

ReadiUse™ TMB Substrate Solution *Optimized for ELISA Assays with HRP Conjugates*

 Catalog number: 11003, 11012
 Unit size: 1 L, 100 ml

Component	Storage	Amount (Cat No. 11003)	Amount (Cat No. 11012)
ReadiUse™ TMB Substrate Solution *Optimized for ELISA Assays with HRP Conjugates*	Refrigerated (2-8 °C), Minimize light exposure	1 bottle (1 L)	1 bottle (100 mL)

OVERVIEW

Horseshoe peroxidase (HRP) and HRP conjugates facilitate the ABTS oxidation in the presence of hydrogen peroxide, turning ABTS into its blue-green oxidized product. ReadiUse™ TMB Substrate Solution is a premixed solution of TMB substrate with hydrogen peroxide. It produces a blue product upon interaction with HRP or HRP conjugates without the addition of hydrogen peroxide. The soluble blue product can be quantitated at 650 nm. Use of a stop solution enhances sensitivity 2-4 fold and the resulting yellow solution can be read at 450 nm. ReadiUse™ TMB Substrate Solution provides a convenient and ultrasensitive quantitative substrate system.

AT A GLANCE

Important

Warm ReadiUse™ TMB Solution to room temperature before use.

Note The reagent is to be used as supplied, no dilution is required.

KEY PARAMETERS

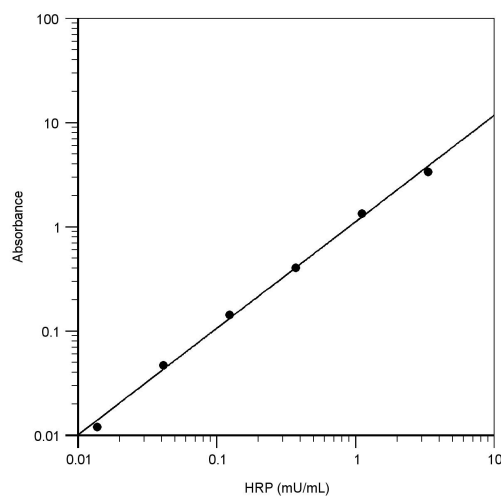
Absorbance microplate reader

Absorbance 650 nm
 Recommended plate Clear bottom

SAMPLE EXPERIMENTAL PROTOCOL

1. Wash the assay plate following the incubation of HRP-labeled reagent.
2. Add 100 µL of ReadiUse™ TMB Solution into each well.
3. Incubate the plate at room temperature for 15 – 30 min or until the desired color develops. *Note:* The incubation time varies depending on the assay conditions.
4. Measure the absorbance signal at 650 nm with an ELISA microplate reader. *Note:* If desired, the reaction can be stopped by adding an equal volume of 2M sulfuric acid to each well. Stopped reaction should be read at 450 nm.

EXAMPLE DATA ANALYSIS AND FIGURES



HRP dose response was measured with ReadiUse™ TMB Substrate Solution in a clear 96-well plate using a SpectraMax microplate reader (Molecular Devices).

Image generated with Quest™ Graph. ©2019 AAT Bioquest

Figure 1. HRP dose response was measured with ReadiUse™ TMB Substrate Solution in a clear 96-well plate using a SpectraMax microplate reader (Molecular Devices).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.