

Product Number:	101476
Lot Number:	104329
Expiration:	12-Feb-2022
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 25 Feb 2021
Signature/Date