

<b>Product Number:</b>	101474
<b>Lot Number:</b>	301712
<b>Expiration:</b>	17-Jan-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 Khilka  
Name

QC Manager  
Title

Sonia Khilka / 19 JAN 2023  
Signature/Date