

Product Number:	103158
Lot Number:	519817
Expiration:	17-Jan-2027
Storage:	2-8°C

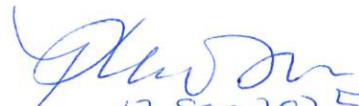
Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.40	7.35 - 7.45	PASS

Review/Approval

Kathleen Cornell
Name

Quality Control
Supervisor
Title


12 Sep 2025
Signature/Date