

Product Number:	101362
Lot Number:	035010
Expiration:	15-Dec-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 06 Jan 2021
Signature/Date