

<b>Product Number:</b>	100452
<b>Lot Number:</b>	120804
<b>Expiration:</b>	26-Jul-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 04 Aug 2021  
Signature/Date