

Product Number:	101353
Lot Number:	426016
Expiration:	16-Mar-2026
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.35 - 7.45	PASS

Review/Approval

*Heidi Caswell*

*Director, QC*

*H/Caswell 11Feb2025*

Name

Title

Signature/Date