

Product Number:	103208
Lot Number:	124513
Expiration:	2-Sep-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.201 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 13 Dec 2021
Signature/Date

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