

Product Number:	101954
Lot Number:	217004
Expiration:	19-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	PASS	7.75 - 7.85

Review/Approval

Matt Smart
Name

QC
Title

M. Smart 24 Jun 2022
Signature/Date