

Product Number:	101954
Lot Number:	507616
Expiration:	17-Sep-2026
Storage:	2-8°C

Laboratory Analysis

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

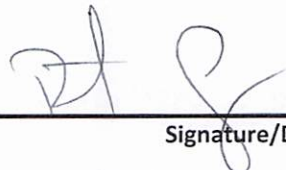
Review/Approval

Robert Somogyi

QC Manager

Name

Title


Signature/Date