

Product Number:	103704
Lot Number:	117113
Expiration:	20-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC Supervisor
Title

Shivani Goel 13 Jul 2021
Signature/Date