

| | |
|------------------------|------------|
| Product Number: | 103611 |
| Lot Number: | 108822 |
| Expiration: | 5-Dec-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-------------------------------|
| Concentration | Multisizer | 2.49x10 ⁹ Beads/mL |

Review/Approval Shivani Goel QC Manager Shivani Goel 14 Apr 2021
 Name Title Signature/Date

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