

| | |
|------------------------|-------------|
| Product Number: | 103527 |
| Lot Number: | 123804 |
| Expiration: | 26-Aug-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|-----------------------|--------------------------|---------------------------------|
| Concentration | Multisizer | 1.18 x 10 ⁹ Beads/mL |

Review/Approval Shivani Goel QC Shivani Goel 15 Sep 2021
 Name Title Signature/Date

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