

Product Number:	101955
Lot Number:	309414
Expiration:	4-Apr-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan      QC Manager      Sonia Khillan 12 MAY 2023  
Name    Title    Signature/Date