

| | |
|-----------------|------------|
| Product Number: | 103612 |
| Lot Number: | 130717 |
| Expiration: | 5-Dec-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|--------------------------------|
| Concentration | Multisizer | 3.23 x10 ⁹ Beads/mL |

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 20Dec2021
Signature/Date

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