

Product Number:	103527
Lot Number:	036603
Expiration:	31-Dec-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.22 x 10 ⁹ Beads/mL

Review/Approval

Junpil Park

QC Supervisor



15 Feb 2021

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.