

Product Number:	101954
Lot Number:	104806
Expiration:	17-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 01 Mar 2021
Signature/Date