

Product Number:	101474
Lot Number:	201215
Expiration:	12-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

SG 18 Jan 2022
Shivani Goel 18 Jan 2022

Signature/Date