

Product Number:	100458
Lot Number:	514005
Expiration:	20-Nov-2026
Storage:	2-8°C

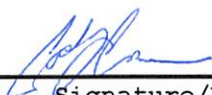
Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Cody Carlson
Name

QC Manager
Title

 29 MAY 2025
Signature/Date