

| | |
|-----------------|------------|
| Product Number: | 103612 |
| Lot Number: | 130720 |
| Expiration: | 5-Jun-2025 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|-----------------------------|
| Concentration | Multisizer | 3.27×10^9 Beads/mL |

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan

17 Oct 2022

Signature/Date