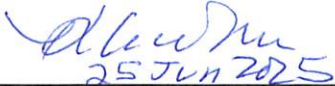


Product Number:	101359
Lot Number:	516201
Expiration:	11-Dec-2026
Storage:	2-8°C

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

RESULT	
pH Reading	7.42

Review/Approval

<u>Kathleen Cornell</u>	<u>Quality Control Supervisor</u>	<u> 25 Jun 2025</u>
Name	Title	Signature/Date