

| | |
|-----------------|------------|
| Product Number: | 103208 |
| Lot Number: | 125016 |
| Expiration: | 7-Sep-2022 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|---------------------------------|
| Concentration | Multisizer | 1.43 x 10 ⁹ Beads/mL |

Review/Approval Shivani Goel QC Shivani Goel 29 NOV 2021
Name Title Signature/Date

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