

Product Number:	103158
Lot Number:	205006
Expiration:	19-Feb-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	PASS

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 01Mar2022  
Signature/Date