

<b>Product Number:</b>	101475
<b>Lot Number:</b>	404416
<b>Expiration:</b>	13-Aug-2025
<b>Storage:</b>	2 - 8°C

Laboratory Analysis

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

Review/Approval

<i>Cody Carlson</i>	QC Manager	 23 FEB 2024
Name	Title	Signature/Date