



CERTIFICATE OF ANALYSIS

TMB Super Sensitive 1-Component HRP Microwell Substrate (SUBS)			
Catalog#	Size	Lot#	Expiry at 2-8°C
3324	Bulk	20L31	01/06/2023
6275	100 mL	NA	NA
6328	500 mL	NA	NA
6329	1 L	NA	NA
6700	5 L	20L32	01/06/2023
6330	10 L	NA	NA

Specifications	Criteria	Lot Data	Results
pH	3.70 ± 0.20	3.67	Pass
Absorbance at 285 nm	0.600 ± 0.150	0.505	Pass
Absorbance at 650 nm	≤ 0.050	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.

Quality Assurance 
 Sally Hed Dahlquist

Date 7/29/2020
 7/29/2020