

Product Number:	103158
Lot Number:	102619
Expiration:	26-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park

Supervisor, QC



04 Mar 2021

Name

Title

Signature/Date