

Product Number:	100376
Lot Number:	316508
Control Number:	316508A
Expiration:	14-Jun-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

*Sonia Khillaw*  
Name

QC Manager  
Title

*Sonia Khillaw 22 JUN 2023*  
Signature/Date