

Product Number:	101362
Lot Number:	506312
Expiration:	4-Sep-2026
Storage:	2-8°C

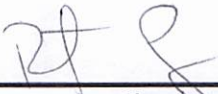
Laboratory Analysis		
Characteristic	pH Target	PASS / FAIL
pH Measurement	7.75 - 7.85	PASS

RESULT	
pH Reading	7.81

Review/Approval

Robert Somogyi

QC Manager

 10 MAR 2025

NameTitleSignature/Date