

CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
---	--	--	--

Catalog#	Size	Lot#	Expiry
3324	Bulk	17N37	09/21/2018
6700	5 L	17N38	09/21/2018

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.70	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.489	Pass
Absorbance at 650 nm	≤ 0.050	0.001	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
 Quality Assurance

Date 06/26/2017
 06/26/2017