

Product Number:	100460
Lot Number:	401708
Expiration:	17-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 08 FEB 2024

Signature/Date