

Product Number:	100453
Lot Number:	509411
Expiration:	4-Oct-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Kathleen Cornell
Name

*Quality Control
Supervisor*
Title

Kathleen Cornell 23 May 2025
Signature/Date