

Product Number:	101353
Lot Number:	105014
Expiration:	19-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Jumpil Park

Name

QC Supervisor

Title

[Signature] 08 Mar 2021

Signature/Date