

| | |
|------------------------|------------|
| Product Number: | 101631 |
| Lot Number: | 028207 |
| Expiration: | 8-Oct-2021 |
| Storage: | 2 - 8°C |

Laboratory Analysis

| Characteristic | Result | pH Target |
|-----------------------|---------------|------------------|
| pH | Pass | 6.8-7.8 |

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 09 Oct 2020
Signature/Date