

Product Number:	103317
Lot Number:	728502
Expiration:	27-Jan-2020
Storage Conditions:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.88×10^9 Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor



29 Oct 2018

Name

Title

Signature/Date

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