

| | |
|-----------------|-------------|
| Product Number: | 101359 |
| Lot Number: | 416317 |
| Expiration: | 11-Dec-2025 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | pH Target | PASS / FAIL |
|----------------|-------------|-------------|
| pH Measurement | 7.35 - 7.45 | PASS |

RESULT

| | |
|------------|------|
| pH Reading | 7.39 |
|------------|------|

Review/Approval

Robert Somogyi

QC Manager

 26 JUN 2024

Name

Title

Signature/Date