

Product Number:	103158
Lot Number:	431612
Expiration:	11-May-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Robert Somogyi

QC Manager

Name

Title

 19 Nov 2024
Signature/Date