

Product Number:	101476
Lot Number:	404609
Expiration:	15-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.98	8.95 - 9.05	PASS

Review/Approval

Cody Carlson
Name

QC Manager
Title

Cody Carlson 23 FEB 2024
Signature/Date