

Product Number:	101353
Lot Number:	402218
Expiration:	22-Jul-2025
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.40	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 14 FEB 2024

Name

Title

Signature/Date