



# CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
Catalog#	Size	Lot#	Expiry at 2-8°C
3324	Bulk	19L3	12/04/2020
6275	100 mL	19L4	12/04/2020
6328	500 mL	NA	NA
6329	1 L	NA	NA
6700	5 L	NA	NA
6330	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
<b>pH</b>	3.70 ± 0.20	3.67	Pass
<b>Absorbance at 285 nm</b>	0.600 ± 0.150	0.471	Pass
<b>Absorbance at 650 nm</b>	≤ 0.050	0.000	Pass
<b>Appearance</b>	Colorless to very light yellow solution	Conforms	Pass
<b>Storage</b>	Refrigerated	Refrigerated	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 7/1/19  
 7/01/19