

Product Number:	101359
Lot Number:	402216
Expiration:	22-Jul-2025
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



Signature/Date

25 JAN 2024