

Product Number:	103704
Lot Number:	119705
Expiration:	16-May-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 23 Jul 2021

Signature/Date