

Product Number:	102908
Lot Number:	404703
Expiration:	16-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 20FEB2024

Signature/Date