

Product Number:	101359
Lot Number:	206616
Expiration:	7-Mar-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 23 Mar 2022

Signature/Date