

Product Number:	100460
Lot Number:	234211
Expiration:	8-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 12/11/23 10FEB2023  
Signature/Date