


Product Number:	101358
Lot Number:	405510
Expiration:	24-Aug-2025
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

ROBERT SOMOGYI QC MANAGER  16 APR 2024

Name

Title

Signature/Date