

Product Number:	103158
Lot Number:	121718
Expiration:	5-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Jampil Park

Name

QC Supervisor

Title



01 Sep 2021

Signature/Date