

Product Number:	101474
Lot Number:	202708
Expiration:	27-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 01Feb2023

Signature/Date