

Product Number:	102912
Lot Number:	207609
Expiration:	17-Mar-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Richa Shrestha  
Name

QC Supervisor  
Title

 21 APR 2022  
Signature/Date