

Product Number:	100460
Lot Number:	416316
Expiration:	11-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager

 24 JUN 2024

Name

Title

Signature/Date