

Product Number:	103847
Lot Number:	223811
Expiration:	26-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC manager

Sonia Khillan / 20 Aug 2023

Name

Title

Signature/Date