

Product Number:	100458
Lot Number:	234304
Expiration:	9-Dec-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan / 22 DEC 2022

Signature/Date