

ImmPRESS® Polymer Detection Systems

Vector® ImmPRESS Micropolymer Systems are carefully produced to ensure consistent, optimal staining, are rigorously tested using a variety of immunohistochemical applications, and are guaranteed to retain activity during prolonged storage. However, background staining may occur, or specific staining may be absent or diminished in test specimens due to factors intrinsic to the tissue or cell samples (i.e. fixation, antigen expression, endogenous tissue components, etc.).

Not all of the causes of off-target staining or weak staining may be obvious. Trying to solve the problem often becomes a time-consuming and frustrating task. We hope this troubleshooting guide helps to identify and correct the most common sources of problems encountered in IHC/ICC staining.

To evaluate background staining, run these deletion controls.

A Substrate Alone <ul style="list-style-type: none"> • Substrate 	B Detection System: Secondary Antibody Enzyme Conjugate <ul style="list-style-type: none"> • Blocking Serum • ImmPRESS Polymer Detection Reagent • Substrate 	C Complete System <ul style="list-style-type: none"> • Blocking Serum • Primary Antibody • ImmPRESS Polymer Detection Reagent • Substrate
<p>Staining</p> <p>No Staining SEE B</p>	<p>Staining</p> <p>No Staining SEE C</p>	<p>Inappropriate Staining</p>
<p>Endogenous enzyme may be developing the substrate.</p> <ul style="list-style-type: none"> • When running this control, use the incubation time and conditions recommended in the substrate instructions. • Note: Color that develops with long exposure to substrate may not be seen within the development time used for specific staining. <p>Block endogenous enzymes appropriately:</p> <p>HRP and/or AP – Incubate slides in BLOXALL® Endogenous Blocking Solution (SP-6000) for 10 minutes to quench both peroxidase (HRP) and/or alkaline phosphatase (AP) activity.</p> <p>Or use the following protocols:</p> <p>HRP – use 0.3% H₂O₂ in methanol for 30 minutes or 3% H₂O₂ in water for 5 minutes.</p> <p>AP – add Levamisole (SP-5000) to the working solution of the AP substrate. If the endogenous enzyme is the intestinal isoform, use BLOXALL Blocking Solution (SP-6000).</p>	<p>ImmPRESS Polymer Detection Reagent can bind to endogenous immunoglobulins if used to detect a primary antibody that is of the same species as the tissue (e.g. mouse primary on mouse tissue).</p> <p>For mouse primary on mouse tissue, use M.O.M.® ImmPRESS Kit (MP-2400).</p> <p>Cross-reactivity may occur between ImmPRESS Polymer Detection Reagent and tissues that are related in species to primary antibody or have tissue components with related amino acid sequences to the primary.</p> <ul style="list-style-type: none"> • Use reagents that have been adsorbed for the tissue species. • Add 2% or more normal serum from tissue species to the secondary antibody diluent, and/or reduce concentration of the secondary antibody. • Use additional blocking agents such as 2% immunohistochemical grade BSA (SP-5050), nonfat dry milk, gelatin, or 0.1% detergent. <p>Washing steps following the Polymer Detection Reagent may be inadequate.</p> <ul style="list-style-type: none"> • Lengthen the wash time and include several changes of wash buffer. <p>Blocking serum from the wrong species was used.</p> <p>Use serum from the same species in which the polymer detection reagent was produced.</p>	<p>Excess primary antibody has been used.</p> <p>Titer the primary antibody concentration. The optimal antibody concentration should produce clean specific staining with no background.</p> <p>The primary antibody may bind non-specifically or cross-react with other tissue epitopes.</p> <ul style="list-style-type: none"> • Add normal serum, BSA, nonfat dry milk, or detergent to buffer used as the primary antibody diluent. Be sure that the antibody diluent has sufficient salt to minimize nonspecific ionic interactions. Generally diluents should contain from 0.15M (0.9%) to 0.6M sodium chloride. • Change source or species of primary. <p>Some commercial diluents for the primary antibody can contribute to the background.</p> <p>Use diluent compatible with the detection system.</p> <p>If the section shows small, amorphous, punctuate staining, the primary antibody may have some denatured precipitated immunoglobulin.</p> <p>Centrifuge primary antibody; use supernatant.</p> <p>Tissue sections dried out during procedure.</p> <p>Be sure to keep tissue sections moist during all steps in the procedure.</p>
<p>Repeat Control A</p>	<p>Repeat Control B</p>	<p>Repeat Control C</p>

For technical assistance

(800) 227-6666 | technical@vectorlabs.com

US HEADQUARTERS – 6737 Mowry Avenue, Newark, CA. 94560, USA | (650) 697-3600
vector@vectorlabs.com | ORDERS: (800) 227-6666

BLOXALL, ImmPRESS, M.O.M., and Vector are trademarks of Vector Laboratories, Inc.

