

Product Number:	101352
Lot Number:	210923
Expiration:	19-Apr-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

MSA 14 Jun 2022  
Signature/Date