

| | |
|-----------------|------------|
| Product Number: | 103847 |
| Lot Number: | 512713 |
| Expiration: | 7-Nov-2026 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target | PASS / FAIL |
|----------------|--------|-------------|-------------|
| pH | 7.41 | 7.35 - 7.45 | PASS |

Review/Approval

Lady Carlson
Name

QC Manager
Title

[Signature] 21 MAY 2025
Signature/Date