

|                 |             |
|-----------------|-------------|
| Product Number: | 101472      |
| Lot Number:     | 204616      |
| Expiration:     | 15-Feb-2023 |
| Storage:        | 2 - 8°C     |

**Laboratory Analysis**

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

Review/Approval

Richa Shrestha

QC Supervisor

Name

Title

 29A702022  
Signature/Date