

Product Number:	104072
Lot Number:	402219
Expiration:	22-Jul-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

 07-MAY-2024

Signature/Date