

<b>Product Number:</b>	104071
<b>Lot Number:</b>	304005
<b>Expiration:</b>	9-Feb-2024
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH	Pass	7.35 - 7.45

Review/Approval

*Pratish Suresh*  
Name

*EL, SUPERVISOR*  
Title

*[Signature]* 22 MAY 2023  
Signature/Date