

Product Number:	102620
Lot Number:	030012
Expiration:	26-Oct-2021
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Junpil Park
Name

Supervisor, Quality Control
Title

 09 Dec 2020
Signature/Date