

Product Number:	101956
Lot Number:	215212
Expiration:	1-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Matt Smart
Name

QC
Title

MS 22 Jun 2022
Signature/Date