

| | |
|---------------------|--------------|
| Product Number: | 101360 |
| Lot Number: | MF-021218-05 |
| Expiration: | 22-Aug-2018 |
| Storage Conditions: | 2-8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-----------------------------|
| Concentration | Multisizer | 2.38×10^9 Beads/mL |

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

Signature/Date

 12 FEB 2018

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