

Product Number:	103208
Lot Number:	102212
Expiration:	22-Jan-2022
Storage:	2±8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.34 x 10 ⁹ Beads/mL

Review/Approval

Junpil Park

QC, Supervisor



Name

Title

Signature/Date

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