

Product Number:	101472
Lot Number:	133916
Expiration:	5-Dec-2022
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 26Jan

Signature/Date *2022*