

Product Number:	103208
Lot Number:	108518
Expiration:	26-Mar-2022
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.24 x 10 ⁹ Beads/mL

Review/Approval

Junpil Park
Name

QC *Supervisor*
Title

[Signature] 19 Apr 2021
Signature/Date

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