

Product Number:	102914
Lot Number:	434010
Expiration:	5-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.96	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager



Name

Title

Signature/Date