

Product Number:	101631
Lot Number:	126909
Expiration:	26-Sep-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.30-7.50

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 18 Oct 20 21
Signature/Date