

Product Number:	101359
Lot Number:	212705
Expiration:	7-May-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 17 May 2022

Signature/Date