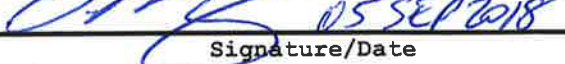


Product Number:	103208
Lot Number:	816902
Expiration:	18-Dec-2018
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.33x10 ⁹ Beads/mL

Review/Approval Sivasankari Murugan QC Supervisor  05 SEP 2018

Name **Title** **Signature/Date**

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