

| | |
|------------------------|-------------|
| Product Number: | 103208 |
| Lot Number: | 005113 |
| Expiration: | 20-Feb-2021 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-------------------------------|
| Concentration | Multisizer | 1.20x10 ⁹ Beads/mL |

Review/Approval

| | | | |
|------------------|---------------------|-------------------|-------------|
| Patrizia Stadler | Assoc. Director, QC | <i>p. Stadler</i> | 04 Feb 2020 |
| Name | Title | Signature/Date | |

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