

|                 |            |
|-----------------|------------|
| Product Number: | 101359     |
| Lot Number:     | 227718     |
| Expiration:     | 4-Oct-2023 |
| Storage:        | 2 - 8°C    |

**Laboratory Analysis**

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

Review/Approval

*Sonia Khillan*

Name

*QC Manager*

Title

*Sonia 12/11/27 OCT 2022*

Signature/Date