

Product Number:	103358
Lot Number:	323313
Expiration:	21-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.05	7.95 - 8.05	PASS

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 30 Aug 2023

Signature/Date