

Product Number:	103358
Lot Number:	501605
Control Lot Number:	501605A
Expiration:	16-Jan-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.97	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager



22 JAN 2025

Name

Title

Signature/Date