



AnaPrep™ IHC Histain Kit-DAB

CATALOG No.: 29700

INTRODUCTION:

IHC Histain kit-DAB is developed for research use only. In general, this kit is used to check the presence of interest targets (including proteins, special modification) in the tissue or cells. With exposure of the antigens, the primary antibody is applied to recognize its target. The result is visualized by the interaction of the substrate with enzyme-linked streptavidin which has high affinity to the biotinylated secondary antibody derived from either mouse or rabbits.

DESCRIPTION:

The reagents in this kit constitute a labeled streptavidin –biotin immunoenzymatic antigen detection system. This technique involves the sequential incubation of the specimen with an unconjugated primary antibody specific to the target antigen, a biotinylated secondary antibody which reacts with the primary antibody, enzyme-labeled streptavidin, and substrate-chromogen.

KIT CONTENTS:

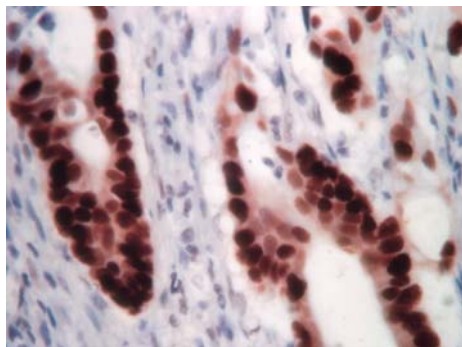
Reagent	Form	Description	Storage
Hydrogen Peroxide Block	Ready to use (15 ml)	To block Endogenous peroxidase	4-8 °C
Protein Block	Ready to use (15 ml)	Serum blocking solution	4-8 °C
Anti-polyvalent	Ready to use (15 ml)	Biotinylated Secondary antibody	4-8 °C Do not freeze
Streptavidin Peroxidase	Ready to use (15 ml)	Streptavidin-enzyme Conjugate with stabilizer	4-8 °C Do not freeze
DAB Chromogen (20x)	1 ml	Carcinogenic, Handle with caution	4-8 °C



DAB Substrate	15 ml	Contains hydrogen peroxide	4-8 °C
---------------	-------	----------------------------	--------

STAINING PROTOCOL

- Deparaffinize and rehydrate Formalin-fixed paraffin-embedded tissue section.
- Add enough drops of **Hydrogen Peroxide Block** to cover the sections. Incubate for 10 minutes, wash 2 times in buffer.
- Perform appropriate pretreatment if required, wash 3 times in buffer.
- Apply **Protein Block** and incubate for 5 minutes at RT to block nonspecific binding. Wash 1 times in buffer.
- Apply primary antibody and incubated according to manufacturer's recommendation.
- Wash 4 times in buffer. Apply **Anti-Polyvalent** and incubate for 10 min at RT. Wash 4 times in buffer.
- Apply **Streptavidin Peroxidase** and incubate for 10 min at RT.
- Rinse 4 times in buffer. Add 50 µl **DAB Chromogen** to 1 ml of **DAB Substrate**. Mix by swirling and apply to tissue. Incubate for 10 min. Rinse 4 times in buffer.
- Add enough drops of Hematoxylin to cover the section, incubate for 1 minute.
- Rinse 7-8 times in tap water. Add Mounting medium to cover this section.
- Read the results under microscope.



Rabbit anti-P53 (CT) (Cat # 29594) stained on colon carcinoma by using the IHC Histain Kit-DAB (Cat# 29700)

STORAGE:

Stored at 4-8 °C

For Research Use Only

