

Product Number:	101355
Lot Number:	031131
Expiration:	6-Nov-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title



08 Mar 2021

Signature/Date