

Product Number:	101358	
Lot Number:	311604	
Control Number:	311604A	
Expiration:	26-Apr-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