

| | |
|---------------------|-------------|
| Product Number: | 103317 |
| Lot Number: | 823908 |
| Expiration: | 27-Jan-2020 |
| Storage Conditions: | 2-8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-----------------------------|
| Concentration | Multisizer | 2.63×10^9 Beads/mL |

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 28 Aug 2018
Signature/Date

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