

Product Number:	103847
Lot Number:	211305
Expiration:	23-Apr-2023
Storage:	2 - 8°C


**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Richa Shrestha

QC Supervisor

 28Apr2022

Name

Title

Signature/Date