

Product Number:	101472
Lot Number:	402615
Expiration:	26-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval

Cody Carlson
Name

QC Manager
Title

 22 FEB 2024
Signature/Date