

Product Number:	101359
Lot Number:	506319
Expiration:	4-Sep-2026
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

Robert Somogyi

QC Manager

Name

Title

 29 APR 2025

Signature/Date