

Product Number:	101985
Lot Number:	820807
Expiration:	27-Jan-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 2/02/2018
Signature/Date

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