

| | |
|------------------------|-------------|
| Product Number: | 103358 |
| Lot Number: | 108516 |
| Expiration: | 26-Mar-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|-----------------------|---------------|------------------|
| pH | Pass | 7.95-8.05 |

Review/Approval

Shivani Goel
Name

Quality Control
Title

Shivani Goel 14 Apr
Signature/Date 2021