

Product Number:	103527
Lot Number:	125611
Expiration:	13-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.29 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 18Jan2022
Signature/Date

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