

Product Number:	103208
Lot Number:	036317
Expiration:	28-Dec-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.12 x 10 ⁹ Beads/mL

Review/Approval

Junpil Park

QC, Supervisor



10 Feb 2021

Name

Title

Signature/Date

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