

| | |
|-----------------|-------------|
| Product Number: | 101474 |
| Lot Number: | 227106 |
| Expiration: | 28-Sep-2023 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-----------|
| pH | Pass | 7.35-7.45 |

Review/Approval

Matt Kaelouris
Name

QC Supervisor
Title

[Signature] 10 OCT 2022
Signature/Date