

Product Number:	102908
Lot Number:	223608
Expiration:	24-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 Khillan / 25AUG2022*

Signature/Date