

<b>Product Number:</b>	101472
<b>Lot Number:</b>	315310
<b>Control Number:</b>	315310A
<b>Expiration:</b>	2-Jun-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result	pH Target	PASS / FAIL
pH	7.39	7.35 - 7.45	PASS

Review/Approval Sonia Khillan QC Manager Sonia Khillan 01SEP2023  
 Name Title Signature/Date