

Product Number:	101359
Lot Number:	300910
Expiration:	9-Jan-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 17 JAN 2023
Signature/Date