

Product Number:	101353
Lot Number:	202210
Expiration:	22-Jan-2023
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel  
Name

Quality Control  
Title

Shwam Goel 15Feb2022  
Signature/Date