

<b>Product Number:</b>	103847
<b>Lot Number:</b>	328304
<b>Expiration:</b>	10-Oct-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

**Review/Approval**

*Sonia Khillan*

**Name**

*QC Manager*

**Title**

*Sonia Khillan 13 Oct 2023*

**Signature/Date**