

Product Number:	101360
Lot Number:	MF-030218-03
Expiration:	22-Aug-2018
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.64×10^9 Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 06 MAR 2018

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