

Product Number:	101474
Lot Number:	234812
Expiration:	14-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 20 DEC 2022
Signature/Date