

Product Number:	100376
Lot Number:	121719
Expiration:	5-Aug-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Jungil Park

Name

QC Supervisor

Title

[Signature] 01 Sep 2021

Signature/Date