

Product Number:	101476
Lot Number:	201804
Expiration:	18-Jan-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 21Jan2022

Signature/Date