



# CERTIFICATE OF ANALYSIS

<b>Antigen-Down HRP Conjugate Stabilizer, 5X</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry at 2-8°C</b>
586	Bulk	20N10	06/2022
6169	25 mL	NA	NA
6102	100 mL	20T20	06/2022
6103	500 mL	21F18	06/2022
6104	1 L	20N11	06/2022
6105	10 L	NA	NA

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
<b>pH at 5X</b>	7.2-7.6	7.41	Pass
<b>Appearance</b>	Clear to light yellow liquid	Conforms	Pass
<b>Conductivity <math>\mu</math>MHOs</b>	55,000-70,000	63,300	Pass
<b>A<sub>280</sub> of 1:100 dilution</b>	0.3-0.6	0.417	Pass
<b>Storage</b>	Refrigerated	Refrigerated	Pass
<b>Documentation</b>	Approved	Approved	Pass

This is an assurance that the specified lot conforms to all requirements of the product specifications. This information is believed to be accurate and is offered in good faith.

Quality Assurance Tracy Murphy  
Tracy Murphy

Date 2/4/2021  
02/04/2021