

Certificate of Analysis

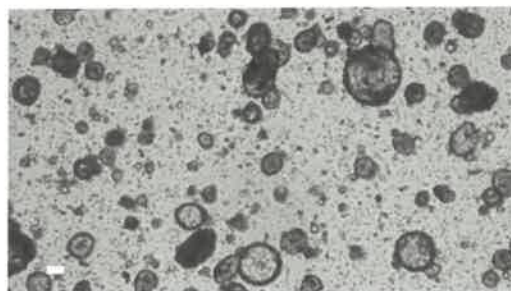
Item number: OES-ISO50-CXP1
 Lot number: CXP1-090
 Line ID: ISO50

CofA Issue: 1

Donor Age/Sex: 75/Male
 Species: Human
 Tissue: Colorectal Cancer
 Volume per vial: ≥0.5 mL
 Date frozen: 31 JUL 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	P31
Average organoids per vial	≥100,000	148,007
Average organoid size post thaw	Report results	44.56 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-ISO50-CXP1
Lot number: CXP1-090
Line ID: ISO50

CofA Issue: 1

Written by:

Name	Role	Signature	Date
Kim Letchford	Senior Scientist		18 SEP 2024

QA approval:

Name	Role	Signature	Date
Ben Jones	QA specialist		18 Sep 2024