

Product Number:	103358
Lot Number:	520306
Expiration:	22-Jul-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.00	7.95 - 8.05	PASS

Review/Approval

Kathleen Cornell

Name

*Quality Control
Supervisor*

Title

[Signature]
04 Aug 2025

Signature/Date