

Product Number:	101986
Lot Number:	822707
Expiration:	15-Feb-2019
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Bead Concentration (beads/mL)	Multisizer	$1.31 \times 10^9$

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 12SEP2018  
Signature/Date

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