

Product Number:	101474
Lot Number:	210902
Expiration:	19-Apr-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 20 Apr 2022

Name

Title

Signature/Date