

Product Number:	101472
Lot Number:	122009
Expiration:	8-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.20-7.60

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 18Oct2021

Signature/Date