

Product Number:	100376
Lot Number:	424714
Expiration:	3-Mar-2026
Storage:	2-8°C



Laboratory Analysis

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

RESULT

pH Reading	7.38
------------	------

Review/Approval

---

Name: ROBERT SOMOGYI      Title: QC MANAGER      Signature/Date: 16 SEP 2024