

Product Number:	101359
Lot Number:	426013
Expiration:	16-Mar-2026
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*
*Director, QC*
*H. Caswell 26 Nov 2024*

---

Name
Title
Signature/Date