

Product Number:	103208
Lot Number:	825610
Expiration:	13-Mar-2019
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.24x10 ⁹ Beads/mL

On behalf of:
Sivasankari Murugan
QC Supervisor

Sr. Director,
Manufacturing
Operations

[Handwritten Signature]
12 Oct 2018

Review/Approval

Bud Drummey		
Name	Title	Signature/Date

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