

Product Number:	103661
Lot Number:	203713
Expiration:	6-Feb-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 08 Mar 2022

Name

Title

Signature/Date