

Product Number:	101472
Lot Number:	514711
Expiration:	27-Nov-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.44	7.35 - 7.45	PASS

Review/Approval

Greg Simone

Name

Quality Events and Training Specialist

Title

[Signature] 09 Oct 2025

Signature/Date