

Product Number:	100460
Lot Number:	306812
Expiration:	9-Mar-2024
Storage:	2-8 °C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan QC manager Sonia 22/6 11 APR 2023
Name Title Signature/Date