

Product Number:	101362
Lot Number:	413608
Expiration:	15-Nov-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS

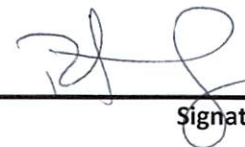
RESULT

pH Reading	7.85
------------	------

Review/Approval

Robert Somogyi

QC Manager



23 MAY 2024

Name

Title

Signature/Date