

Product Number:	100457
Lot Number:	024110
Expiration:	28-Aug-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 16 Feb 2021
Signature/Date