

Product Number:	101761
Lot Number:	830512
Expiration:	1-May-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.54×10^9 Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 09NOV2018
Signature/Date

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