

Product Number:	101474
Lot Number:	312906
Control Number:	312906A
Expiration:	9-May-2024
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

*Sonia Khillan*

Name

*QC Manager*

Title

*Sonia K/L 23MAY2023*

Signature/Date