

VitroView™ LSAB IHC/DAB Kit (Goat IgG) (150 Tests)

Cat. No. VB-6021D

Introduction

Immunohistochemistry (IHC) is a method of detecting the presence of specific proteins in cells of a tissue section by exploiting the principle of antibodies binding specifically to antigens in biological tissues. IHC is widely used in the diagnosis of abnormal cells and basic research to understand the distribution and localization of biomarkers and differentially expressed proteins in different parts of a biological tissue.

LSAB IHC Kit is based on labeled streptavidin-biotin (LSAB) method. This method utilizes a biotinylated secondary antibody that links primary antibodies to streptavidin-peroxidase conjugate. In this method a single primary antibody is subsequently associated with multiple biotin molecules. Therefore, an increase in sensitivity is achieved compared to direct peroxidase-conjugate methods.

Kit Components

1. RTU normal rabbit serum: 15ml.
 2. RTU biotinylated anti-goat secondary antibody: 15ml
 3. RTU streptavidin-HRP: 15ml
 4. DAB stock solution (40x): 0.75 ml
 5. Stable H₂O₂ solution (40x): 0.75ml
 6. DAB enhancer buffer: 30ml
 7. RTU hematoxylin solution: 15ml
- Note: RTU=ready-to-use

Reagents and Materials Required but Not Provided

- Xylene and ethanol
- Distilled or deionized water
- 30% hydrogen peroxide
- 10 mM phosphate-buffered saline (PBS), pH 7.4
- Triton X-100
- Mini PAP Pen
- Primary antibody
- Mounting Media

Storage

Store at 2-8°C.

Application

Immunohistochemistry for detecting a primary antibody made in goat.

Protocol

1. Preparation of Slides

1) Cell Lines

- Grow cultured cells on sterile glass cover slips or slides overnight at 37 °C
- Wash briefly with PBS
- Fix as desired. Possible procedures include:
 - a. 20 minutes with 10% formalin in PBS (keep wet)

- b. 10 minutes with ice cold methanol, allow to air dry
- c. 10 minutes with ice cold acetone, allow to air dry
- Wash in PBS

2) Frozen Sections

- Snap frozen fresh tissues in liquid nitrogen or isopentane pre-cooled in liquid nitrogen, embedded in OCT compound in cryomolds. Store the frozen tissue block at -80°C until ready for sectioning.
- Transfer the frozen tissue block to a cryotome cryostat (e.g. -20°C) prior to sectioning and allow the temperature of the frozen tissue block to equilibrate to the temperature of the cryotome cryostat.
- Section the frozen tissue block into a desired thickness (typically 5-10 µm) using the cryotome.
- Place the tissue sections onto glass slides suitable for immunohistochemistry (e.g. Superfrost).
- Sections can be stored in a sealed slide box at -80°C for later use.
- Before staining, warm slides at room temperature for 30 minutes and fix in ice cold acetone or ice cold methanol for 10 minutes. Air dry for 30 minutes.
- Wash in PBS

3) Paraffin Sections

- Deparaffinize sections in xylene, 3x5min.
- Hydrate with 100% ethanol, 2x2min.
- Hydrate with 95% ethanol, 2x2min.
- Rinse in distilled water.
- Follow procedure for pretreatment as required.

2. Antigen retrieval

Most formalin-fixed tissues require an antigen retrieval step before immunohistochemical staining can proceed. Heat-mediated and enzymatic antigen retrievals are common methods.

- **For Citrate:** Bring slides to a boil in 10 mM sodium citrate buffer, pH 6.0; maintain at a sub-boiling temperature for 10 minutes. Cool slides on bench top for 30 minutes.
- **For EDTA:** Bring slides to a boil in 1 mM EDTA, pH 8.0: follow with 15 minutes at a sub-boiling temperature. No cooling is necessary.
- **For TE:** Bring slides to a boil in 10 mM TE/1 mM EDTA, pH 9.0: then maintain at a sub-boiling temperature for 18 minutes. Cool at room temperature for 30 minutes.
- **For Pepsin:** Digest for 10 minutes at 37°C.

