

Product Number:	101475
Lot Number:	510607
Expiration:	16-Oct-2026
Storage:	2 - 8°C

Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

 24 APR 2025