

Product Number:	102914
Lot Number:	215603
Expiration:	5-Jun-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Matt Smart
Name

QC

Title

MSA 14Jun2022
Signature/Date