

Product Number:	101472
Lot Number:	220206
Expiration:	21-Jul-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan* / 26 JUL 2022  
Signature/Date