

Product Number:	100376
Lot Number:	402514
Expiration:	25-Jul-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.41
------------	------

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 14 FEB 2024

Signature/Date