

Product Number:	102910
Lot Number:	433719
Expiration:	2-Dec-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.03	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

Name

Title

Signature/Date

 05 DEC 2024