

Product Number:	101362
Lot Number:	532911
Expiration:	25-May-2027
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

Kathleen
Cornell

Quality Control
Supervisor

 09 Dec 2025

Name

Title

Signature/Date