

Product Number:	102914
Lot Number:	403710
Expiration:	6-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

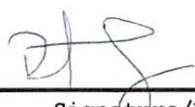
Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 14 FEB 2024

Signature/Date