

Product Number:	101476
Lot Number:	328612
Expiration:	13-Oct-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.95	8.95 - 9.05	PASS

Review/Approval

Sonia Khillan

Senior QC Analyst

Sonia Khillan 20OCT2023

Name

Title

Signature/Date