

Product Number:	103527
Lot Number:	114117
Expiration:	21-May-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.36 x 10 ⁹ Beads/mL

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 27 Jul 2021
Signature/Date

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