

Product Number:	103847
Lot Number:	524605
Expiration:	3-Mar-2027
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Heidi Caswell

Name

Director, QC

Title

Heidi Caswell 08 Sep 2025

Signature/Date