

Product Number:	103611
Lot Number:	031521
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.17×10^9 Beads/mL

Review/Approval Shivani Goel QC Manager *Shivani Goel* 01 Dec 2020
Name Title Signature/Date

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