

Product Number:	102912
Lot Number:	403708
Expiration:	6-Feb-2025
Storage:	2-8°C

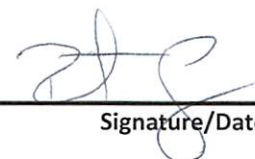
Laboratory Analysis

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