

<b>Product Number:</b>	103172
<b>Lot Number:</b>	314411
<b>Control Number:</b>	314411A
<b>Expiration:</b>	24-May-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

Sonia Khillan  
Name

QC Manager  
Title

Soni Kalla 18 JUL 2023  
Signature/Date