

Product Number:	100458
Lot Number:	104805
Expiration:	17-Feb-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 21 Mar 2021
Signature/Date