

Product Number:	103316
Lot Number:	823907
Expiration:	15-Aug-2020
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.55 x 10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date



[Handwritten Signature] 28 Aug 2018

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