

Product Number:	101474
Lot Number:	425605
Expiration:	12-Sep-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.45
------------	------

Review/Approval

Robert Somogyi

QC Manager

24SEP2024

Name

Title

Signature/Date