

Product Number:	103847
Lot Number:	104921
Expiration:	18-Feb-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Guel  
Name

QC  
Title

Shivani Guel 13Dec2021  
Signature/Date