

| | |
|-----------------|-------------|
| Product Number: | 101631 |
| Lot Number: | 208109 |
| Expiration: | 22-Mar-2023 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-----------|
| pH | Pass | 7.30-7.50 |

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 30 Mar 2022

Name

Title

Signature/Date