

Product Number:	101955
Lot Number:	109215
Expiration:	2-Apr-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Junpil Park
Name

Quality Control
Title

Supervisor  28 Apr 2021
Signature/Date