

Product Number:	103207
Lot Number:	MF-042018-02
Expiration:	30-Jan-2021
Storage Conditions:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$3.23 \times 10^9$ Beads/mL

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 23 APR 2018  
Signature/Date

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