

Product Number:	103360
Lot Number:	223103
Expiration:	19-Aug-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Matt Smart
Name

QC
Title

MA 22 Aug 22
Signature/Date