

## CERTIFICATE OF ANALYSIS

<b>TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)</b>			
<b>Catalog#</b>	<b>Size</b>	<b>Lot#</b>	<b>Expiry</b>
3324	Bulk	18B51	12/04/2020
6700	5 L	18B52	12/04/2020

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.67	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.471	Pass
Absorbance at 650 nm	≤ 0.050	0.000	Pass
Appearance	Colorless to very light yellow	Conforms	Pass
Storage	Refrigerated	Refrigerated	Pass

We hereby certify that this material fully complies with ICT's technical requirements.  
 This information is believed to be accurate and is offered in good faith.

Quality Control                     *Tracy Murphy*                      
 Tracy Murphy

Date                     11/30/2018                      
 11/30/2018