

Product Number:	101956
Lot Number:	110008
Expiration:	10-Apr-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
pH Measurement	PASS

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 20 Apr 2024
Signature/Date