

Product Number:	101353
Lot Number:	105614
Expiration:	25-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Jungil Park
Name

Quality Control Supervisor
Title

[Signature] 08 Mar 2022
Signature/Date