

Product Number:	103847
Lot Number:	106822
Expiration:	9-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 2021