

| | |
|-----------------|-------------|
| Product Number: | 101362 |
| Lot Number: | 512008 |
| Expiration: | 30-Oct-2026 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | pH Target | PASS / FAIL |
|----------------|-------------|-------------|
| pH Measurement | 7.75 - 7.85 | PASS |

RESULT

| | |
|------------|------|
| pH Reading | 7.83 |
|------------|------|

Review/Approval

Heidi Caswell

Director, QC

Heidi Caswell 07JUL2025