

Product Number:	101631
Lot Number:	230418
Expiration:	31-Oct-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.30 - 7.50

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 14 DEC 2022
Signature/Date