

Product Number:	100376
Lot Number:	404318
Expiration:	12-Aug-2025
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.43
------------	------

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 29 MAR 2024

Signature/Date