

Product Number:	101362
Lot Number:	215605
Expiration:	5-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan 21 JUL 2022

Signature/Date