

Product Number:	101474
Lot Number:	429508
Expiration:	21-Oct-2025
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

31 OCT 2024

Name

Title

Signature/Date