

<b>Product Number:</b>	101556
<b>Lot Number:</b>	328317
<b>Expiration:</b>	10-Oct-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH	7.36	7.35 - 7.45	PASS

**Review/Approval**

*Sonia Khillan*

**Name**

*Senior QC Analyst*

**Title**

*Sonia Khillan 06 NOV 2023*

**Signature/Date**