

# CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
Catalog#	Size	Lot#	Expiry
3324	Bulk	15D30	03/21/2018
6275	100 mL	NA	NA
6328	500 mL	NA	NA
6329	1 L	NA	NA
6330	10 L	15D31	03/21/2018

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.68	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.484	Pass
Absorbance at 650 nm	≤ 0.050	0.001	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  
 Tracy Murphy

Date: 12/16/2015  
 12/16/2015