

# CERTIFICATE OF ANALYSIS

## TMB 1-Component HRP Microwell Substrate (SUBT)

Catalog#	Size	Lot#	Expiry
3325	Bulk	15S36	08/18/2019
6276	100 mL	15S37	08/18/2019
6337	1 L	NA	NA
6331	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	3.55 ± 0.20	3.55	Pass
Absorbance at 285 nm	0.450 ± 0.100	0.416	Pass
Absorbance at 650 nm	≤ 0.020	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.

Signature: \_\_\_\_\_

*Tracy Hanson*

Quality Assurance

Date: \_\_\_\_\_

*09/22/2015*

09/22/2015