

Product Number:	101359
Lot Number:	108821
Expiration:	29-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval Shivani Guel QC Manager Shivani Guel 14 Apr 2021
Name Title Signature/Date