

Product Number:	103208
Lot Number:	134927
Expiration:	15-Dec-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.28 x 10 ⁹ Beads/mL

Review/Approval Shivani Goel QC Shivani Goel 17 Feb 2022
Name Title Signature/Date

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