

Product Number:	101353
Lot Number:	218619
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 / 08 AUG 2022

Signature/Date