

Product Number:	101358
Lot Number:	117227
Expiration:	21-Jun-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Gael
Name

QC
Title

Shivani Gael 12 Oct 2021
Signature/Date