

Product Number:	101362
Lot Number:	509803
Expiration:	8-Oct-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.81
------------	------

Review/Approval

Cody Carlson
Name

QC Manager
Title

 09 Apr 2025
Signature/Date