

Product Number:	103172
Lot Number:	205007
Expiration:	19-Feb-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 08 Mar 2022

Name

Title

Signature/Date