



# Single Component TMB Peroxidase EIA Substrate Kit

## Catalog Numbers

**172-1068**

**172-1072**

**172-1073**



# Specifications

---

<b>Description</b>	This one-component substrate develops a deep blue color when reacted with peroxidase labeled conjugates in microplates.
<b>Contents</b>	3,3',5,5'-tetramethylbenzidine in an acidic buffer. Concentration of $H_2O_2$ is 0.01%.
<b>Storage conditions</b>	This substrate solution should be stored at 4 °C.
<b>Shelf life</b>	1 year at 4 °C.
<b>Use</b>	Warm substrate to room temperature before using. Recommended volume is 100 $\mu$ l. Color will develop immediately.
<b>Kinetic assays</b>	The TMB substrate produces a blue color. Read at a wavelength of 655 nm.

## Endpoint assays

After adding a stop solution, the TMB substrate produces a yellow color. Read at a wavelength of 450 nm.

## Stop solutions

Recommended stop solution is an equal volume of 0.18 M sulfuric acid [ $H_2SO_4$ ]. Other commonly used acids (0.3 M phosphoric [ $H_3PO_4$ ], or 0.12 M hydrochloric [HCl]), may also be used but are less efficient in stopping color development.

## Managing reaction rates

To reduce the intensity of the substrate reaction, it is recommended that the conjugate and/or antibodies in the immunoassay be further diluted. **Dilution of substrate is not recommended.**