

Product Number:	101732
Lot Number:	803001
Expiration:	30-Jul-2018
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.27×10^9 Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date

 30 OCT 2018

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