

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101556      |
| <b>Lot Number:</b>     | 423203      |
| <b>Expiration:</b>     | 19-Feb-2026 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| <b>Characteristic</b> | <b>Result</b> | <b>pH Target</b> | <b>PASS / FAIL</b> |
|-----------------------|---------------|------------------|--------------------|
| pH                    | 7.42          | 7.35 - 7.45      | PASS               |

**Review/Approval**

**Robert Somogyi**

**QC Manager**

 23 AUG 2024

**Name**

**Title**

**Signature/Date**