

Certificate of Analysis

Item number: OES-BCO037-CXP1
 Lot number: BCXP1-021
 Line ID: BCO-037

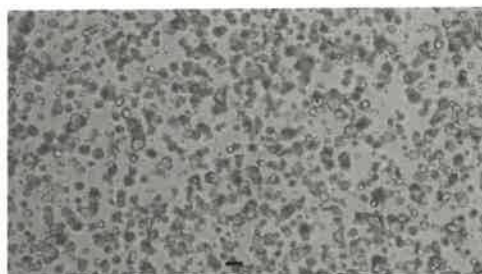
CofA Issue: 2

Donor Age/Sex: Unknown/F
 Species: Human
 Tissue: Breast Cancer
 Volume per vial: ≥0.5 mL
 Date frozen: 19 JUL 2022
 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

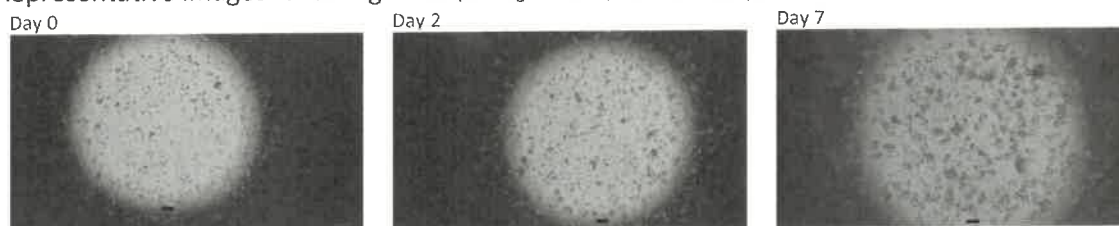
Characteristic	Specification	Result
Passage number	Not applicable	P15
Average organoids per vial	≥250,000	369,399
Average organoid size post thaw	Report results	25.33 μm
Growth after recovery (Visual observation)	Growth observed over 7 days	Growth observed over 7 days
Mycoplasma (PCR based assay)	Not 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 = 100μm)



Certificate of Analysis



Item number: OES-BCO037-CXP1
Lot number: BCXP1-021
Line ID: BCO-037

CofA Issue: 2

Written by:			
Name	Role	Signature	Date
Mairian Thomas	Project Manager		31 MAY 2024.

QA approval:			
Name	Role	Signature	Date
Ben Jones	QA Specialist	CB Jones	31 May 2024