

Product Number:	101353
Lot Number:	231914
Expiration:	15-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Name

Title

Signature/Date

Sonia Kellan *QC Manager* *Sonia Kellan* *30 NOV 2022*