

Product Number:	103172
Lot Number:	213210
Expiration:	12-May-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khilkin
Name

QC

Title

Sonia Khilkin / 08 AUG 2022
Signature/Date