

Product Number:	101474
Lot Number:	531410
Expiration:	10-Nov-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.35 - 7.45	PASS

RESULT

pH Reading	7.35
------------	------

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


19 Nov 2025
Signature/Date