

Product Number:	101472
Lot Number:	117711
Expiration:	26-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.20-7.60

Review/Approval

Shivani Guel

Name

QC

Title

Shivani Guel 23 Aug 2021

Signature/Date