

| | |
|-----------------|------------|
| Product Number: | 103612 |
| Lot Number: | 230419 |
| Expiration: | 8-Dec-2026 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|----------------|-------------------|---------------------------------|
| Concentration | Multisizer | 3.15 x 10 ⁹ Beads/mL |

Review/Approval

Sonia Khilka

Name

QC

Title

Soni Khilka 29 NOV 2023

Signature/Date