

Product Number:	101352
Lot Number:	318108
Control Number:	318108A
Expiration:	30-Jun-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 14 JUL 2023
Name Title Signature/Date