

Product Number:	101353
Lot Number:	031415
Expiration:	9-Nov-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Junpil Park

Name

Quality Control, Supervisor

Title

 05 Feb 2024

Signature/Date