

Product Number:	102912
Lot Number:	320102
Expiration:	20-Jul-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 27 JUL 2023
Signature/Date