

Product Number:	100376
Lot Number:	407412
Expiration:	14-Sep-2025
Storage:	2-8°C

Laboratory Analysis

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

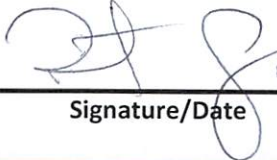
RESULT

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

Review/Approval

ROBERT SOMOSYI

QC MANAGER

 10 APR 2024

Name

Title

Signature/Date