




<b>Product Number:</b>	101631
<b>Lot Number:</b>	325511
<b>Expiration:</b>	12-Sep-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.30 - 7.50	PASS

Review/Approval

 Name	QC  SUPERVISOR Title	 Signature/Date	27 SEP 2023
---	--	---	-------------