

Product Number:	102912
Lot Number:	434017
Expiration:	5-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 14 MAR 2025

Name

Title

Signature/Date