

Product Number:	101782
Lot Number:	032601
Expiration:	21-Nov-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 05May2021
Signature/Date