

| | |
|------------------------|-------------|
| Product Number: | 100376 |
| Lot Number: | 104729 |
| Expiration: | 16-Feb-2022 |
| Storage: | 2-8°C |

Laboratory Analysis

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

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 18 Mar 2021

Signature/Date