

Product Number:	103527
Lot Number:	108825
Expiration:	29-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.20 x 10 ⁹ Beads/mL

Review/Approval

Jungil Park
Name

QC Supervisor
Title

 19 Apr 2021
Signature/Date

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