

Product Number:	100376
Lot Number:	217316
Expiration:	22-Jun-2023
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

Signature/Date