

Product Number:	103358
Lot Number:	421925
Expiration:	6-Aug-2025
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	8.01	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 15 AUG 2024

Name

Title

Signature/Date