

Product Number:	101473
Lot Number:	115231
Expiration:	1-Jun-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 01 Sep 2021
Signature/Date