LIT3035.Rev00

ImmPRESS Polymer Detection Systems

If staining is weak or absent, use these tests.

A Enzyme/Substrate	B Primary Antibody	C Polymer Detection Reagent (Secondary Antibody Enzyme Conjugate)	Notes
<p>For Peroxidase Substrate:</p> <p>Add 1-2 drops of ImmPRESS Reagent to 1 ml peroxidase substrate working solution. Color of solution should change within about 5 seconds.</p> <p>For Any Substrate:</p> <p>Place 1 drop ImmPRESS Reagent on a small piece of nitrocellulose and then immediately dip the nitrocellulose into substrate. A colored spot will develop where the ImmPress Reagent was dotted.</p> <p>If color develops, SEE B. If no color develops, SEE BELOW.</p> <p>Deionized water can contain inhibitors of the peroxidase reaction. Even if the water has very low conductivity, the peroxidase reaction can be severely compromised.</p> <p>Use glass distilled water for the preparation of the substrate solution.</p> <p>Check the pH of the substrate buffer. Buffers of different pH values are recommended for different substrates. Use clean glassware to prepare substrate; traces of chlorine, cleaning solutions, etc. may inhibit the peroxidase reaction.</p> <p>The substrate should be made according to instructions.</p>	<p>Use the primary antibody at the optimal concentration. If activity of the primary is lost over time, a higher concentration may be required to achieve optimal staining. Treatments such as freeze/thawing, especially with monoclonal antibodies, may result in partial or complete inactivation of the antibody. High concentrations of antibodies may also reduce staining.</p> <p>Testing the antibody on a known positive sample may provide information on the activity of the antibody. If the known positive sample is positive, but the test section is negative, SEE NOTES.</p> <p>If the pH of the diluent for the primary antibodies is incorrect, the antibody may not bind well to the antigen.</p> <p>Check the pH of the diluent. Generally TBS or PBS, pH 7.0-8.2, is recommended.</p> <p>If the primary antibody recognizes an antigen in the diluent, it may bind to the antigen in solution rather than on the tissue section. Common diluent additives such as normal serum, fetal bovine serum, or nonfat dry milk may contain significant antigen concentrations that are recognized by the primary antibody.</p> <p>Take care that the diluent for the antibody does not contain the antigen.</p>	<p>ImmPRESS Polymer Reagents are provided ready-to-use (R.T.U.) at an optimal concentration for most applications.</p> <p>Inappropriately high dilutions of secondary antibody enzyme conjugate can result in diminished staining.</p> <p>Generally a 1:200 to 1:500 dilution of our secondary antibody conjugates should give optimal staining..</p> <p>ImmPRESS Polymer Reagents are provided ready-to-use in an optimal diluent for most applications.</p> <p>If the diluent contains any neutralizing antibodies, diminished staining could result. For example, Anti-Mouse IgG should not be diluted in mouse serum. The immunoglobulins in mouse serum will bind the Anti-Mouse IgG and prevent this secondary antibody from binding to the primary antibody.</p> <p>Remove source of neutralizing antibodies.</p> <p>If the secondary antibody is incorrect, no staining will occur. The secondary antibody should be specific for the species in which the primary antibody is made. For example, Anti-Rabbit IgG should be used with primary antibodies made in rabbit.</p> <p>Use correct Polymer Detection Reagent or Secondary Antibody Enzyme Conjugate.</p>	<p>Blocking</p> <p>Some animals from which blocking serum was obtained may have developed antibodies to the antigen in question. If present, the antibodies may bind to the antigen and prevent the primary antibody from binding. Try other blocking proteins such as an immunohistochemical grade of BSA, gelatin, fetal bovine serum, nonfat dry milk, etc., or 1% detergent.</p> <p>Fixation Check</p> <p>Be sure that the method employed for preparing the sample is appropriate to preserve the primary antibody target antigen. Use a high temperature antigen unmasking technique with an appropriate Antigen Unmasking Solution (Citrate-based, H-3300; or Tris-based, H-3301).</p> <p>Counterstain/Mounting</p> <p>Some enzyme reaction products are soluble in alcohol, xylenes or other solvents used for non-aqueous permanent mounting. Be certain that the enzyme reaction product is compatible with the counterstain and mounting medium. A substrate/counterstain compatibility chart is available on our website: vectorlabs.com</p>
	<p style="text-align: center;">Repeat Control C</p>	<p style="text-align: center;">If negative, see notes.</p>	