

|                 |             |
|-----------------|-------------|
| Product Number: | 103847      |
| Lot Number:     | 404319      |
| Expiration:     | 12-Aug-2025 |
| Storage:        | 2 - 8°C     |

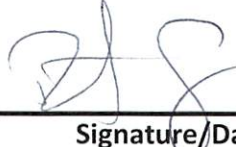
**Laboratory Analysis**

| Characteristic | Result | pH Target   | PASS / FAIL |
|----------------|--------|-------------|-------------|
| pH             | 7.36   | 7.35 - 7.45 | PASS        |

Review/Approval

ROBERT SOMOGYI  
Name

QC MANAGER  
Title

  
Signature/Date  
11 APR 2024