

Product Number:	103158
Lot Number:	110010
Expiration:	10-Apr-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Tumpil Park
Name

QC Supervisor
Title

[Signature] 22 Apr 2021
Signature/Date