

Product Number:	101472
Lot Number:	033713
Expiration:	2-Dec-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 15Jan 2021

Name

Title

Signature/Date