


Product Number:	101475
Lot Number:	313204
Control Number:	313204A
Expiration:	12-May-2024
Storage:	2 - 8°C


**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

  
PRATISH S. MEHTA  
Name

  
QC SUPERVISOR  
Title

  
22 MAY 2023  
Signature/Date