

Product Number:	103358
Lot Number:	412402
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

QC Manager

 15 MAY 2024

Name

Title

Signature/Date