

<b>Product Number:</b>	101358
<b>Lot Number:</b>	220915
<b>Expiration:</b>	28-Jul-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Sonia Khillan      QC Manager      Sonia Khillan / 06 SEP 2022  
Name                                  Title                                  Signature/Date