

Product Number:	103359
Lot Number:	106706
Expiration:	8-Mar-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel
Name

QC Manager
Title

15 Mar 2021
Signature/Date