

Product Number:	103208
Lot Number:	MF-042418-09
Expiration:	24-Oct-2018
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.25x10 <sup>9</sup> Beads/mL

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

 14 Oct 2018  
Signature/Date

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