

Product Number:	101359
Lot Number:	422026
Expiration:	7-Feb-2026
Storage:	2-8°C

Laboratory Analysis		
Characteristic	pH Target	PASS / FAIL
pH Measurement	7.30- 7.50	PASS

RESULT	
pH Reading	7.43

Review/Approval

Heidi Caswell

Director, QC

 15 Apr 2025

Name

Title

Signature/Date