

| | |
|-----------------|------------|
| Product Number: | 101355 |
| Lot Number: | 227908 |
| Expiration: | 6-Oct-2023 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-------------|
| pH Measurement | PASS | 7.35 - 7.45 |

Review/Approval

Matt Koudouris

Name

QC Supervisor

Title

M. Koudouris 10 Oct 2022

Signature/Date