

| | |
|------------------------|-------------|
| Product Number: | 101476 |
| Lot Number: | 121110 |
| Expiration: | 30-Jul-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|-----------------------|---------------|------------------|
| pH | Pass | 8.95-9.05 |

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 03 Aug 2021
Signature/Date