

| | |
|-----------------|-------------|
| Product Number: | 101358 |
| Lot Number: | 510808 |
| Expiration: | 18-Oct-2026 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-------------|
| pH | Pass | 7.35 - 7.45 |

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

 29 APR 2025