

| | |
|------------------------|-------------|
| Product Number: | 101955 |
| Lot Number: | 118021 |
| Expiration: | 29-Jun-2022 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result |
|-----------------------|---------------|
| pH Measurement | Pass |

Review/Approval

Jupil Park

Name

Quality Control

Title

Supervisor  *08 Jul 2021*

Signature/Date