

Product Number:	101360
Lot Number:	MF-021518-02
Expiration:	22-Aug-2018
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.87×10^9 Beads/mL

Review/Approval

Sivasankari Murugan

Name

QC Supervisor

Title

 15 FEB 2018

Signature/Date

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