

<b>Product Number:</b>	103516
<b>Lot Number:</b>	317307A
<b>Expiration:</b>	22-Dec-2024
<b>Storage:</b>	2-8°C

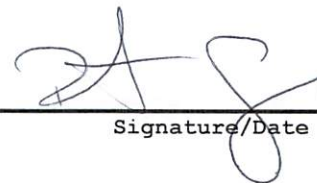
**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
pH Measurement	PASS

Review/Approval

Robert Somogyi

QC Manager



11 JUN 2024

Name

Title

Signature/Date