

Product Number:	100993
Lot Number:	018904
Expiration:	7-Jul-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	5.10x10 ⁶ Beads/mL

Review/Approval Shivani Goel QC Manager Shivani Goel 28 Jul 2020
 Name Title Signature/Date

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