

Product Number:	103847
Lot Number:	233508
Expiration:	1-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khilkin

QC Manager

Soni 12/6/19DEC2022

Name

Title

Signature/Date