

| | |
|------------------------|------------|
| Product Number: | 100458 |
| Lot Number: | 118710 |
| Expiration: | 6-Jul-2022 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result |
|-----------------------|---------------|
| pH Measurement | PASS |

Review/Approval

Jungil Park

Name

QC Supervisor

Title

[Signature] 30 Jul 2021

Signature/Date