

Product Number:	101362
Lot Number:	327614
Expiration:	3-Oct-2024
Storage:	2-8°C

Laboratory Analysis

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

RESULT

pH Reading	7.75
------------	------

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 14/NOV2023
Signature/Date