

Product Number:	103358
Lot Number:	404704
Expiration:	16-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.05	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 20FEB2024

Signature/Date