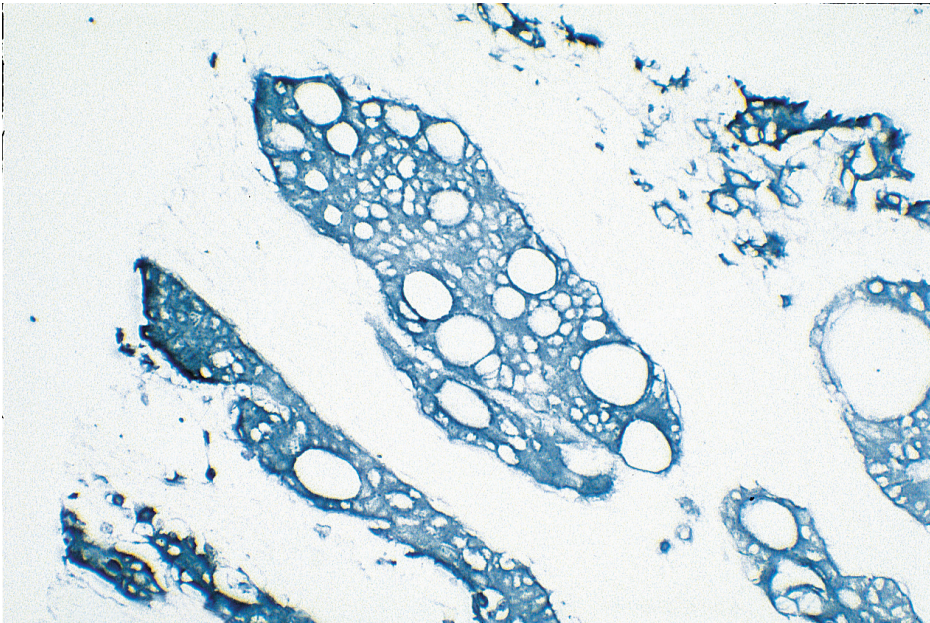




TMB Substrate Kit, Peroxidase (HRP), (3,3', 5,5'-tetramethylbenzidine)

SK-4400

[Product Images](#)



Short Description

TMB (3,3', 5,5' tetramethylbenzidine dihydrochloride) is a sensitive chromogen substrate for peroxidase and a non-carcinogenic substitute for benzidine. TMB can be used for blotting, immunohistochemical, and *in situ* hybridization applications. Slides stained with TMB can be dehydrated and permanently mounted and the soluble blue product, produced in the absence of the stabilizing solution, is useful for ELISA assays.

The TMB HRP Substrate Kit (SK-4400) contains the reagents needed to produce either a precipitating blue reaction product or a soluble blue reaction product. This kit also contains a stabilizing solution which, when added to the working solution, allows the development of a blue precipitate. Kit materials are pre-packaged in convenient dropper bottles.

Features:

- Greater sensitivity than conventional substrates
- Stock solutions supplied in convenient dropper bottles promoting ease of handling
- No wait times for mixing and dissolving powders or tablets
- Sufficient reagents to produce 300 ml of working solution

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, In situ hybridization, Blotting Applications, Elispot, ELISAs
Detection Enzyme(s)	Peroxidase
Substrate Reaction Product	Precipitating Reaction Product
Chromogen Color	Blue
Microscopy	Brightfield, Spectral Imaging
Mounting	Non-Aqueous (Permanent)
Safety Title	Cancer

