

| | |
|------------------------|-------------|
| Product Number: | 103360 |
| Lot Number: | 322806 |
| Expiration: | 16-Aug-2024 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target | PASS / FAIL |
|-----------------------|---------------|------------------|--------------------|
| pH | 8.05 | 7.95 - 8.05 | PASS |

Review/Approval

Sonia Khillan

Name

QC manager

Title

Sonia Khillan 30AUG2023

Signature/Date