

# ImmPRESS® Excel Amplified Polymer Kit

## Peroxidase (Anti-Rabbit IgG)

<b>Cat. No.</b>	<b>MP-7601</b>
<b>Storage</b>	Store reagents in original bottles at 2-8 °C. Do not freeze.
<b>Description</b>	<p>Instructions for Immunohistochemical staining using rabbit primary antibodies.</p> <p>The ImmPRESS Excel Amplified Polymer Kit is an enzymatic, non-biotin amplification system that produces crisp, highly sensitive, specific staining with low background.</p> <p>The reagents in the ImmPRESS Excel Amplified 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.</p> <p>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.</p>

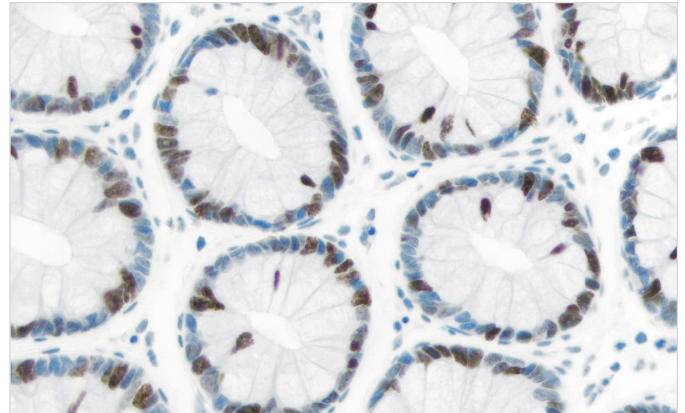
### Kit Components

Product Name	Volume
BLOXALL® Blocking Solution	15 ml or 50 ml
Normal Horse Serum, 2.5%	15 ml or 50 ml
Amplifier Antibody (Goat Anti-Rabbit IgG)	15 ml or 50 ml
ImmPRESS Horse Anti-Goat IgG Polymer Reagent	15 ml or 50 ml
ImmPACT® DAB EqV Reagent 1 (Chromogen)	10 ml or 30 ml
ImmPACT DAB EqV Reagent 2 (Diluent)	10 ml or 30 ml

The ImmPRESS Excel Amplified Polymer Kit (15 ml) will stain approximately 75-150 sections. The ImmPRESS Excel Amplified Polymer Kit (50 ml) will stain approximately 250-500 sections.

### Staining Procedure

- 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.



Colon (FFPE): Ki67 (rm) detected with ImmPRESS Excel Amplified Polymer Kit (Anti-Rabbit IgG) and ImmPACT DAB EqV (brown). Counterstained with Vector Hematoxylin QS (blue).

- 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).
- If quenching of endogenous peroxidase activity is required, incubate the sections in BLOXALL Blocking Solution (SP-6000) 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 15 minutes with Amplifier Antibody.
- Wash in buffer for 5 minutes.
- Incubate for 30 minutes with ImmPRESS 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](http://vectorlabs.com)