

Product Number:	103847
Lot Number:	202310
Expiration:	23-Jan-2023
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 17Feb2022  
Signature/Date