

Product Number:	101355
Lot Number:	501713
Control Lot Number:	501713A
Expiration:	17-Jul-2026
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Heidi Caswell
Name

Director, QC
Title

H/Caswell 11/16/2025
Signature/Date