

## CERTIFICATE OF ANALYSIS

<b>TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)</b>			
Catalog#	Size	Lot#	Expiry
3324	Bulk	18B7	01/30/2021
6700	5 L	18B8	01/30/2021

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.69	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.479	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   
 Kristi Strandberg

Date 11/2/18  
 11/2/2018