Note: Do not use this pretreatment with frozen sections or cultured cells that are not paraffin-embedded.

3. Staining Procedure

- 1) Rinse sections in PBS-Triton X-100 (0.025%) for 2x2min
- 2) **Serum Blocking:** incubate sections with 3-4 drops of RTU normal rabbit serum for 30 minutes to block non-specific binding of immunoglobulin.
- 3) **Primary Antibody:** incubate sections with primary antibody at appropriate dilution in antibody dilution buffer (Cat#: VB-6002) for 1-2 hour at room temperature or overnight at 4 °C. Rinse in PBS.
- 4) **Peroxidase Blocking (optional):** incubate sections in 0.3% hydrogen peroxide in PBS for 10 minutes at room temperature. Rinse in PBS.
- 5) **Secondary Antibody:** incubate sections with 3-4 drops of RTU biotinylated anti-goat secondary antibody for 30 minutes at room temperature.
- 6) Rinse in PBS for 3x2min.
- 7) **Detection:** incubate sections with 3-4 drops of RTU streptavidin-HRP for 30 minutes at room temperature.

- 8) Rinse in PBS for 3x2min.
- 9) **Chromogen/Substrate:** incubate sections with 3 drops of DAB solution for 2-8 minutes. Monitor signal development under a microscope
Note: DAB solution is made by mixture of 20 µl of DAB stock solution and 20 µl of stable H₂O₂ solution with 1ml of DAB enhancer buffer (dark-brown stain).
- 10) Rinse in distilled water 2x2 min
- 11) **Counterstain:** For using Hematoxylin Nuclear Counterstaining Kit (CAT#: VB-6004), incubate sections with 3 drops of RTU hematoxylin solution for 1-2 minutes. Rinse in tape water 2x2 min.
- 12) Dehydrate through 75% ethanol for 2 min, 95% ethanol for 2 min, and 100% ethanol for 2x3min. Clear in xylene for 2x5min.
- 13) Coverslip with mounting medium.

IHC Troubleshooting

1. High background staining

Possible Cause	Solution
Endogenous peroxidase activity was incompletely blocked.	Incubate sections in 0.3% hydrogen peroxide in methanol or PBS for 10-30 minutes at room temperature.
Deparaffinization was incomplete.	Prepare new sections and deparaffinize according to standard laboratory protocol using fresh xylene or xylene substitute.
Inadequate rinsing of slides.	Gently rinse slide with wash buffer bottle and place in wash bath for 5 minutes. Gentle agitation of the wash bath may increase effectiveness.
Over-development of substrate.	Reduce incubation time.
Dehydration of specimen during staining.	Keep section wet.

2. Negative staining on positive slides

Possible Cause	Solution
Steps in the staining protocol were performed in incorrect sequence.	Repeat the procedure.
Primary or secondary antibody incubation steps were omitted.	Repeat the procedure.
Labile antigens were destroyed.	Use fresh cutting slides. Use a paraffin wax with a melting temperature ~55-58°C. Wax used for embedding should not be exceed 60 °C.
Specimen was improperly fixed and/or processed.	Check manufacture's specifications regarding recommended fixative
Specimen dehydrated during staining.	Repeat the procedure by following the manufacture's protocol.

3. Weak staining on all slides

Possible Cause	Solution
Specimen retained excess liquid after rinsing steps.	Remove excess liquid after rinsing steps.
Incubation times were insufficient.	Prolong incubation time.
Substrate prepared improperly.	Check compatibility of buffer ingredients with enzyme and substrate-chromogen reagents. Repeat staining.
9Deparaffinization was incomplete	Prepare new sections and deparaffinize according to

(staining may be accompanied by high background).	standard laboratory protocol using fresh xylene or xylene substitute.
---	---

Warning

DAB is a possible carcinogen. Please take necessary precautions.

Precautions

Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.