

Product Number:	101357
Lot Number:	100703
Expiration:	7-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 14 Jan 2021
Signature/Date