

# Certificate of Analysis

Item number: OES-ISO72-CXP1

Lot number: CXP1-100

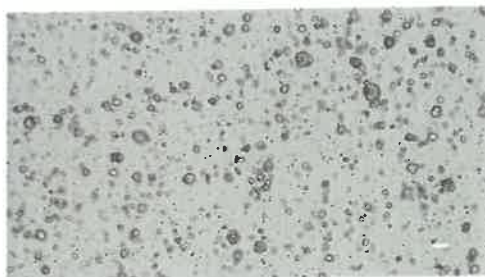
Line ID: ISO72

CofA Issue: 1

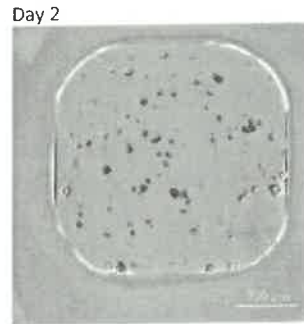
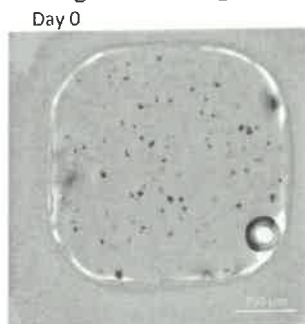
**Donor Age/Sex:** 83/Female  
**Species:** Human  
**Tissue:** Colorectal Cancer  
**Volume per vial:** 0.5 mL  
**Date frozen:** 21 NOV 2024  
**Format:** Organoids stored in freezing medium  
**Storage conditions:** Store at -150°C  
**Usage:** For research use only – organoids are single use only and not to be passaged for repeated use or for other organoid expansion processes

| Characteristic   | Specification  | Result   |
|--|--|--|
| Passage number   | Not applicable   | P33  |
| Average organoids per vial   | ≥ 100,000  | 127,555  |
| Average organoid size post thaw  | Report results   | 32.63 μm   |
| Growth after recovery<br>(Visual observation)                                    | Growth observed over 7 days  | Growth observed over 7 days  |
| Mycoplasma<br>(PCR based assay)  | Not detected   | Not detected   |
| Viral screening<br>HIV-1, HIV-2<br>HBV<br>CMV<br>EBV<br>HPV<br>(PCR based assay) | Not detected<br>Not detected<br>Not detected<br>Not detected<br>Not detected | Not detected<br>Not detected<br>Not detected<br>Not detected<br>Not detected |
| Sterility<br>Bacteria, yeast and fungi<br>(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-ISO72-CXP1

Lot number: CXP1-100

Line ID: ISO72

CofA Issue: 1

## Written by:

| Name           | Role      | Signature | Date        |
|----------------|-----------|-----------|-------------|
| Harman Chagger | Scientist |           | 20 DEC 2024 |

## QA approval:

| Name      | Role          | Signature | Date        |
|-----------|---------------|-----------|-------------|
| Ben Jones | QA Specialist | C B Jones | 20 Dec 2024 |