


Product Number:	103208
Lot Number:	032309
Expiration:	18-Nov-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.16x10 ⁹ Beads/mL

Review/Approval

Junpil Park	QC Supervisor		18 Dec 2020
Name	Title	Signature/Date	

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