

Product Number:	103359
Lot Number:	204201
Expiration:	11-Feb-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.95 - 8.05

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 15Feb2022  
Signature/Date