

<b>Product Number:</b>	101473
<b>Lot Number:</b>	215103
<b>Expiration:</b>	31-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

MSA 03 Jun 2022  
Signature/Date