

Product Number:	103317
Lot Number:	729002
Expiration:	27-Jan-2020
Storage Conditions:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$2.51 \times 10^9$ Beads/mL

Review/Approval

Sivasankari Murugan

Name

QC Supervisor

Title

 17 OCT 2018

Signature/Date

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