

Product Number:	102199
Lot Number:	118305
Expiration:	2-Jul-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 06Oct2021

Signature/Date