

Product Number:	101732
Lot Number:	MF-013018-17
Expiration:	30-Jul-2018
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	$1.25 \times 10^9$ Beads/mL

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 25 MAY 2018  
Signature/Date

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