

Product Number:	102002
Lot Number:	311102
Control Number:	311102A
Expiration:	21-Apr-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 26 APR 2023*  
Signature/Date