

Product Number:	100453
Lot Number:	528907
Expiration:	16-Apr-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	6.20	6.18 - 6.22	PASS

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature]
20 Nov 2025

Signature/Date