

Product Number:	103847
Lot Number:	431107
Expiration:	6-May-2026
Storage:	2 - 8°C

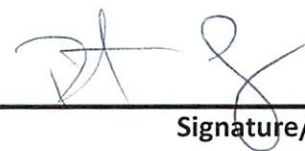
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.37	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi

QC Manager



21 JUN 2024

Name

Title

Signature/Date