

Product Number:	100376
Lot Number:	100805
Expiration:	8-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 02 Feb 2021
Signature/Date