

Product Number:	103612
Lot Number:	102531
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.14×10^9 Beads/mL

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 22 Feb 2021
Signature/Date

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