

| | |
|------------------------|------------|
| Product Number: | 103527 |
| Lot Number: | 015308 |
| Expiration: | 1-Jun-2021 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|---------------------------------|
| Concentration | Multisizer | 1.10 x 10 ⁹ Beads/mL |

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

11 Jun 2020

Name

Title

Signature/Date

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