

Product Number:	102912
Lot Number:	417608
Expiration:	24-Jun-2025
Storage:	2-8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.97	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 27 JUN 2024

Name

Title

Signature/Date