

Product Number:	103612
Lot Number:	031523
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.08×10^9 Beads/mL

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 14 Dec 2020
Signature/Date

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