

| | |
|-----------------|------------|
| Product Number: | 103612 |
| Lot Number: | 015310 |
| Expiration: | 5-Dec-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-----------------------------|
| Concentration | Multisizer | 3.36×10^9 Beads/mL |

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 15 Jun 2020

Name

Title

Signature/Date

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