

Product Number:	104101
Lot Number:	310104
Control Number:	310104A
Expiration:	11-Apr-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan QC Manager Sonia Khillan 05 JUN 2023
Name Title Signature/Date