

Product Number:	100453
Lot Number:	416608
Expiration:	14-Dec-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.21	6.18 - 6.22	PASS

Review/Approval

Cody Carlson  
Name

QC Manager  
Title

[Signature] 18 JUN 2024  
Signature/Date