

## EUROPE

LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00  
• Fax: +32 4 372 75 00 • Toll-free: + 800 666 00 123 • info@eurogentec.com  
• www.eurogentec.com

## NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560  
• Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com  
• www.anaspec.com

## Technical Data Sheet

### iD TMB Substrate ID-SUTMB1-060

Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics.

#### Description

Tetramethyl Benzidine (TMB), a noncarcinogenic analog of benzidine, is routinely used as the key substrate for enzyme immunoassays involving horseradish peroxidase-labelled (HRP-labelled) antibodies and molecular probes. The iD TMB Substrate is a ready-to-use, ultra-sensitive, single-solution substrate for the detection of horseradish peroxidase (HRP) in Western blot and Dot blot assays.

#### Key Features

- Ready-to-use
- High sensitivity and low background
- Reproducible results

#### Storage

Store the kit at 2 - 8°C. It will remain stable for 12 months

#### Content

Components	ID-SUTMB1-060
TMB Solution	60 mL
Enough Reagents for	600 cm <sup>2</sup> of membrane*

\* based on 0.1 ml of the working solution per cm<sup>2</sup>

#### Protocol

This substrate is designed for the development of membranes from Western and Dot blots after the final wash step.

1. Drain off the excess of wash solution from the membrane by holding it vertically with forceps and touching its edge against a tissue.
2. Place the membrane on a clean, flat surface, and cover it with the TMB solution (0.1 ml/cm<sup>2</sup> recommended).
3. Incubate until the desired color is reached.
4. Stop the reaction by rinsing the membrane 3 times for 30 seconds each in 20 ml of deionized water.
5. Drain off the excess water
6. Transfer the membrane to a piece of paper towel.
7. Air-dry the membrane in a dark place.

**For further information please contact our Customer Help Desk:**

#### For Europe:

E-mail: info@eurogentec.com  
Tel: +32 4 372 76 65

#### For USA:

E-mail: service@anaspec.com  
Tel: +1 510 791 9560