

Product Number:	103847	
Lot Number:	308220	
Expiration:	23-Mar-2024	
Storage:	2 - 8°C	

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan 01MAY2023

Name

Title

Signature/Date