

Product Number:	103360
Lot Number:	208302
Expiration:	24-Mar-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95-8.05

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel 28 Mar 2022*  
Signature/Date