

<b>Product Number:</b>	101556
<b>Lot Number:</b>	203209
<b>Expiration:</b>	1-Feb-2023
<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 14 Apr 2022*

Name

Title

Signature/Date