

Product Number:	104071
Lot Number:	414905
Expiration:	28-Nov-2025
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager

 24 FEB 2025

Name

Title

Signature/Date