

Product Number:	101359
Lot Number:	325612
Expiration:	13-Sep-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 Khillar
Name

QC Manager
Title

Sonia Khillar 03 OCT 2023
Signature/Date