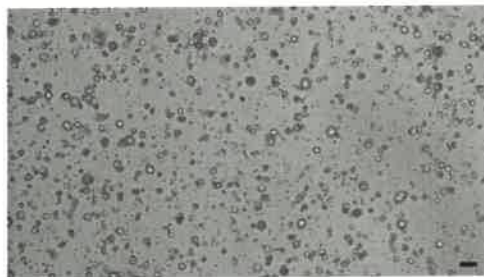


Item number: ISO72-CXP1
Lot number: CXP1-080
Line ID: ISO72
Donor Age/Sex: 83/F
Species: Human
Tissue: Colorectal Cancer
Volume per vial: ≥0.5 mL
Date frozen: 09 JUN 2023
Expiry date: 09 JUN 2024
Format: Organoids stored in freezing medium
Storage conditions: Store at -150°C
Usage: For research use only – 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 | P30 |
| Average organoids per vial | ≥100,000 | 125,072 |
| Average organoid size post thaw | Report results | 33.81µm |
| 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: ISO72-CXP1
Lot number: CXP1-080
Line ID: ISO72

| Written by: | | | |
|----------------|-----------------|-----------|-------------|
| Name | Role | Signature | Date |
| Mairian Thomas | Project Manager | | 19 OCT 2023 |

| QA approval: | | | |
|--------------|---------------|-----------|-------------|
| Name | Role | Signature | Date |
| Ben Jones | QA Specialist | | 19 oct 2023 |