

Product Number:	101473
Lot Number:	308223
Expiration:	23-Mar-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

Senior QC Analyst

Sonia Khillan 30 OCT 2023

Name

Title

Signature/Date