

<b>Product Number:</b>	101472
<b>Lot Number:</b>	122011
<b>Expiration:</b>	8-Aug-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.20-7.60

Review/Approval

Shivam Goel  
Name

QC  
Title

Shivam Goel 16 Nov 2021  
Signature/Date