

Product Number:	102910
Lot Number:	403706
Expiration:	6-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

RS 14 FEB 2024
Signature/Date