

# CERTIFICATE OF ANALYSIS

<b>10X ELISA Wash Buffer (WB1)</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry at room temperature</b>
527	Bulk	17D41	12/2019
6251	25 mL	NA	NA
650	100 mL	NA	NA
651	500 mL	NA	NA
652	1 L	17D42	12/2019
676	10 L	17D43	12/2019

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH at 10X	6.8 – 7.0	6.84	Pass
pH at 1X	7.2 – 7.6	7.42	Pass
Appearance	Clear liquid	Conforms	Pass
Conductivity uMHOs	99,000 - 121,000 uMHO	116,233	Pass
Storage	Room temperature or refrigerated	Room temperature	Pass

Quality Control Tracy Murphy  
 Tracy Murphy

Date 12/18/2017  
 12/18/2017