

Product Number:	101631
Lot Number:	234814
Expiration:	14-Dec-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.30 - 7.50

Review/Approval

Matt Kocouris  
Name

QC Supervisor  
Title

MK 17Jan2023  
Signature/Date