

Product Number:	101761
Lot Number:	819911
Expiration:	18-Jan-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.20x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 17 Oct 2018

Signature/Date

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