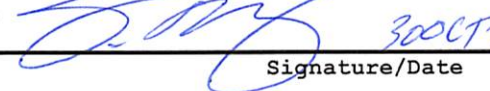


Product Number:	101986
Lot Number:	822709
Expiration:	15-Feb-2019
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Bead Concentration	Multisizer	1.32 x 10 ⁹

Review/Approval	Sivasankari Murugan	QC Supervisor	 30 Oct 2018
	Name	Title	Signature/Date

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