

<b>Product Number:</b>	104101
<b>Lot Number:</b>	408112
<b>Expiration:</b>	21-Sep-2025
<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

Robert Somogyi  
Name

QC Manager  
Title

 27 JUN 2024  
Signature/Date