

Product Number:	101353
Lot Number:	102710
Expiration:	27-Jan-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Junpil Park
Name

Quality Control, Supervisor
Title

 18 Feb 2021
Signature/Date