

<b>Product Number:</b>	101476
<b>Lot Number:</b>	102915
<b>Expiration:</b>	29-Jan-2022
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	8.95-9.05

Review/Approval

Shivani Goel

Manager, QC

Shivani Goel 05 Feb 2021

Name

Title

Signature/Date