

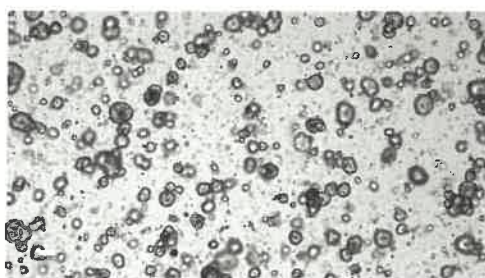
Certificate of Analysis

Item number: OES-ISO49-CXP1
Lot number: CXP1-089
Line ID: ISO49 **CofA Issue: 1**
Donor Age/Sex: 77/Male
Species: Human
Tissue: Colorectal Cancer
Volume per vial: 1 mL
Date frozen: 18 JUN 2024
Format: Organoids stored in freezing medium
Storage conditions: Store at -150°C
Usage: For research use only – organoids are single use only and not to be passaged for repeated use or for other organoid expansion processes

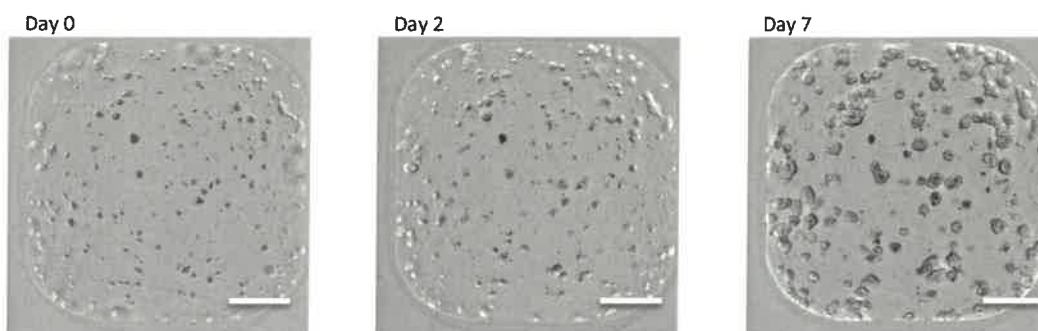
Characteristic	Specification	Result
Passage number	Not applicable	P25
Average organoids per vial	≥100,000	133,686
Average organoid size post thaw	Report results	47.6 μm
Growth after recovery (Visual observation)	Growth observed over 7 days	Growth observed over 7 days
Mycoplasma (PCR based assay)	Not detected	Not detected
Viral screening HIV-1, HIV-2 HBV CMV EBV HPV (PCR based assay)	Not detected Not detected Not detected Not detected Not detected	Not detected Not detected Not detected Not detected Not detected
Sterility Bacteria, yeast and fungi (Direct inoculation)	Not detected	Not detected

Representative image of organoids in culture:

4x magnification
scale bar = 100μm



Representative images following thaw (4x magnification, scale bar = 700μm)



Certificate of Analysis



Item number: OES-ISO49-CXP1
Lot number: CXP1-089
Line ID: ISO49

CofA Issue: 1

Written by:

Name	Role	Signature	Date
GIUSY TORNILLO	SENIOR SCIENTIST	<i>amy z</i>	26 JUL 2024

QA approval:

Name	Role	Signature	Date
Ben Jones	QA specialist	CB Jones	26 Jul 2024