

CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
Catalog#	Size	Lot#	Expiry
3324	Bulk	15L60	09/10/2017
6275	100 mL	NA	NA
6328	500 mL	NA	NA
6329	1 L	NA	NA
6330	10 L	15L61	09/10/2017

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.71	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.512	Pass
Absorbance at 650 nm	≤ 0.050	0.002	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: *Mark Degen*

Date: 7/28/15

Quality Assurance

07/28/2015