



# 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 at 2-8°C</b> |
| 3324                                                                  | Bulk        | 19S20       | 04/30/2021             |
| 6275                                                                  | 100 mL      | NA          | NA                     |
| 6328                                                                  | 500 mL      | NA          | NA                     |
| 6329                                                                  | 1 L         | NA          | NA                     |
| 6700                                                                  | 5 L         | 19S21       | 04/30/2021             |
| 6330                                                                  | 10 L        | NA          | NA                     |

| <b>Specifications</b>       | <b>Criteria</b>                         | <b>Lot Data</b> | <b>Results</b> |
|-----------------------------|-----------------------------------------|-----------------|----------------|
| <b>pH</b>                   | 3.70 ± 0.20                             | 3.67            | Pass           |
| <b>Absorbance at 285 nm</b> | 0.600 ± 0.150                           | 0.484           | Pass           |
| <b>Absorbance at 650 nm</b> | ≤ 0.050                                 | -0.001          | Pass           |
| <b>Appearance</b>           | Colorless to very light yellow solution | Conforms        | Pass           |
| <b>Storage</b>              | 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 Tracy Murphy  
 Tracy Murphy  
 Date 9/23/19  
 9/23/2019

9401 James Avenue S., #155  
 Bloomington, MN 55431 USA  
 TOLL-FREE 1-800-829-3194  
 LOCAL 952-888-8788