

Product Number:	101474
Lot Number:	204908
Expiration:	18-Feb-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 24Feb2022

Signature/Date