

Product Number:	101359	
Lot Number:	031411	
Expiration:	9-Nov-2021	
Storage:	2 - 8°C	

Laboratory Analysis

Characteristic	Result
pH Measurement	Pass

Review/Approval

Shivani Goel
Name

Quality Control Manager
Title

Shivani Goel 19 NOV
Signature/Date 2020