

Product Number:	103359
Lot Number:	218913
Expiration:	8-Jul-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval

*Sonia Khillan*

Name

QC

Title

*Sonia Khillan 12/5/2022*

Signature/Date