

# CERTIFICATE OF ANALYSIS

<b>1X Neptune HRP Conjugate Stabilizer Diluent</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry at 2-8°C</b>
5169	Bulk	17T36	10/2019
6347	100 mL	17T37	10/2019
6348	1 L	17T38	10/2019
6349	10 L	NA	NA

<b>Specifications</b>	<b>Criteria</b>	<b>Lot Data</b>	<b>Results</b>
pH	7.2-7.6	7.35	Pass
Appearance	Clear to slightly cloudy, yellow tinted liquid	Conforms	Pass
Conductivity	Report Results	15,720	Pass
Spec A <sub>280</sub> (1:20)	Report Results	0.561	Pass
Storage	Refrigerated	Refrigerated	Pass
Documentation	Approved	Approved	Pass

Quality Control

*Tracy Murphy*  
 Quality Assurance

Date

*10/31/2017*  
 10/31/2017