

Product Number:	103316
Lot Number:	110613
Expiration:	4-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.55 x10 ⁹ Beads/mL

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 19 Jul 2021
Signature/Date

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