

Product Number:	102620
Lot Number:	104021
Expiration:	9-Feb-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Jumpil Park
Name

QC Supervisor
Title

[Signature] 05 Apr 2021
Signature/Date