

Product Number:	102910
Lot Number:	218102
Expiration:	30-Jun-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 07JUL2022
Signature/Date