

Product Number:	101631
Lot Number:	234816
Expiration:	14-Dec-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.30 - 7.50

Review/Approval

*Sonia Khillan*

Name

*QC Manager*

Title

*Sonia 12/16/23 FEB 2023*

Signature/Date