

Product Number:	100376
Lot Number:	330321
Expiration:	30-Oct-2024
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

Sonia Khillan

Name

Senior QC Analyst

Title

Sonia Khillan 14 NOV 2023

Signature/Date