

Product Number:	100453
Lot Number:	100709
Expiration:	7-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

QC Supervisor
Title


Signature/Date

04 Mar 2021