

Product Number:	101476
Lot Number:	109213
Expiration:	2-Apr-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

 20 Apr 2021
Signature/Date