

Product Number:	100460
Lot Number:	117226
Expiration:	21-Jun-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 17 Aug 2021
Signature/Date