

Product Number:	101987
Lot Number:	825710
Expiration:	14-Mar-2019
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Sivasankari Murugan

QC Supervisor

 10DEC2018

Name

Title

Signature/Date

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