

Product Number:	103208
Lot Number:	122921
Expiration:	17-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30 x 10 ⁹ Beads/mL

Review/Approval Shivani Goel QC Shivani Goel 15 Sep 2021
Name Title Signature/Date

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