

Product Number:	100376
Lot Number:	433720
Expiration:	2-Jun-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

Robert Somogyi

QC Manager



11/25/2024

Name

Title

Signature/Date