

Product Number:	103361
Lot Number:	321905
Expiration:	7-Aug-2024
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 21 Aug 2023
Signature/Date