

| | |
|------------------------|-------------|
| Product Number: | 101631 |
| Lot Number: | 105206 |
| Expiration: | 21-Feb-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

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

Review/Approval

Tunpil Park
Name

QC Supervisor
Title

[Signature] 23 Apr 2021
Signature/Date