

Product Number:	103847
Lot Number:	300606
Expiration:	6-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan / 17 JAN 2023

Signature/Date