

# Echelon Biosciences Inc.

---

## TMB Precipitating

K-TMBP

Support: echelon@echelon-inc.com

### **Description:**

A precipitating chromogen substrate that reacts with horseradish peroxidase (HRP) producing a permanent blue colored spot or band. This detection reagent is best used with an HRP conjugated secondary or tertiary antibody on a dot blot or western blot.

### **Applications:**

For the detection of a protein-lipid interaction on PIP Strips™ and related membrane-based products.

### **Properties:**

Size – 22 mL

Storage – Store at 4 °C. Warm to room temperature prior to use.

Specificity – Horseradish peroxidase, HRP

### **Protocol:**

To be used in conjunction with a protein-lipid overlay assay protocol found on the webpage of each individual product. See related products table for a list of products.

1. Incubate membrane with lipid binding protein, primary antibody and/or HRP conjugated secondary antibody. Wash membrane 3x 5 min with ample amount of wash buffer before adding K-TMBP.
2. After discarding the final wash buffer, add 2 mL TMB Precipitating (K-TMBP) per membrane and incubate with gentle agitation. Protein-lipid interaction will develop within 2-20 minutes with the spots turning blue. Watch for color development closely and avoid overdevelopment. When overdeveloped, blue color may run off the spot causing increased background.
3. Stop the reaction by adding an ample amount of dH<sub>2</sub>O (≥ 15mL).
4. Record an image of the membrane.

**Note:** Do not let the membrane dry at any time during the protein-lipid overlay assay. For Western blots add 1 mL per 10 cm<sup>2</sup> of membrane.

### **Related Products:**

Product	Catalog Number
PIP Strips™	P-6001
Membrane Lipid Strips™	P-6002
SphingoStrips™	S-6000
Custom Lipid Array Strip	P-6111
PIP Arrays™	P-6100
Membrane Lipid Arrays™	P-6003
Sphingo Arrays™	S-6001
PIP Pack	P-P100

---

Technical Data Sheet Rev. 3, 07-17-19 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.com