

| | |
|-----------------|-------------|
| Product Number: | 101355 |
| Lot Number: | 311602 |
| Control Number: | 311602A |
| Expiration: | 26-Apr-2024 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-------------|
| pH Measurement | PASS | 7.35 - 7.45 |

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 10 MAY 2023
Signature/Date