

<b>Product Number:</b>	102620
<b>Lot Number:</b>	331018
<b>Expiration:</b>	6-May-2025
<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

ROBERT SOMOSYI

Name

QC MANAGER

Title

 24 JAN 2024

Signature/Date