

Product Number:	103172
Lot Number:	417211
Expiration:	20-Jun-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager



20 NOV 2024

Name

Title

Signature/Date