

Product Number:	103208
Lot Number:	010716
Expiration:	16-Oct-2020
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.44x10 ⁹ Beads/mL

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 19 May 2020

Name

Title

Signature/Date

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