

<b>Product Number:</b>	103847
<b>Lot Number:</b>	134212
<b>Expiration:</b>	8-Dec-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35-7.45

Review/Approval

*Shivani Goel*

Name

QC

Title

*Shivani Goel* <sup>2022</sup> 26 Jan 2022  
Signature/Date *SG*