

|                 |            |
|-----------------|------------|
| Product Number: | 103360     |
| Lot Number:     | 431102     |
| Expiration:     | 6-Nov-2025 |
| Storage:        | 2-8°C      |

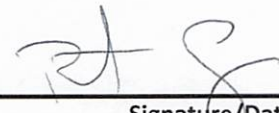
Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager



19NOV2024

Name

Title

Signature/Date