

Product Number:	101474
Lot Number:	526904
Expiration:	26-Sep-2026
Storage:	2-8°C

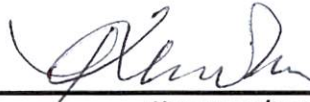
Laboratory Analysis		
Characteristic	pH Target	PASS / FAIL
pH Measurement	7.35 - 7.45	PASS

RESULT	
pH Reading	7.43

Review/Approval

Kathleen Cornell

Quality Control Supervisor

 30 Sep 2025

Name

Title

Signature/Date