

Product Number:	102914
Lot Number:	232003
Expiration:	16-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia/28/16 30NOV2022

Signature/Date