

Product Number:	100376
Lot Number:	501523
Control Lot Number:	501523A
Expiration:	15-Jul-2026
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.39
------------	------

Review/Approval

Robert Somogyi

QC Manager

Robert Somogyi 29 JAN 2025

Name

Title

Signature/Date