

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101476      |
| <b>Lot Number:</b>     | 401608      |
| <b>Expiration:</b>     | 16-Jan-2025 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| <b>Characteristic</b> | <b>Result</b> | <b>pH Target</b> | <b>PASS / FAIL</b> |
|-----------------------|---------------|------------------|--------------------|
| pH                    | 8.95          | 8.95 - 9.05      | PASS               |

Review/Approval

*Cody Carlson*  
Name

*QC Manager*  
Title

*[Signature]* 18JAN2024  
Signature/Date