

Product Number:	103359
Lot Number:	235006
Expiration:	16-Dec-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.95 - 8.05

Review/Approval

Sonia Khilkin

Name

QC Manager

Title

Sonia Khilkin / 03 JAN 2023

Signature/Date