

|                 |             |
|-----------------|-------------|
| Product Number: | 101473      |
| Lot Number:     | 107724      |
| Expiration:     | 18-Mar-2022 |
| Storage:        | 2 - 8°C     |

**Laboratory Analysis**

| Characteristic | Result | pH Target |
|----------------|--------|-----------|
| pH             | Pass   | 7.35-7.45 |

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 24 May 2021

Signature/Date