

Product Number:	103207
Lot Number:	822906
Expiration:	30-Jan-2021
Storage Conditions:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$3.04 \times 10^9$ Beads/mL

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date

 29 AUG 2018

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