

Product Number:	101359
Lot Number:	218620
Expiration:	5-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan / 09 AUG 2020

Signature/Date