

Product Number:	101732
Lot Number:	MF-013018-13
Expiration:	30-Jul-2018

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.31×10^9 Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 20 MAR 2018
Signature/Date

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