

Product Number:	102908
Lot Number:	433704
Expiration:	2-Dec-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

Manager, Quality Control

Name

Title

 09 DEC 2024

Signature/Date