

Product Number:	101955
Lot Number:	511811
Expiration:	28-Oct-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.41	7.30 - 7.50	PASS

Review/Approval

<u>Heidi Crowell</u>	<u>Director, QC</u>	<u>Heidi Crowell 08/29/25</u>
Name	Title	Signature/Date