

Product Number:	104605
Lot Number:	321317
Expiration:	1-May-2025
Storage:	2 - 8°C

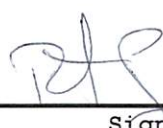
**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 15 FEB 2018  
Signature/Date