

| | |
|------------------------|-------------|
| Product Number: | 103528 |
| Lot Number: | 126313 |
| Expiration: | 20-Sep-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Means of Analysis | Result |
|-----------------------|--------------------------|---------------------------------|
| Concentration | Multisizer | 1.23 x 10 ⁹ Beads/mL |

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 01 Apr 2022
Signature/Date