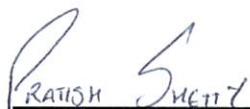


Product Number:	101473
Lot Number:	308218
Expiration:	23-Mar-2024
Storage:	2 - 8°C


**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

  
PRATIK S. JETHI  
Name

QC  
SUPERVISOR  
Title

  
Signature/Date  
20 APR 2023