

Product Number:	100376
Lot Number:	302512
Expiration:	25-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 / 02 FEB 2023

Signature/Date