

Product Number:	101356
Lot Number:	328412
Expiration:	11-Oct-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 17 OCT 2023

Signature/Date