

Certificate of Analysis

Item number: OES-DP41N2-CXP1B
 Lot number: FCXP1-028
 Line ID: DP41N2

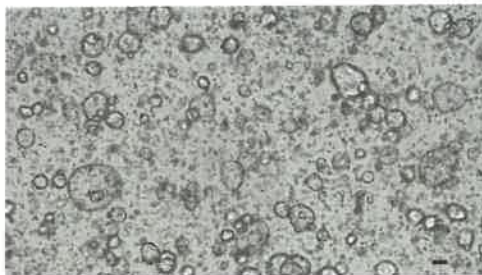
CofA Issue: 1

Donor Age/Sex: 39/Male
 Species: Human
 Tissue: Duodenum
 Volume per vial: ≥1mL
 Date frozen: 06 Sep 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	P23
Average organoids per vial	≥100,000	138,781
Average organoid size post thaw	Report results	38.08 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

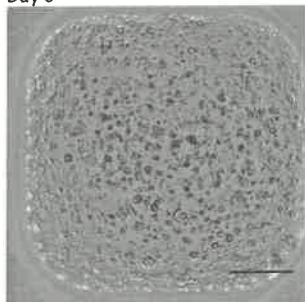
Representative image of organoids in culture:

4x magnification
 scale bar = 100μm

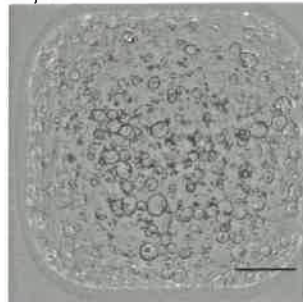


Representative images following thaw (4x magnification, scale bar = 700μm)

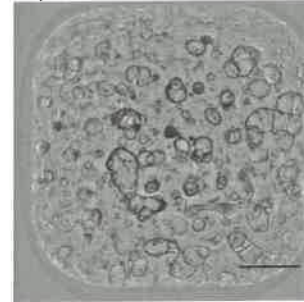
Day 0



Day 2



Day 7



Certificate of Analysis



Item number: OES-DP41N2-CXP1B
Lot number: FCXP1-028
Line ID: DP41N2

CofA Issue: 1

Written by:

Name	Role	Signature	Date
William Newell	Senior Bioprocess Engineer	<i>wnewell</i>	25 Oct 2024

QA approval:

Name	Role	Signature	Date
Ben Jones	QA Specialist	<i>C B Jones</i>	25 Oct 2024