

<b>Product Number:</b>	101472
<b>Lot Number:</b>	024417
<b>Expiration:</b>	31-Aug-2021
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 02NOV2020*

Name

Title

Signature/Date