

Product Number:	101358
Lot Number:	429903
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


Signature/Date

12/01/2024