

Product Number:	101359
Lot Number:	210114
Expiration:	11-Apr-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH Measurement	Pass	7.35 - 7.45

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 13 Apr 2022

Signature/Date