

Product Number:	100458
Lot Number:	402924
Expiration:	29-Jul-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

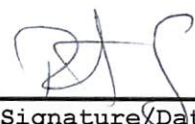
Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

 02 FEB 2024

Signature/Date