

Product Number:	101986
Lot Number:	822708
Expiration:	15-Feb-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.40×10^9

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date

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