

Product Number:	101358
Lot Number:	429909
Expiration:	25-Apr-2026
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager

 17 Dec 2024

Name

Title

Signature/Date