

Product Number:	103612
Lot Number:	031526
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.80×10^9 Beads/mL

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 15 Feb 2021
Signature/Date

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