



3050 Spruce Street
Saint Louis, Missouri 63103 USA
Telephone 800-325-5832 • (314) 771-5765
Fax (314) 286-7828
email: techserv@sial.com
sigma-aldrich.com

Product Information

VIMENTIN

For Immunohistochemical Demonstration of Vimentin in Paraffin-embedded and Frozen Human Tissue Sections

Product Number **IMMH-10**

Store at 2-8 °C

TECHNICAL BULLETIN

Product Description

The introduction of immunohistochemical techniques has ushered a new era of staining into the laboratory based upon sensitive, specific methods.^{1,2} Using antigen-antibody relationships, tissue components previously undetected can be precisely identified.

In the Sigma procedure, an antigen-specific primary antibody is applied to deparaffinized or frozen hydrated tissue sections. Following a brief wash, the section is incubated with a biotinylated secondary antibody. Upon the addition of an ExtrAvidin[®] Peroxidase reagent, a stable avidin-biotin complex is formed with the bound biotinylated antibody. The sites of antibody deposition are visualized by the addition of freshly prepared substrate which contains Hydrogen Peroxide and the chromogen 3-Amino-9-Ethylcarbazole (AEC; an electron donor). The bound peroxidase catalyzes the oxidation of the AEC to form a reddish-brown insoluble precipitate at the antigen sites. Compared to the classic PAP procedures, avidin-biotin techniques are particularly valuable as background staining is virtually eliminated while the specific reaction is amplified.^{3,4}

Vimentin is a 57 kDa molecular weight intermediate filament protein which is expressed in various cells and tissues of mesenchymal origin. Expression of vimentin can frequently be demonstrated in tumors derived from mesenchymal cells.^{5,6}

Materials Provided

Primary Antibody, Vial 1: Mouse Monoclonal Anti-vimentin in buffered saline. Sodium Azide, 0.1%, added as preservative.

Biotinylated Secondary Antibody, Vial 2: Goat Anti-mouse Immunoglobulin in buffered saline. Sodium Azide, 0.1%, added as preservative.

Peroxidase Reagent, Vial 3: ExtrAvidin-conjugated Peroxidase in buffered saline. Preservative added.

Acetate Buffer, Vial 4a: Acetate Buffer, 2.5 mol/L, pH 5.0.

AEC Chromogen, Vial 4b: 3-Amino-9-Ethylcarbazole (AEC) in N,N-Dimethylformamide.

Hydrogen Peroxide, Vial 5: 3% H₂O₂ in deionized water.

Mixing Vial, Vial 6

Materials Required But Not Provided

Blocking Reagent: Normal Goat Serum (Product No. G 9023), 1% (v/v) in buffered saline

Negative Control: Normal Mouse Serum (Product No. M 5905), 5% (v/v) in buffered saline

Phosphate Buffered Saline, pH 7.4 (available as tablets, Product No. P 4417)

Deionized water

Mayer's Hematoxylin Solution (Product Code MHS-1)

Glycerol Gelatin (Product Code GG-1)

Slides, coverslips

Humidity chamber (Product No. H 6644)

Light microscope

Precautions and Disclaimer

Primary Antibody and Biotinylated Secondary Antibody contain Sodium Azide. Due to the Sodium Azide content a material safety data sheet (MSDS) for this product has been sent to the attention of the safety officer of your institution.

AEC chromogen contains 3-Amino-9-Ethylcarbazole, a possible carcinogen. Consult the MSDS for information regarding hazards and safe handling practices.

Preparation Instructions

Tissues fixed in 10% Neutral Buffered Formalin, B-5 Fixative or Bouin's Solution¹² can be used. Cut tissue sections at 4-6 microns. Acetone fixed frozen sections may also be used. For detection of antigens requiring unmasking, digest formalin fixed material with 0.1% Trypsin (Product No. T 7409) or 0.1% Protease (Product No. P 5147) prior to the first step in the procedure.

