoids



Breast Cancer Organoids



Healthy Intestinal Organoids



Get started with confidence

Receive full protocols and comprehensive scientific support with each vial, enabling even new users to begin their research quickly and effectively.



Streamline your research

Simplify your workflow with an integrated HUB Organoid license for proof-of-concept studies, saving time, reducing costs, and eliminating complexity.



Ready for immediate use

Our organoids are delivered in convenient frozen vials containing at least 100,000 organoids, can be stored in your freezer and are ready to use when you need them—no culturing required, just thaw and seed into your microplate.



Achieve reliable results

Our quality-assured, tested and traceable organoids minimize variability, ensuring consistent batch-to-batch results for reliable, reproducible data.



Get started with 3D Ready Organoids today – speak to an expert!

<https://bit.ly/3yNTzL4>

Available organoid lines

Organoid line	Description	Minimum number of organoids/vial	Part Number
3D Ready Breast Cancer	BCO-005, Primary Tissue, Invasive Carcinoma Mixed	250,000	OES-BCO005-CXP1
	BCO-037, PDX Derived, Invasive Ductal Carcinoma	250,000	OES-BCO037-CXP1
	BCO-160, PDX Derived, Invasive Ductal Carcinoma/DCIS	250,000	OES-BCO160-CXP1
3D Ready Colorectal Cancer	ISO38, Sigmoid Colon; Duke's Stage C1; Mut <i>APC</i> , <i>TP53</i> , <i>SMAD4</i>	100,000	OES-ISO38-CXP1
	ISO48, Caecum; Duke's Stage C2; Mut <i>PIK3CA</i> , <i>CTNNB1</i>	100,000	OES-ISO48-CXP1
	ISO49, Lower Sigmoid; Duke's Stage C1; Mut <i>APC</i> , <i>TP53</i> , <i>KRAS</i> (G12D), <i>PIK3CA</i> , <i>FBXW7</i>	100,000	OES-ISO49-CXP1
	ISO50, Rectum; Duke's Stage C2; Mut <i>APC</i> , <i>TP53</i> , <i>KRAS</i> (G12V), <i>SMAD4</i> , <i>SMAD2</i>	100,000	OES-ISO50-CXP1
	ISO57, Upper Rectum; Duke's Stage A; Mut <i>APC</i> , <i>TP53</i> , <i>KRAS</i> (G13D)	100,000	OES-ISO57-CXP1
	ISO68, Sigmoid Colon; Duke's Stage C1; Mut <i>APC</i> , <i>TP53</i> , <i>KRAS</i> (G13D), <i>SMAD4</i>	100,000	OES-ISO68-CXP1
	ISO72, Caecum; Duke's Stage B; Mut <i>APC</i> , <i>KRAS</i> (G12V), <i>ERBB3</i>	100,000	OES-ISO72-CXP1
	ISO75, Caecum; Duke's Stage C1; Mut <i>BRAF</i> , <i>ARID1A</i> , <i>MSH3</i> , <i>RNF43</i>	100,000	OES-ISO75-CXP1
	ISO78, Caecum; Duke's Stage C1; Mut <i>APC</i> , <i>TP53</i> , <i>KRAS</i> (G12D)	100,000	OES-ISO78-CXP1
3D Ready Normal Duodenum from Familial Adenomatous Polyposis (FAP) patient	DP41N2, FAP Normal Duodenum (heterozygous Mut <i>APC</i>)	200,000	OES-DP41N2-CXP1
	DP41N2, FAP Normal Duodenum (heterozygous Mut <i>APC</i>)	100,000	OES-DP41N2-CXP1B

Need a different model?

We also offer **organoid expansion services!** Generate large numbers of customized, high-quality, low-variability organoids in a convenient assay-ready format.



Get started with 3D Ready Organoids today – speak to an expert!

<https://bit.ly/3yNTzL4>

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

