

Product Number:	101359
Lot Number:	423219
Expiration:	19-Feb-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.42
------------	------

Review/Approval

Robert Somogyi

QC Manager



29 AUG 2024

Name

Title

Signature/Date