

Product Number:	101362
Lot Number:	423609
Expiration:	23-Feb-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.78
------------	------

Review/Approval

Robert Somogyi

QC Manager



29 Aug 2024

Name

Title

Signature/Date