

Product Number:	102620
Lot Number:	124105
Expiration:	29-Aug-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 19 Oct 2021

Signature/Date