

Product Number:	100458
Lot Number:	412312
Expiration:	2-Nov-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Robert Somogyi

QC Manager

 17 MAY 2024

Name

Title

Signature/Date