

Product Number:	103207
Lot Number:	MF-041018-11
Expiration:	7-Dec-2019
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	3.21×10^9 Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 12 APR 2018
Signature/Date

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