

Product Number:	101955
Lot Number:	031910
Expiration:	14-Nov-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Junpil Park

Quality Control, Supervisor

 14 Jan 2021

Name

Title

Signature/Date