

ImmPRESS® HRP Horse Anti-Rabbit IgG PLUS Polymer Kit



Peroxidase

Cat. No. MP-7801

Storage Store reagents in original bottles at 2-8 °C. Do not freeze.

Description Instructions for Immunohistochemical staining using mouse or primary antibodies.

The ImmPRESS HRP PLUS Polymer Kit is an enzymatic, non-biotin amplification system that produces crisp, highly sensitive, specific staining with low background.

Reagents in the ImmPRESS PLUS Polymer Kit are ready-to-use—no mixing or titering is necessary to obtain optimal staining. Dilution of the reagents or changes in suggested incubation time may affect performance. The reagents are supplied in convenient dropper bottles. (To remove the drop dispenser tip, press laterally with thumb until tip snaps off).

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 and is especially recommended for use with automated stainers.

Kit Components

| Product Name | Volume |
|--|--------|
| BLOXALL® Endogenous Blocking Solution | 15 ml |
| Normal Horse Serum, 2.5% | 15 ml |
| ImmPRESS HRP Horse Anti-Rabbit IgG Polymer Reagent | 15 ml |
| ImmPACT® DAB EqV Reagent 1 (Chromogen) | 7.5 ml |
| ImmPACT DAB EqV Reagent 2 (Diluent) | 7.5 ml |

The ImmPRESS PLUS Polymer Kit (15 ml) will stain approximately 75-150 sections.

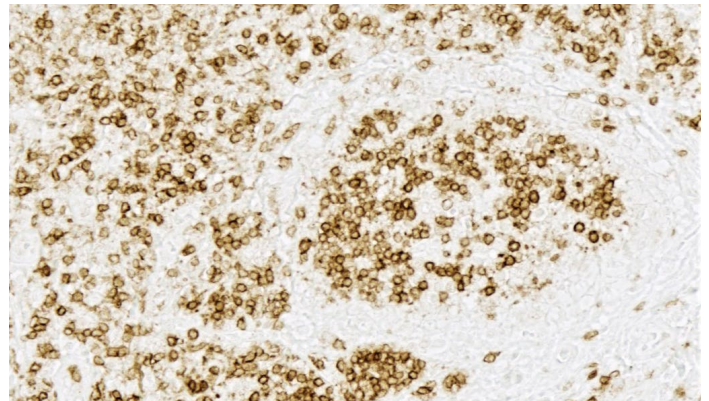
Staining Procedure

1. For paraffin sections, deparaffinize and hydrate tissue sections 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, pH 6.0 (H-3300) or Tris-based, pH 9.0 (H-3301).



Tonsil (FFPE): Stained for CD3 antigen using a rabbit monoclonal antibody and ImmPRESS® HRP Horse Anti-Rabbit IgG Plus Polymer Kit (MP-7801). No counterstain.

- If quenching of endogenous peroxidase activity is required, incubate the sections in BLOXALL Blocking Solution for 10 minutes.
- Wash in buffer for 5 minutes.
- Incubate sections for 20 minutes with Normal Horse Serum, 2.5%.
- Tip off excess serum from sections.
- Incubate with rabbit primary antibody diluted in appropriate antibody diluent solution, such as diluted normal horse serum or BSA.
- Wash in buffer for 5 minutes.
- Incubate for 30 minutes with ImmPRESS HRP Horse Anti-Rabbit IgG Polymer Reagent.
- Wash for 2 x 5 minutes in buffer.
- Combine equal volumes of ImmPACT DAB EqV Reagent 1 with Reagent 2. Mix well.
- Incubate in ImmPACT DAB EqV working solution until desired stain intensity develops, approximately 2-10 minutes.
- Wash for 2 x 5 minutes in buffer.
- Rinse sections in tap water.
- Counterstain (optional), clear and mount.

Detailed product listings, specifications, protocols and additional information is available on our website: vectorlabs.com

For research use only. Not intended for animal or human therapeutic or diagnostic use.

©2022 Vector Laboratories. All rights reserved.

BLOXALL, ImmPACT, ImmPRESS, and Vector are trademarks of Vector Laboratories, Inc.

Tween is a trademark of Atlas Chemical Industries.

MP-7801.LBL-00284.Rev01