

|                 |            |
|-----------------|------------|
| Product Number: | 103359     |
| Lot Number:     | 403707     |
| Expiration:     | 6-Feb-2025 |
| Storage:        | 2 - 8°C    |

**Laboratory Analysis**

| Characteristic | Result | pH Target   | PASS / FAIL |
|----------------|--------|-------------|-------------|
| pH             | 8.02   | 7.95 - 8.05 | PASS        |

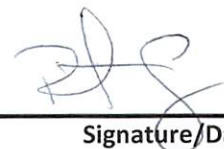
Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title



Signature/Date

14 FEB 2024