

Product Number:	101473
Lot Number:	015436
Expiration:	2-Jun-2021
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 01 Dec 2020

Signature/Date