

Product Number:	101475
Lot Number:	119015
Expiration:	9-Jul-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

Name

QC

Title

Shwam Goel 23 Jul 2021

Signature/Date