

Product Number:	104101
Lot Number:	403209
Expiration:	1-Aug-2025
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

16 FEB 2024

Signature/Date