

Product Number:	101359
Lot Number:	416318
Expiration:	11-Dec-2025
Storage:	2-8°C

Laboratory Analysis

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

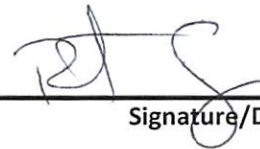
RESULT

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

Review/Approval

Robert Somogyi

QC Manager



09SEP2024

Name

Title

Signature/Date