

Product Number:	101761
Lot Number:	830513
Expiration:	1-May-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.36x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor



Name

Title

Signature/Date

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