

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 102620      |
| <b>Lot Number:</b>     | 030012      |
| <b>Expiration:</b>     | 26-Oct-2021 |
| <b>Storage:</b>        | 2 - 8°C     |

**Laboratory Analysis**

| <b>Characteristic</b> | <b>Result</b> | <b>pH Target</b> |
|-----------------------|---------------|------------------|
| pH                    | Pass          | 7.35-7.45        |

Review/Approval

Junpil Park  
Name

Supervisor, Quality Control  
Title

 09 Dec 2020  
Signature/Date