

Product Number:	101359
Lot Number:	327711
Expiration:	4-Oct-2024
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.38
------------	------

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 11 Oct 2023

Signature/Date