


Product Number:	103317
Lot Number:	729001
Expiration:	27-Jan-2020
Storage Conditions:	2-8°C

**Laboratory Analysis**

Characteristic	Means of Analysis	Result
Concentration	Multisizer	2.41 x 10 <sup>9</sup> Beads/mL

Review/Approval      Sivasankari Murugan      QC Supervisor            26 SEP 2018

---

Name      Title      Signature/Date

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