

Product Number:	103208
Lot Number:	031812
Expiration:	13-Nov-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.17×10^9 Beads/mL

Review/Approval

Junpil Park

QC, Supervisor

 14 Dec 2020

Name

Title

Signature/Date

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