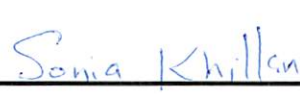

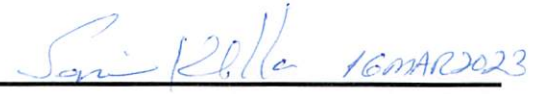


Product Number:	101362
Lot Number:	305413
Expiration:	23-Feb-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Name

Title

Signature/Date