

Certificate of Analysis

Item number: OES-ISO48-CXP1
 Lot number: CXP1-091
 Line ID: ISO48

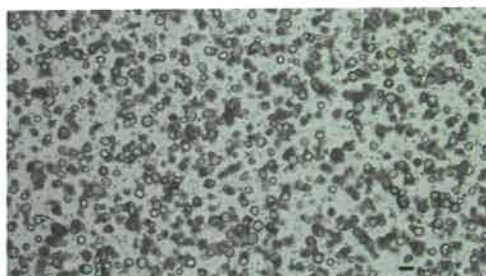
CofA Issue: 1

Donor Age/Sex: 43/Male
 Species: Human
 Tissue: Colorectal Cancer
 Volume per vial: ≥0.5mL
 Date frozen: 04 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	P35
Average organoids per vial	≥100,000	163,713
Average organoid size post thaw	Report results	42.42 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

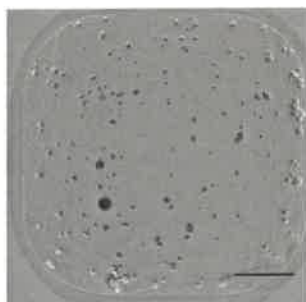
Representative image of organoids in culture:

4x magnification
 scale bar = 100μm

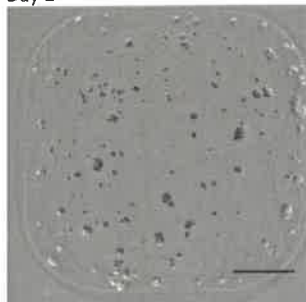


Representative images following thaw (4x magnification, scale bar = 700μm)

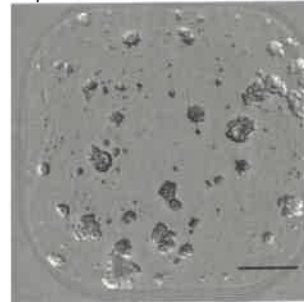
Day 0



Day 2



Day 7



Certificate of Analysis



Item number: OES-ISO48-CXP1
Lot number: CXP1-091
Line ID: ISO48

CofA Issue: 1

Written by:

Name	Role	Signature	Date
William Newell	Senior Bioprocess Engineer	<i>W Newell</i>	15 AUG 2024

QA approval:

Name	Role	Signature	Date
Ben Jones	QA Specialist	<i>CB Jones</i>	15 Aug 2024