

<b>Product Number:</b>	100458
<b>Lot Number:</b>	116408
<b>Expiration:</b>	13-Jun-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
pH Measurement	PASS

Review/Approval

Shwani Goel  
Name

QC  
Title

Shwani Goel 17 Aug 2021  
Signature/Date