


# TMB Stabilized Chromogen

Catalog Number SB01

Pub. No. MAN0018289 Rev. A.0 (30)

## Product description

Volume	25 mL
Product Description	SB01, TMB Stabilized Chromogen, is a one-component horseradish peroxidase substrate containing 3,3',5,5"-tetramethylbenzidine plus H <sub>2</sub> O <sub>2</sub> in a mildly acidic buffer. This solution is ready for use. Dilution of this product prior to use is not recommended.
Applications	This solution is intended for use in enzyme assays which employ horseradish peroxidase. It may be used in assays in which optical density is monitored over several time points (kinetic assays) and in endpoint assays. It is suitable for use in microwell assays, such as ELISA.
Directions for Use	<p>The solution should be colorless or pale blue prior to use. In preparation for use of this product, the volume required should be decanted from the bottle into a clean dispensing trough. This step is necessary to prevent contamination of the contents of the bottle. The decanted volume should be allowed to warm to room temperature (22°C) prior to use. In microwell assays, pipette 100 µL of the warmed solution into each well. In the presence of peroxidase, the solution develops a blue reaction product. When color formation is monitored over several time points (kinetic assays), the optical density of the solution is measured at two wavelengths: 492 nm and 650 nm. When color formation is monitored at an endpoint, an equal volume of an acidic stop solution, such as 1 N HCl or Stop Solution (Cat. No. SS01) is added to the wells and allowed to mix with the solution in the wells. Addition of the stop solution results in the rapid formation of a yellow end product. The optical density of this yellow end product is measured at 450 nm.</p> <p> <b>WARNING!</b> Hazardous materials: TMB Stabilized Chromogen is highly flammable and extremely toxic, if swallowed. In case of accident, call a physician immediately.</p>
Storage	Store at 2–8° C. Protect from light and from exposure to metal ions. It is highly recommended that the volume required for each assay be carefully decanted into a clean dispensing trough, and warmed to room temperature (22°C) prior to use. The unused solution remaining in the trough should be discarded.
Expiration Date	See label.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria  
 For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.