

Imm**PRESS**[™] REAGENT Anti-Goat IgG

PEROXIDASE Cat. No. MP-7405

INSTRUCTIONS FOR IMMUNOHISTOCHEMICAL STAINING USING GOAT PRIMARY ANTIBODIES

Introduction

The ImmPRESS[™] Polymerized Reporter Enzyme Staining System is an enzymatic, non-biotin, one-step detection kit that provides very high sensitivity with very low background staining in immunohistochemical and immunocytochemical applications.

The ImmPRESS[™] Reagent employs a novel, proprietary approach to conjugate horseradish peroxidase micropolymers to our affinitypurified, extensively cross-adsorbed secondary antibodies. This conjugation technology avoids the use of large dextrans or other macromolecules as a backbone and allows a higher density of enzymes per antibody to bind to the target with minimal steric interference. The overall result is an increase in target accessibility, binding specificity, and signal intensity with low background staining.

COMPONENTS

Reagents supplied:

- ImmPRESS™ (Peroxidase) Polymer Anti-Goat IgG Reagent (made in horse, ready-to-use)
- 2.5% Normal Horse Serum for blocking (readyto-use)

The **ImmPRESS**[™] (15 ml) Kit will stain approximately 75-150 sections and the ImmPRESS[™] (50 ml) Kit will stain approximately 250-500 sections.

Unless labeled otherwise, ImmPRESS[™] reagents are designed for laboratory use only.

Storage:

Store $ImmPRESS^{\rm TM}$ Kit Reagents at 2-8 °C (do not freeze).

Reagents not supplied:

- Primary Antibody
- Buffer
- Peroxidase Substrate

Imm**PRESS**[™] DETECTION SYSTEM

The **ImmPRESS**[™] Reagent is ready-to-use — No mixing or titering of the **ImmPRESS**[™] reagent is necessary to obtain optimal immunohistochemical staining. Dilution of this reagent or changes in suggested incubation time may affect performance. The reagent is supplied in a convenient dropper bottle. (To remove the drop dispenser tip, press laterally with thumb until tip snaps off).

The staining procedure should be performed at room temperature (20 - 25 °C). For optimal performance, the **ImmPRESS**[™] Reagent should be equilibrated to room temperature before use.

A number of different wash buffers can be used with the **ImmPRESS**[™] Reagents. One of the most common is 10 mM sodium phosphate, pH 7.5, 0.9% saline (PBS). 0.1% Tween 20 detergent may be added to the wash buffer. Adding 0.1% Tween 20 is especially recommended for use with automated stainers.

STAINING PROCEDURE

1. For paraffin sections, deparaffinize and hydrate through xylenes or other clearing agents and graded alcohol series.

For frozen sections or cell preparations, fix with acetone or an appropriate fixative for the antigen under study, if necessary.

Wash for 5 minutes in tap water.

- 2. If antigen unmasking is required, perform this procedure using a Vector[®] Antigen Unmasking Solution, Citrate-based (H-3300) or High pH-based (H-3301).
- 3. If quenching of endogenous peroxidase activity is required, incubate the slides in **BLOXALL™** Blocking Solution (SP-6000) for 10 minutes. If endogenous peroxidase activity does not present a problem, this step may be omitted. For alternative quenching procedures please see Note 3.
- 4. Wash in buffer for 5 minutes.
- 5. Incubate for 20 minutes with ready-to-use (2.5%) normal horse blocking serum or blocking solution of choice. Blot excess serum from sections.
- 6. Incubate with goat primary antibody diluted in appropriate antibody diluent solution. See Note 4.
- 7. Wash for 5 minutes in buffer.
- 8. Incubate for 30 minutes with ImmPRESS[™] reagent.
- 9. Wash for at least 2x 5 minutes in buffer.
- 10. Incubate in peroxidase substrate solution until desired stain intensity develops. For a list of peroxidase substrates see reverse.
- 11. Rinse in tap water.
- 12. Counterstain, clear and mount.

Vector Laboratories, Inc.

30 Ingold Road • Burlingame, CA 94010 Tel: (650) 697-3600 • Fax: (650) 697-0339 E-mail: vector@vectorlabs.com • www.vectorlabs.com

NOTES:

- 1. ImmPRESS[™] reagents are ideal for multiple antigen labeling applications. A brochure is available with protocols - "Discovery Through Color". Please request a free printed copy or download from our website: www.vectorlabs.com. Additional information on Enzyme Substrate Combinations, Counterstain/Substrate Compatibility, and Relative Substrate Sensitivity is also available on our website.
- 2. Use only freshly prepared buffers. Bacterial contamination which can occur in buffers stored at room temperature may affect the quality of the staining.
- 3. Alternative peroxidase guenching procedures:

For formalin fixed cells and tissues, incubate in 3% H_2O_2 in tap water for 5 minutes or 0.3% H_2O_2 in either methanol or water for 30 minutes.

For frozen tissue or cell preparations, use 0.3% H₂O₂ in 0.3% normal serum in PBS for 5 minutes, or 0.3% H₂O₂ in methanol for 30 minutes or use other published methods. (eq. Andrew, S.M., Jasani, B., Histochem J. 1987, 19, 426-30).

- 4. To avoid adsorption of the antibody to the plastic or glass container in which the final dilution is made, the primary antibody may be diluted in buffers containing diluted (2.5%) normal serum. Alternatively, 0.1% immunohistochemical grade Bovine Serum Albumin (SP-5050) can be used. Other grades of BSA can contain undesired impurities.
- 5. For thicker sections, longer incubation times may be required for optimal staining.
- 6. To prevent sections from detaching from the glass, slides can be treated with VECTABOND[™] Reagent (SP-1800), a non-protein tissue section adhesive.

