

Product Number:	100376
Lot Number:	506219
Control Lot Number:	506219A
Expiration:	3-Sep-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

Robert Somogyi
Name

QC Manager
Title

 05 MAR 2025
Signature/Date