

Product Number:	103358
Lot Number:	433705
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

 09 Dec 2024

Name

Title

Signature/Date