

<b>Product Number:</b>	101359
<b>Lot Number:</b>	433807
<b>Expiration:</b>	3-Jun-2026
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH Measurement	7.30 - 7.50	PASS

**RESULT**

<b>pH Reading</b>	7.40
-------------------	------

**Review/Approval**

Robert Somogyi

QC Manager

 11/25/2024

Name

Title

Signature/Date