

| | |
|-----------------|-------------|
| Product Number: | 102912 |
| Lot Number: | 207611 |
| Expiration: | 17-Mar-2023 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|----------------|--------|-------------|
| pH | Pass | 7.95 - 8.05 |

Review/Approval

Matt Smart
Name

QC
Title

MRA 10/19/2022
Signature/Date