

Product Number:	101358
Lot Number:	100818
Expiration:	8-Jan-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 05 Feb 2021
Signature/Date