

Product Number:	100376
Lot Number:	510511
Control Lot Number:	510511A
Expiration:	15-Oct-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

Name

Title

Signature/Date

 18 APR 2025