

<b>Product Number:</b>	101353
<b>Lot Number:</b>	412313
<b>Expiration:</b>	2-Nov-2025
<b>Storage:</b>	2 - 8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.43	7.35 - 7.45	PASS

Review/Approval

ROBERT SOMOGYI

QC MANAGER

*[Signature]* 15 MAY 2024

Name

Title

Signature/Date