

Product Number:	101474
Lot Number:	102912
Expiration:	29-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