

Product Number:	102252
Lot Number:	106717
Expiration:	8-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 15 Nov 2021

Signature/Date