

| | |
|------------------------|-------------|
| Product Number: | 101357 |
| Lot Number: | 210809 |
| Expiration: | 18-Apr-2023 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|-----------------------|---------------|------------------|
| pH Measurement | PASS | 6.18 - 6.22 |

• Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 20 Apr 2022
Signature/Date