

Product Number:	103207
Lot Number:	819702
Expiration:	30-Jan-2021
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.25x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date

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