

Product Number:	101352
Lot Number:	220310
Expiration:	22-Jul-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Killian  
Name

QC  
Title

Sonia Killian / 26 Jul 2022  
Signature/Date