

Product Number:	102910
Lot Number:	223612
Expiration:	24-Aug-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 129 AUG 2022

Signature/Date