

Product Number:	101476
Lot Number:	501517
Expiration:	15-Jan-2026
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.95	8.95 - 9.05	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 22 JAN 2025

Name

Title

Signature/Date