

Product Number:	103359
Lot Number:	412404
Expiration:	3-May-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.03	7.95 - 8.05	PASS

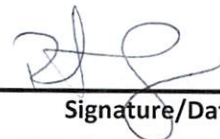
Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title



15 MAY 2024

Signature/Date