

# Certificate of Analysis

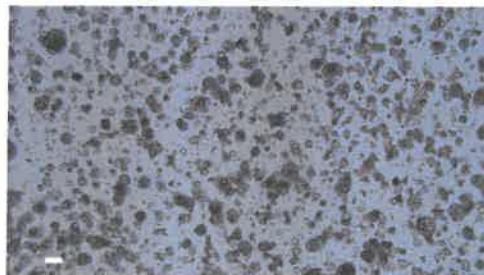
**Item number:** OES-BCO160-CXP1  
**Lot number:** BCXP1-025  
**Line ID:** BCO160

**CofA Issue: 1**

**Donor Age/Sex:** Unknown/Female  
**Species:** Human  
**Tissue:** Breast Cancer  
**Volume per vial:** ≥0.5 mL  
**Date frozen:** 17 SEP2024  
**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

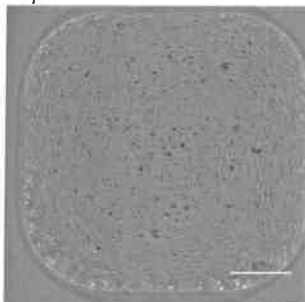
Characteristic	Specification	Result
<b>Passage number</b>	Not applicable	P11
<b>Average organoids per vial</b>	≥250,000	483,795
<b>Average organoid size post thaw</b>	Report results	28.91 μm
<b>Growth after recovery</b> (Visual observation)	Growth observed over 7 days	Growth observed over 7 days
<b>Mycoplasma</b> (PCR based assay)	Not detected	Not detected
<b>Viral screening</b> 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
<b>Sterility</b> Bacteria, yeast and fungi (Direct inoculation)	Not detected	Not detected

Representative image of organoids in culture:  
 4x magnification  
 scale bar = 100μm

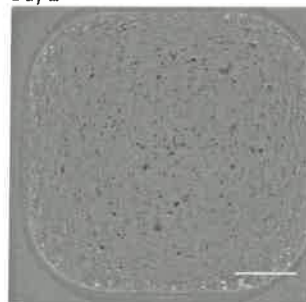


Representative images following thaw (4x magnification, scale bar = 700μm)

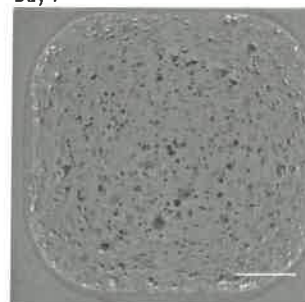
Day 0



Day 2



Day 7



# Certificate of Analysis



Item number: OES-BCO160-CXP1  
Lot number: BCXP1-025  
Line ID: BCO160

CofA Issue: 1

Written by:			
Name	Role	Signature	Date
Carly Bunston	Scientist		25 Oct 2024

QA approval:			
Name	Role	Signature	Date
Ben Jones	QA specialist		25 Oct 2024