

Product Number:	101359
Lot Number:	330621
Expiration:	2-Nov-2024
Storage:	2-8°C

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

RESULT	
pH Reading	7.41

Review/Approval

Sonia Khillan

Senior QC Analyst

Sonia 12/1/23 14 Nov 2023

Name

Title

Signature/Date