

Product Number:	100458
Lot Number:	404321
Expiration:	12-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Binh Vu

QC Supervisor

[Signature] 29 APR 2024

Name

Title

Signature/Date