

Product Number:	103847
Lot Number:	427010
Expiration:	26-Mar-2026
Storage:	2 - 8°C

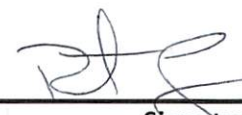
**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi

QC Manager

 09 Oct 2024

Name

Title

Signature/Date