

Product Number:	101362
Lot Number:	502410
Control Lot Number:	502410A
Expiration:	24-Jul-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.79
------------	------

Review/Approval

Heidi Caswell

Name

Director, QC

Title

H. Caswell 11 Feb 2025

Signature/Date