

Product Number:	101359
Lot Number:	523823
Expiration:	26-Feb-2027
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.41
------------	------

Review/Approval

Heidi Caswell
Name

Director, QC
Title

Heidi Caswell 08 Sep 2025
Signature/Date