

Product Number:	103847
Lot Number:	303309
Expiration:	2-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan / 27 FEB 2023

Name

Title

Signature/Date