

| | |
|------------------------|-------------|
| Product Number: | 103847 |
| Lot Number: | 226219 |
| Expiration: | 19-Sep-2023 |
| Storage: | 2 - 8°C |

Laboratory Analysis

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

Review/Approval

Sonia Khillan QC Manager Sonia Khillan / 13 DEC 2022

Name

Title

Signature/Date