

| | |
|------------------------|-------------|
| Product Number: | 100458 |
| Lot Number: | 134609 |
| Expiration: | 12-Dec-2022 |
| Storage: | 2-8°C |

Laboratory Analysis

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

Review/Approval Shivani Goel QC Shivani Goel 20Dec2021
Name Title Signature/Date