Note: Since tissue sections have a tendency to fall off during immunohistochemical procedures, Poly-L-Lysine (Product No. P 8920) may be used as a tissue adhesive.

CONTROLS

For the correct interpretation of the staining results it is necessary to run a positive control tissue section known to contain the antigen in question and a negative control test section incubated with the negative control reagent.

Note: All rinses are with Phosphate Buffered Saline (PBS), pH 7.4. Following incubations slides should be washed gently with PBS from a wash bottle, avoid a direct jet of water which may wash off or loosen sections. Carefully wipe each slide free of excess fluid before the application of the next reagent. Avoid touching the tissue section. Be certain to apply enough drops of the reagent to cover the sections. DO NOT allow the tissue sections to dry out at any time during the procedure. It is recommended the incubations be performed in a humidity chamber. All incubations are at room temperature unless otherwise specified. Wash steps can include placing slides in a PBS bath for 2 minutes.

Procedure

1. Deparaffinize and hydrate sections to water.
2. Quench endogenous peroxide with 2 drops 3% Hydrogen Peroxide (Vial 5) for 5 minutes. Wash and wipe slides. Incubate with Blocking Reagent for 10 minutes. Wipe off excess reagent but do not wash slides.
3. Apply 2 drops Primary Antibody (Vial 1) or Negative Control and incubate 60 minutes. Wash and wipe slides.
4. Apply 2 drops Biotinylated Secondary Antibody (Vial 2) and incubate 20 minutes. Wash and wipe slides.
5. Apply 2 drops Peroxidase Reagent (Vial 3) and incubate 20 minutes. Wash and wipe slides.
6. Prepare Substrate Reagent in Mixing Vial (Vial 6)
In order add:
4 ml deionized water
2 drops Acetate Buffer (Vial 4A)
1 drop AEC Chromogen (Vial 4B)
1 drop 3% Hydrogen Peroxide (Vial 5)
7. Apply 2 drops Substrate Reagent incubate up to 10 minutes. Check slide microscopically for adequate chromogen development.
8. When sufficient staining has been achieved rinse slides in deionized water for 5 minutes. Wipe off excess.
9. Counterstain with Mayer's Hematoxylin for 2 minutes.
10. Rinse in gently running tap water to "blue" the hematoxylin.
11. Apply glycerol gelatin or other aqueous mounting media and carefully cover with coverslip.

Nuclei will be blue, while the cytoplasm of positive cells will be rose-red to brownish-red.

References

1. Taylor, C. R., Immunoperoxidase Techniques. Practical and Theoretical Aspects, Arch. Pathol. Lab. Med., **102**, 113 (1978).
2. Guesdon, J. L., et al., The use of avidin-biotin interaction in immunoenzymatic techniques, J. Histochem. Cytochem., **27**, 1131 (1979).
3. Naritoku, W. Y., and Taylor, C. R., A comparative study of the use of monoclonal antibodies using three different immunohistochemical methods: an evaluation of monoclonal and polyclonal antibodies against human prostatic acid phosphatase, J. Histochem. Cytochem., **30**, 253 (1982).
4. Hsu, S. M., et al., Use of avidin-biotin peroxidase complex (ABC) in immunoperoxidase techniques: A comparison between ABC and unlabeled antibody (PAP) procedures, J. Histochem. Cytochem., **29**, 577 (1981).
5. Lazarides, E., Intermediate Filaments: A chemically heterogenous developmentally regulated class of proteins., Annu. Rev. Biochem., **51**, 219 (1982).
6. Denk, H., et al., Proteins of intermediate filaments. An immunohistochemical and biochemical approach to the classification of soft tissue tumors. AM. J. Pathol., **110**, 193 (1983).
7. Culling, C. F., et al., The effect of various fixatives and trypsin digestion upon the staining of routine paraffin embedded sections by the peroxidase-anti-peroxidase and immunofluorescent technique, J. Histotechnology, **3**, 10 (1980).

JWM/VLE 10/02

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.