

Product Number:	101358
Lot Number:	214005
Expiration:	20-May-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Matt Smart

Name

QC

Title

MSA 24 May 2022

Signature/Date