

Product Number:	101362
Lot Number:	300608
Expiration:	6-Jan-2024
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sania Khillan

QC Manager

Sania Khillan / 17 JAN 2023

Name

Title

Signature/Date