

# CERTIFICATE OF ANALYSIS

<b>5X HRP Conjugate Stock Stabilizer Diluent (5X CS6)</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry at 2-8°C</b>
52	Bulk	18B46	11/2020
6173	25 mL	NA	NA
667	100 mL	18B47	11/2020
668	500 mL	18B48	11/2020
669	1 L	18B49	11/2020
670	10 L	NA	NA

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH at 5X	7.2-7.6	7.41	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity uMHOs	60,000-80,000 uMHO	63,400	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Tracy Murphy

Date

11/28/2018  
 11/28/2018