

Product Number:	103360
Lot Number:	221704
Expiration:	5-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Kella / 10AUG2022
Signature/Date