

<b>Product Number:</b>	104969
<b>Lot Number:</b>	422020
<b>Expiration:</b>	7-Feb-2026
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Robert Somogyi

QC Manager

*Robert Somogyi* 25 MAR 2025

Name

Title

Signature/Date