

Product Number:	101631
Lot Number:	404415
Expiration:	13-Feb-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.30 - 7.50	PASS

Review/Approval

ROBERT SONOGYI

QC MANAGER

 21 FEB 2024

Name

Title

Signature/Date