

Product Number:	103208
Lot Number:	829015
Expiration:	17-Apr-2019
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.30x10 ⁹ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 30 OCT 2018
Signature/Date

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