

Product Number:	100460
Lot Number:	100824
Expiration:	8-Jan-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 11 Mar 2021

Signature/Date