

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101474      |
| <b>Lot Number:</b>     | 214901      |
| <b>Expiration:</b>     | 29-May-2023 |
| <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

Matt Smart  
Name

QC  
Title

M. Smart 06 Jun 2022  
Signature/Date