

Product Number:	101732
Lot Number:	MF-010818-14
Expiration:	8-Jul-2018

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.32x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 30 Jan 2018
Signature/Date

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