

Product Number:	101359
Lot Number:	416319
Expiration:	11-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.30 - 7.50	PASS

RESULT

pH Reading	7.41
------------	------

Review/Approval

Robert Somogyi

QC Manager

 24 SEP 2024

Name

Title

Signature/Date