

Product Number:	102002
Lot Number:	234907
Expiration:	15-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan

QC manager

Jan 12/16 / 22 DEC 2022

Name

Title

Signature/Date