

Product Number:	102914
Lot Number:	102510
Expiration:	25-Jan-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Junpil Park  
Name

Supervisor, Quality Control  
Title

 02 Feb 2021  
Signature/Date