

<b>Product Number:</b>	101353
<b>Lot Number:</b>	502213
<b>Expiration:</b>	22-Jul-2026
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.38	7.35 - 7.45	PASS

Review/Approval

Robert Somogyi

QC Manager

 06 MAR 2025

Name

Title

Signature/Date