

CERTIFICATE OF ANALYSIS

TMB 1-Component HRP Microwell Substrate (SUBT)			
Catalog#	Size	Lot#	Expiry
3325	Bulk	17R37	08/23/2020
6276	100 mL	17R38	08/23/2020
6337	1 L	17R39	08/23/2020
6331	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	3.55 ± 0.20	3.63	Pass
Absorbance at 285 nm	0.450 ± 0.100	0.439	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.
 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: 
 Kristi Strandberg

Date: 03/22/17
 03/22/2017