

| | |
|-----------------|--------------|
| Product Number: | 101761 |
| Lot Number: | MF-072717-10 |
| Expiration: | 16-Oct-2018 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-------------------------------|
| Concentration | Multisizer | 1.38x10 ⁹ Beads/mL |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date

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