

Product Number:	101362
Lot Number:	312407
Control Number:	312407A
Expiration:	4-May-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Sonia Khillan 12 MAY 2023  
Signature/Date