

Product Number:	103360
Lot Number:	403709
Expiration:	6-Feb-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.98	7.95 - 8.05	PASS

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 14 FEB 2024

Signature/Date