

Product Number:	103360
Lot Number:	106701
Expiration:	8-Mar-2022
Storage:	2 - 8°C

**Laboratory Analysis**

Characteristic	Result
pH	Pass

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 16Mar 2021  
Signature/Date