

CERTIFICATE OF ANALYSIS

TMB 1-Component HRP Microwell Substrate (SUBT)			
Catalog#	Size	Lot#	Expiry
3325	Bulk	18J9	11/17/2021
6276	100 mL	18J10	11/17/2021
6337	1 L	18J11	11/17/2021
6331	10 L	NA	NA


Specifications	Criteria	Lot Data	Results
pH	3.55 ± 0.20	3.64	Pass
Absorbance at 285 nm	0.450 ± 0.100	0.460	Pass
Absorbance at 650 nm	≤ 0.020	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:


 01/09/2018