

Product Number:	101631
Lot Number:	429902
Expiration:	25-Oct-2025
Storage:	2 - 8°C

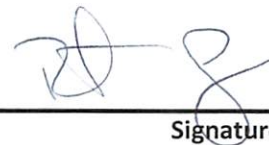
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.33	7.30 - 7.50	PASS

Review/Approval

Robert Somogyi

QC Manager



12 NOV 2024

Name

Title

Signature/Date