

Product Number:	103359
Lot Number:	104327
Expiration:	12-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Junpil Park
Name

Supervisor, Quality Control
Title

 5 Feb 2021
Signature/Date