

| | |
|------------------------|-------------|
| Product Number: | 101985 |
| Lot Number: | 820806 |
| Expiration: | 27-Jan-2019 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|-----------------------|--------------------------|-------------------------------|
| Concentration | Multisizer | 1.22x10 ⁹ Beads/mL |


Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 22 Aug 2018

Signature/Date

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