

<b>Product Number:</b>	103847
<b>Lot Number:</b>	409216
<b>Expiration:</b>	1-Oct-2025
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH	7.41	7.35 - 7.45	PASS

**Review/Approval**

Robert Somogyi

QC Manager

 18 JUL 2024

**Name**

**Title**

**Signature/Date**