

Product Number:	101352
Lot Number:	323415
Expiration:	22-Aug-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.35 - 7.45	PASS

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 01SEP2023

Signature/Date