Peroxidase Substrates:

A variety of chromogens can be used to localize peroxidase in tissue or cell preparations. All Vector Laboratories' substrates are supplied in convenient, easy to use dropper bottles. Vector Laboratories offers conventional as well as proprietary substrates producing the colors listed.

Note: A chart of the Relative Sensitivity of Substrates in Immunohistochemistry and further description of substrate properties is available on our website: www.vectorlabs.com

ImmPACT [™] DAB <i>EqV</i> (Brown)	400 ml	SK-4103
ImmPACT [™] DAB (Brown)	120 ml	SK-4105
ImmPACT [™] AEC (Red)*	120 ml	SK-4205
ImmPACT [™] AMEC Red (Red)*	120 ml	SK-4285
ImmPACT [™] VIP (Purple)	120 ml	SK-4605
ImmPACT [™] SG (Blue-Gray)	120 ml	SK-4705
ImmPACT [™] NovaRED [™] (Red)	120 ml	SK-4805
DAB/Ni Substrate (Brown or Gray-Blac	:k) 1 Kit	SK-4100
AEC Substrate (Red)*	1 Kit	SK-4200
TMB Substrate (Blue)	1 Kit	SK-4400
Vector [®] VIP Substrate (Purple)	1 Kit	SK-4600
Vector [®] SG Substrate (Blue-Gray)	1 Kit	SK-4700
Vector [®] NovaRED [™] Substrate (Red)	1 Kit	SK-4800

* AEC. ImmPACT[™] AEC and ImmPACT[™] AMEC Red must be mounted in aqueous mounting media. All other substrates may be dehydrated, cleared, and permanently mounted.

Kits contain sufficient reagent to prepare approximately 300 ml of working solution.

Imm**PRESS**[™] Reagents:

Amplified Peroxidase:

ImmPRESS[™] Excel Reagent Anti-Rabbit IgG 15 ml MP-7601 (made in horse)

ImmPRESS[™] Excel Reagent Anti-Mouse IgG 15 ml MP-7602 (made in horse)

Peroxidase:

ImmPRESS™ Reagent Anti-Rabbit IgG	15 ml MP-7401
(made in horse)	50 ml MP-7401
ImmPRESS™ Reagent Anti-Rabbit IgG	15 ml MP-7451
(made in goat)	50 ml MP-7451
ImmPRESS [™] Reagent Anti-Mouse IgG	15 ml MP-7402
(made in horse)	50 ml MP-7402
ImmPRESS [™] Reagent Anti-Mouse IgG	15 ml MP-7452
(made in goat)	50 ml MP-7452
ImmPRESS [™] Reagent Anti-Mouse IgG (rat adsorbed, made in horse)	15 ml MP-7422
ImmPRESS [™] Reagent Anti-Goat IgG	15 ml MP-7405
(made in horse)	50 ml MP-7405
ImmPRESS™ Reagent Anti-Rat IgG (made in goat)	50 ml MP-7404

ImmPRESS [™] Reagent Anti-Rat IgG (mouse adsorbed, made in goat)	15 ml MP-7444
ImmPRESS [™] Universal Reagent Anti-Mouse/Rabbit IgG (made in horse)	15 ml MP-7500 50 ml MP-7500
Vector [®] M.O.M. [™] ImmPRESS [™] Peroxidase Kit (made in horse)	1 Kit MP-2400
<u>Alkaline Phosphatase</u> : ImmPRESS [™] -AP Anti-Rabbit IgG (made in horse)	15 ml MP-5401 50 ml MP-5401
ImmPRESS [™] -AP Anti-Mouse IgG (made in horse)	15 ml MP-5402 50 ml MP-5402

Related Reagents:

Antigen Unmasking Solution (dilutes to 25 liters)			
Citrate-based	250 ml	H-3300	
High pH	250 ml	H-3301	
BLOXALL [™] Blocking Solution	100 ml	SP-6000	
ImmEdge™ Hydrophobic Barrier Pen	2-pen set	H-4000	
ImmPrint [™] Histology Pen	5-pen set	H-6100	
Vectabond [™] Reagent (dilutes to 350 ml)	7 ml	SP-1800	
VectaMount [™] Mounting Medium	60 ml	H-5000	
VectaMount [™] AQ Mounting Medium	60 ml	H-5501	
Vector [®] Hematoxylin	500 ml	H-3401	
Vector [®] Hematoxylin QS	100 ml	H-3404	
Vector [®] Methyl Green	500 ml	H-3402	
Vector [®] Nuclear Fast Red	500 ml	H-3403	

Heat-treated, ultrafiltered normal serum:

<u>near treated, attainterea normal seram</u> .		
Normal Goat Serum	20 ml	S-1000
2.5% Normal Goat Serum	50 ml	S-1012
Normal Horse Serum	20 ml	S-2000
2.5% Normal Horse Serum	50 ml	S-2012
Normal Chicken Serum	20 ml	S-3000
Normal Swine Serum	20 ml	S-4000
Normal Rabbit Serum	20 ml	S-5000
Control Antibodies:		
Rabbit IgG Control Antibody	5 mg	I-1000
Mouse IgG Control Antibody	1 mg	I-2000
Rat IgG Control Antibody	1 mg	I-4000
Goat IgG Control Antibody	5 mg	I-5000

Trademarks: Vector, VECTABOND, BLOXALL, ImmPRESS, ImmEdge, ImmPACT, ImmPrint, VectaMount, NovaRED, and M.O.M. are trademarks of Vector Laboratories.

Detailed product listings, specifications and protocols are available on our website. A complete catalog listing is also available upon request.

Visit our website: www.vectorlabs.com