

Product Number:	103361
Lot Number:	412408
Expiration:	3-May-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	8.04	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 16 MAY 2024

Name

Title

Signature/Date