

Product Number:	101359
Lot Number:	408515
Expiration:	25-Sep-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.40
------------	------

Review/Approval

Cody Carlson

QC Manager

[Signature] 29/MAR/2024

Name

Title

Signature/Date