

Product Number:	101362
Lot Number:	323616
Expiration:	24-Aug-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.79	7.75 - 7.85	PASS

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 30 Aug 2023

Signature/Date