

Product Number:	101362
Lot Number:	434502
Expiration:	10-Jun-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS


RESULT

pH Reading	7.84
------------	------

Review/Approval

Robert Somogyi

QC Manager



17 Dec 2024

Name

Title

Signature/Date