

Product Number:	103359
Lot Number:	511401
Expiration:	24-Apr-2026
Storage:	2 - 8°C

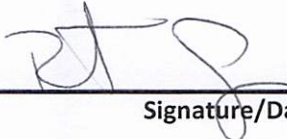
Laboratory Analysis

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

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 25 APR 2025
Signature/Date