

Product Number:	101474
Lot Number:	521712
Expiration:	5-Aug-2026
Storage:	2-8°C

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


RESULT	
pH Reading	7.39

Review/Approval

Kathleen Cornell

Quality Control Supervisor

NameTitle

 18 Aug 2025

Signature/Date