

<b>Product Number:</b>	101475
<b>Lot Number:</b>	107808
<b>Expiration:</b>	19-Mar-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Gael  
Name

QC Manager  
Title

Shivani Gael 30Mar 2021  
Signature/Date