

Product Number:	100458
Lot Number:	414211
Expiration:	21-Nov-2025
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

*Thanh Vu* QC Supervisor

Name

Title

*[Signature]* 20JUN2024

Signature/Date