

Product Number:	103847
Lot Number:	110809
Expiration:	18-Apr-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 16Nov2021  
Signature/Date