

Product Number:	101631
Lot Number:	105205
Expiration:	21-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.30-7.50

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel

Signature/Date

24 MAR 2021