

<b>Product Number:</b>	102910
<b>Lot Number:</b>	417701
<b>Expiration:</b>	25-Jun-2025
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH Measurement	7.95	7.95 - 8.05	PASS

Review/Approval

Robert Somogyi

QC Manager

 01 Jun 2024

Name

Title

Signature/Date