

Product Number:	101473
Lot Number:	230115
Expiration:	28-Oct-2023
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 / 19 DEC 2022
Signature/Date