

Product Number:	103847
Lot Number:	101604
Expiration:	16-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

 27Jan2021
Signature/Date