

Product Number:	103208
Lot Number:	MF-042418-08
Expiration:	24-Oct-2018
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.26x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 14 Jun 2018
Signature/Date

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