

<b>Product Number:</b>	101355
<b>Lot Number:</b>	323413
<b>Expiration:</b>	22-Aug-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>	<b>PASS / FAIL</b>
pH	7.41	7.35 - 7.45	PASS

Review/Approval

*Sonia Killam*  
Name

*QC Manager*  
Title

*Sonia Killam 01SEP2023*  
Signature/Date