

Product Number:	101987
Lot Number:	817009
Expiration:	19-Dec-2018
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	1.29 x 10 ⁹

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 17 Oct 2018

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