

# CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
Catalog#	Size	Lot#	Expiry
3324	Bulk	17G45	01/18/2020
6275	100 mL	17G46	01/18/2020
6328	500 mL	NA	NA
6329	1 L	17G47	01/18/2020
6330	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.63	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.501	Pass
Absorbance at 650 nm	≤ 0.050	0.000	Pass
Appearance	Colorless to very light yellow solution	Conforms	Pass
Storage	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.  
 Product is for research use only. Not for use in diagnostic procedures.

Quality Assurance:   
 Tracy Murphy

Date: 08/30/2017  
 08/30/2017