

Product Number:	103612
Lot Number:	130716
Expiration:	5-Dec-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.22 x10 ⁹ Beads/mL

Review/Approval Shivani Gnel QC Shivani Gnel 20Dec2021
Name Title Signature/Date

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