

Product Number:	101362
Lot Number:	022413
Expiration:	11-Aug-2021
Storage:	2-8°C


Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 05 Feb 2021
Signature/Date