

<b>Product Number:</b>	100376
<b>Lot Number:</b>	308609
<b>Expiration:</b>	27-Mar-2024
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>	<b>pH Target</b>
pH Measurement	PASS	7.35 - 7.45

Review/Approval

Sonia Khillan  
Name

QC Manager  
Title

Sonia Khillan 06 APR 2023  
Signature/Date