

Product Number:	101987
Lot Number:	MF-030118-06
Expiration:	1-Sep-2018
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beans/mL)	Multisizer	1.28 x 10 ⁹

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 14 MAY 2018
Signature/Date

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