

Product Number:	103361
Lot Number:	232002
Expiration:	16-Nov-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Name

Title

Signature/Date

Sonia Khilkin

QC manager

Sonia/20/6 30 Nov 2022