

Product Number:	101474
Lot Number:	124003
Expiration:	28-Aug-2022
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Tunpil Park
Name

QC Supervisor
Title

[Signature] 01 Sep 2021
Signature/Date