

Product Number:	101901
Lot Number:	107716
Expiration:	18-Mar-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
Signature/Date
21 Mar
2021