

Product Number:	101631
Lot Number:	034201
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

 27 Jan 2021
Signature/Date