

Product Number:	101631
Lot Number:	034204
Expiration:	7-Dec-2021
Storage:	2 - 8°C


Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 27Jan2021
Signature/Date