

Product Number:	102908
Lot Number:	412401
Expiration:	3-May-2025
Storage:	2 - 8°C

**Laboratory Analysis**

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

Review/Approval


Robert Somogyi

QC Manager

Name

Title

Signature/Date

 15 MAY 2024