

Product Number:	102908
Lot Number:	320006
Expiration:	19-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