

Product Number:	102103
Lot Number:	128012
Expiration:	7-Oct-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 21Jan2022

Signature/Date