

Product Number:	100376
Lot Number:	108323
Expiration:	24-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Jumpil Park

Name

QC

Supervisor

Title

[Signature] 28 Apr 2021

Signature/Date