

Product Number:	102910
Lot Number:	211307
Expiration:	23-Apr-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

*[Signature]* 26 Apr 2022

Signature/Date