

Product Number:	101353
Lot Number:	303216
Expiration:	1-Feb-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC Manager

Soni 12/16 13FEB2023

Name

Title

Signature/Date