

Product Number:	103847
Lot Number:	104914
Expiration:	18-Feb-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 12Mar2021
Signature/Date