

Product Number:	103704
Lot Number:	227802
Expiration:	5-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 / 02 Nov 2022
Signature/Date