

Product Number:	101352
Lot Number:	521010
Expiration:	29-Jan-2027
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell
Name

*Quality Control
Supervisor*
Title

[Signature]
19 Aug 2025
Signature/Date