

Product Number:	101355
Lot Number:	408806
Expiration:	28-Sep-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.36	7.35 - 7.45	PASS

Review/Approval

Binh Vu
Name

QC supervisor
Title

[Signature] 10 APR 2024
Signature/Date