

<b>Product Number:</b>	101631
<b>Lot Number:</b>	416218
<b>Expiration:</b>	10-Jun-2025
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.30 - 7.50	PASS

Review/Approval

Robert Somogyi  
Name

QC Manager  
Title

 09 SEP 2024  
Signature/Date