

Product Number:	101954
Lot Number:	408504
Expiration:	25-Sep-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.84	7.75 - 7.85	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 27 MAR 2024

Name

Title

Signature/Date